



# Konzentrations-Messsysteme MicroPolar (++) LB 566

Betriebsanleitung Geräteanleitung 41986BA1

Rev. Nr.: 08, 07/2023

Die gelieferten Geräte dürfen nur vom Service der Firma Berthold Technologies oder durch von der Firma Berthold Technologies autorisierte Techniker instand gehalten werden!

Im Störungsfall wenden Sie sich bitte an unseren zentralen Kundendienst (Kontaktadresse siehe unten). The units supplied should not be repaired by anyone other than Berthold Technologies Service engineers or technicians by Berthold Technologies.

In case of operation trouble, please address to our central service department (address see below).

Die komplette Betriebsanleitung besteht aus der Geräte- und Softwareanleitung.

#### Die Geräteanleitung beinhaltet:

- Mechanische Komponenten
- > Montage
- > Elektrische Installation
- > Technische Daten
- > Elektrische und mechanische Zeichnungen

#### Die Softwareanleitung beinhaltet:

- > Bedienung der Auswerteeinheit
- > Parameter-Beschreibung
- Grundeinstellung
- Kalibrierung
- > Fehlermeldungen

#### Der vorliegende Teil beinhaltet die Geräteanleitung.

Änderungen vorbehalten

BERTHOLD TECHNOLOGIES GmbH & Co. KG	
Calmbacher Str. 22 · D-75323 Bad Wildbad	
Zentrale:	Service:
Tel. +49 7081 177 0	Tel. +49 7081 177 111
Fax +49 7081 177 100	Fax +49 7081 177 339
industry@Berthold.com	Service@Berthold.com
www.Berthold.com	

# **Inhaltsverzeichnis**

		Seite
Kapitel 1. Sich	nerheitshinweise	7
1.1	Kennzeichnung und Warnhinweise	7
1.2	Generelle Hinweise	8
1.5	Algemente Sichemens- und Wahnninweise	5
Kapitel 2. Allg	jemeines	11
2.1	Bestimmungsgemäßer Gebrauch	11
2.2	Frequenzzulassung	12
2.3	Begriffserklärung	14
Kapitel 3. Sys	tembeschreibung	17
3.1	Messprinzip	17
3.2	Messwertberechnung	18
3.3	lemperaturkompensation	19
3.5	Mechanische Komponenten	20
	3.5.1 Die Auswerteeinheiten	23
	3.5.2 Die FlowCell	25
	3.5.3 Die Behältersonde	28
3.6	Messanordnung am Rohr	34
3.7	Messanordnung am Behälter	35
Kapitel 4. Inb	etriebnahme	36
4.1	Transport	36
4.2	Inbetriebnahme FlowCell	37
	4.2.1 Installation der FlowCell	37
	4.2.3 Anschluss der HF-Kabel	40 41
4.3	Inbetriebnahme Behältersonde	43
	4.3.1 Installation der Behältersonden	43
	4.3.2 Installation der AWE	45
4.4	Anschluss der AWE	43
	4.4.1 Anschlussbelegung der Steckerleiste	48
	4.4.2 Digitale Ausgänge, Relais	50
Kapitel 5. Wa	rtung	53
5.1	Allgemeines	53
5.2	verschleibtelle Gerätereinigung	54
5.4	Batterie	57
5.5	Sicherung austauschen	58

Kapitel 6. Tee	chnische Daten	60
6.	1 Technische Daten AWE	60
6.	2 Technische Daten Sensoren	63
6.	3 Technische Daten HF-Kabel	69
6.	4 Format der Datenausgabe RS232 und RS485	71
Kapitel 7. Zei	rtifikate	73
7.	1 EG-Konformitätserklärung	73
7.	2 Frequenzzulassung	76
Kapitel 8. Te	chnische Zeichnungen	85
8.	1 Maßbilder der AWE-Wandgehäuse	85
8.	2 Elektrischer Anschlussplan	87
8.	3 Maßbilder FlowCell	88
	8.3.1 Typ LB 5660-102-00x, FlowCell DN 50 Flansch, FDA	88
	8.3.2 Typ LB 5660-202-00x, FlowCell DN 65 Flansch, FDA	89
	8.3.3 Typ LB 5660-402-00x, FlowCell DN 80 Fiditsch, FDA	90
	8.3.4 Typ LB 5660-502-00x, FlowCell DN 100 Flansch, FDA	92
	8.3.6 Typ LB 5660-602-00x, FlowCell DN 150 Flansch, FDA	93
	8.3.7 Typ LB 5660-112-00x, FlowCell DN 50 G-BS/M	94
	8.3.8 Typ LB 5660-212-00x, FlowCell DN 65 G-BS/M	95
	8.3.9 Typ LB 5660-312-00x, FlowCell DN 80 G-BS/M	96
	8.3.10 Typ LB 5660-412-00x, FlowCell DN 100 G-BS/M	97
	8.3.11 Typ LB 5660-512-00x, FlowCell DN 125 G-BS/M	98
	8.3.12 Typ LB 5660-612-00x, FlowCell DN 150 G-BS/M	100
	8.3.13 Type LB 5660-132-00X DN 50	101
	8.3.14 Type LB 5660-332-00X DN 80	101
	8 3 16 Type LB 5660-432-00X DN 100	102
	8.3.17 Type LB 5660-532-00X DN 125	104
	8.3.18 Type LB 5660-632-00X DN 150	105
8.4	4 Maßbilder Behältersonden	106
	8.4.1 Typ LB 5650-01	106
	8.4.2 Typ LB 5650-02	107
	8.4.3 Typ LB 5650-03	108
	8.4.4 Typ LB 5650-04	109
	8.4.5 Typ LB 5650-05	110
	8.4.0 Typ LD 3030-09 8.4.7 Einbausituation in Pohrloitungon	117
8	5. Maßhilder Snülsonden	113
0.	8.5.1 Tvp I B 5651-01	113
	8.5.2 Tvp LB 5651-02	114
	8.5.3 Typ LB 5651-03	115
	8.5.4 Typ LB 5651-04	116
	8.5.5 Typ LB 5651-05	117
	8.5.6 Einbausituation in Rohrleitungen	118
8.	6 Montageblätter für LB 5650 (Behältersonde)	119
8.	/ Montageblatter für LB 5651 (Spülsonde)	121



# Kapitel 1. Sicherheitshinweise

### 1.1 Kennzeichnung und Warnhinweise

Der Begriff Berthold Technologies steht in dieser Betriebsanleitung stellvertretend für die Firma Berthold Technologies GmbH & Co.KG.

Zur Vermeidung von Personen- und Sachschäden beachten Sie unbedingt die in dieser Betriebsanleitung gegebenen Warn- und Sicherheitshinweise. Sie sind mit GEFAHR, WARNUNG, VORSICHT oder HINWEIS besonders gekennzeichnet.



# 

Enthält Anwendertipps und andere besonders nützliche Informationen.

Bedeutung weiterer, in dieser Dokumentation verwendeter Symbole

Warnhinweis: Kein Eingriff, nichts verändern

Gebot: Spannungsfrei schalten



Gebot: Sicherheitsschuhe tragen

## **1.2 Generelle Hinweise**

In dieser Betriebsanleitung sind die wichtigsten Sicherheitsmaßnahmen zusammengefasst. Sie ergänzt die entsprechenden Vorschriften, zu deren Studium das verantwortliche Personal *verpflichtet* ist.

Beachten Sie unbedingt:

- > die nationalen Sicherheits- und Unfallverhütungsvorschriften
- > die nationalen Montage- und Errichtungsvorschriften
- > die allgemein anerkannten Regeln der Technik
- die Angaben zu Transport, Montage, Betrieb, Wartung, Instandhaltung
- > die Sicherheitshinweise und Angaben in dieser Betriebsanleitung
- sowie die beiliegenden technischen Zeichnungen und Verdrahtungspläne
- die Kennwerte, Grenzwerte und die Angaben f
  ür die Betriebsund Umgebungsbedingungen auf den Typenschildern und in den Datenblättern
- > die Hinweisschilder auf den Geräten
- > die landesspezifischen Zulassungsbestimmungen





#### 

Ist Flüssigkeit in das Innere des Geräts gelangt, so ist die Stromzufuhr zu unterbrechen. Das Gerät muss durch eine autorisierte Servicestelle kontrolliert und gereinigt werden.

	Elektrische Gefahren:
	Während der Installation und für Servicearbeiten ist die Stromver- sorgung abzuschalten, damit Berührungen mit spannungsführen- den Teilen verhindert werden.
	Vor jedem Öffnen des Gerätes ist die Stromzufuhr zu unterbrechen. Arbeiten am geöffneten und unter Spannung stehenden Gerät sind verboten.
HINWEIS	Achtung! Mögliche Gefährdung, Sachschäden! Betrifft den Gerätetyp:
	Bei Anschluss der 24 V DC Hilfsenergie, müssen die + und – Pole richtig angeschlossen werden. Es besteht kein Verpolungsschutz!
HINWEIS	Ersatzsicherungen müssen die vom Gerätehersteller vorgegebenen Werte aufweisen. Ein Kurzschließen oder Manipulieren ist nicht gestattet.
	<b>WICHTIG</b> Das LB 566 und alle Zusatzgeräte müssen geerdet am Netz ange- schlossen sein.
	<b>i WICHTIG</b> Das Konzentrations-Messgerät I B 566 darf ausschließlich von sach-
	kundigen Personen installiert, gewartet und repariert werden.
Sachkundige Personen	Sachkundig sind Personen dann, wenn sie durch ihre fachliche Aus- bildung ausreichende Kenntnisse auf dem geforderten Gebiet besit- zen und mit den einschlägigen nationalen Arbeitsschutzvorschriften, Unfallverhütungsvorschriften, Richtlinien und anerkannten Regeln der Technik vertraut sind. Sie müssen in der Lage sein, die Ergeb- nisse ihrer Arbeit sicher beurteilen zu können und sie müssen mit dem Inhalt dieser Betriebsanleitung vertraut sein.

# Kapitel 2. Allgemeines

# 2.1 Bestimmungsgemäßer Gebrauch

	Das MicroPolar LB 566 wurde als Konzentrations-Messgerät entwi- ckelt und darf nur für diesen Zweck eingesetzt werden. Werden die Geräte auf eine Weise verwendet, die nicht in der vorliegenden Be- triebsanleitung beschrieben wird, so ist der Schutz der Geräte be- einträchtigt und der Garantieanspruch geht verloren.
	Berthold Technologies haftet bzw. garantiert lediglich, dass die Ge- räte seinen veröffentlichten Spezifikationen entsprechen. Das LB 566 darf nur in unbeschädigtem, trockenem und sauberem Zu- stand eingebaut werden. Umbauten und Veränderungen an den Systemkomponenten sind nicht gestattet.
	Das LB 566 ist nicht als "sicherheitsgerichtete Messung" qualifiziert.
Normenkonformität	Die Normen und Richtlinien, denen das LB 566 entspricht, sind in dieser Geräteanleitung in <i>Kapitel 2.2 Frequenzzulassung</i> und <i>Kapitel 7.1 EG-Konformitätserklärung</i> aufgeführt.
Druckgeräterichtlinie	Die FlowCell und die Behältersonde wurden als Druckgeräte nach Artikel 4 Absatz 3 der Richtlinie 2014/68/EU eingestuft. Vor der Ver- wendung ist zu prüfen, ob der Einsatzfall dieser Einstufung ent- spricht. Insbesondere ist die Medienverträglichkeit der Fluid berüh- renden Teile zu prüfen.
Schutzart	Der Schutzgrad des LB 566 nach IEC 60529 beträgt maximal IP 65.
Warnung	Bestimmungswidrig und zu verhindern ist:
vor Fehlgebrauch	Die Verwendung unter anderen als den durch den Hersteller in seinen technischen Unterlagen, Datenblättern, Betriebs- und Montageanleitungen und in anderen spezifischen Vorgaben ge- nannten Bedingungen und Voraussetzungen.
	Die Verwendung nach Instandsetzung durch Personen, die nicht von Berthold Technologies autorisiert wurden.
	Die Verwendung in beschädigtem oder korrodiertem Zustand.
	Der Betrieb mit geöffnetem oder mit unzureichend verschlosse- nem Deckel.
	Der Betrieb mit unzureichend festgezogenen Adaptern und Ka- belverschraubungen.
	Der Betrieb ohne die vom Hersteller vorgesehenen Sicherheits- vorkehrungen.
	<ul> <li>Bestehende Sicherheitseinrichtungen zu manipulieren oder zu umgehen.</li> </ul>
Autorisierte Personen	Autorisierte Personen sind Personen, die entweder aufgrund gesetz- licher Vorschriften für die entsprechende Tätigkeit vorgesehen sind oder durch Berthold Technologies für bestimmte Tätigkeiten zuge- lassen wurden.

## 2.2 Frequenzzulassung

Das Gerät erfüllt die Vorschriften der FCC<sup>1</sup>, Part 15 und die der kanadischen Industrie nach lizenzfreien RSS-Standard(s). Der Betrieb unterliegt folgenden zwei Bedingungen:

- (1) Dieses Gerät darf keine schädlichen Störungen verursachen und
- (2) Dieses Gerät muss alle empfangenen Störungen akzeptieren, einschließlich solcher, die einen unerwünschten Betrieb verursachen können.

BERTHOLD TECHNOLOGIES MicroPolar (Brix) FCC ID: R9ZFCC01X01 IC: 4777A-IC01X01

BERTHOLD TECHNOLOGIES MicroPolar (Brix) ++ FCC ID: R9ZFCC01X12 IC: 4777A-IC01X12

Dieses Gerät wurde getestet und entspricht den Grenzwerten für ein digitales Gerät der Klasse B gemäß Teil 15 der FCC-Bestimmungen. Diese Grenzwerte sind so ausgelegt, dass sie einen angemessenen Schutz gegen schädliche Störungen in einer Wohnanlage bieten. Dieses Gerät erzeugt und verwendet Hochfrequenzenergie und kann diese ausstrahlen, und wenn es nicht entsprechend den Anweisungen installiert und verwendet wird, kann es zu schädlichen Störungen der Funkkommunikation kommen. Es besteht jedoch keine Garantie dafür, dass bei bestimmte Installationen keine Störungen auftreten. Wenn dieses Gerät schädliche Störungen des Radio- oder Fernsehempfangs verursacht, die durch Ein- und Ausschalten des Gerätes festgestellt werden können, wird der Benutzer aufgefordert, die Störung durch eine oder mehrere der folgenden Maßnahmen zu beheben:

- Empfangsantenne neu ausrichten oder verlegen.
- Vergrößerung des Abstands zwischen Gerät und Empfänger.
- Schließen Sie das Gerät an eine Steckdose an, die nicht mit dem Stromkreis des Empfängers verbunden ist.
- Wenden Sie sich an den Händler oder einen erfahrenen Radio-/Fernsehtechniker.

Änderungen oder Modifikationen an diesem Gerät, die nicht ausdrücklich von Berthold Technologies genehmigt wurden, können

## Commission

FCC<sup>1</sup> und IC<sup>2</sup> Zulassungsschilder

 $<sup>^{1}\ \</sup>mathrm{FCC}\ \ldots$  Federal Communications Commission

<sup>&</sup>lt;sup>2</sup> IC Industry Canada

zum Erlöschen der FCC-Zulassung zum Betrieb dieses Gerätes führen.

Das MicroPolar und MicroPolar ++ sind konform nach der R&TTE-Richtline 1999/5/EG und erfüllt somit alle Anforderungen für derartige Hochfrequenzgeräte. Die Geräte tragen zur Kennzeichnung der Konformität nach dem CE Zeichen die Nr. 0682 der Zertifizierungsstelle. Das Zertifikat befindet sich unter *Kapitel 7.3 Frequenzzulassung*.

# (€ 0682

## WICHTIG

Die Frequenzzulassungen und Konformitäten sind nur gültig in Kombination mit den Behältersonden und der FlowCell.

Zur Einhaltung der Konformität nach ETSI (European Telecommunications Standards Institute) darf das LB 566 ++ in Kombination mit Horn- und Spiralantennen nur mit einem 10 dB Dämpfungsglied im Sendepfad (MTX) betrieben werden.

## 🚺 WICHTIG

Das LB 566 wurde in Übereinstimmung mit den Sicherheitsanforderungen für Mikrowellengeräte hergestellt. Bestehen besondere gesetzliche Bestimmungen zur Verwendung von Mikrowellen, so ist es Aufgabe des Benutzers, diese einzuhalten.



## 

Eine Veränderung der Frequenz oder anderweitige Manipulation an dem Mikrowellengerät führt zum Verlust der Frequenzzulassung und kann zu strafrechtlichen Folgen führen.

Die Mikrowellenmodule besitzen keine austauschfähigen Komponenten und dürfen nicht geöffnet werden.

## 2.3 Verwendungszweck

Mit dem LB 566 können der Wasser- bzw. Feuchtegehalt und die Konzentration von nahezu beliebigen Materialien bestimmt werden. Dafür stehen folgende Sensor und AWE Varianten zur Verfügung:

- Die Behältersonden sind für den Einbau in Behältern und Rohrleitungen mit Nennweite ≥ 200 mm konstruiert. Die Sonde wird so angebracht, dass die beiden Messstäbe (Sender und Empfänger) in das Messgut eintauchen.
- 2. Bei der Messzelle handelt es sich um eine Rohrsonde, die direkt in das bestehende Rohrleitungssystem oder in einem Bypass eingebracht wird.

Die AWE steht in zwei Varianten zur Verfügung. Das Standardmodel MicroPolar und die High Dynamik Version MicroPolar ++. Das Micro-Polar ++ erfordert eine Mikrowellensignaldämpfung von mindestens 40 dB. Bei geringerer Dämpfung muss das Standardmodel MicroPolar angewandt werden.

Die Konzentrationsmessgeräte MicroPolar und MicroPolar ++ sendet beim Betrieb elektromagnetische Strahlung im Frequenzbereich von 2.4-2.5 GHz aus (Bereichseinschränkungen, je nach Länderbestimmung). Die Mikrowellen sind für Mensch und Umwelt ungefährlich (Leistungsabstrahlung < 10 mW). Die Mikrowellen treten dabei gerichtet aus dem Mikrowellenfenster heraus, das Produkt wird dabei nicht durch die Mikrowellen verändert.

Um eine einwandfreie Funktion des Messgerätes zu gewährleisten ist unbedingt auf folgendes zu achten:



- Das zu messende Material darf nur begrenzt elektrisch leitfähig sein.
- In dem Produkt dürfen sich keine Gaseinschlüsse befinden, bzw. müssen bei Rohrleitungsmessung mit genügend großem Druck komprimiert werden.
- Die Ionen-Konzentration, z.B. Salzgehalt muss annähernd konstant sein.

# 2.4 Begriffserklärung

AWE	Auswerteeinheit
Dämpfung	Schwächung der Mikrowellensignale, Mikrowellen-Messeffekt
Fabrikeinstellung	Siehe Werkseinstellung
HF-Kabel	Hochfrequenz-Kabel
FlowCell	Rohrsonde zur einfachen Integration in das bestehende Rohrlei- tungssystem
Mikrowellen	Bezeichnung für elektromagnetische Wellen in einem bestimm- ten Frequenzbereich
Phase	Phase oder auch Phasenverschiebung, Mikrowellen-Messeffekt
Quad Kabel	Zusammenfassung von vier gleichlangen HF-Kabeln in einem Wellrohr.
Spülsonde	Behältersonde mit Spülvorrichtung
Softkeys	Software zugeordneten Tasten
ТК	Temperaturkompensation
Werkseinstellung	In der Werkseinstellung sind alle Parameter mit Standard-Wer- ten voreingestellt. In den meisten Fällen ist damit die Kalibrie- rung des Gerätes wesentlich erleichtert. Trotz Werkseinstellung muss <b>immer</b> eine Anpassung durchgeführt werden.

Kapitel 3. Systembeschreibung

## 3.1 Messprinzip

Die Mikrowellen durchstrahlen das Messgut und werden in ihrer Ausbreitungsgeschwindigkeit verlangsamt (= Phasenverschiebung) und in ihrer Intensität gedämpft (= Dämpfung). Die Abb. 3-1 veranschaulicht das Messprinzip: Im Vergleich zu einem Referenzsignal erfahren die Mikrowellen beim Durchgang durch das Messprodukt eine Verringerung ihrer Ausbreitungsgeschwindigkeit (Phasenverschiebung) und ihrer Intensität (Dämpfung).



*Abb. 3-1: Prinzipdarstellung: Veränderung der Mikrowelle durch das Messgut* 

Voraussetzung dafür ist, dass das Messgut dielektrische Eigenschaften aufweist. Im Allgemeinen ist dies das Wasser, welches ein stark ausgeprägtes Dielektrikum darstellt. Die Wasser- bzw. Trockenmassekonzentration kann deshalb durch die Messung der Phasenverschiebung und/oder Dämpfung ermittelt werden.

Die im Messgut nachzuweisende Konzentration hängt in guter Näherung linear von der auftretenden Phasenverschiebung und der Dämpfung ab. Aus diesem Grund kann mit einer linearen Kalibrierung die Konzentration bzw. der Trockensubstanzgehalt des Messguts gemessen werden (siehe *Kapitel 3.2 Messwertberechnung*).

## 3.2 Messwertberechnung

Die Mikrowellenmesseffekte Phase und Dämpfung werden nach einer automatischen Plausibilitätsanalyse kalibriert.

Bei der Kalibrierung werden der Phase und oder der Dämpfung ein Konzentrationswert (oder Dichtewert) durch Probenahme zugewiesen. Die Kalibrierung verläuft automatisch und die Probeaufnahme wird durch die AWE unterstützt.

Welcher der Parameter (Phase, Dämpfung oder beide) zur Kalibrierung benutzt werden, hängt von der Größe und Störbeeinflussung des Messeffekts ab. So z.B. ist die Dämpfung deutlich empfindlicher auf elektrolytischer Leitfähigkeit (Salzgehalt).

In vielen Fällen wird die reine Phasenmessung empfohlen und der Messwert wie folgt berechnet:

#### Messwert = $A \cdot Phase + C$ Gl. 3-1

wobei gilt:

MesswertKonzentration / Feuchte / TrockenmasseA, CKoeffizienten der jeweiligen Kalibrierfunktion

Das LB 566 erlaubt die Kalibrierung, Anzeige und Ausgabe von zwei Konzentrationen Kon1 und Kon2. Die Kalibrierkoeffizienten müssen für die Konzentrationen 1 und 2 getrennt eingegeben werden. Näheres zu diesem Thema finden Sie in der Software Beschreibung.

Einschränkungen.

- Schwach gebundenes Wasser kann in Abhängigkeit der Stärke der Bindung nachgewiesen werden. Somit kann der Messeffekt von der auftretenden Korngrößenverteilung und der chemischen Zusammensetzung des Messgutes abhängen, wenn hierdurch die Bindung des Wassers an den Feststoff verändert wird.
- Zu durchstrahlende Wandungen aus Kunststoff, Gummi oder Isolierstoffen mit möglichst niedriger Dielektrizitätszahl haben kaum Einfluss auf den Messeffekt, und werden konstant einkalibriert.
- Die Messung von Eis und Kristallwasser ist nicht möglich, da hier die Wassermoleküle nicht frei rotieren können (Eis und Kristallwasser sind trocken).

Leitfähige Materialien wie z. B. Graphit oder Koks können nicht mit Mikrowellen durchstrahlt werden. Ebenso ist eine Durchstrahlung von Metallwandungen nicht möglich.

Kompensation

Neben dem Wassergehalt haben auch u.a. die Produkttemperatur, Produktdichte und eine schwankende Materialbelegung (schwankender Mikrowellendurchstrahlungsweg) einen Einfluss auf die Phasen und Dämpfung. Dieser Einfluss muss bei der Kalibrierung kompensiert werden.

## 3.3 Temperaturkompensation

Eine Temperaturkompensation (TK) ist bei schwankender Produkttemperatur erforderlich. Generell empfiehlt es sich eine Temperaturkompensation vorzusehen, d.h. ein Temperatursignal (0/4...20 mA oder PT100) an die AWE anzuschließen und gegebenenfalls die Kompensation in der AWE zu aktivieren. Die AWE ist dafür so ausgelegt, dass die erforderlichen TK's selbstständig berechnet werden können. Ab welcher Temperaturschwankung eine TK zwingend erforderlich ist, hängt von dem Produkt und von dem Wassergehalt ab. In erster Näherung sollte  $\pm$  2 °C als Schwankungsgrenze angesetzt werden.

Wird beispielsweise die Produkttemperatur über den PT100-Eingang erfasst, so wird die Gl. 3-1 wie folgt erweitert:

#### Messwert = $A \cdot Phase + D \cdot T_{mess} + C$ Gl. 3-2

wobei gilt:

Messwert	Konzentration / Feuchte / Trockenmasse
A, D, C	Koeffizienten der Kalibrierfunktion
T <sub>mess</sub>	Produkttemperatur

Der Umgang mit der Temperaturkompensation wird ausführlich in dem Softwareteil beschrieben.

## 3.4 Durchsatz-Berechnung und Ausgabe

Das LB 566 bietet bei Rohrleitungsapplikationen die Möglichkeit den Durchsatz (Massenstrom) zu berechnen und u. a. über Stromausgang auszugeben.

Die Berechnung erfolgt aus dem Mikrowellenmesswert, korreliert dieser mit der Produktdichte, so kann mit Zusatzinformationen der Durchsatz berechnet werden. Die benötigten Zusatzinformationen sind der Rohrleitungsinnendurchmesser/Querschnitt und die Produktgeschwindigkeit. Die Produktgeschwindigkeit muss dabei über Stromeingang eingespeist werden.

Näheres hierzu finden Sie im *Softwareteil unter Kapitel 4.2.12 Massenstrom*.



## 3.5 Mechanische Komponenten

Das Messsystem besteht aus einer AWE, einer Sonde/Antennenpaar und einem Satz spezieller Hochfrequenzkabel (kurz HF-Kabel). Die AWE steht in zwei Varianten zur Verfügung: Das Standardmodel MicroPolar LB 566 und die High Dynamik Version MicroPolar ++ LB 566, siehe Abb. 3-4 und Abb. 3-5.



Abb. 3-4: AWE MicroPolar LB 566

Abb. 3-5: AWE MicroPolar ++ LB 566



Die Sonden/Antennenpaare gibt es in verschiedenen Varianten, als Rohrleitungs- und Behältersonde mit oder ohne Spülvorrichtung (siehe Abb. 3-6 bis 3-7).



*Abb. 3-6: Von links: Behältersonde LB 5650 und LB 5651 mit Spülvorrichtung* 



Abb. 3-7: FlowCell LB 5660-102-00x Nennweite 50mm mit V-Flansch

Die FlowCell steht in den Rohrnennweiten 50 -150 mm (50, 65, 80, 100, 125 und 150) zur Verfügung. Die Anschlüsse sind in folgenden Varianten ausgeführt:

- Hygiene Milchrohrverschraubung DIN 11853-1
- V-Flansch EN 1092-1/11

#### 3.5.1 Die Auswerteeinheiten

Die Auswerteeinheiten bestehen aus dem Auswerterechner mit Mikrowelleneinheit. In der Mikrowelleneinheit werden die Mikrowellen erzeugt, empfangen und ausgewertet. Im Auswerterechner erfolgen die Signalverarbeitung und die Kommunikation. Für die einfache Bedienung des Messgerätes sind ein Display, 4 Softkeys und eine alphanumerische Tastatur vorgesehen. Den Softkeys werden über das Display verschiedene Funktionen zugeordnet.

Unterschiede zwischen MicroPolar ++ und MicroPolar

Die AWE MicroPolar ++ hat ergänzend zum Standardmodel ein zusätzliches HF-Verstärker-Modul, wodurch das Wandgehäuse größer ist (Außenmaße siehe *Kapitel 6.1 Technische Daten AWE*). Ansonsten unterscheiden sich die AWE's nur noch in Ihrer Anwendung.

MicroPolar ++ Mit der High Dynamik Version MicroPolar ++ sind höhere Produktdämpfungen erlaubt. Damit können größere Messwege durchstrahlt werden, z.B. eine FlowCell größerer Nennweite eingesetzt werden. Die Anwendungen der beiden AWE's werden durch die Produktdämpfung vorgegeben. Bis zu einer Dämpfung von 50 dB wird MicroPolar und darüber hinaus MicroPolar ++ eingesetzt. Das Micro-Polar ++ benötigt generell eine Dämpfung von 40 dB. Wird diese unterschritten, so weist die Software durch eine Fehlermeldung darauf hin.

Auf der Geräteunterseite steht eine RS232-Schnittstelle zur Verfügung.



*Abb. 3-8: Frontansicht der AWE* 



#### Interpretation der LED's

Auf der Gerätefront sind fünf LED's angebracht, die den jeweiligen Gerätezustand übersichtlich anzeigen.

- Run
- Error
- Signal 1

*Abb. 3-10: LED's auf der Frontseite der AWE* 

- Signal 2
- Comm

LED	Funktion
Run	<u>Leuchtet</u> : Gerät im Messmodus <u>Blinkt + ERROR LED aus</u> : Gerät im Warn-, Halte-, Pausen- oder Beladungsunterschreitungs-Zu- stand. Eine Displaymeldung mit Fehlercode gibt die Ursache an ( <i>siehe Softwareteil Kapitel 11.</i> <i>Fehlerlisten und Gerätezustände</i> ).
Error	<u>Leuchtet</u> : Gerät im Fehlerzustand. Eine Display- meldung mit Fehlercode gibt die Ursache an ( <i>siehe</i> <i>Softwareteil Kapitel 11. Fehlerlisten und Gerätezu-</i> <i>stände</i> ). Erlischt nach Quittierung oder Fehlerbehebung
Signal 1	Anzeige entsprechend der ausgewählten Funktion von Relais 1, mögliche Funktionen: Fehler, kein Produkt, Alarm min., Alarm max., Messung angehalten, Beladungsunterschreitung
Signal 2	Anzeige entsprechend der ausgewählten Funktion von Relais 2, mögliche Funktionen: Fehler, kein Produkt, Alarm min., Alarm max., Messung angehalten, Beladungsunterschreitung
Comm	Kommunikation aktiv. z.B. über RS232 und RS485

#### Anschlussklemmleiste

Die elektrischen Anschlüsse des LB 566 befinden sich auf einer Steckerleiste im Wandgehäuse. Die Anschlussleiste ist von vorne zugänglich, indem man den Deckel des Gehäuses öffnet. Dort befinden sich außerdem der Testschalter zur Netzunterbrechung und die Sicherungen. Die Hochfrequenzanschlüsse sind außen am Gehäuse. Alle übrigen Elemente, insbesondere der spannungsführenden Elemente (auf dem Motherboard) sind mit einer Schutzabdeckung versehen.

#### 3.5.2 FlowCell

Die FlowCell gibt es in den Nennweiten von 50 bis 150 mm (siehe Abb. 3-11). Für die Anschlüsse stehen die Varianten V-Flansch EN 1092-1/11 oder Hygiene-Milchrohrverschraubung DIN 11853-1 zur Verfügung. Die technischen Daten sind im *Kapitel 6.2 Technische Daten Sensoren* aufgeführt.



*Abb. 3-11: Varianten FlowCell* 

- A: mit V-Flansch EN 1092-1/11
- B: mit Hygiene-Milchrohrverschraubung DIN 11853-1
- C: mit Schweissflansch

Die FlowCell besteht aus einem stabilen Edelstahlkörper. Der Mikrowellen-Sender und -Empfänger sind fest an der Rohraußenseite verschweißt. Das gesamte Produktrohr erfüllt die speziellen Anforderungen für den Einsatz in Lebensmittel.

Es ragen keine Gegenstände in das Rohr hinein (wie z.B. Messfühler). Die FlowCell kann, je nach Variante, mittels V-Flansch, Milchrohrverschraubung oder Schweissflansch in die Rohrleitung eingebaut werden. Für die Variante mit V-Flansch sind ASA-Adapterflansche als Zubehör erhältlich.

Die FlowCell verfügt über zwei HF-Anschlüsse, über denen die Mikrowellensignale eingespeist und ausgegeben werden. Dabei kann der Ein- und Ausgang beliebig zugeordnet werden (M-Tx, M-Rx). Die Mikrowellensignale durchstrahlen das Produkt über den gesamten Rohrleitungsquerschnitt.

Für sämtliche Varianten gibt es folgendes Zubehör:

- 1. Rohranlege-PT100 oder Inline PT100
- 2. Leitfähigkeitsmessgerät
- 3. Probenahmeventil (Kombination mit 1. und 2. möglich)

#### Übersicht Zubehör (siehe auch Kap. 6.2):

Leitfähigkeitsmessgerät





Inline-PT100







Rohranlege-PT100

Probenahmeventil

Kombination Leitfähigkeitsmessgerät mit Probenahmeventil

Kombination Inline-PT100 mit Probenahmeventil

#### 3.5.3 Behältersonde

Die Behältersonde gibt es in den Varianten mit und ohne Spülvorrichtung (siehe Abb. 3-12). Die technischen Daten sind im *Kapitel* 6.2 Technische Daten Sensoren aufgeführt.



*Abb. 3-12: Varianten Behältersonde* 

A: Hochfrequenzanschlüsse

B: Prozessanschluss, Flansche unterschiedlicher Größen

#### Behältersonde vom Typ LB 5650 und Typ LB 5651

Die Behältersonde ist speziell für die Konzentrationsmessung in Behältern entwickelt worden. Die beiden Messstäbe tauchen dabei in das Messgut hinein. Die Mikrowellen werden aus einem Stabende abgestrahlt und von dem zweiten Stabende empfangen, dabei treten die Mikrowellen nur in Richtung des gegenüberliegenden Stabendes aus. Diese Richtcharakteristik der Sonde minimiert den störenden Einfluss von Metallteilen im Umfeld der Sonde und ermöglicht den Einbau auch bei wenig Platzreserve. So kann z.B. die Konzentration eines Zuckersudes kontinuierlich gemessen werden, um den geeigneten Impfzeitpunkt zu finden.

Die Kunststoffkappen der Messstäbe werden den speziellen Anforderungen für den Einsatz in Lebensmitteln gerecht.

Es stehen zwei unterschiedliche Sondentypen zur Verfügung:

- > Der Standardtyp ist die Behältersonde ohne Spülvorrichtung
- für Prozesse in denen vermehrt Ablagerungen z.B. durch Inkrustationen zu erwarten sind, wird die Spülsonde eingesetzt. Die Spülvorrichtung verhindert eine Ablagerung an den Mikrowellenaustrittsfenstern.



Die Strömungsrichtung des Messguts muss für den in Abb. 3-13 dargestellten Fall senkrecht zur Zeichnungsebene verlaufen. Damit ist gewährleistet, dass sich bei guter Durchmischung repräsentatives Messgut zwischen den Messstäben befindet.



*Abb. 3-13: Behältersonde LB 5650* 

HINWEIS

PT100

Nur die Behältersonde LB 5650 ist mit einem PT100-Anschluss ausgestattet und kann mit der AWE über ein 4-adriges Kabel verbunden werden. Der Anschlussplan für den PT100 wird in *Kapitel 4.6.1 Anschlussbelegung der Steckerleiste* beschrieben. Um die Inkrustationsgefahr in unmittelbarer Nähe der Messstäbe zu minimieren, ist die Spülsonde nicht mit einem PT100 ausgestattet.

Achtung, mögliche Sachschäden! Die Deckelschrauben auf der Frontseite der Behältersonden dürfen nicht gelöst werden, siehe Abb. 3-14.



HINWEIS

#### Behältersonde vom Typ LB 5651 mit Spülvorrichtung

Die Spülsonde LB 5661 ist für Prozesse entwickelt worden, bei denen Ablagerungen z.B. durch Inkrustationen an der Sonde zu erwarten sind.

Die Spülsonde weist zwei Spülkanäle auf, welche die Messstäbe frei von Inkrustationen halten, damit die Mikrowellen direkt mit dem Messgut in Berührung kommen. Alle produktberührenden Teile der Sonde werden den speziellen Anforderungen für den Einsatz in Lebensmittel gerecht. Abb. 3-15 zeigt den Sondenaufbau.

Achtung, Kommunikationssignal gestört! mögliche Sachschäden! Achten Sie beim Einbau der Sonde mit Spülung darauf, dass der HF-Stecker immer trocken ist.



Abb. 3-15: Spülsonde LB 5651

2 x 3/8" Spülanschlüsse, Innengewinde



Die Spülschlitzbreite ist für beide Sondenstäbe gleich ausgelegt und in Abb. 3-16 angegeben.



*Abb. 3-16: Stabspitze mit Spülrohr* 

#### 3.5.4 Das Hochfrequenzkabel

Die Hochfrequenzkabel (kurz HF-Kabel) dienen der Übertragung der Mikrowellensignale.

Die HF-Kabel ändern ihre Leitfähigkeit (für Mikrowellen) mit der Temperatur und würden deshalb bei Schwankungen der Umgebungstemperatur einen Messfehler erzeugen. Dieser Fehler wird durch die automatische Kabelkompensation ausgeglichen. Dabei werden die Einflüsse der Umgebungstemperatur auf die Messkabel mittels des Referenzkabels kompensiert. Zu diesem Zweck wird die Summe der Referenzkabel genauso lang gewählt, wie die Summe der Messkabel.

Es stehen zwei verschiedene HF-Kabeltypen unterschiedlicher Längen zur Verfügung.

**Variante 1:** Das sogenannte HF-Kabel quad: Es steht aus vier einzelnen HF-Kabeln gleicher Länge, deren Enden mit je einem HF-Stecker (N-Typ) abgeschlossen sind. Verfügbare Kabellängen: 2, 4, 6, 8 und 10 m (siehe Abb. 3-17).

Die Referenzleitung des HF-Kabels quad wird auf der Sondenseite mittels N-Verbinder kurzgeschlossen (siehe Abb. 3-18).



Abb. 3-17: HF-Kabel quad



*Abb. 3-18: HF-Kabel quad sondenseitig: Die Enden der Referenzleitung R-Rx und R-Tx sind mit einem N-Verbinder kurzgeschlossen.* 

**Variante 2:** Es besteht aus einem einzelnen HF-Kabel, dessen Enden mit einem HF-Stecker (N-Typ) versehen sind. Verfügbare Längen: 2, 2.5, 3, 3.5 und 4 m (siehe Abb. 3-19).



## 3.6 Messanordnung am Rohr

Die Auswerteeinheit wird in unmittelbarer Nähe zur FlowCell installiert, damit die HF-Kabel zwischen Auswerteeinheit und Sonde möglichst kurz sind. Je kürzer die Kabelverbindung, desto besser ist die Stabilität der Messung. Das HF-Kabel muss jedoch mindestens 2 m lang sein. Die Standardlänge ist 2 m und die maximale Länge der HF-Kabel beträgt 10 m.

Die FlowCell wird in das bestehende Rohrleitungssystem oder in einem Bypass integriert. Die Orientierung der FlowCell kann sowohl senkrecht als auch waagerecht sein. Um jedoch mögliche sedimentäre Ablagerungen zu vermeiden, wird der senkrechte Einbau in eine Steigleitung bevorzugt (siehe Abb. 3-20).

Der Einbau sollte möglichst nahe einer Probeentnahmestelle sein, damit die Probenahme repräsentativ für die Kalibrierung ist.

Für eine Produkttemperaturkompensation sollte ein repräsentatives Temperatursignal (Stromsignal oder PT100) an die Auswerteeinheit angeschlossen werden.



## 3.7 Messanordnung am Behälter

Die Auswerteeinheit wird in unmittelbarer Nähe zur Behältersonde installiert, damit die HF-Kabel zwischen Auswerteeinheit und Sonde möglichst kurz sind. Je kürzer die Kabelverbindung, desto besser ist die Stabilität der Messung. Das HF-Kabel muss jedoch mindestens 2 m lang sein. Die Standardlänge ist 2 m und die maximale Länge der HF-Kabel 10 m.

Der Einbau sollte möglichst nahe einer Probeentnahmestelle sein, damit die Probenahme repräsentativ für die Kalibrierung ist. Für eine evtl. erforderliche Produkttemperaturkompensation sollte ein repräsentatives Temperatursignal (Stromsignal oder PT100) an die Auswerteeinheit angeschlossen werden.

Als Beispiel ist in Abb. 3-21 die Messanordnung an einem Prozessbehälter dargestellt. Die Sonde wird durch Flanschankopplung an der Behälterwand so befestigt, dass die beiden Messstäbe in das Messgut getaucht sind.



*Abb. 3-21: Beispiel Messanordnung am Prozessbehälter* 

# Kapitel 4. Inbetriebnahme

#### 4.1 Transport

 

 HINWEIS
 Achtung, mögliche Sachschäden! Beim Transport können Systemteile beschädigt werden.

 Alle Komponenten in der Originalverpackung transportieren. Teile vor Erschütterungen schützen. Insbesondere die Messstäbe der Behältersonden müssen vor mechanischen Stößen geschützt werden!

 Nach dem Auspacken sind alle Teile laut Packliste auf Vollständigkeit und Beschädigung zu kontrollieren und falls erforderlich zu reinigen. Bei Schäden sofort das Transportunternehmen und den Hersteller verständigen.

Das Gewicht der Systemkomponenten kann je nach Ausführung über 25 Kg betragen. Tragen Sie daher Sicherheitsschuhe.

MicroPolar (++) LB 566
🕄 WARNUNG

## 4.2 Inbetriebnahme FlowCell

## 4.2.1 Installation der FlowCell

Verletzungsgefahr durch austretendes Betriebsmedium

Die FlowCell muss richtig montiert werden.

Insbesondere muss das richtige Anzugsmoment für die Flansch-Schrauben beachtet werden.

Voraussetzungen für die hygienische Installation der FlowCell:

- > Die Einbaulage muss eine Selbstentleerung gewährleisten.
- Das Gerät wurde für Cleaning in Place (CIP)-Anwendungen entwickelt und darf zur Reinigung nicht zerlegt werden.
- Verwenden Sie keine Reinigungsgeräte, die kratzend oder scheuernd sind, um eine Beschädigung der Produktkontaktflächen zu vermeiden.
- Verwenden Sie keine aggressiven Reinigungsmittel oder Chemikalien, die die Produktkontaktfläche angreifen können.
- Der Betreiber ist verpflichtet, die FlowCell vor dem Kontakt mit Lebensmitteln ordnungsgemäß zu reinigen.
- Um die Anforderungen für die EHEDG-Zertifizierung zu erfüllen, muss der Sensor mit Prozessanschlüssen gemäß dem EHEDG-Positionspapier "Easy Cleanable Pipe Couplings and Process Connections" (www.ehedg.org) verbunden werden.

Beachten Sie folgende Punkte bei der Montage der FlowCell:

- Die FlowCell ist an einer geeigneten Stelle in das Rohrleitungssystem integriert. Vorzugsweise sollte unmittelbar nach der FlowCell eine Materialentnahme für die Kalibrierung möglich sein.
- Die FlowCell sollte möglichst in einer senkrechten Steigleitung eingebaut werden. Es muss sichergestellt sein, dass sich an den Rohrwandungen keine Materialablagerungen bilden können und im Produkt keine Blasenbildung auftritt.

Bei horizontaler Anordnung ist auf die Entleerbarkeit der Rohrleitung zu achten, dazu wird die FlowCell entsprechend der Abb. 4-2 eingebaut.

Vor und nach der FlowCell sollte eine gerade Rohrstrecke von min. 300 mm gleicher Nennweite vorhanden sein, um ein möglichst homogenes Strömungsprofil zu gewährleisten und um evtl. auftretende Mikrowellenreflexionen in der Rohrleitung zu vermeiden. Diese Rohrbereiche müssen ebenfalls frei von Einbauten sein.

- In dem Produkt sollten möglichst keine Gaseinschlüsse vorhanden sein. Sollten sich diese nicht vermeiden lassen, so ist ein Rohrleitungsdruck von min. 4 bar erforderlich, um den Einfluss der Gasblasen vernachlässigen zu können. Beachten Sie den max. zulässigen Betriebsdruck, siehe Kapitel 6.2 Technische Daten Sensoren.
- Die Hochfrequenzkabel sollten vorzugsweise von unten an die FlowCell angeschlossen werden, um nachfließendes Wasser nicht auf die Anschlussbuchsen zu führen.
- Die Mess- und Referenzkabel sollten soweit wie möglichst den gleichen Weg verfolgen, um gleicher Temperatur ausgesetzt zu sein und nicht mit warmen Rohrleitungen in Berührung kommen. Am besten verlegt man die HF-Kabel durch ein gemeinsames Schutzrohr. Bei dem HF-Kabel quad übernimmt diese Funktion in guter Näherung bereits das Wellrohr.







## 4.2.2 Installation der AWE

Beachten Sie folgende Punkte bei der Montage der AWE:

- Die AWE muss entsprechend der HF-Kabellänge in der Nähe der Mikrowellensonde positioniert werden.
- Die AWE muss vibrationsarm aufgestellt werden. In einigen Fällen empfiehlt es sich, die AWE an einem von dem Rohrleitungssystem getrennten Ständer anzubringen.
- Sehen Sie bei der Geräteinstallation eine Trennvorrichtung vor, mit der Sie das Gerät einfach und schnell von der Netzversorgung trennen können.
- Sehen Sie eine automatische Trennvorrichtung (Leitungsschutzschalter) vor, die das Gerät im Fehlerfall in 0,03 Sekunden vom Netz trennt. Die Trennvorrichtung muss auf den Leitungsquerschnitt der Zuleitung angepasst sein, mindestens aber für 1 A Dauerstrom ausgelegt sein.
- Wenn Sie die AWE an einer warmen/heißen Wand montieren müssen, so verwenden sie eine Distanzschiene, um die Wärmestrahlung und –leitung zu minimieren. Siehe Abb. 4-3.
- Bei Aufstellen der AWE im Freien muss die AWE vor direkten Sonnstrahlen und Regen geschützt werden, etwa durch ein genügend großes Schutzdach.



*Abb. 4-3: Ansicht von oben: Montage der AWE an einer heißen Behälterwand* 

## 4.2.3 Anschluss der HF-Kabel

Die FlowCell wird mit der AWE über die HF-Kabel verbunden. Dabei gibt es zwei verschiedene Möglichkeiten:

Variante 1:	$1 \times HF$ -Kabel quad und $1 \times N$ -Verbinder
Variante 2:	2 x HF-Kabel (als Messkabel)
	1 x HF-Kabel (als Referenzkabel)

Die Entscheidung der Kabelvariante trifft Berthold Technologies bei der Projektierung. Sie ist Abhängig von der Applikation und gewünschten Entfernung zwischen AWE und FlowCell.

Voraussetzung für eine einwandfreie Messung ist die korrekte Kabelverlegung! Bitte beachten Sie dafür folgende Hinweise:

	<b>TIPP</b> Das Kabel auf seiner gesamten Länge (Wellrohr- und Einzelkabel- abschnitt nach Splitting), darf nicht mit heißen Leitungen in Berüh- rung kommen, z.B. direkter Kontakt mit der Apparatewand (uniso- liert). Nur so ist sichergestellt, dass alle Einzelkabel denselben Um- gebungsbedingungen ausgesetzt sind und damit die Kom-pensation der Kabeldrift einwandfrei arbeiten kann.
HINWEIS	Das HF-Kabel niemals knicken! Der Biegeradius sollte 100 mm nicht unterschreiten. Fixieren Sie das HF-Kabel mit Kabelbindern oder an- deren geeigneten Mitteln, so dass das Kabel nicht mehr bewegt wer- den kann!
HINWEIS	Achten Sie bei der Montage darauf, dass das Ende des Wellrohres auf der Sondenseite nach unten gebogen ist. Auf diese Weise wird das Eindringen von Flüssigkeiten in das Wellrohr verhindert.
Für Variante 1	Das HF-Kabel quad und die HF-Anschlüsse an der AWE sind beschrif- tet. Verbinden Sie gemäß der Abb. 4-4 die FlowCell mit der AWE, dabei werden nur gleichlautende Beschriftungen zusammengefügt. Die beiden Anschlüsse an der FlowCell sind nicht beschriftet, die Zu- ordnung der beiden Kabelstecker M-Tx und M-Rx ist beliebig. Die Kabelstecker R-Tx und R-Mx werden mit dem N-Verbinder verbun- den (kurzgeschlossen).
Für Variante 2	Verbinden Sie gemäß der Abb. 4-5 die FlowCell mit der AWE, dabei beachten Sie, dass das Referenzkabel (Ringleitung) an R-Tx und R-Mx angeschlossen ist.
	Das Referenzkabel muss so lang sein wie die Summe der beiden Messkabel.



Drehen Sie alle Schraubverbindungen der HF-Kabel handfest an (2 Nm = 0,2 kg/m) ! Vor dem Festziehen die Kabel vorsichtig per Hand anschrauben **-Vorsicht! Verschraubung verkantet leicht.** 

Die Schraubverbindungen gelegentlich auf festen Sitz überprüfen. Bei vibrierenden Installationen könnte sich sonst die Schraubverbindung lockern und damit Messungenauigkeiten oder Korrosionen der Anschlüsse verursachen.

Während die Kabel nicht angeschlossen sind, müssen die Koaxialbuchsen mit Kunststoffkappen und die Kabelstecker durch geeignete Maßnahmen gegen Feuchte und Verschmutzung geschützt werden. 

## 4.3 Inbetriebnahme Behältersonde

## 4.3.1 Installation der Behältersonden

Verletzungsgefahr durch austretendes Betriebsmedium

Die Behältersonde muss richtig montiert werden.

Insbesondere muss das richtige Anzugsmoment für die Flansch-Schrauben beachtet werden.

Beachten Sie folgende Punkte bei der Montage der Sonden:

- Der Einbauort ist so zu wählen, dass eine gute Durchmischung, ein homogenes Produkt und keine Blasenbildung an der Sonde gewährleistet sind. Ein Probenhahn sollte in nächster Nähe für repräsentative Probeentnahmen zur Verfügung stehen.
- Die Sonde muss so am Behälter angeflanscht werden, dass das Messgut zwischen den beiden Messstäben hindurch fließt. D. h. die Gabel (beide Messstäbe) muss senkrecht zum Materialstrom angeordnet sein.
- Der Abstand zwischen der Messstabspitzen und jeglicher metallisierten Wandung (Heizstäbe, Rührer, Behälterwand) sollte mindestens 60 mm betragen.
- Für den Sondeneinbau sind folgende Einbaulochmaße im Anschlussflansch erforderlich:

Flansch	Minimales Einbaulochmaß $\varnothing$ (mm)
DN 65 / PN 6	$102\pm0.3$
sonst	$104\pm0.5$

- Weitere Einbaumaße, wie die erforderliche Einbautiefe, befinden sich in Kapitel 8.6 und 8.7 Montageblätter.
- Verwenden Sie die zugehörige Flachdichtung (Standard Zubehör). Diese gleicht geringe Oberflächentoleranzen in dem Anschlussflansch gut aus.

Installation am Prozessbehälter

Die Abb. 3-26 zeigt die Position der Behältersonde am Behälter. Die Position gilt gleichermaßen für die Spülsonde. Beachten Sie die Hinweise unter *Kapitel 3.9 Messanordnung am Behälter*.

Hilfreich für die Montage ist das Montageblatt unter *Kapitel 8.6 und 8.7*. Auf diesem sind die Montageinformationen übersichtlich zusammengefasst.

#### Demontage bei festgeklebtem Flansch

Prozessbedingt kann es vorkommen, dass der Sondenflansch am Prozessflansch festklebt. Entfernen Sie in diesem Fall bei der Demontage zunächst die beiden Ringschrauben. Drehen Sie dann in deren Bohrungen zwei längere Schrauben (M8) bis zum Prozessflansch ein. Wenn Sie die Schrauben nun weiterdrehen, drücken Sie die Sonde vom Prozessflansch weg.

### Installation in der Rohrleitung

Die Behältersonden können mit Hilfe eines Adapterflanschs in Rohrleitungen mit einer Nennweite  $\geq$  200 mm eingesetzt werden. Beachten Sie dabei die Position und die Orientierung der Behältersonde, siehe dazu die technischen Zeichnungen unter *Kapitel 8.4.7 und 8.5.6 Einbausituation in Rohrleitungen*.

### Anschließen der Spülleitungen

Die Behältersonde mit Spülvorrichtung besitzt zwei Spülanschlüsse mit G3/8" Innengewinde (DIN ISO 228-1). Die Spülanschlüsse sind ausschließlich an diesem Gewinde abzudichten. Eine Abdichtung zum Sondendeckel, etwa mit Silikon ist unzulässig.

Wird der Spülanschluss an der Behältersonde nicht angeschlossen, muss er durch einen Blindstopfen verschlossen werden.

### Spülparameter (nur für Spülsonde)

Für die Spüleinstellungen, wie z.B. Spülintervall und Spüldauer ist der Grad der Ablagerung bzw. Inkrustation maßgeblich. Die Spülparameter müssen auf das Produkt und auf den Prozess abgestimmt werden.

Folgende produkt- und prozessunabhängige Spülparameter müssen zwingend erforderlich eingehalten werden:

Spüllösung	Wasser, Kondensat
Temperatur der	Maximal 120 °C
Spüllösung	
Druck	≥ 3 bar, maximal 8 bar
Anschlüsse	2 x G3/8" Innengewinde (DIN ISO 228-1)
Zuleitung	≥ 1/2 Zoll

**Allgemein gilt:** Die Spülanschlüsse können gleichzeitig oder zeitlich versetzt gespült werden. Die Spülparameter gelten für jeden Spülanschluss.





Bei der Spüldauer sind eventuelle Trägheiten des Systems, z.B. Ventilöffnungen zu berücksichtigen. Die Spülleitungen müssen gut wärmeisoliert sein, um eine anfänglich kältere Spüllösung zu vermeiden.

Wassermengen

Bei einem Spüldruck von 5 bar beträgt die Wassermenge pro Spülanschluss ca. 0.8 Liter pro Sekunde.

#### 4.3.2 Installation der AWE

Installation der AWE wie unter Kapitel 4.2.2 beschrieben.

### 4.3.3 Anschluss der HF-Kabel

Die Behältersonde wird mit der AWE über das HF-Kabel quad verbunden.

Voraussetzung für eine einwandfreie Messung ist die korrekte Kabelverlegung! Bitte beachten Sie dafür folgende Hinweise:



Das Kabel auf seiner gesamten Länge (Wellrohr- und Einzelkabelabschnitt nach Splitting), darf nicht mit heißen Leitungen in Berührung kommen, z.B. direkter Kontakt mit der Apparatewand (unisoliert). Nur so ist sichergestellt, dass alle Einzelkabel denselben Umgebungsbedingungen ausgesetzt sind und damit die Kom-pensation der Kabeldrift einwandfrei arbeiten kann.

#### HINWEIS

Das HF-Kabel niemals knicken! Der Biegeradius sollte 100 mm nicht unterschreiten. Fixieren Sie das HF-Kabel mit Kabelbindern oder anderen geeigneten Mitteln, so dass das Kabel nicht mehr bewegt werden kann!

Das HF-Kabel, die HF-Anschlüsse an der AWE und Sonde sind beschriftet. Verbinden Sie gemäß Abb. 4-6 die Sonde mit der AWE, dabei werden nur gleichlautende Beschriftungen zusammengefügt.



Auswerteeinheit, Unterseite



Drehen Sie alle Schraubverbindungen der HF-Kabel handfest an (2 Nm = 0,2 kg/m) ! Vor dem Festziehen die Kabel vorsichtig per Hand anschrauben - **Vorsicht! Verschraubung verkantet leicht.** 

Die Schraubverbindungen gelegentlich auf festen Sitz überprüfen. Bei vibrierenden Installationen könnte sich sonst die Schraubverbindung lockern und damit Messungenauigkeiten oder Korrosionen der Anschlüsse verursachen.

Während die Kabel nicht angeschlossen sind, müssen die Koaxialbuchsen sofort mit Kunststoffkappen und die Kabelstecker durch geeignete Maßnahmen gegen Feuchte und Verschmutzung geschützt werden.

## 4.4 Anschluss der AWE

	Elektrische Gefahren:
	Während der Installation und für Servicearbeiten ist die Stromver- sorgung abzuschalten, damit Berührungen mit spannungsführen- den Teilen verhindert werden. Vor jedem Öffnen des Gerätes ist die Stromzufuhr zu unterbre- chen. Arbeiten am geöffneten und unter Spannung stehenden Ge- rät sind verboten
HINWEIS	Achtung! Mögliche Gefährdung, Sachschäden! Betrifft den Gerätetyp: LB 566-12 MicroPolar ++ (IdNr. 51833-02)
	Bei Anschluss der 24 V DC Hilfsenergie, müssen die + und – Pole richtig angeschlossen werden. Es besteht kein Verpolungsschutz!
	· · · · · · · · · · · _ · · _ ~ _ · _ ~ _ ~
	Der Leitungsquerschnitt für den Netzanschluss muss min. 1.0 mm <sup>2</sup> betragen.
	<ul> <li>Der Leitungsquerschnitt für den Netzanschluss muss min. 1.0 mm<sup>2</sup> betragen.</li> <li>Schließen Sie alle gewünschten Ein- und Ausgangssignale ent- sprechend der nachfolgenden Seiten an die Klemmleiste an. Be- nutzen Sie dabei die vorgesehenen M-Durchführungen, damit der Schutzgrad erhalten bleibt.</li> </ul>
	<ul> <li>Der Leitungsquerschnitt für den Netzanschluss muss min. 1.0 mm<sup>2</sup> betragen.</li> <li>Schließen Sie alle gewünschten Ein- und Ausgangssignale entsprechend der nachfolgenden Seiten an die Klemmleiste an. Benutzen Sie dabei die vorgesehenen M-Durchführungen, damit der Schutzgrad erhalten bleibt.</li> <li>Überprüfen Sie, ob die auf dem Typenschild zugelassene Spannung mit Ihrer Versorgungsspannung übereinstimmt.</li> </ul>
	<ul> <li>Der Leitungsquerschnitt für den Netzanschluss muss min. 1.0 mm<sup>2</sup> betragen.</li> <li>Schließen Sie alle gewünschten Ein- und Ausgangssignale entsprechend der nachfolgenden Seiten an die Klemmleiste an. Benutzen Sie dabei die vorgesehenen M-Durchführungen, damit der Schutzgrad erhalten bleibt.</li> <li>Überprüfen Sie, ob die auf dem Typenschild zugelassene Spannung mit Ihrer Versorgungsspannung übereinstimmt.</li> <li>Schließen Sie das spannungsfrei geschaltete Netzkabel an die Klemmen 3(L1), 2(N) und 1(PE) an.</li> </ul>
	<ul> <li>Der Leitungsquerschnitt für den Netzanschluss muss min. 1.0 mm<sup>2</sup> betragen.</li> <li>Schließen Sie alle gewünschten Ein- und Ausgangssignale entsprechend der nachfolgenden Seiten an die Klemmleiste an. Benutzen Sie dabei die vorgesehenen M-Durchführungen, damit der Schutzgrad erhalten bleibt.</li> <li>Überprüfen Sie, ob die auf dem Typenschild zugelassene Spannung mit Ihrer Versorgungsspannung übereinstimmt.</li> <li>Schließen Sie das spannungsfrei geschaltete Netzkabel an die Klemmen 3(L1), 2(N) und 1(PE) an.</li> <li>Überprüfen Sie, ob sich der Testschalter (Netzunterbrechung) in Position "EIN" befindet (siehe Abb. 5-1).</li> </ul>

### 4.4.1 Anschlussbelegung der Steckerleiste

Auf der Steckerleiste der AWE sind folgende Anschlüsse zu finden:



Abb. 4-7: Anschlussplan LB 566

Netzanschluss: Klemmen 3 (L1, +), 2 (N, -) und 1 (PE, 🛞)

Für MicroPolar, je nach Geräteausführung, siehe Typenschild auf der Gehäuse Außenwand.

- 1.) 100...240 V AC, 45...65 Hz
- 2.) 24 V DC: 18...36 V
  - 24 V AC: -20%, +5%, 40...440 Hz

Für MicroPolar ++, je nach Geräteausführung, siehe Typenschild auf der Gehäuse Außenwand.

- 1.) 100...240 V AC, 45...65 Hz
- 2.) 24 V DC: 18...36 V, kein Verpolungsschutz

#### Stromeingang Nr. 1 (Klemmen 20+ und 8-), isoliert Stromeingang Nr. 2 (Klemmen 22+ und 10-), nicht isoliert

Eingabe als 0/4 - 20 mA Signal. Z.B. zur Temperaturkompensation oder Referenzsignalaufzeichnung.

#### Stromausgang Nr. 1 (Klemmen 27+ und 15-), isoliert

Ausgabe als 4 - 20 mA Signal. Ausgabemöglichkeiten: Konzentrationen (1 / 2), Stromeingangssignale (1 / 2), PT100 Signal, Massenstrom

#### Stromausgang Nr. 2 (Klemmen 19+ und 7-), isoliert

Ausgabe als 0/4 - 20 mA Signal. Ausgabemöglichkeiten entsprechend Stromeingang Nr. 1.

#### PT100 (Klemmen 23+ und 11-)

Anschluss zur Temperaturmessung.

() Klemmen-Nr.

Im Inbetriebnahmefall der Behältersonde schließen Sie das 4-adrige Kabel des PT100 an die Steckerleiste der AWE an (siehe Abb. 4-8). Verbinden Sie das andere Ende des Kabels mit dem entsprechenden Stecker an der Behältersonde (siehe Abb. 4-6).



*Abb. 4-8: PT100 Anschluss Behältersonde* 



Das PT100-Kabel hat im Lieferzustand eine Länge von 10 Metern. Es wird empfohlen, das Kabel auf die benötigte Länge zu kürzen.

Fassen Sie die beiden Aderpaare entsprechend der Abbildung 4-8 zusammen.

Nach dem Anschluss des PT100-Kabels muss der mitgelieferte Ferritkern um das Kabel gelegt werden. Die Position hierfür ist innerhalb des Gehäuses, möglichst nahe an der Kabeldurchführung.



#### Digitaleingang 1: DI1 (Klemmen 24+ und 12-)

Nur für potentialfreie Kontakte! Konfigurationsmöglichkeiten:

- Keine Funktion
- Messung: Start (geschlossen) und Stopp (offen)

#### Digitaleingang 2: DI2 (Klemmen 25+ und 13-)

Nur für potentialfreie Kontakte! Konfigurationsmöglichkeiten:

- Keine Funktion
- > Mittelwert: Halten (geschlossen) und weiter mitteln (offen)
- Produktauswahl: Produkt 1 (offen) und Produkt 2 (geschlossen)

Abb. 4-9: Ferritkern

### Digitaleingang 3: DI3 (Klemmen 26+ und 14-)

Nur für potentialfreie Kontakte! Konfigurationsmöglichkeiten:

- Keine Funktion
- Probeaufnahme starten, offen: keine Aktion, geschlossen: einmalige Aufnahme startet
- Produktauswahl

### Relais 1: (Klemmen 4, 5 und 6) und Relais 2: (Klemmen 16, 17 und 18)

Wechselkontakte (SPDT), isoliert, Konfigurationsauswahl:

- keine Funktion
- Fehlermeldung
- Messung anhalten
- Grenzwert min. und max.
- Kein Produkt
- > Beladungsgrenze unterschritten

#### RS485 Schnittstelle (Klemmen 21 (RS1) und 9 (RS2)) und RS232 Schnittstelle (an Geräteunterseite)

Serielle Datenschnittstelle zur Ausgabe der Live-Daten (alle Messdaten für jeden Sweep (Messzyklus)), des Protokolls und Datenlogs. Datenformat: Datenübertragungsrate 38400 Baud, 8 Datenbits, 1 Stoppbit, keine Parität, kein Handshake

## 4.4.2 Digitale Ausgänge, Relais

Den Status der Messung wird über zwei Relais ausgegeben:

- Fehler
- Alarm (Alarm min. und max.)
- ➢ Kein Produkt

Unter Menüpunkt Plausibilität, lässt sich für die Pausenerkennung (z.B. Prozesspause, kein Produkt vorhanden) eine min. Dämpfung angeben, bei Unterschreitung wird über Relais "kein Produkt" signalisiert und der Stromausgang fällt auf 0 bzw. 4 mA ab.

- Messung angehalten
- Beladungsunterschreitung

Der jeweilige Schaltzustand wird auch über LED's auf der Frontplatte signalisiert (LED's: Signal 1 und 2).



Relais Nr.	Fehler, Alarm, kein Pro- dukt, Messung angehal- ten, Beladungsunterschreitung, stromloser Zustand	Normal
1	4 0	4 0 5 0 com 6 0
2	16 17 18 	16 0

Die Relais mit Wechselkontakten können wahlweise als Schließer, Klemmen 4 & 5 (offen bei Fehler, Alarm ...) oder als Öffner, Klemmen 5 & 6 (geschlossen bei Fehler, Alarm ...) betrieben werden.



## Kapitel 5. Wartung

## 5.1 Allgemeines

Sollte eine Fehlfunktion des Messsystems vorliegen, muss nicht unbedingt ein Gerätedefekt vorliegen. Oft führen Fehlbedienungen, ungeeignete Installationen oder Unregelmäßigkeiten im Messprodukt zu Messfehlern. Falls doch einmal eine Fehlfunktion vorliegt, erleichtert das Messgerät die Fehlerbeseitigung durch Fehlermeldungen in der Anzeige, womit Bedienfehler und elektronische Defekte signalisiert werden.

Defekte Module der Auswerteeinheit können normalerweise nicht repariert werden, sondern sind auszutauschen. Das Mikrowellen-Modul ist mit einer Abschirmhaube fest verschraubt und darf nicht geöffnet werden.

Für den Fall der Geräteentsorgung, wenden Sie sich bitte an den Service von Berthold und beantragen einen Recyclingpass.

## 5.2 Verschleißteile

Die AWE besitzt keine Verschleißteile und Komponenten, die besonderer Wartung bedürfen.

Die Kunststoffkappen der Messstäbe der Behältersonden und die PEEK Mikrowellenfenster der FlowCell können je nach Messgut mit der Zeit einen Abrieb erfahren. Ein geringer bis mittlerer Abrieb beeinflusst die Messung vernachlässigbar gering bzw. wird durch die Kalibrierung ausgeglichen. Überprüfen Sie daher in einem Intervall von etwa 2 Jahren die Verschleißteile. Bei starkem Verschleiß können die Kunststoffkappen der Messstäbe und die Mikrowellenfenster der FlowCell ausgetauscht werden.

#### Austausch Kunstoffkappen Behältersonde

- 1 Schrauben Sie die beiden Kunststoffkappen von den Messstäben ab und entfernen Sie die vier Dichtungsringe (siehe Abb. 5-1, gelber Pfeil).
- 2 Reinigen Sie die vier Nuten der Dichtungsringe mit einem fusselfreien, ggf. angefeuchteten Tuch.
- **3** Legen Sie pro Messstab zwei neue Dichtungsringe in die Nuten.
- 4 Schrauben Sie die beiden neuen Kunststoffkappen auf.





Abb. 5-1: Kunststoffkappen der Messstäbe tauschen

Kappenset für Behältersonde	
IdNr. 66049-S	2 Stück PEEK Kunststoffkappen mit 4 Dichtungsringen

#### Austausch Mikrowellenfenster FlowCell

Gehen Sie wie folgt vor:

- 1 Lösen Sie die Verschlussklammer (siehe Abb. 5-2, Pos. 1).
- 2 Entfernen Sie die Antenne (siehe Abb. 5-2, Pos. 2), das Mikrowellenfenster (siehe Abb. 5-2, Pos. 3) und die O-Ringe.
- 3 Befestigen Sie das neue Mikrowellenfenster, die neuen O-Ringe und die Antenne wieder mit der Verschlussklammer an der Flow-Cell, entsprechend der Darstellung in Abb. 5-3. Ab einem Fettgehalt von 8 % muss anstelle des EPDM-O-Rings ein HNBR-O-Ring verwendet werden.



Abb. 5-2: Antenne und Mikrowellenfenster

Mikrowellenfenster

Abb. 5-3: Montage



*Abb. 5-4: Montage Mikrowellenfenster mit PEEK* 

HINWEIS

HINWEIS

Es wird empfohlen, immer den gesamten Satz O-Ringe auszutauschen.

Ab einem Fettgehalt von 8 % muss anstelle des EPDM-O-Rings ein HNBR-O-Ring verwendet werden.

Mikrowellenfenster-Set für FlowCell	
IdNr. 66624-S	2 Stück PEEK Mikrowellenfenster mit 6 Dichtungs-O-Ringen
IdNr. 66625-S	2 Stück PEEK Mikrowellenfenster GF30 glasfaserverstärkt mit 6 Dichtungs-O-Ringen
IdNr. 75514-S	2 Stück PEEK EHEDG Mikrowellenfenster mit 8 Dichtungs-O-Ringen

## 5.3 Gerätereinigung

Reinigen Sie alle Systemkomponenten ausschließlich mit einem angefeuchteten Tuch ohne chemische Reinigungsmittel. Die produktberührenden Teile (im Normalbetrieb) können mit warmen Wasser gereinigt werden, dabei sind die Temperaturgrenzen siehe *Kapitel 6.2 Technische Daten Sensoren* zu berücksichtigen.



## 5.4 Batterie

Falls das Messgerät LB 566 längere Zeit ohne Netzversorgung (Stromausfall oder vom Netz getrennt) ist, wird die Systemuhr über die Lithium-Batterie auf dem Motherboard versorgt.

Ist die Batteriespannung nicht mehr ausreichend, wird nach einem Wiedereinschalten der AWE die Fehlermeldung CODE 14 "Batteriespannung" angezeigt. Nach Quittieren der Fehlermeldung arbeitet das Gerät einwandfrei weiter, jedoch sollte das Datum und die Uhrzeit überprüft und gegebenenfalls korrigiert werden. Messdaten, die über eine der seriellen Schnittstellen ausgegeben werden, können durch fehlerhafte Datums- und Uhrzeit-Angaben unbrauchbar werden. Wir empfehlen die Batterie umgehend zu tauschen.

Die Lebensdauer der Batterie, selbst bei Dauerbeanspruchung, beträgt ca. 8 Jahre. Ein Austausch der Batterien darf nur bei spannungsfreiem Gerät vorgenommen werden.

Batterietyp: 3 Volt Lithium Cell (Knopfzelle), Type CR2032

## 5.5 Sicherung austauschen

Die Netzsicherungen des LB 566 befinden sich im Wandgehäuse. Die Sicherungen dürfen nur im spannungslosen Zustand getauscht werden und müssen dem angegebenen Wert entsprechen.

Sicherungen nur mit korrektem Wert verwenden, siehe Kapitel 6.1

HINWEIS

Ersatzsicherungen müssen die vom Gerätehersteller vorgegebenen Werte aufweisen. Ein Kurzschließen oder manipulieren ist nicht gestattet.



Abb. 5-4: Ansicht bei geöffneten Wandgehäuse MicroPolar



# Kapitel 6. Technische Daten

Allgemeine Spezifikationen	
Verfahren	Mikrowellen-Transmissionsmessung
Arbeitsfrequenz	2.4 – 2.5 GHz (ISM-Band), je nach Landes- bestimmung
Sendeleistung	MicroPolar: < 0.1 mW (< - 10 dBm) MicroPolar ++: < 10 mW (< 10 dBm) Alles koaxiale Leitungsleistung
Applikationen	Konzentrations-, Feuchtemessung in Behäl- tern und Rohrleitungen

## 6.1 Technische Daten AWE

Auswerteeinheit	
Gehäuse	Wandgehäuse aus Edelstahl, siehe Maßbild unter Kapitel 8 MicroPolar: HxBxT: 300 x 323 x 140 mm MicroPolar++: HxBxT: 400 x 338 x 170 mm
Schutzart	IP 65
Gewicht	MicroPolar: ca. 6.5 kg MicroPolar ++: ca. 8.0 kg
Umgebungsbedin- gungen im Betrieb	-20 +60 °C (253333 K), keine Betau- ung Relative Luftfeuchte: max. 80 % Höhenlage: max. 2000 m
Umgebungsbedin- gung bei Lagerung	-20 +70 °C (253343 K), keine Betau- ung, Relative Luftfeuchte: max. 80 %
Erreichbare Genauigkeit	
Anzeige	Dot-Matrix LC Display, 114 mm x 64 mm, 240 x 128 Pixel, mit Hintergrundbeleuch- tung, automatische Kontrasteinstellung
Tastatur	Freizugängliche Folientastatur, lichtstabil und witterungsfest: alphanumerische Tasta- tur und 4 Softkeys (software zugeordneten Tasten)



Hilfsenergie	<ul> <li>Für MicroPolar, je nach Geräteausführung:</li> <li>1.) 100 240 V AC, 45 65 Hz</li> <li>2.) 24 V DC: 1836 V; 24 V AC: -20%, +5%, 40 440 Hz</li> <li>Für MicroPolar++, je nach Geräteausführung:</li> <li>1.) 100240 V AC, 4565 Hz</li> <li>2.) 24 V DC: 1836 V, kein Verpolungsschutz</li> </ul>
Leistungsaufnahme	Für MicroPolar: max. 30 VA (AC/DC), je nach Konfiguration Für MicroPolar ++: max. (48/60) VA (AC/DC), je nach Konfigura- tion
Sicherungen	Für MicroPolar: 2 x 2.0 A / Tr. Für MicroPolar ++: 2 x 2.0 A / Tr. bei 100 240 V AC oder 2 x 6.3 A / Tr. bei 24 V DC
Batterietyp	3 V Lithium Knopfzelle, Typ CR2032
Messwert	z.B. Konzentration, Trockengehalt
Ein- und Ausgänge	
Leiterquerschnitt	min. 1.0 mm <sup>2</sup> (Netzanschluss)
Kabeleinführung	2 x M20x1.5 für Kabel 514 mm (je nach Einsatz) 4 x M16x1.5 für Kabel 58 mm (je nach Ein- satz)
Sensoranschluss	Ein- und Ausgänge für Mess- und Referenzkanal, 50 $\Omega$ N-Buchse
HF-Kabel	Kabellängen: 2, 4, 6 und 10 m; 50 $\Omega$ ; beidseitig mit 4 N-Stecker
Stromeingang	2 x Stromeingang 0/420 mA, Bürde 50 $\Omega$ , 1x isoliert, 1x Gerätemasse z.B. für Temperaturkompensation
Stromausgang	Stromausgang 1: 420 mA, Bürde max. 800 $\Omega$ , isoliert Stromausgang 2: 0/420 mA, Bürde max. 800 $\Omega$ , isoliert z.B. zur Messwert- oder Temperaturausgabe
PT100 Anschluss	Messbereich: -50 +200 °C (223 473 K); Messtoleranz: < 0.4 °C

Digitaler Eingang	3 x digitale Eingänge (DI13), für potential- freie Kontakte (nicht an einer Spannungsver- sorgung anschließen). <u>Konfigurationsmöglichkeiten:</u> DI1: Keine, Messung Start/Stopp DI2: Keine, Messung halten, Produktauswahl DI3: Keine, Probeaufnahme, Produktauswahl
	<ul> <li><u>Funktionsbeschreibung:</u></li> <li>1. Messung (Start/Stopp), <u>offen:</u> Messung gestoppt, <u>geschlossen:</u> Messung gestartet bzw. Messung läuft</li> <li>2. Messung anhalten, <u>offen:</u> Messung läuft, geschlossen: Messung angehalten, d h</li> </ul>
	Mittelwerte und Stromausgang sind einge- froren
	3. Produktauswahl über ein DI: offen: Produkt 1 (P1), <u>geschlossen:</u> P2
	Produktauswahl über zwei DI's: <u>DI2 &amp; DI3 offen</u> : P1 <u>DI2 geschlossen &amp; DI3 offen</u> : P2 <u>DI2 offen &amp; DI3 geschlossen</u> : P3 DI2 & DI3 geschlossen; P4
	4. Probeaufnahme: <u>offen</u> : keine Aktionen, geschlossen: einmalige Aufnahme startet
Relaisausgänge	2 x Relais (SPDT), isoliert <u>Konfigurationsmöglichkeiten:</u> - Sammelstörmeldung - Messung anhalten - Grenzwert (min. und max.) - kein Produkt - Beladungsunterschreitung
	Belastbarkeit:AC:max. 400VADC:max. 90WAC / DC:max. 250V, max. 2A, induktionsfrei≥ 150V:Spannung muss erdbezogen sein
	Das an den Relaisausgang verwendete Kabel muss einer Netzanschlussleitung entsprechen.
	Einschränkungen bei 24 V AC/DC (DC: 18 36 V; AC: 24 V +5 %, -20 %) Netzversor- gung, wenn der Schutzleiter nicht an die Klemme 1 (PE) angeschlossen ist: AC: max. 50 V DC: max. 70 V
Serielle	RS232 auf der Geräteunterseite,
Schnittstellen	RS485 über Klemmenleiste Datenformat: 38400 Bd, kein Handshake, 8 Datenbits, 1 Stoppbit, keine Parität

FlowCell	
Anwendung	Mikrowellen-FlowCell verschiedener Nennweiten und Flansche zur Konzentrationsmessung an Rohrleitungen
Material	Inline-Gehäuse aus Edelstahl 1.4404 poliert (AISI 316L)
	Mikrowellenfenster aus PEEK
	Produktberührende Dichtung aus EPDM
Prozessankopplung	Zwei Varianten:
	1. Hygiene Milchrohrverschraubung DIN 11853-1 EHEDG-Zulassung
	2. Flansch nach EN 1092-1/11 (V-Flansch) FDA-zugelassene Materialien
	Optionale Adapter für die V-Flansch Variante mit ASA-Flansch
Prozessdruck	Bis 16 bar (relativ)
Temperaturbereich	Produkttemperatur: 10130 °C (283403 K), kurzzeitig bis 140°C (413 K)
	Umgebungstemperatur: -2060 °C (253333 K) Lagertemperatur: 1080 °C (283353 K)
Anschlüsse	2 x HF-Anschlüsse: N-Buchse, 50 $\Omega$
	Für HF-Kabel mit max. 10 m Länge
Varianten	Rohrnennweiten von 50 150 mm
Маве	Siehe Maßzeichnungen unter Kapitel 8.

## 6.2 Technische Daten Sensoren

Bezeichnung	IdNr.	Nennweite [mm]	Flansch	Druck [bar]
LB 5660-102-00x	66744-001	50	DN 50 / PN 16	
LB 5660-202-00x	66744-002	65	DN 65 / PN 16	
LB 5660-302-00x	66744-003	80	DN 80 / PN 16	
LB 5660-402-00x	66744-004	100	DN 100 / PN 16	
LB 5660-502-00x	66744-005	125	DN 125 / PN 16	
LB 5660-602-00x	66744-006	150	DN 150 / PN 16	
LB 5660-402-200 FlowCell Vfl. FDA Eintauchkappe	66744-031	100	DN 100	16
LB 5660-502-200 FlowCell Vfl. FDA Eintauchkappe	66744-032	125	DN 125	
LB 5660-602-200 FlowCell Vfl. FDA Eintauchkappe	66744-033	150	DN 150	

### Übersicht FlowCell mit V-Flansch

## Übersicht FlowCell mit Milchrohrverschraubung

Bezeichnung	IdNr.	Nennweite [mm]	Druck [bar]
LB 5660-112-00x	66744-013	50	
LB 5660-212-00x	66744-014	65	
LB 5660-312-00x	66744-015	80	
LB 5660-412-00x	66744-016	100	
LB 5660-512-00x	66744-017	125	
LB 5660-612-00x	66744-018	150	
LB 5660-412-200 G-BS-M Eintauchkappe (hygie- nisch)	66744-034	100	16
LB 5660-512-200 G-BS-M Eintauchkappe (hygie- nisch)	66744-035	125	
LB 5660-612-200 G-BS-M Eintauchkappe (hygie- nisch)	66744-036	150	



Bezeichnung	IdNr.	Nennweite [mm]	Druck [bar]
LB 5660-132-00X	66744-025	50	
LB 5660-232-00X	66744-026	65	
LB 5660-332-00X	66744-027	80	
LB 5660-432-00X	66744-028	100	
LB 5660-532-00X	66744-029	125	
LB 5660-632-00X	66744-030	150	16
LB 5660-432-200 Eintauchkappe (hygienisch)	66744-037	100	
LB 5660-532-200 Eintauchkappe (hygienisch)	66744-038	125	
LB 5660-532-200 Eintauchkappe (hygienisch)	66744-039	150	

### Übersicht FlowCell mit Schweißenden

#### Übersicht Mikrowellenfenster-Set

Die Mikrowellenfenster der FlowCell sind für Anwendungen mit hohem Abrieb auch in einer glasfaserverstärkten Ausführung erhältlich. Diese ist nicht für den Lebensmittelbereich zugelassen.

IdNr.	Beschreibung
66624-S	2 Stück PEEK Mikrowellenfenster mit 6 Dichtungs-O-Ringen
66625-S	2 Stück PEEK Mikrowellenfenster GF30 glasfaserverstärkt mit 6 Dichtungs-O-Ringen
75514-S	2 Stück PEEK EHEDG Mikrowellenfenster mit 6 dichtenden O-Ringen

### HINWEIS

Sonden mit PEEK EHEDG Mikrowellenfenstern sind erst ab einer Nennweite von DN100 einsetzbar.

#### Übersicht ASA-Flansch-Adapter

Bezeichnung	IdNr.	
ASA-Flansch-Adapter-Set für Flowcell 50	62324	
ASA-Flansch-Adapter-Set für Flowcell 65	62319	
ASA-Flansch-Adapter-Set für Flowcell 80	62328	
ASA-Flansch-Adapter-Set für Flowcell 100	62331	
ASA-Flansch-Adapter-Set für Flowcell 150 62335		
Das Set besteht aus zwei Adaptern, Schrauben und zwei Dichtungen.		

#### Übersicht Inline-Gehäuse, FDA

für Temperatur-, Leitfähigkeitssensor oder Probenahmeventil

Rozeichpung	VFL	G-BS/M
bezeichnung	IdNr.	IdNr.
Inline-Gehäuse für Flowcell 50	67078	67084
Inline-Gehäuse für Flowcell 65	67079	67085
Inline-Gehäuse für Flowcell 80	67080	67086
Inline-Gehäuse für Flowcell 100	67081	67087
Inline-Gehäuse für Flowcell 125	67082	67088
Inline-Gehäuse für Flowcell 150	67083	67089

#### Übersicht Oberflächen-Temperaturfühler

Selbstklebender PT100 Temperaturfühler mit Befestigungsmaterial, Anschlusskabel 10 m, 4-Leiter (lose Enden)

Temperaturbereich: -50 - +200 °C

Bezeichnung	IdNr.
Selbstklebender Temperaturfühler für DN 50	66655
Selbstklebender Temperaturfühler für DN 65	66656
Selbstklebender Temperaturfühler für DN 80	66657
Selbstklebender Temperaturfühler für DN 100	66658
Selbstklebender Temperaturfühler für DN 125	66659
Selbstklebender Temperaturfühler für DN 150	66660

### Übersicht Sensoren

Bezeichnung	IdNr.
Leitfähigkeitssensor hygienisch, Clamp-Flansch	66693
Induktives Leitfähigkeitsmessgerät für flüssige Medien in hygienischen Anwendungen Messbereich: 0-999 mS/cm Prozessanschluss: Clamp-Flansch Betriebsdruck: max. 16 bar Hilfsspannung: 18-36 V DC max. 190 mA Ausgang: 4-20 mA	
Temperatursensor EHEDG, Clamp-Flansch	66694
PT100 Temperatursensor für hygienische An- wendungen Messbereich: -50 - +250 °C Länge 20 mm, Durchmesser 4 mm Prozessanschluss: Clamp-Flansch Mit Anschlusskabel 10 m (lose Enden)	

#### Übersicht Probenahmeventil und Zubehör

Bezeichnung	IdNr.
Probenahmeventil aseptisch, Clamp-Flansch	66738
Aseptisches Inline Probenahmeventil Edelstahl 1.4404 (AISI 316L), Faltenbalg PTFE Auslaufstutzen S-DN 10 ohne Spülanschluss Prozessanschluss: Clamp-Flansch	
Clamp-Blindflansch	66737
Clamp-Blindflansch für Inline-Gehäuse DN 50-150 Edelstahl 1.4306 (AISI 304L)	
Clamp-Klammer	66736
1 Stück Clamp-Klammer für Inline-Gehäuse DN 50-150 Edelstahl 1.4306 (AISI 304L)	

Behältersonden		
Anwendung	Behältersonden mit und ohne Spülvorrichtung zur Konzentrationsbestimmung in Prozessbehältern und in Rohrleitungen mit NW $\geq$ 200 mm.	
Material	Kunststoffkappen, Edelstahl PT100 Anschlusskabel: Silicon / Teflon	
Prozessankopplung	Flansch nach DIN EN 1092 Typ 05 DN65 / PN6, DN 80, 100, 150 / PN16; ASA Flansch 2.5" / 150 PSI (weiter auf Anfrage)	
Prozessdruck	Bis 16 bar, je nach Ausführung	
Temperaturbereich	Produkttemperatur: 10120 °C (283393 K) Umgebungstemperatur: -2060 °C (253333 K) Lagertemperatur: 1080 °C (283353 K)	
Anschlüsse	4 x HF-Anschlüsse: N-Buchse, 50 $\Omega$ Für HF-Kabel mit max. 10 m Länge	
Маве	Siehe Maßzeichnungen unter Kapitel 8.	
Zubehör Dichtungs	sscheibe	
Material	Klingersil C-4400	
Dicke	3 mm	
Kappenset für Behältersonde		
IdNr. 66049-S	2 Stück PEEK Kunststoffkappen mit 4 Dichtungsringen	

## Übersicht Behältersonden und Dichtungsscheiben

Bezeichnung	IdNr.	Flansch	Druck [bar]	IdNr. Dichtungen
LB 5650-01	65464-01	DN 65 / PN6	6	32175
LB 5650-02	65464-02	DN 80 / PN16	16	33717
LB 5650-03	65464-03	DN 100 / PN16	16	46661
LB 5650-04	65464-04	DN 150 / PN16	16	46664
LB 5650-05	65464-05	ASA 2.5" / 150 PSI	16	46665
LB 5650-09	65464-09	ASA 3" / 150 PSI	16	50659
LB 5651-01	65937-01	DN 65 / PN6	6	32175
LB 5651-02	65937-02	DN 80 / PN16	16	33717
LB 5651-03	65937-03	DN 100 / PN16	16	46661
LB 5651-04	65937-04	DN 150 / PN16	16	46664
LB 5651-05	65937-05	ASA 2.5" / 150 PSI	16	46665



## 6.3 Technische Daten HF-Kabel

HF-Kabel Quad			
Material	Wellrohr: Polyamid (PA6)		
	Kabelmantel: Polyethylen (PE)		
Schutzart	IP 66		
Temperatur	Bei Betrieb: -30 +70 °C Bei Installation: -20 +70 °C		

Kabellänge [m]	IdentNr.
2	43431
4	43432
6	43433
8	43434
10	43435

HF-Kabel Quad, hygienisch			
Material	Wellrohr: Polyamid (PA6)		
	Kabelmantel: Polyethylen (PE)		
Schutzart	IP 66		
Temperatur	Bei Betrieb: -30 +70 °C Bei Installation: -20 +70 °C		

Kabellänge [m]	IdentNr.
2	67048
4	67049

HF-Kabel (Festmantelkabel)			
Material	Kabelmantel: Polyethylen (PE)		
Schutzart	IP 68 im aufgeschraubten Zustand		
Temperatur	Bei Betrieb: -40 +85 °C Bei Installation: -40 +85 °C		
Dämpfungsbe- lag	ca. 0.3 dB/m		



Kabellänge [m]	IdentNr.
2.0	11476
2.5	11477
3.0	11478
3.5	11479
4.0	11480

#### Format der Datenausgabe RS232 und RS485 6.4

### Titelzeile

 $Date \cdot Time \rightarrow State \rightarrow Status \rightarrow Synchronizer \rightarrow Product \rightarrow Att \rightarrow Phi \rightarrow R2 \rightarrow Tint \rightarrow IN1 \rightarrow IN2 \rightarrow PT100 \rightarrow PT100$  $C \rightarrow Cm \rightarrow C2 \rightarrow C2m \rightarrow MF1 \rightarrow MF2$ 

¶

### Folgezeilen

 $01.01.2005 \cdot 00: 00: 00 \rightarrow 0 \rightarrow 0 \rightarrow 0 \rightarrow 1 \rightarrow 0.43 \rightarrow 5.30 \rightarrow 0.07 \rightarrow 0.0 \rightarrow 0$ 

1 2 3 4 5 6 7 8 9 10 11 12

 $75.36 {\rightarrow} 75.00 {\rightarrow} 0.00 {\rightarrow} 0.00 {\rightarrow} 0.000 {\rightarrow} 0.000 \P$ 

13 14 15 16 17 18

Spalte Nr.	Beschreibung	Format
1	Datum und Uhrzeit	TT.MM.JJ·HH:MM:SS
2	State	4 –Stellig, HEX
3	Status: Information über die Qualität der letzten Messung	0 : Messung OK < 0 : Fehler
4	Produktsynchronisation	5: nicht aktiv 1: noch unsynchron 0: alle Werte synchron -1: Fehler -2: Zeit zu klein für syn. -3: Geschw. außer Ber.
5	Produktnummer	X (1 bis 4)
6	Dämpfung [dB]	X.XX
7	Phase [°/GHz]	X.XX
8	Streuung der Phasenregression	X.XX
9	Gerätetemperatur [Temperatureinheit]	X.X
10	Stromeingang 1 [Einheit des Stromeingangs]	X.X
11	Stromeingang 2 [Einheit des Stromeingangs]	X.X
12	PT100 Temperatur [Temperatureinheit]	X.X
	[] mit Auswahl der Einheit g/cm <sup>3</sup>	
13	Konzentration 1 live	X.XX [X.XXXX]
14	Konzentration 1 gemittelt	X.XX [X.XXXX]
15	Konzentration 2 live	X.XX [X.XXXX]
16	Konzentration 2 gemittelt	X.XX [X.XXXX]
17	Massenstrom für Konzentration 1	X.XXX
18	Massenstrom für Konzentration 2	X.XXX

### Sonderzeichen

" $\rightarrow$ " Tabulation

"¶" Carriage return + Line feed "·" Leerzeichen
### Kapitel 7. Zertifikate

#### 7.1 EG-Konformitätserklärung



BERTHOLD TECHNOLOGIES GmbH & Co.KG

Calmbacher Sm 22 25522 Bad westbad, Grenwoy

Planter (54) 2021 (277-0) Las (54) 2091 (277-196) Selegi (54) (Seld) (Charao) Www.Bart (Self) (Sharao) Www.Bart (Self) (Sharao)

### EG – Konformitätserklärung

Hiermit erklären wir, daß die Bauart des(r) nachfolgend bezeichneten Gerätes / Systems / Anlage in der von uns in den Verkehr gebrachten Ausführung den unten genannten einschlägigen EG-Richtlinien entspricht.

Durch nicht mit uns abgestimmte Änderungen oder nicht bestimmungsgemäßen Gebrauch verliert diese Erklärung ihre Güftigkeit.

Bezeichnung: Konzentrations-Messsystem Micro-Polar 1

Тур: LB 566-XX

Einschlägige EG-Richtlinien:

89/336/EWG (Elektromagnetische Verträglichkeit) geändert. 91/263/EWG, 92/31/EWG, 93/68/EWG, 93/97/EWG

73/23/EWG (Niederspannungsnchtlinie) geändert: 93/68/EWG

Zur Beurteilung des Erzeugnisses wurden folgende Normen herangezogen:

EN 55011.1998 + A1:1999 + A2:2002 EN 61010-1:2002-08 EN 61006-6-2:2001 EN 61000-4-2:1995 + A1:1998 + A2:2001 EN 61000-4-3:2002 + A1:2002 EN 61000-4-4:1995 + A1:2001 EN 61000-4-6:1996 + A1:2001 EN 61000-4-11:1994 + A1:2001

Diese Erklärung wird verantwortlich für den Hersteller:

BERTHOLD TECHNOLOGIES GmbH & Co. KG Calmbacher Strasse 22 D-75323 Bad Wildbad

abgegeben durch

Speld Dr. J. Bnggmann Entwicklungsleker Prozessmesslechnik

Prozessmessiechnik

Bad Wildbad, den 04.05.2004

Providence of the antibal sector of the sector of the sector of the Personal in terms of the sector of the first of the basic sector of the the sector of the sector of

H. Tetzlaff Pachiorati CE Zerilitzarung

(1) We (Constant) (M) (CONSTANT), and (MACONSTANT) of the formation of the Constant of the Million of the Constant of the Macon Constant of the Constant of the Constant of the Decomposition of the Constant of the Constant of the Decomposition of the Constant of the C

Deg Ber Back – 2018 v forward – 3 en 28 v v v j (28 v v v for 12 mer 2 for each e by Swill war out stress for park to the volument of the stress of the stre





BERTHOLD TECHNOLOGIES GmbH & Co.KG

Calmbacher Str. 22 75323 Bad Wildbad, Sermany

Phoise +49 7081 (77-0 Fax +49 7081 177-100 info@Berthold.com www.Berthold.com

### EG-Konformitätserklärung

Hiermit erklären wir, dass die Bauart des(r) nachfolgend bezeichneten Geräte / Systems / Anlage In der von uns in den Verkehr gebrachten Ausführung den unten genannten einschlägigen EG-Richtlinien entspricht.

Durch nicht mit uns abgestimmte Änderungen oder nicht bestimmungsgemäßen Gebrauch verliert diese Erklärung ihre Gültigkeit.

Produktbezeichnung:

Konzentrations-Messsysteme Micro-Polar ++

Typenbezeichnung: LB 566-XX

	Richtlinie	und Änderungen	angewendete Normen		
EMV	2004/108/EG		EN 55011	1998 +A1:1999 +A2:2002	
			EN 61326-1 EN 61000-4-2	2006-05	
				+A1:1998 +A2:2001	
			EN 61000-4-3	2006-12	
			EN 61000-4-4	2004	
			EN 61000-4-5	1995 +A1:2001	
			EN 61000-4-6	1996 +A1:2001	
			EN 61000-4-11	1994-08 +A1:2001-02	
			Namur NE21	2004	
Niederspg.	73/23/EWG	93/68/EWG	EN 61010 Teil 1	2002-08	
R&TTE	1999/5/EG		ETSI EN 300 440-1 ETSI EN 300 440-2	2007-08 2007-08	

Diese Erklärung wird verantwortlich für den Hersteller

BERTHOLD TECHNOLOGIES GmbH & Co. KG Calmbacher Str. 22, D-75323 Bad Wildbad

aktyalisiert durch/ ١.

Dr. Wilfried Reuter Technischer Geschäftsführer Bad Wildbad, den 06.06.2008

Registergenisht / Court of Pegistration Persönlich halter de Gelet schaftenn / Elity Jobie Associates Resistergevisint / Court of Registration Gaschsterkerveg / Waragemeet Beinste vorsikzender / Chasman of the Board Ostria-Ne / WAT Reg. Ne Devisite Stouernammer / German Tax Ne WEE-Reg. No.

Stuttgert HkA-330001 BERTHOLD TECHNOLIXSJES Verwaltungs-GmtH Stuttgert HRB-331520 Anno 3, Diterhotev, Dr. Wiltingd Senter Dr. Tritz Berthold DE91302051: 49038/08038 DE99468500 DE99468500 D299468500

Dresdner Pank 751.03 Pforzheim Konto/Account No. 6511120 (SLZ 666 600 19) //SWIFT-BIC DREG DREF 660 IBAN: DE05 6668 0013 0651 1120 33 Svarkasse PF-CW 75323 Bad Wildbad Konto/Account No. 80 450 03 (BLZ 666 500 S5) //SWIFT-BIC PZHSDEE6 IBAN: DE37 6665 0005 00.36 3450 03

	Date of issue: 1 August 2023 Valid unvil: 31 December 2024
C	EHEDG hereby declares that the product
	microwave sensor FlowCell, type LB5660 with planar and immersion cap windows made of PEEK and EPDM O-ring for ball housings diameter 68 mm
	from
	Berthold Technologies GmbH & Co.KG, Calmbacher Straße 22, 75323 Bad Wildbad, Germany
	has/have been evaluated for compliance and meets/meet the current criteria for Hygienic Equipment Design of the EHEDG
	Certificate No. EHEDG-C2300049
	Signed President EHEDG
	Signed EHEDG Certification Officer
	EHEDG Karspeldreef 8 1101 CJ Amsterdam Netherlands ©EHEDG

### 7.2 Frequenzzulassung

### **CETECOM ICT Services GmbH**

EC Identification number 0682

authorized by the German Government



to act as Notified Body in accordance with the R&TTE Directive 1999/S/EC of 09. March 1999.

#### CERTIFICATE EXPERT OPINION

Registration-No.:	E814059R-EO
Certificate Holder:	Berthold Technologies GmbH & Co KG Calmbacher Strasse 22
	D-75323 Bad Wildbad
Product Designation:	LB 465-xx, LB 466-xx, LB 565-xx, LB 566-xx
Product Description:	Short Range Devices
Product Manufacturer:	Berthold Technologies GmbH & Co KG Calmbacher Strasse 22
	D-75323 Bad Wildbad

Essential requirements	Specifications / Standards	Submitted documents	Result
EMC (R&TTE, Article 3.1b)	EN 55011:1998+A1:1999 (class A) EN 61000-6-2:2001	Test Report	conform
Radio spectrum (R&TTE, Article 3.2)	EN 300 440-1 V1.3.1 (2001-09) EN 300 440-2 V1.1.1 (2001-09)	Test Report	Conform

Marking: The product shall be signed with CE and our notified body number as shown right hand.

# <€ 0682

The scope of this evaluation relates to the submitted documents oply 7 The certificate is only valid in con unction with the following mimber of annexes Number of annexes: 1

Number of annexes: Saarbrücken, 24.06.2004

Place, Date of Issue

. • Dí annexes.

Signed by Ernst flussinge Nonfied Body



CETECOM ICT Services GnibH, Unterlärkheimer Straße 6-10, D-66117 Saupbrucken, Germany http://www.cetecom.de

55

### **CETECOM ICT Services GmbH**



### **CERTIFICATE OF CONFORMITY**

Registration-No :	E814059R-CC	Number of annexes:
Certificate Holder:	Berthold Technologies GmbH & Co KG Calmbacher Strasse 22	
	D-75323 Bad Wildbad	
Product Designation:	LB 465-xx, LB 466-xx, LB 565-xx, LB 566	- <i>ž</i> z
Product Description:	Short Range Devices (humidity sensor)	
Product Manufacturer:	Berthold Technologies GmbH & Co KG Calmbacher Strasse 22	
	D-75323 Bad Wildbad	

Specifications and test reports:			
Specification	Test report no. & date	Name of test laboratory	Notes
EN 55011:1998+A1(1999 (class A)	2003-731-1182-REN dated Sept. 9, 2003	ELMAC GmbH	conform
EN 61000-6-2:2001			
EN 300 440-1 V1.3.1 (2001-09)	2-3389-01-01/03 dated May 14, 2004	CETECOM ICT	conform
EN 300 440-2 V1.1.1 (2001-09)			

Statement

This equipment fulfils the requirements or parts thereof in the above mentionend specifications.

CETECOM ICT Services is authorized to act as Notified Body in accordance with the R&TTE Directive 1999/5/EC of 09. March 1999

Signed by Ernst Hussinger



Saarbrücken, 24.06.2004 -Place, Date of Issue

Notified Body

CETECOM ICT Services GmbH, Untertürkheimer Stieße 6-10, D-66117 Saarbrücken, Germany

### **CETECOM ICT Services GmbH**



### **CERTIFICATE OF CONFORMITY**

Registration-No.:	E815580V-CC	Number of annexes:
Certificate Holder:	BERTHOLD TECHNOLOGIE: Calmbacher Strasse 22 75323 Bad Wildbad Germany	S GmbH & Co KG
Product Designation:	Micro-Polar ++ LB 566-1x Micro-Polar Brix ++ LB 565-1x	
Product Description:	Short Range Devices (humidity sensor)	
Product Manufacturer:	BERTHOLD TECHNOLOGIES Calmbacher Strasse 22 75323 Bad Wildbad Germany	S GmbH & Co KG

Specifications and test reports:				
Specification	Test report no. & date	Name of test laboratory	Notes	
EN 300 440-1 V1.4.1 (2007-08)	2-4837-01-02/07	CETECOM ICT	conform	
EN 300 440-2 V1.2.1 (2007-08)	dated February 07, 2008			

Statement

Saarbrücken, 2008-12-05 Place, Date of Issue

This equipment fulfils the requirements or parts thereof in the above mentionend specifications.

CETECOM ICT Services is authorized to act as Notified Body in accordance with the R&TTE Directive 1999/5/EC of 09, March 1999

ñ d Signed by Ernst Hussinger CHE BENAMN Notified Body

CETECOM 1CT Services GmbH, Untertürkheimer Straße 6-10, D-66117 Saarbrücken, Germany



#### **CETECOM ICT Services GmbH**

**EC Identification Number 0682** 

authorized by the German Government

Bundesnetzagentur

BNetzA-bS-02/51-52

to act as Notified Body in accordance with the R&TTE Directive 1999/5/EC of 9. March 1999.

#### **EXPERT OPINION**

Registration-No.:	E815580V-EO
Certificate Holder:	BERTHOLD TECHNOLOGIES GmbH & Co KG Calmbacher Strasse 22 75323 Bad Wildbad Germany
Product Designation:	Micro-Polar ++ LB 566-1x Micro-Polar Brix ++ LB 565-1x
Product Description:	Short Range Devices
Product Manufacturer:	BERTHOLD TECHNOLOGIES GmbH & Co KG Calmbacher Strasse 22 75323 Bad Wildbad Germany

Essential requirements	Specifications / Standards	Submitted documents	Result
Radio spectrum	EN 300 440-1 V1.4.1 (2007-08)	Test Report	conform
(R&TTE, Article 3.2)	EN 300 440-2 V1.2.1 (2007-08)		

Marking:	The product shall be sig number as shown right	ned with CE and our notified body hand.	(6	0682
The scope of the The certificate Number of a Saarbrücken, Place, Date of	his evaluation relates to the sub is only valid in conjunction w annexes: <u>1</u> 2008-12-05 Issue	mitted documents only ith the following number of annexes.		Struke BERN NOTIE

CETECOM ICT Services GmbH, Untertürkheimer Straße 6-10, D-66117 Saarbrücken, Germany http://www.cetecom-ict.de

cal Acceptance Certificate German www.ctcadvanced.com CTC advanced GmbH (formerly CETECOM ICT Services GmbH) 66117 **Jntertuerkheimer Str.** 

СТС	advanced member of RWTÜV group		
Certificate Holder:	BERTHOLD TECHNOLOGIES GMBH & CO. KG CALMBACHER STR. 22 75323 BAD WILDBAD Germany	<b>(†</b> )	Bundesnetzagentur
	Germany		BNetzA-CAB-03/22-51
ISED Certification Number:	4777A-IC01X01		authorized by the German Government to act as CAB
CTC Registration Number:	1865		in accordance with the MRA EU Canada of 1st November 1998.
OATS Facility ID Number:	3462C-1		
OATS Facility:	CTC advanced GmbH		
	66117 Saarbruecken		
	Germany		

Phone: +49 681 598-0

Fax:

+49 681 598-8775

Email: info@ctcadvanced.com

**Product Description:** 

Saarbrücken

Concentration, Dry Mass, Water, Density measuring system

Certification of equipment means only that the equipment has met the requirements of the above-noted specification. Licence applications, where applicable to use certified equipment, are acted on accordingly by the ISED issuing office and will depend on the existing radio environment, service and location of operation. This certificate is issued on condition that the holder complies and will continue to comply with the requirements and procedures issued by ISED. The equipment for which this certificate is issued shall not be manufactured, imported, distributed, leased, offered for sale or sold unless the equipment complies with the applicable technical specifications and procedures issued by ISED.

La certification du matériel signifie seulement que le matériel a satisfait aux exigences de la norme indiquée ci dessus. Les demandes de licences nécessaires pour l'utilisation du matériel certifié sont traitées en conséquence par le bureau de délivrance d'ISDE et dépendent des conditions radio ambiantes, du service et de l'emplacement d'exploitation. Le présent certificat est délivré à la condition que le titulaire satisfasse et continue de satisfaire aux exigences et aux procédures d'ISDE. Le matériel à l'égard duquel le présent certificat est délivré ne doit pas être fabrique, importé, distribué, loué, mis en vente ou vendu à moins d'être conforme aux procédures et aux spécifications techniques applicables publiées par ISDE.

I hereby attest that the subject equipment was tested and found in compliance with the above-noted specification. J'atteste par la présente que le matériel a fait l'objet d'essai et jugé conforme à la spécification ci-dessus.

Stich CERTIFICATION BOD + GABID NO DEOOD CTC advanced GmbH cn=Gerald Schmidt, o=CTC advanced GmbH, ou=SCT-141216, email=gerald.schmidt@ctcadvan ced.com, c=DE 2017.09.08 10:20:11 +02'00' Certification Body







# тсв

GRANT OF EQUIPMENT AUTHORIZATION

Certification Issued Under the Authority of the Federal Communications Commission

By:

CTC advanced GmbH (former CETECOM ICT Services Gmb Untertuerkheimer Strasse 6-10 66117 Saarbruecken, Germany

Date of Grant: 09/08/2017

тсв

Application Dated: 09/08/2017

Berthold Technologies Calmbacher Str. 22 75323 Bad Wildbad Germany Bad Wildbad, 75323 Germany

Attention: Dirk Moermann , Dr.

#### NOT TRANSFERABLE

EQUIPMENT AUTHORIZATION is hereby issued to the named GRANTEE, and is VALID ONLY for the equipment identified hereon for use under the Commission's Rules and Regulations listed below.



## TCB

GRANT OF EQUIPMENT AUTHORIZATION Certification Issued Under the Authority of the Federal Communications Commission

By:

CTC advanced GmbH (former CETECOM ICT Services Gmb Untertuerkheimer Strasse 6-10 66117 Saarbruecken, Germany

Date of Grant: 09/08/2017 Application Dated: 09/08/2017

TCB

Berthold Technologies Calmbacher Str. 22 75323 Bad Wildbad Germany Bad Wildbad, 75323 Germany

Attention: Dirk Moermann , Dr.

#### NOT TRANSFERABLE

EQUIPMENT AUTHORIZATION is hereby issued to the named GRANTEE, and is VALID ONLY for the equipment identified hereon for use under the Commission's Rules and Regulations listed below.



## Kapitel 8. Technische Zeichnungen



### 8.1 Maßbilder der AWE-Wandgehäuse





BERTHOLD

### 8.2 Elektrischer Anschlussplan





### 8.3 Maßbilder FlowCell

#### 8.3.1 Typ LB 5660-102-00x, FlowCell DN 50 Flansch, FDA







#### 8.3.2 Typ LB 5660-202-00x, FlowCell DN 65 Flansch, FDA



#### 8.3.3 Typ LB 5660-302-00x, FlowCell DN 80 Flansch, FDA



#### 8.3.4 Typ LB 5660-402-00x, FlowCell DN 100 Flansch, FDA





#### 8.3.5 Typ LB 5660-502-00x, FlowCell DN 125 Flansch, FDA



#### 8.3.6 Typ LB 5660-602-00x, FlowCell DN 150 Flansch, FDA

# 5 5 LB 5660-112-00X FlowCell DN 50 G-BS/M 3.10 С 66744 M13 Va fielab GLOHTABE R co. 228 ۲ 囙 (0. 164 O Kanten-2 Form- u oleranz-Z768-m ĺ۵ nkelmasse,F en ohne To DIN ISD 2 Б Dateiname Wodell: 66744-013-WI3 Dateiname Zeichnung: 66744-013-WI3 nenearad nattiab then traiqosi sabwa prummituta edulitinta endo trab drunnta sa so Diserettin di antervesti printa antervetti di antervetti antervetti antervetti antervetti antervetti antervetti

#### 8.3.7 Typ LB 5660-112-00x, FlowCell DN 50 G-BS/M





#### 8.3.8 Typ LB 5660-212-00x, FlowCell DN 65 G-BS/M



#### 8.3.9 Typ LB 5660-312-00x, FlowCell DN 80 G-BS/M



#### 8.3.10 Typ LB 5660-412-00x, FlowCell DN 100 G-BS/M



#### 8.3.11 Typ LB 5660-512-00x, FlowCell DN 125 G-BS/M



#### 8.3.12 Typ LB 5660-612-00x, FlowCell DN 150 G-BS/M



#### 8.3.13 Typ LB 5660-132-00X DN 50









#### 8.3.15 Typ LB 5660-332-00X DN 80



#### 8.3.16 Typ LB 5660-432-00X DN 100





#### 8.3.17 Typ LB 5660-532-00X DN 125









### 8.4 Maßbilder Behältersonden

#### 8.4.1 Typ LB 5650-01



#### 8.4.2 Typ LB 5650-02



#### 8.4.3 Typ LB 5650-03


# 8.4.4 Typ LB 5650-04



# 8.4.5 Typ LB 5650-05



### 8.4.6 Typ LB 5650-09





# 8.4.7 Einbausituation in Rohrleitungen

# 8.5 Maßbilder Spülsonden

# 8.5.1 Typ LB 5651-01





# 8.5.2 Typ LB 5651-02



# 8.5.3 Typ LB 5651-03





# 8.5.4 Typ LB 5651-04



# 8.5.5 Typ LB 5651-05





# 8.5.6 Einbausituation in Rohrleitungen

BERTHOLD

# 8.6 Montageblätter für LB 5650 (Behältersonde)





# 8.7 Montageblätter für LB 5651 (Spülsonde)





BERTHOLD

# Indexverzeichnis

### A

Abstand zu Metall · 43 Adapterflansch · 44 Anschlussflansch · 43 Anschlussplan · 87 ASA-Flansch-Adapter · 66 Auswerteeinheit · 23 AWE · 15

### B

Batterie · 57 Behältersonde · 28

### С

CE Zeichen · 13 Clamp-Stutzen · 26

### D

Datenformat RS232 · 71 Datenübertragungsrate · 50 Digitale Ausgänge · 50 Distanzschiene · 40 Durchsatz-Berechnung · 20

### E

EG-Konformitätserklärung · 73 Einbausituation in Rohrleitungen · 112, 118 Einbautiefe · 43 Einschränkungen · 18

### F

Flachdichtung · 43 Format RS232 · 71 Frequenzzulassung · 13, 76

### G

Gaseinschlüsse · 14, 38 Genauigkeit · 60 Geräteentsorgung · 53 Gerätereinigung · 56

### Η

HF-Kabel · 32 HF-Kabel Quad · 32 Hochfrequenzkabel · 32

### Ι

Inkrustation · 30 Installation AWE, Behältersonde · 45 Installation AWE, Messzelle · 40 Installation der Behältersonden · 43 Installation der Messzelle · 37

### K

Kabelverlegung · 41, 45 Kompensation · 19 Komponenten · 21

### L

Lebensmitteltauglich · 28 LED's · 24, 50 Leistungsabstrahlung · 14 Leitfähige Materialien · 19

### М

Messanordnung am Behälter · 35 Messanordnung am Rohr · 34 Messprinzip · 17 Messwertberechnung · 18 Messzelle · 25 Montageblätter · 119

### N

Netzanschluss · 47 Netzsicherungen · 58

### R

Recyclingpass · 53 Relais · 50 Rohrleitungsdruck · 38 RS232 Schnittstelle · 50

### S

Salzgehalt · 14 Sendeleistung · 60 Sicherheitshinweise · 7 Sicherungen · 58 Spülparameter · 44 Spülsonde · 15, 30 Steckerleiste · 48 Steigleitung · 37 Symbole · 7

### Т

Technische Daten · 60



Technische Daten HF-Kabel · 69 Technische Daten Sensoren · 63 Technische Zeichnungen · 85 Temperaturkompensation · 19 Transport · 36

### U

Übersicht Behältersonden · 68 Übersicht Dichtungsscheiben · 68

### V

Verschleißteile · 54

### W

Warnhinweise · 7 Wartung · 53 Werkseinstellung · 15

### Ζ

Zertifikate · 73

# Notizen



		 	 	 		 	 	 		 			 	 	 	 _			
<u> </u>		 	 	 		 	 	 		 		 	 	 	 	 			
		 				 		 	_		_	 _	 _						
<u> </u>		 	 	 		 	 	 		 		 	 	 	 	 _		_	
<u> </u>	 	 	 	 		 	 	 		 		 	 	 	 	 			
<u> </u>		 <u> </u>			-			 		 		 	 						
													 				_		



# Notizen








# Konzentrations-Messsysteme MicroPolar (++) LB 566

Betriebsanleitung Softwareanleitung 41986BA1

Rev. Nr.: 07, 03/2018



Die gelieferten Geräte dürfen nur vom Service der Firma Berthold oder durch von der Firma Berthold autorisierte Techniker instand gehalten werden!

Im Störungsfall wenden Sie sich bitte an unseren zentralen Kundendienst (Kontaktadresse siehe unten). The units supplied should not be repaired by anyone other than Berthold Service engineers or technicians by Berthold.

In case of operation trouble, please address to our central service department (address see below).

Die komplette Betriebsanleitung besteht aus der Geräte- und Softwareanleitung.

### Die Geräteanleitung beinhaltet die

- Komponentenbeschreibung
- Montageanleitung
- > Elektrische Installationsbeschreibung
- Technische Daten
- > Zertifikate
- Maßzeichnungen

### Die Softwareanleitung beinhaltet die Beschreibung der

- Bedienung
- Softwarefunktionen
- > Kalibrierung
- > Fehlermeldungen

### Der vorliegende Teil beinhaltet die Softwareanleitung.

Änderungen vorbehalten.

 BERTHOLD TECHNOLOGIES GmbH & Co. KG

 Calmbacher Str. 22 · D-75323 Bad Wildbad

 Zentrale:
 Service:

 Tel. +49 7081 177 0
 Tel. +49 7081 177 111

 Fax +49 7081 177 100
 Fax +79 7081 177 339

 industry@Berthold.com
 Service@Berthold.com

 www.Berthold.com
 Service@Berthold.com

# **Inhaltsverzeichnis**

		Seite
Kapitel 1. Sic	cherheitshinweise	7
Kapitel 2. Ko	mmunikation mit MicroPolar	9
Kapitel 3. Lei	itfaden zur Inbetriebnahme	11
Kapitel 4. So	ftwarefunktionen	13
	4.1 Hinweise zur Menü-Struktur	13
	4 2 Menüstruktur	14
	4.2.1 Startmenü	16
	4.2.2 Diagnose	16
	4.2.3 Einstellungen	18
	4.2.4 Zugriffsebene	19
	4.2.5 Sprache	19
	4.2.6 Konfiguration	20
	4.2.7 Allgemeines	21
	4.2.8 Messung	21
	4.2.9 Plausibilität	22
	4.2.10 Phasenmessung	23
	4.2.11 Offline Modus	23
	4.2.12 Massenstrom	24
	4.2.13 Kalibrierung	25
	4.2.14 Systemabgleich	25
	4.2.15 Kalibriere Konzentration	26
	4.2.16 Probentabelle	27
	4.2.17 Erweiterte Probedaten	28
	4.2.18 Extras	28
	4.2.19 Kalibrierung	29
	4.2.20 EIN- / Ausgange	30
	4.2.21 Stromausgang	31
	4.2.22 Stromausgang 1	31
	4.2.25 Stromaingang	23 22
	4.2.24 Stromoingang 1	22 23
	4.2.25 Stromeingang 2	33
	4 2 27 PT100	34
	4.2.27 Digitalausgang	34
	4.2.29 Digitaleingang	35
	4.2.30 Service	37
	4.3 Trendanzeige	39
Kapitel 5. Ko	nfiguration	41
I	5.1 Konfigurationseinstellungen	41
	5.1.1 Allgemeines	41
	5.1.2 Messung	42
	5.1.3 Plausibilität	42
	5.1.4 Mikrowelle	43
	5.1.5 Einheit	43
	5.1.6 Marker	44
	5.2 Start-Kalibrierkoeffizienten	45
	5.3 Plausibilitäten konfigurieren	46

# BERTHOLD ...

Kapitel 6. Kalibrierung	47
6.1 Systemabgleich	47
6.1.1 Überprüfung der Referenz	werte 49
6.2 Probenanme 6.2 1 Eingabe der Laborwerte	49
6.3 Kalibrierung	52
Kapitel 7. Inbetriebnahme MicroPolar +	+ 53
Kapitel 8. Kalibrierung und Extras	55
8.1 Plausibilität Verhältnis Phi/D 8.1.1 Verhältnis Phi/D 8.1.2 Prozessaufnahme	konfigurieren 55 55 56
8.2 Anpassung der Kalibrierung	58 Destabulla
8.3 Ausgabe des Indetriednanme 8.4 Kalibrierung	Protokolis 60
8.4.1 Kalibrierung mit zwei Kon	zentrationen 64
8.4.2 Kalibrierung mit Splitwert	68
8.5 Typische Kalibrierkoeffiziente	n/Startwerte 70
Kapitel 9. Passwort	71
9.1 Passwort vergessen	71
Kapitel 10. Fehlerlisten und Gerätezustä	inde 73
10.1 Stromausfall	73
10.2 Temperaturfehler	73
10.3 Hardwarefehler	74
10.4 Sensorrenier	/5
10.6 Fehler Messbereich	75
10.7 Fehler Zusatzmessung	77
10.8 Fehler Bereich Analogausga	ng 77
10.9 Watchdog-Fehler	77
10.10 Systemfehler	78
10.11 Dichterenier 10.12 Eingabefehler	/8 78
10.13 Gerätezustände	79
Kapitel 11. Inbetriebnahme-Protokoll	81
11.1 Beispiel Inbetriebnahme-Pro	otokoll 84
11.2 Probentabelle	89



# Kapitel 1. Sicherheitshinweise

Berücksichtigen Sie alle Sicherheitshinweise im *Hardwareteil*, insbesondere die unter *Kapitel 1. Sicherheitshinweise*.

HINWEIS

### Parametereinstellungen

Es dürfen keine Änderungen an den Parametereinstellungen vorgenommen werden ohne genaue Kenntnis dieser Betriebsanleitung mit allen Hinweisen und genauer Kenntnis über das Verhalten eines angeschlossenen Reglers und die möglichen Einflüsse auf den zu steuernden Betriebsprozess.

# Kapitel 2. Kommunikation mit MicroPolar

Die Kommunikation mit MicroPolar und MicroPolar ++ erfolgt über 4 Softkeytasten. Entsprechend der Position im Menü ändert sich die Funktion der einzelnen Tasten. Werte und Texte werden über eine alphanumerische Tastatur eingegeben. Der Gerätezustand wird mittels 5 LED's angezeigt.



Nutzen Sie die Hilfetaste ? in der Displayfußzeile, sie liefert nützliche Informationen.



# Kapitel 3. Leitfaden zur Inbetriebnahme

Zur Inbetriebnahme führen Sie der Reihe nach die unten aufgeführten Schritte durch.

Bei Inbetriebnahme der High Dynamik Version MicroPolar ++ beachten Sie zunächst das *Kapitel 8. Inbetriebnahme Micro-Polar* ++.

Das *Kapitel 4. Softwarefunktionen* beschreibt alle Softwarefunktionen und dient unter anderem als Nachschlagewerk.

### 1. Schritt

Konfigurieren Sie je nach Erfordernis die analogen Eingänge: Stromeingänge 1, 2 und den PT100. Siehe dazu *Kapitel 4.2.22 Ein- / Ausgänge*.



Alle analogen Ein- und Ausgänge sind bereits ab Werk abgeglichen. Daher sind bei der Inbetriebnahme keine Abgleichsarbeiten erforderlich.

### 2. Schritt

Überprüfen und passen Sie die Software-Parameter der Applikation an. Einige Parameter sind bereits ab Werk voreingestellt. Führen Sie dazu das *Kapitel 5. Konfiguration* aus.

### 3. Schritt

Führen Sie die Kalibrierung mit Probenahme durch, siehe *Ka- pitel 6. Kalibrierung*.

Eine Temperaturkompensation ist nur dann erforderlich, wenn die Phase/Dämpfung von der Produkttemperatur deutlich beeinflusst wird, dieses ist abhängig vom Produkt und Wassergehalt.

### 4. Schritt

Konfigurieren Sie je nach Erfordernis die Stromausgänge, digitalen Ein- und Ausgänge.

# Kapitel 4. Softwarefunktionen

# 4.1 Hinweise zur Menü-Struktur

Die Menü-Struktur auf den nächsten Seiten gibt einen Überblick über alle Funktionen des LB 566. Anhand der darauf angegebenen **Seiten-Nummern**, können die Funktionen des dargestellten Fensters nachgeschlagen werden.

Der Wechsel von Ebene "**Nur lesen**" auf "**Benutzer Modus**" ist nur mit Passwort möglich. Die "**Service-Ebene**" ist aufgrund von Zulassungsbestimmungen verschlossen.



Die Auswerteeinheit bzw. die Software beinhaltet Funktionen, welche für das Messsystem LB 566 (++) und deren Messanwendung nicht relevant sind.

Es handelt sich dabei um folgende Funktionen:

- Beladungskompensation

- Synchronisation der Messeinrichtungen

In der Betriebsanleitung werden diese Funktionen und die zugehörigen Menüs nicht beschrieben.



# 4.2 Menüstruktur











### 4.2.1 Startmenü

1   -	LB 566	07.05 -	13:25								
Live Anzeige Diagnose Einstellungen											
Zugr Spra	iffsebene che										
RUN											

### Live Anzeige:

Anwahl schaltet zur Live Anzeige

#### **Diagnose:**

Dieser Menüpunkt verzweigt zum Datenlogger, Fehlerspeicher, zu Geräteinformationen und Inbetriebnahme-Protokoll-Ausgabe.

#### Einstellungen:

Hier werden alle notwendigen Eingaben zur Funktion des Messsystems vorgenommen.

#### Zugriffsebene:

Auswahl der Zugriffsebene. Durch Passwörter geschützte Bereiche können hier freigegeben werden.

### Sprache:

Auswahl der Dialogsprache.

### 4.2.2 Diagnose



### Datenlog:

Der Datenlog zeichnet die Daten entsprechend dem Inhalt der seriellen Datenausgabe RS232 und RS485 (siehe *Hardwareteil Kapitel 6.6*) auf.

Sämtliche Messdaten einer Messung (Sweep) werden über die Mittelungszeit (siehe unten) gemittelt und gespeichert. Diese Zeit ergibt sich aus der ausgewählten Aufzeichnungszeit. Der Inhalt des Datenlogs kann über die Live Anzeige grafisch aufgerufen werden, siehe *Kapitel 4.3 Trendanzeige*. Eine Ausgabe als Text-Datei über RS232 und RS485 sowie über das Memory Tool (optionales Zubehör) ist möglich.

• Logart deaktivieren

einzeln kontinuierlich

stopp bei Fehler

- Logzeit Aufzeichnungszeitraum
- 15 Minuten bis 3 Tage
- Log Neustart Löscht den Datenlog und startet mit obiger Einstellung
- Mittelungszeit Ergibt sich aus der Logzeit
- Log drucken Ausdruck der Tabellen, Ausgabe über RS232 und RS485, Format siehe Hardwareteil Kapitel 6.6

### Wechsel der Datenlog-Einstellungen:

Bei Wechsel der Logart von beliebig auf "einzeln" wird der Datenlog gelöscht und mit der aktuellen Einstellung neu gestartet.

Bei Wechsel aller übrigen Logarten und Logzeiten wird der Datenlog nicht gelöscht und mit den neuen Einstellungen weiter aufgezeichnet.

### Verhalten bei gestoppter Messung:

Wird während des Datenlogs die Messung eine zeitlang gestoppt, so wird bei Logart "einzeln" die Messpause als Logzeit interpretiert. Bei allen übrigen Logarten addiert sich die Messpause zur Logzeit.

### Fehlerlog:

Anzeige der aufgezeichneten Fehler. Gespeichert werden die letzen 20 Fehlermeldungen mit Datum und Uhrzeit.

#### Info:

•

- Messstelle :...
- Gerätetyp : LB 566
  - Lieferant : Berthold Technologies
- Hersteller : Berthold Technologies
- Geräte Nr. : ...
- Fertigungs Nr. : ...-...
- Software Ver. : V...
- SW Rev.datum : ...

#### Protokoll drucken:

Ausgabe des Inbetriebnahme-Protokolls über RS232 und RS485.

Format, Inhalt und Beispiel siehe Kapitel 11. Inbetriebnahme-Protokoll.

RTHOLE

# 4.2.3 Einstellungen



### **Konfiguration:**

Einstellung von

- allgemeinen Daten
- messspezifische Daten
- Plausibilitätsdaten
- Mikrowellendaten
- Einheiten
- Marker

#### **Kalibrierung:**

- Systemabgleich
- Konzentration Kalibrierung
- Extras

### Ein-/Ausgänge:

- Stromausgänge
- Stromeingänge
- PT100
- Digitalausgänge
- Digitaleingänge

### Service:

- Fabrikeinstellung
- General Reset
- Memory Tool (Bedienung des Memory Tools, optionales Zubehör)
- Datenausgabe (über RS232 und RS485, Dateninhalte können gewählt werden)

#### **Produkt:**

Auswahl des Produktes (1–4), mit Umschaltung werden die produktspezifischen Daten geladen: Aus-, Eingänge und Kalibrierung.

Beim erstmaligen Aufruf der Produkte 2 bis 4, werden alle Einstellungen und Inhalte (wie z.B. Systemabgleich, Probentabelle, Datenlog und Kalibrierung) des aktuellen Produktes in das neue Produkt kopiert.

#### Passwort ändern:

Das Passwort für die Zugriffsebenen Benutzer Modus kann hier geändert werden.

Weitere Informationen siehe Kapitel 9. Passwort.
## 4.2.4 Zugriffsebene



#### Nur lesen:

Hiermit kann das Messsystem gegen Zugriff Unberechtigter geschützt werden. Ein Verlassen dieser Ebene ist nur mit Passwort möglich. Das Messgerät kann nicht gestartet und nicht gestoppt werden. Es kann nur vom Hauptmenü zur Diagnose und Zugriffsebene gewechselt werden.

#### **Benutzer Modus:**

- Der Benutzer Modus ist der Standard Modus und ermöglicht den Zugriff auf alle benutzerrelevanten Parameter.
- Von "Nur lesen" aus, wird ein Passwort abgefragt.
- Das Passwort ist veränderbar.

#### Service:

• Diese Ebene ist nur dem Servicepersonal vorbehalten.

## 4.2.5 Sprache



#### Sprache:

Auswahl der Dialogsprache



## 4.2.6 Konfiguration



#### **Allgemeines:**

• Eingabe von Datum, Uhrzeit und Messstelle

#### Messung:

- Messmodus (batch/kontinuierlich)
- Startmodus (Tastatur/extern)
- Mittelung (Anzahl der Messwerte zur Mittelwertbildung)
- Mittelwert Reset ( ja/nein )
- Stromausgang bei Stopp

Weiteres unter Kapitel 4.2.8 Messung

#### Plausibilität:

- Die Prozessgrenzen definieren den zulässigen Bereich, in der die aktuelle Konzentration liegen muss.
- Die Phasenmessung unterliegt einer Plausibilitätsanalyse, welche hier einstellbar ist.
- Aktivierung und Schaltwert der Pausenfunktion

Weiteres unter Kapitel 4.2.9 Plausibilität

#### Mikrowelle:

Kabel (Eingabe der Kabellängen für Messzweig und Referenzzweig). Z.B. für 4 m HF-Kabel quad muss 8 m für beide Längen eingegeben werden.

#### Massenstrom:

Wird in einer Rohrleitungsapplikation die Dichte gemessen (Einheit der Konzentration =  $g/m^3$  ausgewählt), so besteht die Möglichkeit den Massenstrom (Durchsatz) über Stromausgang 1 / 2 (in Tonnen pro Stunde) auszugeben. Details zum Massenstrom siehe *Kapitel 4.2.12 Massenstrom*.

#### Einheiten:

Gemäß Konfiguration sind verschiedene Dimensionen für die Konzentrationen, Stromeingänge und Temperatur wählbar.

Für die Konzentration (1 und 2) sind wählbar: keine, spezifisch, %, %TS, °Bx, g/L, g/cm<sup>3</sup>, °Be

Für Stromeingang 1 sind wählbar: keine, spezifisch, °C, °F, g/cm<sup>3</sup>, kg, t/h

Für Stromeingang 2 sind wählbar: keine, spezifisch, °C, °F, cm, m/s

Für den PT100-Eingang sind wählbar: keine, °C, °F

#### Marker:

Markerwert und Bezeichnung bis 5 Zeichen werden hier eingegeben. Die Darstellung erfolgt in der Live Anzeige und bezieht sich auf das Bardiagramm. Um den Marker zu deaktivieren, wählen Sie einen Markerwert außerhalb der Diagrammgrenzen bzw. der Stromausgangsgrenzen.

## 4.2.7 Allgemeines



#### Datum:

• Eingabe des aktuellen Datums

#### **Uhrzeit:**

• Eingabe der aktuellen Uhrzeit

#### Messstelle:

• Der Name der Messstelle kann hier eingegeben werden. Die Messstelle (max. 8 Zeichen) wird in der Kopfzeile im Display angezeigt.

### 4.2.8 Messung



#### Messmodus:

Auswahl Kontinuierlich oder Batch. Im Batchbetrieb wird ein Mittelwert zwischen Start und Stopp gebildet. Im Kontinuierlichen Modus wird eine gleitende Mittelwertsbildung entsprechend der eingestellten Mittelungszahl durchgeführt.

#### Startmodus:

Das Messgerät kann über externe Klemmen (Digitaleingang) oder via Tastatur gestartet und gestoppt werden.

#### Mittelung:

Hier erfolgt die Eingabe der Mittelungszahl. Sie gibt an, über wie viele Messungen der Konzentrationswert gemittelt wird (gleitende Mittelung). Betrifft nur den Messmodus Kontinuierlich.

#### **Mittelwert Reset:**

Mittelwert wird zurückgesetzt (ja/nein). Bezieht sich auf Batch und Kontinuierlich.

#### Stromausgang bei Stopp:

Auswahl "0/4 mA" oder "Wert halten". Die Auswahl legt fest, wie sich die Stromausgänge bei gestoppter Messung verhalten. Betrifft nur den Messmodus Kontinuierlich.

## 4.2.9 Plausibilität



#### Prozessgrenzen:

Eingabe einer zulässigen Messbereichsüberschreitung. Läuft die Konzentration aus dem Bereich heraus, so wird der Konzentrationsmittelwert eingefroren und eine Fehlermeldung angezeigt (Fehlerzustand). Die Prozessgrenzen sind unabhängig von den Stromausgangsgrenzen.

#### **Phasenmessung:**

Die Phase wird einer Plausibilitätsanalyse unterzogen. Näheres siehe Kapitel 4.2.10 Phasenmessung.

#### **Offline Modus:**

Kann aktiviert oder deaktiviert werden. Schaltgröße ist die Dämpfung, wird die eingegebene min. Dämpfung unterschritten, so geht die AWE in den Offline Modus:

- Stromausgang fällt auf die untere Stromausgangsgrenze (0/4 mA)
- Displaymeldung
- RUN-LED blinkt

Details zur Offline-Funktion siehe Kapitel 4.2.11 Offline Modus.

## 4.2.10 Phasenmessung



#### Sigma max.:

Hier wird die max. Streuung (Sigma) aus der Regression Phase über Frequenz eingestellt. Die Streuung liegt im normalen Messbetrieb zwischen 0 bis

100.

Default: Sigma = 100.

Mit Sigma =0 wird die Plausibilität ausgeschaltet.

#### Verhältnis Phi/D:

Ein weiteres Plausibilitätskriterium ist das Verhältnis aus Phase und Dämpfung. Dabei wird ein fester Verhältniswert Phi/D eingegeben. Ist dieser nicht bekannt, so muss er einmalig aufgenommen werden, siehe nachfolgende Beschreibung "Auto-Einstellung". Die genaue Vorgehensweise wird in *Kapitel 8.1* beschrieben.

#### Auto-Einstellung:

Die automatische Verhältnisaufnahme Phi/D wird hier einund ausgeschaltet. Details zur Funktion, siehe *Kapitel 8.1.* 

### 4.2.11 Offline Modus

#### **Beschreibung:**

Der Offline Modus ist ein Softwarefeature mit der ein bestimmter Produktzustand (z.B. kein Produkt vorhanden) erkannt werden kann. Unterschreitet die gemessene Dämpfung eine definierte minimale Dämpfung, so geht die AWE in den Offline Modus. Der Gerätezustand für diesen Modus wird in *Kapitel 10.4 Gerätezustände* beschrieben.



#### Aktiviert:

Hier wird der Offline Modus ein- und ausgeschaltet.

#### Dämpfung min:

Eingabe der minimalen Dämpfung, bei Unterschreitung geht die Messung in den Offline Modus und bei Überschreitung wird er wieder verlassen.



### 4.2.12 Massenstrom

Das Menü MASSENSTROM erscheint nur unter zwei Voraussetzungen:

- 1. Für die Konzentration ist die Dichteeinheit g/cm^3 ausgewählt.
- 2. Stromeingang 2 ist aktiviert und die Einheit für Stromeingang 2 ist auf m/s angewählt.

Für diesen Fall kann aus dem Dichtemesswert mit Angabe der Rohrquerschnittsfläche oder Rohrinnendurchmesser der Massenstrom (Durchsatz) in Tonnen pro Stunde angezeigt und über Stromausgang ausgegeben werden.

1   -   Massenstrom   07.05–13:25			
Massens Querschi	tr.berec	hn.	aktiviert 0.79cm^2
Rohrinne	ndurchr	nesser	1.0cm

#### Massenstromberechnung:

Hier wird die Berechnung aktiviert oder deaktiviert.

#### Querschnittsfläche:

Hier wird die Querschnittsfläche der Rohrleitung eingegeben.

#### **Rohrinnendurchmesser:**

Hier wird der Rohrinnendurchmesser eingegeben.

Bemerkung: Es ist nur eine Eingabe erforderlich, entweder die Querschnittsfläche oder der Rohrinnendurchmesser, die andere wird automatisch berechnet.

## **I** WICHTIG

Die Massenstromberechnung erfolgt nur mit variierender Produktgeschwindigkeit, welche über Stromeingang 2 eingespeist werden muss.

## 4.2.13 Kalibrierung



#### Systemabgleich:

Der Systemabgleich wird hier gestartet. Details siehe *Kapitel 4.2.14 Systemabgleich*.

#### **Kalibriere Konz:**

Verzweigt zum Kalibriermenü von Konzentration 1

#### Kalibriere Konz 2:

Verzweigt zum Kalibriermenü von Konzentration 2. Die zweite Konzentration wird nur dann angezeigt, wenn unter Menü | EXTRAS | PROZESSART | eine zweite Konzentration angewählt ist.

#### Extras:

Hier werden die Tara Werte, Anzahl der Sweeps bei Probeaufnahme, die Prozessart und der Splitwert eingestellt. Details siehe *Kapitel 4.2.18 Extras*.

### 4.2.14 Systemabgleich



#### Abgleich:

Hier wird der Systemabgleich gestartet. Mit dem Systemabgleich werden Phase und Dämpfung auf null gesetzt, wodurch z.B. alle Kabelparameter berücksichtigt werden. Dieser Abgleich bildet zugleich die Referenz für die Messung.

Der Systemabgleich (= Referenzmessung) muss einmalig durchgeführt werden.

#### **Referenzwerte:**

Nach Abschluss der Referenzmessung können die Referenzwerte für Phase, Dämpfung, Steigung und Sigma ausgelesen werden.

#### **Chart Phi:**

Zeigt die Phase über der Frequenz an.

#### **Chart Dämpfung:**

Zeigt die Dämpfung über der Frequenz an.

Durch einen Systemabgleich wird der Datenlog (siehe *Kapitel* 4.2.2 Diagnose) nicht gelöscht.



## 4.2.15 Kalibriere Konzentration



#### Probentabelle:

Anzeige aller aufgenommenen Proben und Eingabe der Laborwerte.

#### **Kalibrierung:**

Hier erfolgt die:

- die Auswahl der Kalibrierparameter, Temperatur- und Beladungskompensation
- automatische Berechnung der Kalibrierkoeffizienten
- Anzeige der Kalibrierkoeffizienten

Details siehe Kapitel 4.2.19 Kalibrierung.

#### Anpassung:

Eine nachträgliche Korrektur der Anzeige ist durch Eingabe eines Faktors und Offsets möglich. Die Berechnung erfolgt nach folgender Formel:

#### Gl. 4-1:

Korrigiert e Anzeige = Anzeige · Faktor + Offset

#### **Ergebnis:**

Darstellung der Kalibrierkennlinie, Anzeige der Korrelation und Anzeige der Koeffizienten.

#### 4.2.16 **Probentabelle**

1   1/1   Probe Nr.	1   07.05 – 13:25
Weiter	
Aktiv	Ja
Messwert	65.50 %
Laborwert	0.00 %
Erweitert	
Löschen	▲▼√

In der Kopfzeile wird folgendes beschrieben (von links): •

- Produkt-Nr.
- Aktuelle Tabellenposition / Gesamtzahl der Einträge
- Probe-Nr. der aktuellen Tabellenposition
- Datum und Uhrzeit der Probenahme

Bis zu 30 Probeeinträge sind möglich. Die Probe kann durch die Probe-Nr. oder durch Datum/Uhrzeit dem Laborwert zugeordnet werden. Die Probe-Nr. wird fortlaufend vergeben. Wird eine Probe gelöscht, so wird die Probe-Nr. kein zweites Mal vergeben. Es stehen bis zu 999 Probe-Nr. zur Verfügung. Erst mit Vergabe aller Nr. ist eine doppelte Nr.-Vergabe möglich, eine Displaymeldung weißt gegebenenfalls darauf hin.

#### Weiter:

Schaltet zur nächsten Probe.

#### Aktiv:

Auswahlmöglichkeit ob diese Probe in der Kalibrierung berücksichtigt werden soll.

#### Messwert:

Anzeige des Messwertes, berechnet mit den aktuellen Koeffizienten.

#### Laborwert:

Eingabeposition für den Laborwert

#### **Erweitert:**

Schaltet zur nächsten Datenseite

#### Löschen:

Durch kurzes drücken der Softkey Taste kann der angezeigte Probeneintrag gelöscht werden. Durch längeres drücken können alle Probeneinträge gelöscht werden.

## 4.2.17 Erweiterte Probedaten



Stromeingang 1:

Editierbare Anzeige des ersten Kompensationseingangs.

#### Stromeingang 2:

Editierbare Anzeige des zweiten Kompensationseingangs.

#### PT100:

Editierbare Anzeige des PT100 Eingangs.

#### Phi(fm):

Nicht editierbare Anzeige der gemessenen Phase.

#### Dämpfung:

Nicht editierbare Anzeige der gemessenen Dämpfung.

### 4.2.18 Extras



#### **Tara Werte:**

Eingabemöglichkeit von Tara-Werte für Phase und Dämpfung. Die Tara-Werte werden der Phase bzw. der Dämpfung vor der Kalibrierung zugerechnet. Die Berechnung erfolgt wie folgt:

Gl. 4-2 und 4-3

Phase = Phase<sub>mess</sub> - Phi Tara

Dämpfung = Dämpfung<sub>mess</sub> - Dämpfung Tara

#### Anzahl Kalibrier Sweeps:

Frei einstellbare Anzahl der Einzelmessungen [Sweeps], über die ein Kalibrierpunkt (bei der automatischen Probenaufnahme) gemittelt wird.

#### **Prozessart:**

Auswahl der Betriebsart, möglich ist:

- eine Konzentration [1 Messbereich]
- zwei Konzentrationen [2 Messbereiche]
- Split Konzentration [1 Messbereich mit Schaltpunkt (Split Wert) für Koeffizientenumschaltung]

#### Split Wert:

Wertmäßige Einstellung des Schaltpunktes

### 4.2.19 Kalibrierung

Die Kalibrierung erfolgt nach folgender Formel:

GI. 4-4

Messwert =  $A \cdot Phase + B \cdot D$ ämpfung + C + D  $\cdot PT100$ + E  $\cdot Eing1 + F \cdot Eing2 + G \cdot Beladung$ 

wobei gilt:	
Messwert	Konzentration / Feuchte / Trockenmasse /
	Dichte
А	Phasen-Koeffizient
В	Dämpfungs-Koeffizient
С	Offset
D	Kompensationskoeffizient für PT100-Eingang
E	Kompensationskoeffizient für Stromeingang 1
F	Kompensationskoeffizient für Stromeingang 2
G	Kompensationskoeffizient für die Beladung

Die Koeffizienten können manuell eingegeben oder aus den Einträgen der Probetabelle automatisch berechnet werden.



Hier wird die Kalibrierung entsprechend den eingestellten Parametern gestartet und die Koeffizienten automatisch aus den Einträgen der Probetabelle berechnet.

#### Kal. Basis:

Auswahl der Mikrowellensignale, welche für die Kalibrierung berücksichtigt werden. Einstellbar sind:

- Phase
- Dämpfung
- Phase und Dämpfung

Default: Phase



RTHOLD

#### Komp.-Eingang:

Hier können die zur Kompensation benötigten analogen Eingänge (PT100, Stromeingang 1 und 2) ausgewählt werden. Entsprechend den aktivierten analogen Eingängen, steht maximal folgendes zur Auswählbar:

- Kein
- Eing1
- Eing1 +Eing2
- Eing1 + PT100
- Eing1 + Eing2 + PT100
- Eing2
- Eing2 + PT100
- PT100

#### Koeffizienten:

Hier können alle Koeffizienten auch direkt eingegeben werden, z.B. Startkoeffizienten.

Die automatisch berechneten Koeffizienten werden ebenfalls hier abgelegt. Nicht verwendete Koeffizienten werden auf null gesetzt.

### 4.2.20 Ein- / Ausgänge



#### Stromausgang:

In der zu selektierenden Ebene können beide Stromausgänge abgeglichen, zugeordnet und eingestellt werden.

#### Stromeingang:

Aktivierungsebene der Stromeingänge, Abgleich und Anzeige des aktuellen Stromsignals.

#### PT100:

Hier kann ein angeschlossener PT100 aktiviert und abgeglichen werden. Anzeige des aktuellen Temperatursignals.

#### Digitalausgang:

Zuordnung der Relais 1 und 2 und Testfunktion.

#### Digitaleingang:

Statuskontrolle und Zuordnung der Digitaleingänge.

## 4.2.21 Stromausgang



#### 

Bei laufender Messung kann durch die Aktivierung eines nicht belegten oder nicht justierten Stromeingangs ein Fehler verursacht werden.

## 4.2.22 Stromausgang 1

1   -  Stromausg. 1  0	7.05 –13:25	
Zuordnung	keine	
4 mA	60.00	
20 mA	95.00	
Test / Abgleich		
Fehlerstrom		
Strom Belad.unters	schr.	

#### Zuordnung:

Dem Ausgangsstrom können folgende Signale zugeordnet werden:

- keine
- Konzentration
- Konzentration 2 (wenn aktiv)
- Stromeingang 1 oder 2 (wenn aktiv)
- PT100 (wenn aktiv)

#### 4 mA:

Anzeigewert, der dem 4 mA Wert zugeordnet wird.

#### 20 mA:

Anzeigewert, der dem 20 mA Wert zugeordnet wird.



Stromausgang 1 nur 4 – 20mA möglich

Wird die Stromausgangsgrenze überschritten, so geht die Messung in den Warnzustand, siehe *Kapitel 10.4 Geräte-zustände*.

#### Test/Abgleich:

Stromtest, Abgleich und Anzeige des aktuellen Stroms.



Bei Testfunktion sollte die Messung gestoppt sein.

RTHOLD

Zur Überprüfung der Stromschleife und eventuell angeschlossener Fernanzeigen kann über die Testfunktion ein Strom von 4 bis 20 mA eingestellt werden. Wird die Testfunktion verlassen schaltet das System automatisch auf den aktuellen Strom zurück.

#### Fehlerstrom:

Geht die Messung in den Fehlerzustand, so wird über Stromausgang ein Fehlerstrom ausgegeben, dieser kann hier eingestellt werden.

- 22 mA
- 3.5mA
- Halten
- Wert (wählbar)

## 4.2.23 Stromausgang 2

1   -  Stromaus	sg. 2  07.05–13:25	
Zuordnung	keine	
0/4 mA	60.00	
20 mA	95.00	
Bereich	4-20 mA	
Test / Abgleich		
Fehlerstrom		
Strom Belac	l.unterschr.	

#### Alle Funktionen wie bei Stromausgang 1

# TIPP

Stromausgang 2 wahlweise von 0/4 bis 20 mA einstellbar.

#### **Bereich:**

Änderung des Stromausgangs

- 0 20 mA
- 4 20 mA

## 4.2.24 Stromeingang



#### Stromeingang 1:

Bei Anwahl wird zum Aktivierungs- und Abgleichmenü gewechselt.

#### Stromeingang 2:

Wie zuvor beschrieben.

## 4.2.25 Stromeingang 1



#### Zustand:

Durch Auswahl ja/nein wird der Stromeingang aktiviert oder deaktiviert.

#### **Bereich:**

Änderung des Stromausgangs

- 0 20 mA
- 4 20 mA

#### 0/4 mA:

Anzeigewert, der dem 0/4 mA Wert zugeordnet wird.



#### 20 mA:

Anzeigewert, der dem 20 mA Wert zugeordnet wird.

#### Abgleich:

Bei Anwahl des Abgleichs folgen Sie den Anweisungen auf dem Display.

#### Live Strom:

Anzeige des aktuellen Stromsignals.

### 4.2.26 Stromeingang 2

Einstellungen entsprechen Stromeingang 1.

#### 4.2.27 PT100



#### **Aktiviert:**

Bei angeschlossenem PT100 muss der Eingang erst aktiviert werden.

### 💶 WICHTIG

Bei laufender Messung kann durch die Aktivierung eines nicht belegten oder nicht justierten PT100-Eingangs ein Fehler verursacht werden.

#### PT100 Abgleich:

Hierzu benötigen Sie einen 100 Ohm und einen 138.5 Ohm Widerstand. Folgen Sie den Anweisungen auf dem Display.

#### PT100 Live:

Anzeige der aktuellen Temperatur.

Wird eingestellt und aktiviert wie Eingang 1.

### 4.2.28 Digitalausgang

Das Messgerät verfügt über zwei Relais. Relais 1 ist mit der LED Signal 1 und Relais 2 mit LED Signal 2 verknüpft.





#### Relais 1:

Dem Relais 1 können unterschiedliche Funktionen zugeordnet werden.

- Keine
  - Fehler
- Halt
- Kein Produkt
- Alarm min
- Alarm max
- Beladungsunterschreitung

Funktion	Beschreibung
Keine	Relais- und LED Funktion ausgeschaltet
Fehler	Bei Fehler wird Relais und LED gesetzt
Halt	Wenn Haltfunktion aktiviert, wird Relais und LED gesetzt
Offline	Befindet sich die AWE im Offline Modus, so wird dies über das Relais und der LED sig- nalisiert
Alarm min.	Bei unterschreiten des einzustellenden Grenzwertes schaltet das Relais
Alarm max.	Bei Überschreiten des einzustellenden Grenzwertes schaltet das Relais

#### Relais 2:

Zuordnungen wie oben beschrieben möglich.

#### Test:

Der Schaltzustand der Relais kann hier gesetzt und an den zugehörigen Klemmen überprüft werden.

### 4.2.29 Digitaleingang



#### Status:

Zeigt den Zustand der Eingangsbeschaltung an

Das Messgerät verfügt über 3 digitale Eingänge, denen unterschiedliche Funktionen zugeordnet werden können.

auf/zu

#### **DI 1 Funktion:**

DI 1 können nachstehende Funktionen zugeordnet werden

- Keine
- Start (Externer Start)

THOLD

#### **DI 2 Funktion:**

- DI 2 können nachstehende Funktionen zugeordnet werden
- Keine
- Halt (Mittelwertbildung wird angehalten)
- Produkt (Externe Produktauswahl)

#### **DI 3 Funktion:**

- Zuordnungen für DI 3
- Keine
- Probe (Externe Steuerung der Probennahme)
- Produkt (Externe Produktauswahl)

Zur externen Startfunktion muss im Menüfenster **Messung** die Startfunktion auf **extern** gesetzt werden.

*Halt* bedeutet, dass die Mittelwertbildung angehalten wird, die Messung jedoch weiterläuft.

**Probe** bedeutet, dass durch schließen des Kontaktes die Probenahme gestartet wird.

**Produkt** bedeutet, dass durch schließen des Kontaktes in ein anderes Produkt gewechselt wird (Produkt 1 bis 4).



Bei erstmaliger Produktanwahl (Produkt 2 bis 4) werden alle Einstellungen und Inhalte des aktuellen Produkts in das neue Produkt kopiert. Dazu gehören:

- Konfigurationsdaten
- Systemabgleich
- Kalibrierdaten (inklusive Probentabelle)
- Ein-/ Ausgangsdefinitionen

Für die Umschaltung aller 4 Produkte, muss DI 3 ebenfalls auf Produkt eingestellt werden. Die Klemmenbelegung entnehmen Sie bitte der nachfolgenden Tabelle.

Klemmen	DI 2	DI 3
	13 / 25	14 / 26
Produkt 1	offen	offen
Produkt 2	geschlossen	offen
Produkt 3	offen	geschlossen
Produkt 4	geschlossen	geschlossen

## 4.2.30 Service



#### Fabrikeinstellung und General Reset

Siehe Tabelle auf der nächsten Seite.

#### **Memory Tool:**

Bezieht sich auf die Kommunikation mit dem externen Speichermodul (Memory Tool, optionales Zubehör). Die Datenübertragung erfolgt über den 9-poligen SubD-Stecker auf der Geräteunterseite.

- Parameter sichern: Alle Betriebsparameter werden für alle Produkte auf den Memory Tool gespeichert.
- Parameter laden: Alle Betriebsparameter werden vom Memory Tool in die AWE geladen. Damit werden alle Betriebsparameter in der AWE gelöscht.
- Datenlog sichern: Der Datenlog wird auf dem Memory Tool gespeichert.
- Protokoll sichern: Das Inbetriebnahme-Protokoll wird auf dem Memory Tool gespeichert.

HINWEIS

Während der Kommunikation mit dem Memory Tool ist der Konzentrationsmittelwert eingefroren! Damit ist auch der Messwert über Stromausgang eingefroren!

#### Datenausgabe:

Sämtliche Messwerte werden für jede Messung über die seriellen Datenschnittstellen RS232 und RS485 ausgegeben. Die Ausgabe kann wie folgt eingestellt werden:

- Keine (Deaktivierung)
- Zeile (Datenübertragung, siehe *Hardwareteil Kapitel* 6.4)
- Tabelle (Mikrowellendaten für jeden Frequenzpunkt)
- Zeile und Tabelle

Default ist "Zeile".



	Fabrikeinstellung	General Reset
Sprachauswahl	unverändert	unverändert
Zugriffsebene	unverändert	default: Benutzer Modus
Messung	wird gestoppt	wird gestoppt
Passwort	unverändert	default: PASS1
Produktauswahl	unverändert	alle Produkte gelöscht
Fehlerlog	nicht gelöscht	gelöscht
Datenlog	nicht gelöscht, Einstel- lungen default	gelöscht, Einstellungen default
Systemabgleich	nicht gelöscht	gelöscht
Kabellänge	unverändert	default
Probentabelle	nicht gelöscht	gelöscht
Messstellen-Bezeichnung	default	default
Alle Parameter unter Menü: Messung Plausibilität Marker Einheit	default	default
Kalibrier-Koeffizienten	default	default
Alle Einstellungen unter den analogen und digitalen Ein- und Ausgängen	default	default
Justage der analogen Ein- und Ausgänge	unverändert	unverändert
Bemerkung:	wirkt nur auf das aktu- elle Produkt	wirkt für alle Produkte (P1 bis P4)
-		

\*Default: Default Werte, siehe Kapitel 11.1 Beispiel Inbetriebnahme-Protokoll

## 4.3 Trendanzeige



Mit der **ZOOM** Taste kann der eingerahmte Messwert als Großanzeige angezeigt werden.



Durch längeres drücken der **ZOOM** Taste wird der großangezeigte Messwert als Trend über das gesamte Display dargestellt.

Die Trenddarstellung entspricht dem Inhalt des Datenlogs. Der Datenlog muss für die Trenddarstellung aktiviert sein.

HINWEIS

Für die Zeit des Trendaufbaus wird der Messwert bzw. der Stromausgang eingefroren!



## Kapitel 5. Konfiguration

Vor der Kalibrierung müssen die erforderlichen Analogeingänge aktiviert und konfiguriert werden sowie die Konfigurationsparameter überprüft und gegebenenfalls korrigiert werden.

Ohne Aktivierung der erforderlichen Eingänge werden einige Menüs nicht angezeigt und eine ordentliche Konfiguration und Kalibrierung ist unter Umständen nicht möglich. Die Stromausgänge, digitalen Ein- und Ausgänge können auch noch nach der Kalibrierung aktiviert und konfiguriert werden.

Das Messsystem verfügt über zwei unabhängige potentialfreie Stromausgänge.

## 5.1 Konfigurationseinstellungen

1 | - |Einstellungen | 07.05 – 13:25 Konfiguration Kalibrierung Ein-/Ausgänge Service Produkt Passwort ändern zeige über | EINSTELLUNGEN |

Ausgehend vom Hauptmenü erreichen Sie nebenstehende An-

KONFIGURATION

## 5.1.1 Allgemeines



> ALLGEMEINES







#### **Beispiel:**

Es wird der entsprechende Eintrag selektiert, geändert und durch Bestätigung abgespeichert.

> DATUM

Mit **LÖSCHEN** wird der Eintrag gelöscht und anschließend das neue Datum eingegeben. Das geänderte Datum wird mit vestätigt und gespeichert.



Den Doppelpunkt für die Uhrzeiteingabe (z.B. 13:25) rufen Sie mit der Taste [ . ] auf

#### 5.1.2 Messung



Die Einstellungen auf dieser Seite müssen überprüft und den Messbedingungen entsprechend angepasst werden.

Die Mittelung über 20 Messungen ist in der Regel eine gute Wahl.

## 5.1.3 Plausibilität



Die **Prozessgrenzen** müssen angepasst werden. Erlauben Sie eine Messbereichsüberschreitung von  $\pm$  5 % absolut. Beispiel: Der Messbereich beträgt 65-95 %TS. So geben Sie als Prozessgrenzen 60-100 %TS ein.

Die Prozessgrenzen sind unabhängig von den Stromausgangsgrenzen.

Die Einstellungen der **Phasenmessung** siehe *Kapitel 5.3 Plausibilitäten konfigurieren*.

Der **Offline Modus** ist werkseitig ausgeschaltet. Nehmen Sie die Funktion bei Bedarf erst in Betrieb, wenn das Messsystem fehlerfrei in Betrieb genommen wurde. Lesen Sie dazu die Beschreibung unter *Kapitel 4.2.11 Offline Modus*.

## 5.1.4 Mikrowelle



KABEL



Entsprechen die werksseitig eingestellten Kabellängen nicht den tatsächlichen Geometrieverhältnissen, so sind die Werte zu korrigieren.

Beispiel: Für ein 4 m langes HF-Kabel Quad geben Sie 8 m für die Ref- und Messkabellänge ein. Der Eingabewert entspricht immer dem doppelten der Quad Kabellänge.

## 5.1.5 Einheit



Die Einheiten sollten auf die gewünschte Dimension eingestellt werden.

➢ EINHEIT

1  -   Einheiten	07.05 – 13:25
Konz	%
Konz 2	%
Stromeingang 1	keine
Stromeingang 2	keine
Temp. PT100	°C

Die Einheiten der Konzentrationen (Konz 1 und 2) und die der aktivierten analogen Eingänge können ausgewählt werden.

KONZ / KONZ 2



Die Dimensionen können für beide Konzentrationen unterschiedlich eingestellt werden.

≻ %





## 5.1.6 Marker



Es kann ein Marker von max. 5 Zeichen eingegeben werden, der im Live Display den eingestellten Wert kennzeichnet.

> MARKER

## 5.2 Start-Kalibrierkoeffizienten



- |Kalibrierung | 07.05 –13:25

 $\mathbf{\nabla}$ 

Phi

kein

Start Kalibrierung

Komp.-Eingang

Koeffizienten

Kal. Basis Beladungskomp. Ausgehend vom Hauptmenü erreichen Sie nebenstehende Anzeige über: ZURÜCK | EINSTELLUNGEN | KALIBRIERUNG | KALIBRIERE

➢ KALIBRIERUNG

KONZ

#### ➢ KOEFFIZIENTEN

Für die default Einstellung berechnet sich die Konzentration wie folgt:

#### Messwert = $A \cdot Phase + C$

Gl. 5-1

Wobei gilt: A, C: Kalibrierkoeffizienten



Überprüfen Sie die Koeffizienten A und C und korrigieren Sie wenn nötig wie folgt:

C = Mittlerer Messbereichswert (Konzentrationswert)

A = 0

Alle nicht benötigten Koeffizienten werden automatisch auf null gestellt.

Hinweis: Mit diesen Kalibrierkoeffizienten sind während der Inbetriebnahme der Konzentrationsmittelwert und damit auch der Stromausgang eingefroren.

## 5.3 Plausibilitäten konfigurieren

Ausgehend vom Hauptmenü erreichen Sie nebenstehende Anzeige über

| EINSTELLUNGEN | KONFIGURATION | PLAUSIBILITÄT | PHA-SENMESSUNG |

Nebenstehende Anzeige zeigt die Default Einstellungen, diese müssen gegebenenfalls eingegeben werden.

Die **Streuung (Sigma)** liegt im normalen Messbetrieb zwischen 0 bis 100. Daher ist das Sigma<sub>max</sub> = 100 für die meisten Applikationen eine gute Wahl. Sollten Sie während der Messung jedoch laufend Sigma-Werte bis zu 300 erhalten, so kann der Grenzwert bis auf 500 erhöht werden. Bei noch höheren Sigma-Werten, liegt meist eine Störung vor. Wie z.B. ständige Luftblasen, die beseitigt werden müssen.

Das **Verhältnis Phi/D** ist für die meisten Applikationen nicht bekannt, daher muss es einmalig bestimmt werden. Hierfür wird die Auto-Einstellung benötigt, die ansonsten immer auf AUS stehen muss.

Die genaue Vorgehensweise ist unter Kapitel 9.1 beschrieben.





# Kapitel 6. Kalibrierung

Hinweis: Das Messgerät muss die normale Betriebstemperatur erreicht haben (ca. 45 min. Aufwärmzeit).

Es muss sichergestellt sein, dass die Messzelle ganz mit Produkt gefüllt ist, bzw. die Behältersonde ganz mit Produkt umgeben ist.

Die Messeinrichtung muss bei normalem Durchsatz bzw. Materialstrom und dem üblichen Material unter tatsächlichen Betriebsbedingungen laufen.

Voraussetzung: Das **Kapitel 5. Konfiguration** wurde durchgeführt.

## 6.1 Systemabgleich

1   -   LB 566   07.05 – 13:25	Ausgehend vom Hauptmenü:
Live Anzeige Diagnose Einstellungen Zugriffsebene Benutzer M. Sprache Deutsch RUN I I	➢ EINSTELLUNGEN
1   -  Einstellungen   07.05 – 13:25 Konfiguration Kalibrierung Ein-/Ausgänge Service Produkt Passwort ändern	≻ KALIBRIERUNG
1   -   Kalibrierung   07.05 – 13:25 Systemabgleich Kalibriere Konz Extras	➢ SYSTEMABGLEICH

MicroPolar (++) LB 566









Bestätigung



Der Systemabgleich wird nun durchgeführt.

1   -  Systemabgleich 07.05 – 13:25
Abgeglichen!
OK

Mit **OK** bestätigen und mit 3mal **A** zurück zum Hauptmenü.

## 6.1.1 Überprüfung der Referenzwerte



Ausgehend vom Hauptmenü erreichen Sie nebenstehende Anzeige über | EINSTELLUNGEN | KALIBRIERUNG | SYSTEMAB-GLEICH |

➢ REFERENZWERTE

Grenzwerte wichtiger Parameter für die Referenzmessung:

1   -   Referenzwerte  07.05 – 13:25		
Phi(fm) Dämpfung Steigung Sigma	125.00 °/GHz 22.5 dB 120.55 °/GHz 0.00	

Mikrowellen-Durchstrahlung)	
Dämpfung: < 45 dB bei Standard AWE MicroPolar < 65 dB bei High Dynamik AWE MicroPolar+ (Kabel richtig angeschlossen, Produkt durch strahlbar)	·+ -

## 6.2 Probenahme





Im Falle einer Temperaturkompensation muss die Produkttemperatur über eines der analogen Eingänge eingespeist und zusätzlich der entsprechende Eingang aktiviert werden. Ohne dies wird bei den Probenahmen am Gerät die Produkttemperatur nicht gespeichert.

Falls sich das Messsystem noch nicht im Messmodus befindet, so starten Sie jetzt die Messung.

Das Messsystem wird jetzt mit *RUN* gestartet.

Nach Bestätigung der Sicherheitsabfrage 🗸 geht das Gerät in den Messmodus.





Das nebenstehende Display wird nach **RUN** angezeigt.

**Hinweis:** Mit **PROBE** wird die Aufnahme der Rohdaten gestartet. Synchron zur Datenaufnahme muss die Laborprobe genommen und gekennzeichnet werden. Eine Auswertung kann später erfolgen, sofern das Produkt sich dadurch nicht verändert.



Probenahme läuft.....

Mit Taste X kann die Probenahme jederzeit gestoppt werden.



Lief die Probenahme ohne Zwischenfall ab, so wird durch Drücken der Taste die Probe in der Tabelle gespeichert und die Messung läuft weiter.

Der vorab beschriebene Vorgang muss für jede weitere Probe wiederholt werden.

Die Feuchte/Konzentration der Proben sollte über den gesamten Messebereich verteilt sein. Bei zusätzlicher Temperaturkompensation sollte die Temperatur der Proben über den gesamten Temperaturbereich verteilt sein.

In Abhängigkeit von der eingestellten Kalibrierung ergibt sich die min. erforderliche Probenanzahl. Ist die Probenanzahl zu gering, so folgt nach versuchter Kalibrierung eine Fehlermeldung.

Für eine grobe Berechnung der Kalibrierkoeffizienten reichen etwa 6 Proben aus, wenn die Konzentration sich um mindestens 5 % unterscheidet. Für eine Fein-Kalibrierung sowie mit Temperaturkompensation sind mindestens 15 Proben erforderlich.

## 6.2.1 Eingabe der Laborwerte



#### Kapitel 6. Kalibrierung







1   1/1   Probe Nr.1 Laborwert	07.05 –13:25
0.00	%
ZURÜCK ? LO	ÖSCHEN √

PROBENTABELLE

LABORWERT

Vorgabewert mit LÖSCHEN löschen, neuen Wert eingeben und mit 🗸 bestätigen

1   1/1   Probe Nr.1	07.05 –13:25
Weiter	
Aktiv	Ja
Messwert	65.50 %
Laborwert 72,40 %	
Erweitert	
<ul> <li>↓ LÖSCHEN</li> </ul>	

➢ WEITER

und oben beschriebenen Vorgang mit der nächsten Probe wiederholen.

Nach Eingabe der letzten Probe durch Drücken der **T**aste zurück zum Kalibriermenü. (Kurzer Tastendruck eine Seite zurück, längerer Tastendruck sofort zurück zum Kalibriermenü)

## 6.3 Kalibrierung

Gehen Sie vor, wie unter Kapitel 8.4 beschrieben.

## Kapitel 7. Inbetriebnahme MicroPolar ++

Der Systemabgleich und die Kalibrierung erfolgen beim Micro-Polar ++ genauso wie beim Standardmodel MicroPolar. Jedoch gilt es zu beachten, dass die ++ Einheit eine minimale Dämpfung von 40 dB über den gesamten Konzentrationsbereich und beim Systemabgleich benötigt. Bei Unterschreitung ist die Messung nicht mehr eindeutig.

Die gesamte Dämpfung errechnet sich wie folgt:

 $dB_{gesamt} = dB_{Abgleich} + dB_{Live} + 0.4 x Messkabellänge$  **GI. 7-1** 

Wobei gilt:	
dBgesamt:	Gesamte Dämpfung
dB <sub>Abgleich</sub> :	Dämpfung bei Systemabgleich
dB <sub>Live</sub> :	Aktuelle Dämpfung im Messbetrieb
Messkabellänge:	z.B. 4 m HF-Kabel quad ergibt 8 m
-	Messkabellänge (Hin- und Rückweg)

Die Auswerteeinheit überwacht die gesamte Dämpfung automatisch und meldet eine Unterschreitung durch eine Fehlermeldung (Fehler Nr. 55).

#### Abhilfe bei Dämpfungsunterschreitung:

Wird die erforderliche Gesamtdämpfung unterschritten, so besteht die Möglichkeit ein 10 dB Festdämpfungsglied in den Sendezweig zu montieren (siehe Abbildung 7-1). Bei deutlicher Unterschreitung jedoch sollte das Standardmodel Micro-Polar eingesetzt werden.



*Abb. 7-1: Montage des 10 dB Dämpfungsgliedes*
# Kapitel 8. Kalibrierung und Extras

## 8.1 Plausibilität Verhältnis Phi/D konfigurieren





Ausgehend vom Hauptmenü erreichen Sie nebenstehende Anzeige über | EINSTELLUNGEN | KONFIGURATION | PLAUSIBI-LITÄT | PHASENMESSUNG |

Überprüfen bzw. korrigieren Sie entsprechend der nebenstehenden Einstellungen.

➢ VERHÄLTNIS PHI/D

Ist der Verhältniswert bekannt, so geben Sie ihn hier ein, anderenfalls gehen Sie zu *Kapitel 8.1.1 Verhältnis Phi/D*.

Vorgabewert mit LÖSCHEN löschen, neuen Wert eingeben und mit √ bestätigen

### 8.1.1 Verhältnis Phi/D

Ist der Verhältniswert nicht bekannt so gibt es folgende Möglichkeiten:

- 1. Führen Sie eine Prozessaufnahme durch, siehe dazu *Kapitel 8.1.2 Prozessaufnahme*. Voraussetzung hierfür ist, dass während der Aufnahme der Prozess über den gesamten Konzentrationsbereich gefahren wird.
- Ist die Prozessaufnahme aktuell nicht möglich oder ist der Konzentrationsbereich ohnehin klein (< ± 5%), so geben Sie als Startwert Phi/D = 1 ein.

Eine nachträgliche Anpassung oder Prozessaufnahme ist möglich.



#### 8.1.2 Prozessaufnahme

Voraussetzung: Die Kapitel **5. Konfiguration 6.1 Systemabgleich** wurden durchgeführt.

Die Prozessaufnahme dient zur Bestimmung des Verhältnisses aus Phase und Dämpfung (Phi/D), einem Parameter aus der Plausibilitätsanalyse zur korrekten Bestimmung der Phase.

Ist das Verhältnis bereits aus andern Messungen bekannt, so kann es direkt unter Menü PLAUSIBILITÄT eingegeben werden (siehe *Kapitel 4.2.10 Phasenmessung*), die Prozessaufnahme ist in diesem Fall nicht erforderlich.

## **İ** WICHTIG

Die Aufnahme erfolgt automatisch, Sie muss lediglich ge-startet und wieder gestoppt werden. Während der Aufnahme müssen die folgenden Punkte unbedingt beachtet werden.

- Stoppen Sie nicht die Messung
- Ändern Sie die Konzentration nicht sprunghaft (max. 1 %)
- Durchlaufen Sie möglichst den gesamten Messbereich

Ausgehend vom Hauptmenü erreichen Sie nebenstehende Anzeige über

| EINSTELLUNGEN | KONFIGURATION | PLAUSIBILITÄT |

PHASENMESSUNG

AUTO-EINSTELLUNG





BERTHOLD



1 | - |Phasenmessung|07.05 –13:25 Sigma max. 100.00 Verhältnis Phi/D 5.32 Auto- Einstellung EIN ≻ EIN

#### Anhalten der Aufnahme:

Sie haben die Möglichkeit die Aufnahme anzuhalten, indem Sie die Aufnahme ausschalten. Die Aufnahme wird eingefroren und beginnt erst wieder mit erneutem einschalten.

#### Neue Aufnahme starten:

Voraussetzung: die Aufnahme ist ausgeschaltet. Stoppen und Starten Sie die Messung bevor Sie eine neue Aufnahme starten. Die Ergebnisse älterer Aufnahmen sind damit gelöscht.

#### Aufnahme beenden:

Nach der Aufnahme schalten Sie die automatische Aufnahme einfach aus, dass aufgenommene Phi/D wird automatisch abgespeichert.

Führen Sie jetzt, während die Prozessaufnahme läuft, die Probeaufnahmen durch.

Vergessen Sie nicht die Prozessaufnahme auf gleicher Weise wie oben beschrieben wieder auszuschalten!



## 8.2 Anpassung der Kalibrierung

Es kann nachträglich ein Korrekturfaktor und ein Offset eingegeben werden, um eine nachträgliche Anpassung der Kalibrierung zu erreichen (Feinkalibrierung).

Nachfolgend ein Beispiel für eine Offset-Anpassung.

Das nebenstehende Display wird nach **RUN** angezeigt.

Die Displayanzeige wird nun mit dem Analysewert der Laborprobe verglichen. Die Differenz muss vorzeichenrichtig als Offset eingegeben werden.

Berechnung:

Analysewert – Display = Offset

Gl. 8-1



1 | - | Live Anzeige | 07.05 - 13:25

ZURÜCK PROBE ▲▼ ZOOM

M. Konz.

Mittlere Konzentr.

65.50 %

Akt. Konz. 64.35%

Mit ZURÜCK zum Hauptmenü.

> EINSTELLUNGEN

1   -  Einstellungen   07.05 – 13:25
Konfiguration
Kalibrierung
Ein-/Ausgänge
Service
Produkt
Passwort ändern



> KALIBRIERUNG

KALIBRIERE KONZ





> ANPASSUNG

1   -   Anpassung	07.05 –13:25
Faktor	1.00000
Offset	0.000
	7 8

OFFSET

Berechnungsformalismus siehe Kapitel 4.2.16 Kalibriere Konzentration.



Berechneten Offsetwert eingeben, mit 🔨 Taste bestätigen und mit 4-mal Hometaste 🗹 zum Hauptdisplay zurück.



Mit

LIVE ANZEIGE

zurück zur Anzeige.



Der Anzeigewert sollte nun dem Istwert entsprechen.



### 8.3 Ausgabe des Inbetriebnahme-Protokolls



Ausgehend vom Hauptmenü erreichen Sie nebenstehende Anzeige über | DIAGNOSE |

PROTOKOLL DRÜCKEN



Mit 🗸 Taste wird das Protokoll über RS232 und RS485 ausgegeben. Mit 🗙 gehen Sie ohne Ausgabe eine Seite zurück.

Das Inbetriebnahme-Protokoll beinhaltet alle Parameter, Daten des Systemabgleichs, Kalibrierdaten und Einträge der Probentabelle.

Weitere Informationen zum Inbetriebnahme-Protokoll siehe Kapitel 11. Inbetriebnahme-Protokoll.

### 8.4 Kalibrierung

Bei der Kalibrierung gibt es mehrere Einstellmöglichkeiten, Details siehe *Kapitel 4.2.20 Kalibrierung*. Im Weiteren werden die Standardeinstellungen angezeigt, welche in der Regel die beste Wahl sind.

Voraussetzung:	
Die Kapitel	5. Konfiguration
	6.1 Systemabgleich
	6.2 Probenahme
wurden durchae	eführt.



Ausgehend vom Hauptmenü erreichen Sie nebenstehende Anzeige über | EINSTELLUNGEN | KALIBRIERUNG | KALIBRIERE KONZ |

KALIBRIERUNG

KALIBRIER BASIS

- 1 | | Kalibrierung | 07.05 –13:25 Kal. Basis Phase Dämpfung Beide

PHASE (Phasenmessung)

#### Standard für alle Applikationen: Phase



#### Gewünschte Kompensation einstellen:

1   -   Kalibi	rierung	07.05	5 – 13:25
Start K Kal. Ba Beladu	alibrieru sis ngskon	Ing I 1p.	Phase
Komp	Eingan	g.	Kein
Koeffiz	ienten		

Für die Kompensation, wie z.B. **Temperaturkompensation**, gehen sie wie folgt vor:

KOMP.-EINGANG

Hier können die zur Kompensation (z.B. Temperaturkompensation) benötigten analogen Eingänge (PT100, Stromeingang 1 und 2) ausgewählt werden. Auswählbar sind:

- Kein
- Eing1
- Eing1 + Eing2
- Eing1 + PT100
- Eing1 + Eing2 + PT100
- Eing2
- Eing2 + PT100
- PT100

Wenn keine Kompensation erforderlich ist, dann wird hier "kein" ausgewählt.

1 | - | Kalibrierung | 07.05 – 13:25 Start Kalibrierung Kal. Basis Phase Beladungskomp. Komp.-Eingang kein Koeffizienten Starten Sie nach Einstellung der Kompensation die automatische Berechnung der Kalibrierkoeffizienten.

START KALIBRIERUNG



1 | - | Kalibrierung | 07.05 –13:25 Kalibriert! OK Mit **V** Taste wird die Kalibrierung gestartet mit **X** gehen Sie ohne Kalibrierung eine Seite zurück.

OK übernimmt die Kalibrierung und wechselt zur nächsten Anzeige.

Bei der Berechnung der neuen Koeffizienten werden der Faktor auf 1 und der Offset auf 0 zurückgesetzt.

BERTHOLD



Die nebenstehende Kurve zeigt die Darstellung Anzeige über Labor.

|--|



Ausgabe der Korrelation zwischen Anzeige und Labor.

OK



Nach Bestätigung erscheint wieder das Kalibriermenü, von wo aus Sie mit 4-mal d zum Hauptmenü gelangen und die Messung wieder gestartet werden kann.

### 8.4.1 Kalibrierung mit zwei Konzentrationen

1 | - | Kalibrierung | 07.05 – 13:25 Systemabgleich Kalibriere Konz Extras

1 | - | Extras | 07.05 –13:25 Tara Werte Anz. Kal. Sweeps 20 Prozessart 1 Konz.

1 | - | Extras |07.05 –13:25 Prozessart 1 Konz 2 Konz Split Konz ZURÜCK ? ▲▼ √



1   -  Kalibriere Konz1 07.05 –13:25
Probentabelle
Kalibrierung
Anpassung
Ergebnis

> EXTRAS

Display über

PROZESSART

> 2 KONZ

Mit der ✓ Taste wird die selektierte Prozessart übernommen und mit 1-mal ▲ Taste erreichen Sie nachstehendes Display.

Die Kalibrierung für zwei Konzentrationen beginnt mit der Än-

Ausgehend vom Hauptmenü erreichen Sie nebenstehendes

derung der Prozessart wie unten beschrieben.

| EINSTELLUNGEN | KALIBRIERUNG |

> KALIBRIERE KONZ (entspricht der Konzentration 1)

> PROBENTABELLE

1   1/4   Probe Nr.	1   07.05 – 13:25
Weiter	
Aktiv	Ja
Messwert	65.50 %
Laborwert	0.00 %
Erweitert	
<ul> <li>LÖSCHEN</li> </ul>	$\land$

Für beide Kalibrierungen gibt es nur eine Probentabelle. Zu allen Proben, die für die Kalibrierung der Konzentration1 verwendet werden sollen, müssen die Laborwerte eingegeben werden. Alle anderen Proben müssen deaktiviert werden (aktiv: ja/nein).

LABORWERT

1   1/4   Probe Nr. 1   07.05 –	13:25
Laborwert	
60.40	%
ZURÜCK ? LÖSCHEN	$\checkmark$

Vorgabewert mit Löschen löschen, neuen Wert eingeben und mit √ bestätigen

1  1 /4   Probe Nr.1	07.05 –13:25
Weiter	
Aktiv	Ja
Messwert	65.50 %
Laborwert	60.40 %
Erweitert	
◄ LÖSCHEN	$\land$

> WEITER

zur nächsten Probe

1  2 /4   Probe Nr.2	07.05 –13:25
Weiter	
Aktiv	Ja
Messwert	74.35 %
Laborwert	67.80 %
Erweitert	
▲ LÖSCHEN	

Probe deaktivieren

1   2/4   Probe Nr. 2   07.05 – 13:25 Aktiv
Nein
Ja
ZURÜCK ? LÖSCHEN √

> NEIN



1  2 /4   Probe Nr.2	2   07.05 –13:25
Weiter	
Aktiv	Nein
Messwert	74.35 %
Laborwert	67.80 %
Erweitert	
<ul> <li>LÖSCHEN</li> </ul>	$\land$



1   -   Kalibrierung  0	7.05–13:25
Start Kalibr.	
Kal. Basis	Phase
Beladungskomp.	
KompEingang kein	
Koeffizienten	

1   -   Kalibrierung   07.05 – 13:25
Jetzt kalibrieren?
X



Es muss sichergestellt sein, dass alle Proben bearbeitet wurden und nur die für diese Kalibrierung relevanten Proben aktiv sind.

Mit < zurück zur Kalibrierseite

START KALIBRIERUNG

Mit 🗸 Taste wird die Kalibrierung gestartet mit X gehen Sie ohne Kalibrierung eine Seite zurück.

**OK** übernimmt die Kalibrierung und wechselt zur nächsten Anzeige.





Mit 2-mal d gelangen Sie zwei Seiten zurück.



➢ KALIBRIERE KONZENTRATION 2

1   -  Kalibriere Konz2 07.05 –13	25	
Probentabelle		
Kalibrierung		
Anpassung		
Ergebnis		

wiederholen, wobei in der Probentabelle alle Proben wieder aktiviert sind. Es müssen nun alle Proben, die für Konzentration 2 nicht herangezogen werden sollen, deaktiviert werden.

Die Schritte wie oben beschrieben sind für Konzentration 2 zu

➢ PROBENTABELLE



### 8.4.2 Kalibrierung mit Splitwert

Bei dieser Art der Kalibrierung werden in einem Messbereich zwei Kennlinien (Konzentrationen) zusammengefasst, deren Schnittpunkt den Splitwert festlegt.

Die Konz 1 für den unteren und Konz 2 für den oberen Messbereich können nur gemeinsam über Stromausgang herausgegeben werden.

EXTRAS



1 | - | Kalibrierung | 07.05 – 13:25

Systemabgleich Kalibriere Konz

Extras





> PROZESSART

SPLIT KONZ

Mit der Taste wird die selektierte Prozessart übernommen und mit 1-mal Taste erreichen Sie nachstehende Anzeige. Der angezeigte Splitwert ist werkseitig voreingestellt, muss aber der jeweiligen Applikation angepasst werden.

Die Probeaufnahmen sollten so gewählt werden, dass die letzte Probe der unteren Konzentration möglichst dicht bei der ersten Probe der oberen Konzentration liegt. Im Idealfall ist die letzte Probe der Anfangskonzentration die erste Probe der Endkonzentration.



Die Probeaufnahmen erfolgen kontinuierlich über den gesamten Messbereich mit nebenstehender Anzeige. Siehe *Kapitel 6.3 Probenahme.* 

Nach Abschluss der Probenahme werden während der Eingabe der Laborwerte die einzelnen Proben, in Abhängigkeit des eingestellten Splitwertes, aktiviert oder deaktiviert. Alle Proben kleiner oder gleich dem Splitwert werden dem unteren Konzentrationsbereich zugeordnet und alle Proben oberhalb dem oberen Konzentrationsbereich.

Die Zuordnung der Proben wird automatisch durchgeführt, etwa nach Setzen des Splitwerts oder nach Eingabe der Laborwerte (z.B. nach erneuter Probenahme). Die Zuordnung richtet sich nach dem Splitwert und dem Laborwert.

# **WICHTIG**

Mit der Splitwerteingabe können Proben, die zuvor deaktiviert waren, wieder durch die automatische Zuordnung aktiviert werden! In solchen Fällen sollten deaktivierte Proben besser gelöscht oder nach Splitwerteingabe nochmals deaktiviert werden!

Der einzustellende Splitwert muss dem Schnittpunkt der beiden Kalibrierkennlinien entsprechen. Dieser wird nach der Kalibrierung automatisch (in Grenzen) korrigiert.

> SPLITWERT

Eingabe des Splitwertes und mit  $\checkmark$  bestätigen.

Mit Hometaste A zurück zur Kalibrierseite.

	Π		
Spl	it Wert		75.00 %
An: Pro	z. Kal. Sı zessart	weeps S	20 plit Konz
Tar	a Werte		
7   -	Exila	5   07	.05-15.25

1   -   Extras   07.05–13	:25
Split Wert	
75.00	%
ZURÜCK ? LÖSCHEN	$\checkmark$





Kalibrierung Anpassung Ergebnis

 $\sim$ 

KALIBRIERE KONZ

KALIBRIERUNG

Die untere Konzentration wird nun kalibriert. Anschließend KONZ2 anwählen und Kalibriervorgang wiederholen. Zurück zum Hauptmenü und die Messung starten.

### 8.5 Typische Kalibrierkoeffizienten/Startwerte

C: Konzentrationswert bei Systemabgleich

#### Für Applikationen mit Behältersonde oder Messzelle Nennweite DN 50 sowie ohne Temperaturkompensation

- A = 0.19 für die Bestimmung der Konzentration als Trockensubstanz
- A = + 0.19 für die Bestimmung der Konzentration als Feuchte

THOLD

## Kapitel 9. Passwort

Mittels Passwort kann das Messsystem gegen Zugriff Unberechtigter geschützt werden.

Die Zugriffsebenen stellen sich wie folgt dar.

#### Nur lesen

Das Messgerät kann nicht gestartet und nicht gestoppt werden. Es kann nur vom Anzeigedisplay zur Diagnose und zur Zugriffsebene geschaltet werden.

#### **Benutzer Modus**

Der Benutzer Modus ist der Standard Modus und ermöglicht den Zugriff auf alle benutzerrelevanten Parameter.

#### Service

Die Serviceebene ist dem Kundendienst vorbehalten.

Der Zugriffwechsel von "Nur lesen" auf "Benutzer Modus" ist nur durch Passwort möglich. Dieses Passwort ist zum Zeitpunkt der Auslieferung

#### PASS1

Das Passwort kann unter Menü | EINSTELLUNGEN | PASS-WORT ÄNDERN | geändert werden.

#### 9.1 Passwort vergessen

Das Gerät befindet sich im "Nur lesen" Modus und der Benutzer hat das Passwort vergessen. Um einen "Reset" der Benutzerebene durchzuführen, muss man folgendermaßen vorgehen:

Gerät ausschalten.

Gerät einschalten, in dem Moment, wenn alle 5 LED's beim Hochfahren aufleuchten, die 0 (Null) drücken und 8 Sekunden gedrückt halten.

Gerät fährt im "Benutzer Modus" hoch. Sie können jetzt ein neues Passwort eingeben.

### 🚺 WICHTIG

Prüfen Sie Ihren Prozess, bevor Sie das Gerät ausschalten. Die Stromausgänge z.B. fallen auf 0 mA.

# Kapitel 10. Fehlerlisten und Gerätezustände

Die LED's signalisieren den Gerätezustand. Nachdem die Fehler behoben sind, geht die Messung in den Zustand vor Fehlereintritt. Ein Quittieren ist nicht erforderlich.

### 10.1 Stromausfall

Code	Fehler	Mögliche Ursache/ Behebung
10	24 V Betriebsspannung	<i>Bitte an den Service von Berthold Technologies wenden.</i>
11	9 V Betriebsspannung	<i>Bitte an den Service von Berthold Technologies wenden.</i>
12	5 V Betriebsspannung	Bitte an den Service von Berthold Technologies wenden.
13	3 V Betriebsspannung	Bitte an den Service von Berthold Technologies wenden.
14	Batteriespannung	Batterie ist bald leer, um- gehend austauschen, siehe Hardwareteil Kapitel 5.4

### 10.2 Temperaturfehler

Code	Fehler	Mögliche Ursache/ Behebung
20	Achtung: Umgebungstemperatur zu hoch!	Betriebstemperatur der Auswerteeinheit prüfen, zulässiger Bereich: -20 bis 60 °C
21	<i>HF Temperatur außer Bereich</i>	Betriebstemperatur der Auswerteeinheit prüfen, zulässiger Bereich: -20 bis 60 °C



## 10.3 Hardwarefehler

Code	Fehler	Mögliche Ursache/ Behebung
30	Programmspeicher fehlerhaft	<i>Bitte an den Service von Berthold Technologies wenden.</i>
31	<i>Datenspeicher fehler- haft</i>	<i>Bitte an den Service von Berthold Technologies wenden.</i>
32	<i>Parameterspeicher fehlerhaft</i>	<i>Kompatibilitätsüberprü- fung nach Software- Download: Es muss ein General Reset durchge- führt werden.</i>
33	I2C-Bus Kommunika- tion	<i>Bitte an den Service von Berthold Technologies wenden.</i>
34	D/A Wandlung fehler- haft	<i>Bitte an den Service von Berthold Technologies wenden.</i>
35	LCD Kontrasteinstel- lung	Bitte an den Service von Berthold Technologies wenden.
36	LCD Kontroller	Bitte an den Service von Berthold Technologies wenden.
37	Tastatur defekt	<i>Bitte an den Service von Berthold Technologies wenden.</i>
38	HF Kommunikation ge- stört	<i>Bitte an den Service von Berthold Technologies wenden.</i>
39	<i>HF Hardware fehler- haft</i>	Gestörte Kabelverbindung zwischen Motherboard und HF-Modul. Stecker auf dem Motherboard prü- fen. Achtung! Zuvor die AWE vom Netz trennen!
40	I/O Kommunikation gestört	<i>Bitte an den Service von Berthold Technologies wenden.</i>
41	I/O Modul fehlerhaft	<i>Bitte an den Service von Berthold Technologies wenden.</i>
42	HF Board Startup Feh- ler	Bitte an den Service von Berthold Technologies wenden.

## 10.4 Sensorfehler

Code	Fehler	Mögliche Ursache/ Behebung
50	Sigma der Phase ist zu groß	Die gemessene Phase überschreitet den zulässi- gen Grenzwert der Streu- ung.
53	Kein Produkt	<i>Die Auswerteeinheit befin- det sich im Offline-Zu- stand (kein Produkt vor- handen).</i>
54	Kein Systemabgleich vorhanden	Der Systemabgleich wurde noch nicht durch- geführt.
55	Einfügedämpfung ist unterschritten	Siehe Softwareteil Kapi- tel 8.

# 10.5 Fehler Bereich Analogeingang

Code	Fehler	Mögliche Ursache/ Behebung
60	<i>Stromeingang 1 außer Bereich</i>	Der aktivierte Stromein- gang wurde noch nicht abgeglichen oder ist nicht belegt.
61	<i>Stromeingang 2 außer Bereich</i>	Der aktivierte Stromein- gang wurde noch nicht abgeglichen oder ist nicht belegt.
62	<i>PT100 Temperatur au- ßer Bereich</i>	Der aktivierte PT100 Ein- gang wurde noch nicht abgeglichen oder ist noch nicht belegt.



## **10.6 Fehler Messbereich**

Code	e Fehler Mögliche Ursache/ Behebung	
70	<i>Konzentration außer Bereich</i>	<i>Die Konzentration liegt außerhalb der Prozess- grenzen.</i>
71	Konzentration 2 außer Bereich	Die Konzentration 2 liegt außerhalb der Prozess- grenzen.
72	Beladungswert1 klei- ner Mindestbelegung	<i>Unterschreitung der Min- destbeladung für Kon- zentration 1</i>
73	Beladungswert2 klei- ner Mindestbelegung	<i>Unterschreitung der Min- destbeladung für Kon- zentration 1</i>
74	Beladungskompensa- tion inaktiv. Stromein- gang obere&untere Grenze ungültig	<i>Stromeingang liegt außer- halb des Bereichs.</i>
75	Synchronisationszeit zu lang	Einstellungen zur Syn- chronisation überprüfen, siehe Kapitel 7.3 im Hard- wareteil.
76	<i>Synchronisation: Ge- schw. naußer Bereich</i>	Einstellungen zur Syn- chronisation überprüfen, siehe Kapitel 7.3 im Hard- wareteil.
77	<i>Warte auf synchrone Werte</i>	<i>Die Messung ist noch nicht synchronisiert, bitte abwarten.</i>

### 10.7 Fehler Zusatzmessung

Code	Fehler	Mögliche Ursache/ Behebung
78	Massenstrom 1 Be- rechn. inaktiv. Pro- duktgeschw. ist ungül- tig	<i>Bitte an den Service von Berthold Technologies wenden.</i>
79	Massenstrom 1 Be- rechn. inaktiv. Pro- duktdichte ist ungültig	Bitte an den Service von Berthold Technologies wenden.
80	Massenstrom 2 Be- rechn. inaktiv. Pro- duktgeschw. ist ungül- tig	<i>Bitte an den Service von Berthold Technologies wenden.</i>
81	Massenstrom 2 Be- rechn. inaktiv. Pro- duktdichte ist ungültig	Bitte an den Service von Berthold Technologies wenden.

## 10.8 Fehler Bereich Analogausgang

Code	Fehler	Mögliche Ursache/ Behebung
90	<i>Stromausgang 1 außer Bereich</i>	<i>Der aus der Konzentration abgebildete Strom liegt außerhalb des Strombe- reichs.</i>
91	<i>Stromausgang 2 außer Bereich</i>	Der aus der Konzentration abgebildete Strom liegt außerhalb des Strombe- reichs.

## 10.9 Watchdog-Fehler

Code	Fehler	Mögliche Ursache/ Behebung
92	Watchdog	Bitte an den Service von Berthold Technologies
		wenden.



## **10.10** Systemfehler

Code	Fehler	Mögliche Ursache/ Behebung
120	Datum/Uhrzeit nicht eingestellt	Bitte das Datum und die Uhrzeit eingeben.

### **10.11** Dichtefehler

Code	Fehler	Mögliche Ursache/ Behebung
150	<i>Dichte Berechnung: FG (Radiometrie) außer Bereich</i>	Den Messwert der radio- metrische FG überprüfen
151	<i>Dichte Berechnung: Höhensignal außer Be- reich</i>	Den Messwert des Schichtdickensensors überprüfen

# 10.12 Eingabefehler

Fehler	Ursache	
Wert zu groß	Eingabewert ist zu groß	
Wert zu klein	Eingabewert ist zu klein	
Tabelle ist leer	Probentabelle wurde angewählt ohne	
	vorherige Probenaufnahme	
Chart Daten	Während der Kalibrierung hat das	
fehlerhaft	Gerät fehlerhafte Chartdaten ermittelt.	
Keine Chartda-	Die berechneten Diagrammdaten wurden	
ten vorhanden	gelöscht oder Kalibrierung wurde nicht	
	beendet.	
Probentabelle	Es wurde versucht mehr als 30 Proben	
voll	aufzunehmen.	

### 10.13 Gerätezustände

#### Fehlerzustand:

Dieser Zustand tritt u. a. unter den Fehlercodes 50 bis 56, 60 bis 62 und 70 bis 71 auf (siehe obige Tabelle). Die AWE zeigt folgendes Verhalten: <u>LED's</u>: RUN blinkt, ERROR leuchtet, Signal 1

	und 2 je nach Konfiguration.
Stromausgänge:	Fehlerstrom, wie ausgewählt
Display:	Fehlermeldung mit Fehlercodeangabe

#### Warnzustand:

Dieser Zustand tritt u. a. unter den Fehlercodes 14, 21, 90 und 91 auf (siehe obige Tabellen). Die AWE zeigt folgendes Verhalten: LED's: RUN blinkt, ERROR aus, Signal 1 und 2

	kein Zusammenhang.
Stromausgänge:	live
Display:	Fehlermeldung mit Fehlercodeangabe

#### Haltezustand:

Messung wurde über Digitaleingang angehalten. Die AWE zeigt folgendes Verhalten:

Der gemittelte Konzentrationswert wird eingefroren. Die Messung läuft jedoch weiter, sodass eine Messstörung den Fehlerzustand auch aus dem Haltezustand auslösen kann.

<u>LED's</u> :	RUN blinkt, ERROR aus, Signal 1 und 2
	je nach Konfiguration.
Stromausgänge:	eingefroren
Display:	keine Displaymeldung

#### Offline-Zustand:

Die AWE befindet sich im Offline-Zustand, es ist z.B. kein Produkt vorhanden. Die AWE zeigt folgendes Verhalten: LED's: RUN blinkt, ERROR aus, Signal 1 und 2

	je nach Konfiguration.
<u>Stromausgänge</u> :	unterer Stromausgangswert 0/4 mA
Display:	Fehlermeldung mit Code-Nr. 053 An-
	gabe



# Kapitel 11. Inbetriebnahme-Protokoll

Das Protokoll kann über RS232 und RS485 ausgegeben werden. Die Ausgabe erfolgt unter Menü | DIAGNOSE | PROTO-KOLL DRÜCKEN |.

Die serielle Schnittstelle RS232 und RS485 besitzt folgende Anschlusseinstellungen:

Datenübertragungsrate 38400 Bd, 8 Datenbits, Keine Parität, 1 Stoppbit

Das Protokoll wird über Terminalprogramm in eine TXT-Datei gespeichert. Für die Darstellung (z.B. mit Excel<sup>®</sup>) muss folgendes Datenformat berücksichtigt werden.

Trennzeichen: Tabulatoren Dezimaltrennzeichen: . 1000er Trennzeichen: ,

Die nachfolgende Codeliste dient zur Interpretation des Inbetriebnahme-Protokolls, siehe Beispiel eines Protokolls unter *Kapitel 12.1*.

Parameter	Code-	Information
	Nr.	
Log type		Logart:
	0	Deaktiviert
	1	Einzeln
	2	Kontinuierlich
	3	Stopp bei Fehler
Log time		Logzeit:
-	0	15 Minuten
	1	1 Stunde
	2	4 Stunden
	3	8 Stunden
	4	1 Tag
	5	3 Tage
Measuring		Messmodus:
mode	0	Kontinuierlich
	1	Batch
Start mode		Startmodus (Start/Stopp):
	0	Tastatur
	1	Extern

Parameter	Code- Nr.	Information
Calibration		Erw. Kal. Eingangsauswahl:
input	0	Kein
selection	1	Eing. 1
	2	Eing. 1 + Eing. 2
	3	Eing. $1 + PT100$
	4	Eing. 1 + Eing. 2 + PT100
	5	Eing. 2
	6	Eing. 2 + PT100
	7	PT100
Calibration		Kalibrierungsbasis:
variable	0	Phase
	1	Dämpfung
	2	Phase und Dämpfung
Massflow		Durchsatzberechnung:
calculation	0	Aus
mode	1	Ein
Measure		Prozessart:
configuration	0	1 Konzentration
	1	2 Konzentrationen
	2	Split Konzentration
AO Assign		Zuordnung Stromausgang:
Code	0	Keine
	1	Konzentration
	2	Konzentration 2
	3	Stromeingang 1
	4	Stromeingang 2
	5	PT100
AO Alarm se-		Fehlerstrom Stromausgang:
lect code	0	22 mA
	1	3.5 mA
	2	Halten
_	3	Wert
Range		Messbereich Stromausgang:
selection	0	0 20 mA
	1	4 20 mA
AI Range		Messbereich Stromeingang:
selection	0	0 20 mA
	1	4 20 mA

	F;	2	7	-1-	-/	С	)/		$\overline{}$	)		
L		ì	c	H	N	$\tilde{}$	L	-0	G	1	E	s

Parameter	Code-	Information
	Nr.	Z tanduar Champingang 2
AI ENAD-		Zustand von Stromeingang 2
Iea[2]		Eveltien der Disitelen Aussänne.
DO Function		Funktion der Digitalen Ausgange:
	0	Keine
	1	Fehler
	2	Halt
	3	Kein Produkt
	4	Alarm min.
	5	Alarm max.
DO Assign-		Digitale Ausgang: der min./max.
ment		Alarm ist folgendem zugewiesen:
	0	Konzentration
	1	Konzentration 2
	2	Stromeingang 1
	3	Stromeingang 2
	4	PT100
DI Function		Funktion der Digitalen Eingänge:
selection	0	Keine
	1	Start/Stopp
	2	Halt
	3	Probenahme
	4	Produktauswahl
Printout		Form der Datenausgabe:
mode	0	Keine
	1	Zeile
	2	Tabelle
	3	Zeile + <u>Tabelle</u>
Access level		Zugriffsebene:
	0	Nur lesen
	1	Benutzer Modus
	2	Service
Language		Sprachauswahl:
	0	Englisch
	1	Deutsch
	2	Französisch



# **11.1** Beispiel Inbetriebnahme-Protokoll

Menü:	Start of Setup:	Inbetriebnal	nme-Pr	otokol	Interpretation:					
			_			(* nur Service relevant)				
Produkt	Entry	Product1	Pro- duct2	Prod. 3	Prod. 4					
Datenlog	Log type :	0				Logart: siehe Codeliste				
Ū	Log time :	2				Logzeit: siehe Codeliste				
	Number of errors :	0				Anzahl der Einträge im Fehlerlog				
	NTC temperature :	45.3 °C				*				
	max. NTC temperature :	46.7 °C				*				
	9V power supply :	8.94 V				*				
Info	Tag :					Messstelle				
	Device type :	LB 566				Gerätetyn				
	Unique device ID number	761				Ceratetyp				
	Serial number :	1204067000								
	Final assembly number :	4294907000								
	Software version	2 00								
	Software release date	02.02.2016				Software Revisionsdatum				
	Actual date :	10.02.2016				Datum der Aufnahme				
	Actual time :	12.15				Uhrzeit der Aufnahme				
Messung	Measuring mode :	0				Messmodus: siehe Codeliste				
0	Start mode :	0				Startmodus: siehe Codeliste				
	Filter damping value[2] :	20				Mittelungszahl Stromausgang				
	Filter damping value[3] :	20				Mittelungszahl bei Probenahme				
	Reset average :	FALSE				Mittelwert Reset: Ja/Nein				
Plausibilität	Lower limit :	0.00				Min. Prozessgrenze				
	Upper limit :	100.00				Max. Prozessgrenze				
	Raw data average value	5				*				
	Max. phase sigma :	100				Sigma max.				
	Correlation Phi/Att :	1.00				Verhältnis Phase/Dämpfung				
	Auto-set mode :	FALSE				Auto-Einstellung: Ein/Aus				
	Offline Mode detection :	FALSE				Offline-Funktion: Ein/Aus				
	Minimum attenuation :	-15 dB				Grenzwert für den Offline Modus				
Mikrowelle	Ref. cable length :	8.00 m				Referenzkabellänge				
	Meas. cable length :	8.00 m				Messkabellänge				
	Wave band selection :	1				*				
	Start frequency :	2				*				
	Frequency step :	1				*				
	Nbr of freq. points :	5				*				
Marker	Marker name :	Mark1				Markername für Konzentration				
	Marker value :	75.00 %				Markerwert für Konzentration				
	Marker name[2] :	Mark2				Markername für Konzentration 2				
Quatara -h	Marker value[2] :	75.00 %				Markerwert für Konzentration 2				
system-ab- gleich	Nbr of sweeps for reference :	1				*				
AWE Typ	HF amplifier mode :	0				*				
	Minimal insertion loss :	40.00 dB				*				



			-
Kalibriere	Calibration input selection :	0	Erw. Kal. Eingangsauswahl: siehe Codeliste
tration	Calibration mode	0	*
	Calibration variable :	0	Kalibrierungsbasis: siehe Codeliste
	Phase coefficients :	0	Phasen-Koeffizient A
	Attenuation coefficients :	0	Dämpf. Koeffizient B
	Constant coefficient :	0	Konstante C
	d coefficient	0	KompKoeffizient für PT100 Eingang
	e coefficient	0	KompKoeffizient für Stromeingang 1
	f coefficient	0	KompKoeffizient für Stromeingang 2
	a coefficient	0	KompKoeffizient für die Beladung
	Adjust factor :	1	Faktor
	Adjust offset :	0	Offset
	Massflow calculation mode:	0	Durchsatzberechnung: siehe Codeliste
	Loading comp Selection	0	*
	Loading comp. Lower limit:	0	*
Synchron-			
isation	Synchronizer mode:	0	*
	Current input1 distance to uWave	0	*
Kalibriere	Current input1 distance to uWave	0	*
Konzen-	Calibration input selection :	0	Erw. Kal. Eingangsauswahl: siehe Codeliste
tration 2	Lower limit :	0.00 %	Min. Prozessgrenze (Plausibilität für Konz 2)
	Upper limit :	100.00 %	Max. Prozessgrenze (Plausibilität für Konz 2)
	Calibration mode :	0	*
	Calibration variable :	0	Kalibrierungsbasis: siehe Codeliste
	Phase coefficients :	0	Phasen-Koeffizient A
	Attenuation coefficients :	0	Dämpf. Koeffizient B
	Constant coefficient :	0	Konstante C
	d coefficient	0	KompKoeffizient für PT100 Eingang
	e coefficient	0	KompKoeffizient für Stromeingang 1
	f coefficient	0	KompKoeffizient für Stromeingang 2
	g coefficient	0	KompKoeffizient die Beladung
	Adjust factor :	1	Faktor
	Adjust offset :	0	Offset
	Massflow calculation mode:	0	Durchsatzberechnung: siehe Codeliste
	Loading comp. Selection:	0	*
	Loading comp. Lower limit:	0	*
Extras	Tara Phase (°/GHz) :	0.00 °/GHz	
	Tara Attenuation (dB) :	0.00 dB	
	Measure configuration :	0	Prozessart: siehe Codeliste
	Range split value :	75	Splitwert
Stromaus-	AO Assign code :	0	Zuordnung: siehe Codeliste
gang i	AO Upper range value :	95	Obere Grenze
	AO Lower range value :	60	Untere Grenze
	AO Alarm select code :	2	Fehlerstrom: siehe Codeliste
	AO Error current value :	22.00 mA	 Fehlerstrom-Wert
Stromaus-	AO Assign code[2] :	0	 Zuordnung: siehe Codeliste
yang z	AO Upper range value[2] :	95	Obere Grenze
	AO Lower range value[2] :	60	Untere Grenze
	Range selection[2] :	1	Bereich
	AO Alarm select code[2] :	2	Fehlerstrom: siehe Codeliste
	AO Error current value[2] :	22.00 mA	Fehlerstrom-Wert
	<u> </u>		

### Kapitel 11. Inbetriebnahme-Protokoll



Stromein-	AI Enabled :	0	Deaktiviert: 0 Aktiviert: 1
55	AI Range selection :	1	Bereich: siehe Codeliste
	Al Upper range value :	100	Obere Grenze
	AI Lower range value :	0	Untere Grenze
	Analog input filter constant :	5	*
Stromein-	AI Enabled[2] :	0	Deaktiviert: 0 Aktiviert: 1
gang z	AI Range selection[2] :	1	Bereich: siehe Codeliste
	AI Upper range value[2] :	100	Obere Grenze
	AI Lower range value[2] :	0	Untere Grenze
	Analog input filter constant :	5	*
PT100 Ein- gang	AI Enabled[3] :	0	Deaktiviert: 0 Aktiviert: 1
Relais 1	DO Function :	1	Funktion: siehe Codeliste
	DO Assignment :	0	Zuweisung: siehe Codeliste
	DO Threshold :	0.00%	*
	DO Hysteresis :	5.00%	*
Relais 2	DO Function[2] :	2	Funktion: siehe Codeliste
	DO Assignment[2] :	0	Zuweisung: siehe Codeliste
	DO Threshold[2] :	0.00%	*
	DO Hysteresis[2] :	5.00%	*
Digital-ein- gänge	DI Function selection :	0	Funktion Digitaleingang 1: siehe Codeliste
33-	DI Function selection[2] :	0	Funktion Digitaleingang 2: siehe Codeliste
	DI Function selection[3] :	0	Funktion Digitaleingang 3: siehe Codeliste
	Printout mode :	1	Datenausgabe: siehe Codeliste
	Access level :	2	Zugriffsebene: siehe Codeliste
	Language :	1	Sprache: siehe Codeliste
	End of Setup		Ende

Start of Reference Data						Daten des Systemabgleichs:								
Product 1:														
Mean Atten.:		46.8509 dB												
Phase at fm:		42.6285 deg	/GHz											
Phase slope:		380.984 deg	/GHz											
Phase sigma:		0.24575												
Frequency[GHz]		Phase[Deg]	Transformierte Phase[Deg]	Atten.[dB]										
	2.42	35.64	35.64	21.98										
	2.43	361.81	361.81	21.95										
	2.44	689.04	689.04	22.07										
	2.45	1014.44	1014.44	22.36										
	2.46	1339.01	1339.01	22.37										
Start of Sample Data:							Probe	ent	abelle:					
Product 1: Sample Data for	Concenti	ration 1:												
Sample:	Active:	Kon.(%):	Lab.(%):	AIN1:	AIN2:		Temp. (°C):		Phi. (°/GHz):	Att.(dB):				
1 17.08 - 12:37	TRUE	85	40	0		0		0	-0.35	0.02				
2 17.08 - 12:37	TRUE	80	35	0		0		0	30.33	5.08				
3 17.08 - 12:45	TRUE	70	25	0		0		0	59.02	18.98				
Correlation factor between														
lab and meas values.		1												
End of Sample Data		1												
Do not use following data!														

## **11.2** Probentabelle

Nr.	Aktiv	Messwert	Laborwert	Stromeing. 1	Stromeing. 2	PT100	Phi(fm)	Dämpfung	
1									
2									
3									
4									
5									
6									
7									
8									
9									
10									
11									
12									
13									
14									
15									
16									
17									
18									
19									
20									

# Notizen



-		 	-			 		 	 	 	_	 	 	 	 				
<u> </u>		 		 		 	 	 	 	 	_	_							
		 				 	 		 					 				_	
		 							 				 			 		_	
		 				 			 				 		 	 		_	
		 							 				 	 		 		_	
<u> </u>		 					 	 	 			 	 	 	 	 			
<u> </u>		 		 		 	 	 	 	 			 	 	 	 		_	
BERTHOLD ...

## Notizen



	 			 	 		 	 	 	 			 	 		 	 -	_	
	 			 	 		 	 	 	 			 	 		 	 _	_	
								 	 										_
		_					 	 	 _	 	_				_	 			_
<u> </u>				 	 		 	 	 	 						 			_
				 	 		 	 	 _	 	 			_		 		_	_
<u> </u>	 			 	 	 	 	 	 	 			 	 		 	 		
																			_
									_		_				_				
									 										_
		_	_				 				_				_				
							 		 										_
		_									_			_	_				