

# DuoXpert LB 473 Grenzstand

## Technische Information



# Inhaltsverzeichnis

1.	<b>Informationen zur 2-Leiter Technologie.....</b>	<b>2</b>
2.	<b>Auswerteeinheit.....</b>	<b>3</b>
2.1.	Installationsvarianten Wandgehäuse.....	3
2.2.	Einbauvarianten 19" Baugruppenträger.....	4
2.3.	Anschlussplan Anschlussplatine Master.....	5
2.4.	Belegung Klemmenblock AWE.....	7
3.	<b>Wandgehäuse .....</b>	<b>8</b>
4.	<b>19" Baugruppenträger.....</b>	<b>9</b>
5.	<b>AWE .....</b>	<b>10</b>
6.	<b>Nummernschlüssel LB 47x .....</b>	<b>13</b>
7.	<b>Konformitätserklärung LB 47x.....</b>	<b>14</b>
8.	<b>NRTL Zertifikat US/CAN Wandgehäuse .....</b>	<b>16</b>
9.	<b>NRTL Zertifikat US/CAN DuoXpert LB 47x.....</b>	<b>19</b>
10.	<b>Zubehör.....</b>	<b>21</b>

# 1. Informationen zur 2-Leiter Technologie

Das DuoSeries / DuoXpert Messsystem besteht aus einem Detektor mit Szintillatortechnologie sowie einer separaten Auswerteeinheit zur Anzeige und Bedienung.

Die moderne Auswerteeinheit verfügt über ein 3,5" Touchscreen, eine starke Dual Core CPU und verschiedenen Bedien-Optionen. Erweiterte Funktionen zur Selbstdiagnose und Überwachung sorgen zudem für höchste funktionale Sicherheit der Messung im Betrieb. Darüber hinaus können die Betreiber die Datenlog Funktionen für eine detaillierte Prozessanalyse nutzen und so zum Beispiel Trends entwickeln oder Prozessänderungen nachvollziehen.

## Hochentwickeltes Messsystem in 2-Leiter Technologie

- Echte 2-Leiter Technik, nur 2 Adern im Feld
- Verbesserte Diagnosefunktionen und Selbstüberwachung
- Einfache, intuitive Bedienung über Touchscreen
- Kompatibel zu allen 2-Leiter Detektoren LB 4700, LB 44xx und LB 54xx

## 2. Auswerteeinheit

Die Module können entweder in Wandgehäusen oder 19"-Baugruppenträgern eingebaut und kann je nach Bedarf unterschiedlich bestückt werden. Zum elektrischen Anschluss werden die rückwärtigen Klemmenblöcke verwendet.

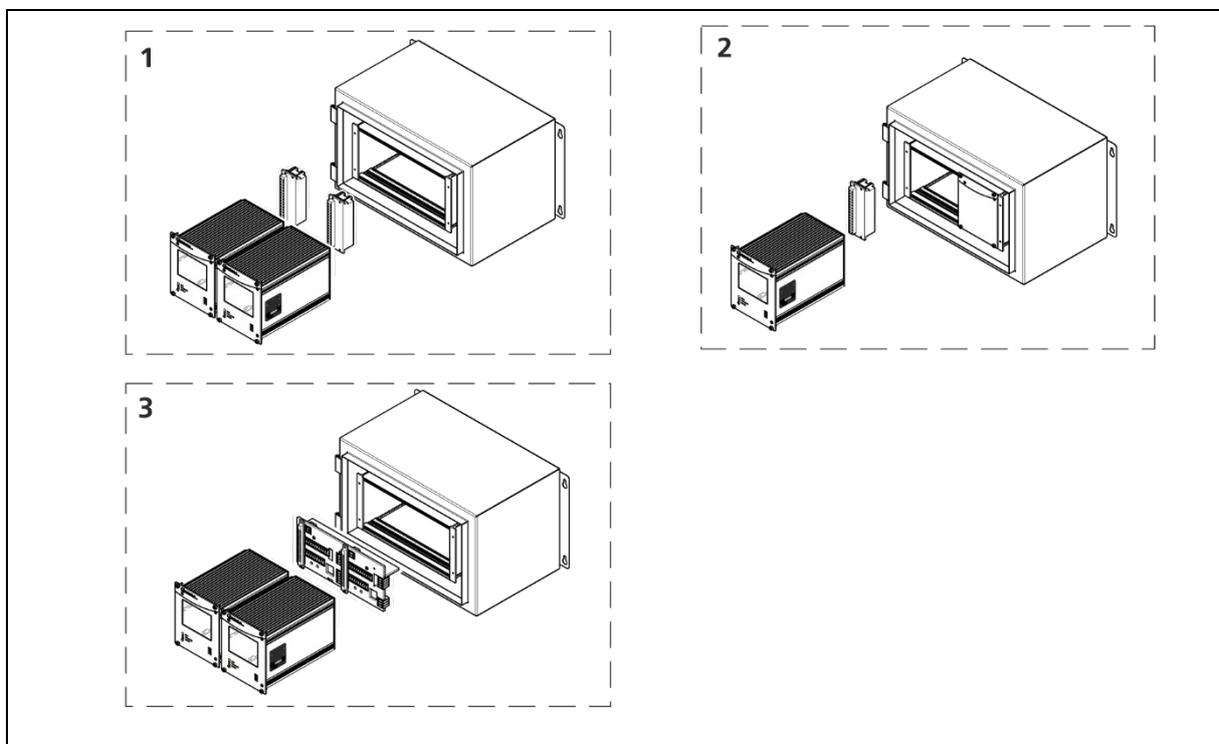
### **⚠️ WARNUNG**



#### Lebensgefahr durch Explosion!

- ▶ Diese Ausführung der Auswerteeinheit ist nicht ex-geschützt ausgeführt und darf nicht in explosionsgefährdete Bereiche verwendet werden.
- ▶ Beachten Sie für Ausführungen der Auswerteeinheit für den eigensicheren Betrieb die mitgeltenden Dokumente, siehe Betriebsanleitung Kapitel 1.1 Mitgeltende Dokumente.

### 2.1. Installationsvarianten Wandgehäuse

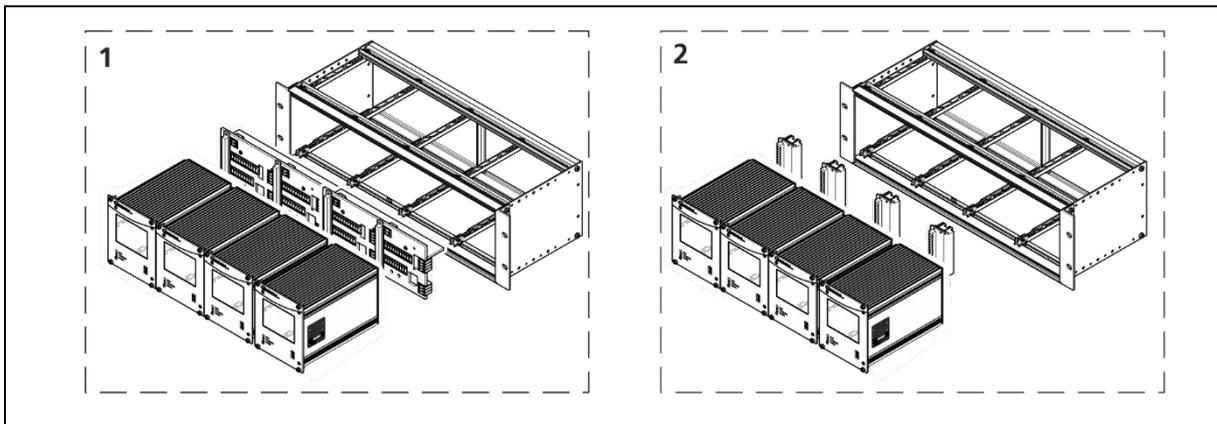


TI-Abb. 1 Installationsvarianten Wandgehäuse

Pos.	Komponenten	Anschluss
1	2 Masters	2 Klemmenblöcke
2	1 Master	1 Klemmenblock
3	2 Master	1 Anschlussplatine Master/Master <sup>1</sup>

<sup>1</sup> NTRL Zertifikat US/CAN

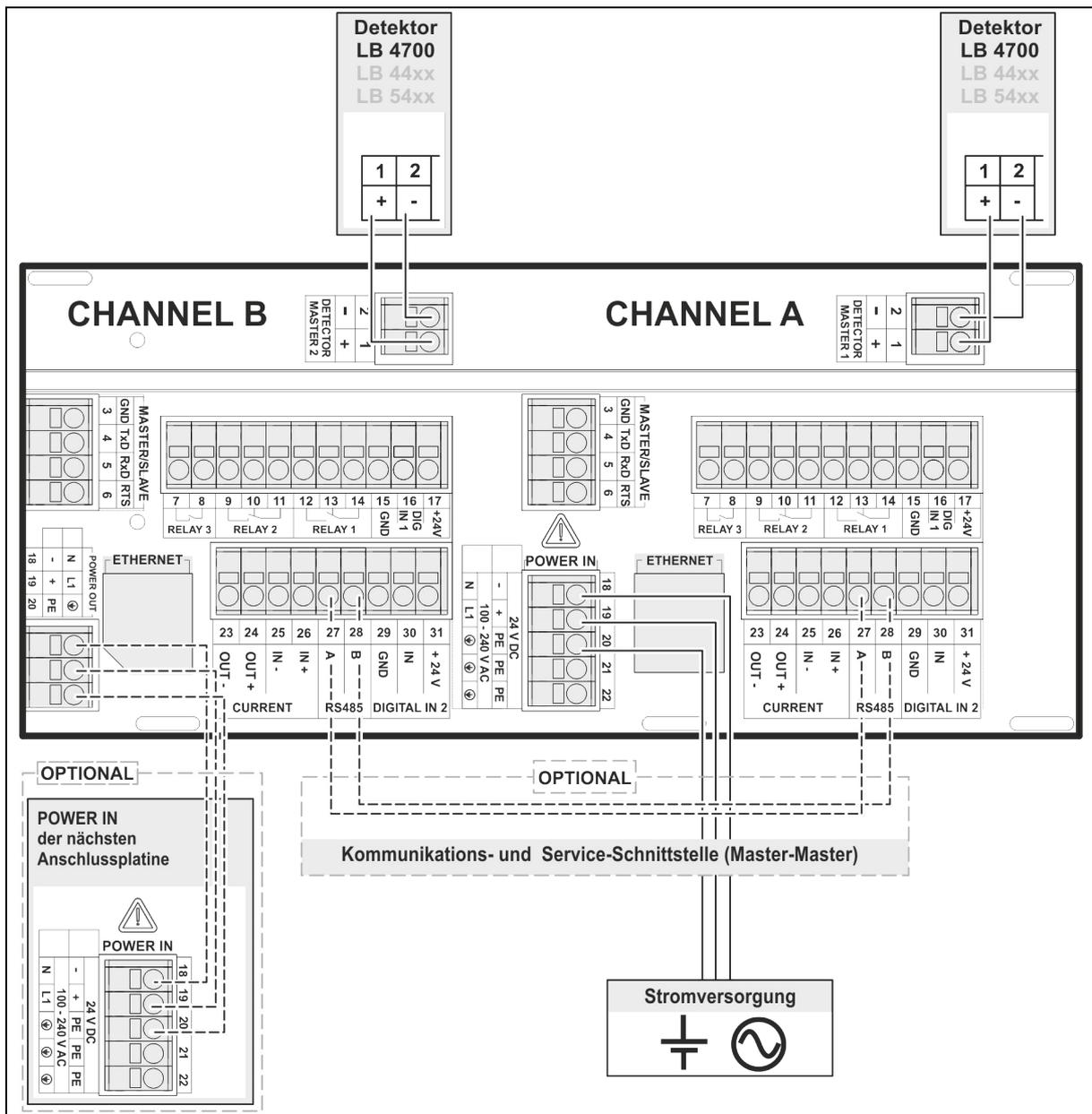
## 2.2. Einbauvarianten 19" Baugruppenträger



TI-Abb. 2 Installationsvarianten 19" Baugruppenträger

Pos.	Komponenten	Anschluss
1	4 Masters	2 Anschlussplatine Master
2	4 Masters	4 Klemmenblöcke

### 2.3. Anschlussplan Anschlussplatine Master



TI-Abb. 3 Anschlussplan Anschlussplatine Master

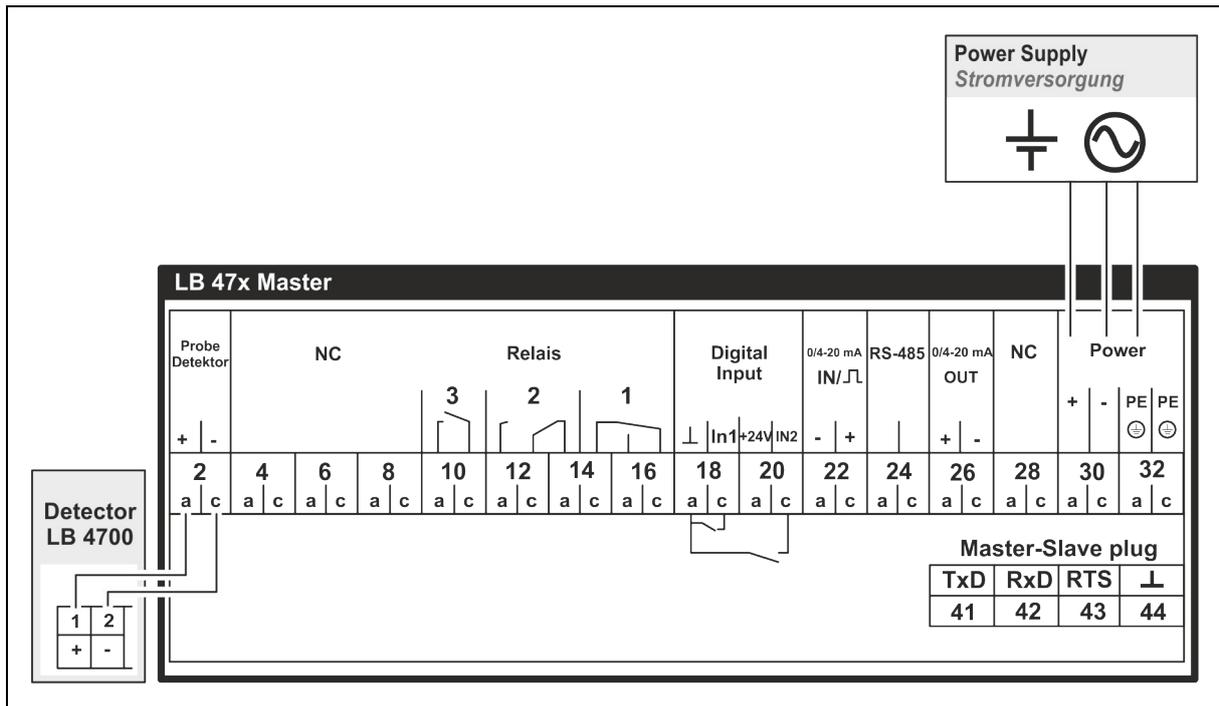
**WICHTIG**

In einem 19" Baugruppenträger für 4 Master (Mat. Nr. 59484), ist neben der oben gezeigten Anschlussplatine für Channel A/B, eine weitere mit Channel C/D enthalten. Die Klemmenbelegung von Channel C/D ist identisch mit der von Channel A/B.

**Klemmen Anschlussplatine Master/Master**

#	Anschluss	Funktion
1	DETECTOR MASTER +	Verbindung Detektor LB 4700 / LB 44xx / LB 54xx
2	DETECTOR MASTER -	
3	MASTER/SLAVE GND	Anschluss von weiteren Slave-Einheiten
4	MASTER/SLAVE TxD	
5	MASTER/SLAVE RxD	
6	MASTER/SLAVE RTS	
7	RELAIS 3 NC	DIGITAL OUT
8	RELAIS 3 COM	
9	RELAIS 2 NC	DIGITAL OUT
10	RELAIS 2 NO	
11	RELAIS 2 COM	
12	RELAIS 1 NC	Fehler DIGITAL OUT
13	RELAIS 1 NO	
14	RELAIS 1 COM	
15	DIGITAL IN 1 GND	GND
16	DIGITAL IN 1 IN	Logic Input
17	+ 24 V (GND → 15)	24 V out (max. 200 mA)
18	POWER DC 24 V – / AC N	24 V DC / 100-240 V AC
19	POWER DC 24 V + / AC L1	
20	PE	
21	PE	
22	PE	
23	CURRENT OUT –	4 mA ... 20 mA
24	CURRENT OUT +	
25	CURRENT IN –	Keine Verwendung bei LB 473
26	CURRENT IN +	
27	RS 485 A	Kommunikations- und Serviceschnittstelle (Master-Master)
28	RS 485 B	
29	DIGITAL IN 2 GND	GND
30	DIGITAL IN 2 IN	Logic Input
31	+ 24 V (GND → 29)	24 V out (max. 200 mA)

## 2.4. Belegung Klemmenblock AWE



TI-Abb. 4 Belegung Klemmenblock Master AWE

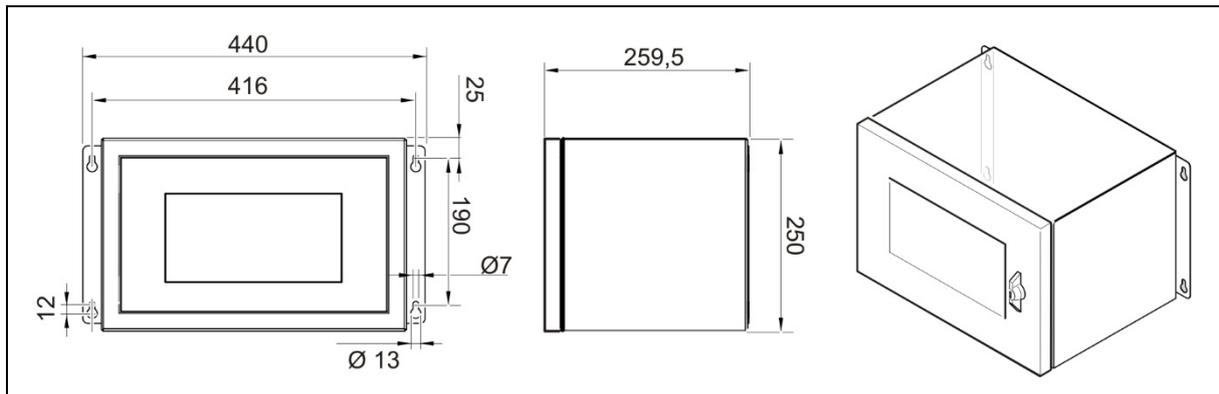
### Klemmenblock

Signal	Pin	Pin	Signal
DETECTOR GND	C - 2	A - 2	DETECTOR +
n.a. *	C - 4	A - 4	n.a. *
n.a. *	C - 6	A - 6	n.a. *
n.a. *	C - 8	A - 8	n.a. *
RELAIS 3 COM	C - 10	A - 10	RELAIS 3 NO
RELAIS 2 COM	C - 12	A - 12	RELAIS 2 NO
RELAIS 1 NC	C - 14	A - 14	RELAIS 2 NC
RELAIS 1 COM	C - 16	A - 16	RELAIS 1 NO
DIGITAL IN 1	C - 18	A - 18	DIGITAL IN GND
DIGITAL IN 2	C - 20	A - 20	+ 24 V (GND → A-18)
CURRENT IN +	C - 22	A - 22	CURRENT IN -
RS 485 B	C - 24	A - 24	RS 485 A
CURRENT OUT -	C - 26	A - 26	CURRENT OUT +
n.a. *	C - 28	A - 28	n.a. *
Netz N AC, DC 24 V (-)	C - 30	A - 30	Netz L1 AC, DC 24 V (+)
PE**	C - 32	A - 32	PE**

\* nicht belegt

\*\* Schutzleiter

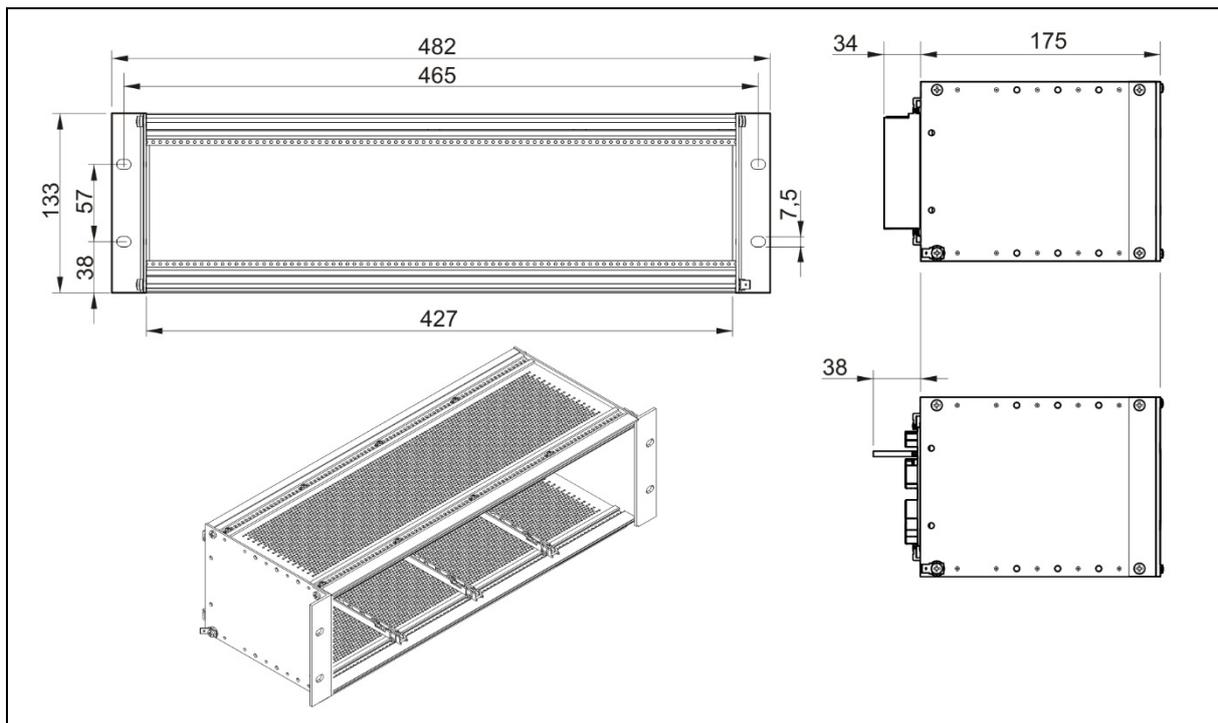
### 3. Wandgehäuse



TI-Abb. 5 Abmessungen Wandgehäuse

Technische Daten	
Max. Bestückung	2 Master
Gewicht (ohne Module)	8,8 kg
Schutzart	IP65
Betriebstemperatur	-20°C ... +40°C
Lagertemperatur	-25°C ... 80°C
Allgemeine Umgebungsbedingungen	Überspannungskategorie: II Verschmutzungsgrad: 2 Höhenlage: bis zu 2000 m Rel. Luftfeuchtigkeit: 93% oder weniger
Oberfläche, Farbe	pulverbeschichtet, grau
Kabeleinführung	8 x M16, 2 x M32

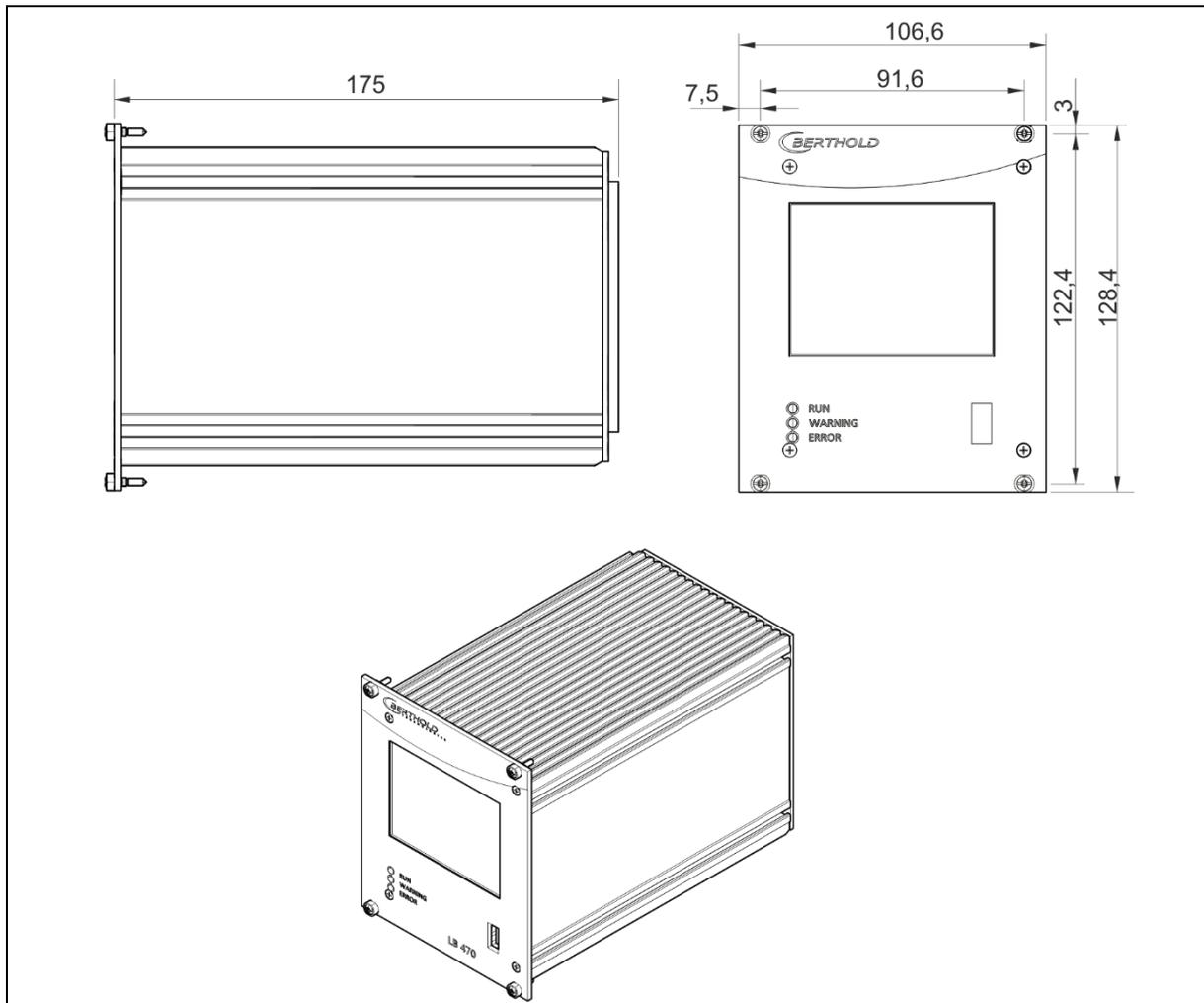
## 4. 19" Baugruppenträger



TI-Abb. 6 Abmessungen 19" Baugruppenträger

Technische Daten	
Normeinheiten	3HE/84TE/5T
Max. Bestückung	4 Master
Gewicht (ohne AWEs)	1,4 kg
Gewicht Klemmenblock	220 g
Betriebstemperatur	-20°C ... +50°C, nicht kondensierend
Lagerungstemperatur	-25°C ... +80°C
Schutzart	IP20

## 5. AWE



TI-Abb. 7 Abmessungen AWE

Technische Daten	
Gewicht	1200 g
Betriebstemperatur	-20°C ... +50°C nicht kondensierend. Direkte Sonneneinstrahlung ist zu vermeiden. Für eine ungehinderte Luftzirkulation ist zu sorgen.
Lagerungstemperatur	-30°C ... +80°C
Allgemeine Umgebungsbedingungen	Überspannungskategorie: II Verschmutzungsgrad: 2 Höhenlage: bis zu 2000 m Rel. Luftfeuchtigkeit: 93% oder weniger
Schutzart	IP20 <sup>1</sup>
Anschlüsse	- USB-Port zum Anschluss von USB-Speichermedium - Master/Slave Buchse (4-polig) und Stecker - RJ45-Buchse für Ethernet (an Rückwand) - 32 polige Stiftleiste nach DIN 19465 Baureihe C

Display	<ul style="list-style-type: none"> <li>- graphisches LCD-Display</li> <li>- 320 x 240 Punkte, 262.000 Farben</li> <li>- Dimmbare LED Hintergrundbeleuchtung</li> <li>- Touchscreen</li> </ul>
Rechnerkern	<ul style="list-style-type: none"> <li>- Prozessor: Dual Core DSP/ARM Controller</li> <li>- Taktfrequenz: 300 MHz intern (20 MHz externer Quarz)</li> <li>- ROM: 512 KByte</li> <li>- RAM: 64 MByte xt. SDRAM, 128 KByte int. shared RAM</li> <li>- FLASH: 8 MByte extern seriell</li> </ul>

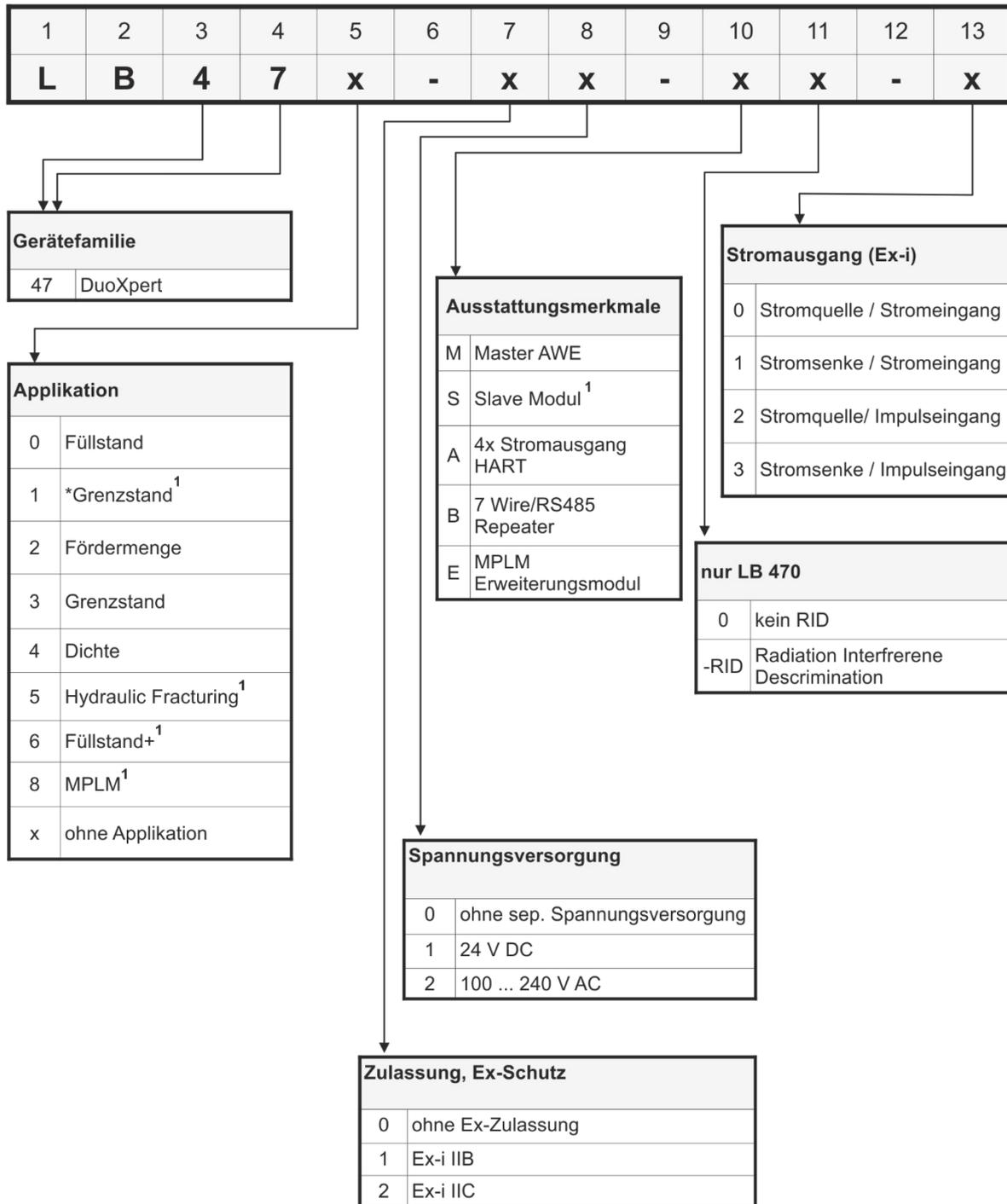
### Stromversorgung

Spannung	100 ... 240 V AC 50/60 Hz (wide range input) +/- 10% 21 ... 32 V DC (24 V DC power input)
Leistungsaufnahme	22 VA, 15 W
Sicherungen	Intern, 2x 250 V, 1A verzögert, 5x20 mm, 1500 A Ausschaltstrom IEC 60127-2, 1x 250 V TR5 T80 mA (Ø 8,5 mm)

<sup>1</sup>Für den Innenbereich konzipiert. Einsatz im Außenbereich nur bei Einbau in Wandgehäuse möglich.

Schnittstellen	
Stromausgang	<p>4-20mA (nach Namur-Empfehlung NE 006 und NE 043) intern von Stromquelle auf Stromsenke umschaltbar. Dip-Schalter auf der Elektronik-Platine in der Auswerteeinheit. Standardeinstellung ist aktiver Stromausgang.</p> <p>Dauerhaft kurzschlussfest und potentialgetrennt (500 V). Innenwiderstand ca. 105 Ohm max. Bürde bei Betrieb als Stromquelle: 850 Ohm. Interne Überwachung des Schleifenstroms und zusätzliche Fehlersignalisierung durch Hardware bei Erkennung eines Fehlerzustands.</p>
Digitale Ausgänge	<p>3 Relais, <math>U_{\max} = 33 \text{ V AC}_{\text{eff}}, 46 \text{ V DC}; I_{\max} = 1 \text{ A}</math></p> <p>Funktionen: Relais 1: SPDT zur Fehlersignalisierung Relais 2: SPDT über Software zuweisbar Relais 3: SPST über Software zuweisbar</p>
Digitale Eingänge	<p>2x gemeinsam potentialgetrennt (500 V), Schalter zwischen DigIn und GND, <math>U_{\text{outmax}}</math> ca. 24 V Funktion über Software konfigurierbar</p>
Externe Versorgung	<p>Ausgangsspannung: 24 V DC Ausgangsstrom: max. 150 mA</p>
RS485	<p>für Prüf-und Testzwecke. Nicht potentialgetrennt von Hauptelektronik und USB-Anschluss potentialgetrennt von restlichen I/Os (500 V)</p>
USB Anschluss	<p>1x USB 2.0 Typ A (Host) über Frontplatte zum Anschluss einer ext. Maus, Tastatur oder Speichermedium <math>U_{\text{out}} = 5 \text{ V}, I_{\text{outmax}} = 0,5 \text{ A}</math></p>
Ethernet	<p>RJ45-Buchse über Rückwand, 10 Mbit, DHCP unterstützt, max. 3 m Vorgesehen für Wartungszwecke. Nicht vorgesehen für den Dauerbetrieb.</p>

## 6. Nummernschlüssel LB 47x



TI-Abb. 8 Nummernschlüssel

<sup>1</sup> Nur verfügbar als Standard-Version

\* Andere Hardware

# 7. Konformitätserklärung LB 47x



BERTHOLD TECHNOLOGIES GmbH & Co. KG  
 Calmbacher Straße 22  
 75323 Bad Wildbad, Germany  
 Phone: +49 7081 177-0  
 Fax: +49 7081 177-100  
 info@Berthold.com  
 www.Berthold.com

**EG-Konformitätserklärung (ORIGINAL)**

Dok.Nr.: CE20028-1

Hiermit erklären wir in alleiniger Verantwortung, dass die Bauart des(r) nachfolgend bezeichneten Geräte / Systems / Anlage / Maschine in der von uns in den Verkehr gebrachten Ausführung den unten genannten einschlägigen Harmonisierungsvorschriften der EU entsprechen.

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

Produktbezeichnung: **radiometrisches Auswertesystem DuoXpert**

Typenbezeichnung / Modell: **LB 47x**

	<b>Richtlinie (Fundstelle)</b>	<b>angewendete Normen und weitere Spezifikationen</b>
NSR	2014/35/EU	EN 61010-1 2010
RoHS	2011/65/EG	
EMV	2014/30/EU	EN 61326-1 2013 EN 61000-4-2 EN 61000-4-3 EN 61000-4-4 EN 61000-4-5 EN 61000-4-6 EN 61000-4-11 EN 61000-3-2 Namur NE21 2012

Diese Erklärung wird verantwortlich für den Hersteller

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

abgegeben durch

Dr. Jürgen Briegleb  
 Leiter Entwicklung  
 Bad Wildbad, den 1. September 2015

Registergericht / Court of Registration Persönlich haftende Gesellschafterin / Fully liable Associates Registergericht / Court of Registration Geschäftsleitung / Management USt.-Id.-Nr. / VAT Reg. No. Deutsche Steuernummer / German Tax No. WEEE-Reg. No.	Stuttgart HRB 330991 BERTHOLD TECHNOLOGIES Verwaltungs-GmbH Stuttgart HRB 331520 Herr: Knäuff, Dr. Dirk Mörmann DE813050511 49038/08038 DE89468690
Sparkasse PF-CW 75323 Bad Wildbad Volksbank 75119 Pforzheim Commerzbank 75105 Pforzheim	Konto/Account No. 8 045 033 (BLZ 655 500 85) Konto/Account No. 957 064 (BLZ 655 900 301) Konto/Account No. 6 511 120 (BLZ 655 800 13)
	SWIFT-BIC PFHS3333 SWIFT-BIC VBP12166 SWIFT-BIC DRES3333
	IBAN: DE37 6565 0085 0038 0450 03 IBAN: DE85 6565 0003 0000 9570 04 IBAN: DE05 6565 0013 0651 1120 00



BERTHOLD TECHNOLOGIES GmbH & Co. KG  
 Calmbacher Straße 22  
 75323 Bad Wildbad, Germany  
 Phone +49 7081 177-0  
 Fax +49 7081 177-100  
 info@berthold.com  
 www.berthold.com

**EG-Declaration of Conformity (ORIGINAL)**

File.No.: CE20028-2

We, hereby declare under our sole responsibility that the design of the following products / systems / units / machines brought into circulation by us comply with the relevant harmonized rules of the EU.

This declaration loses its validity should modifications or unsuitable and improper use take place without our authorisation.

Product name: **radiometric evaluation system  
DuoXpert**

Type / model: **LB 47x**

	<b>directive</b>	<b>applied standards</b>
LVD	2014/35/EU	EN 61010-1 2010
RoHS	2011/65/EG	
EMC	2014/30/EU	EN 61326-1 2013 EN 61000-4-2 EN 61000-4-3 EN 61000-4-4 EN 61000-4-5 EN 61000-4-6 EN 61000-4-11 EN 61000-3-2 Namur NE21 2012

This declaration is issued by the manufacturer

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

released by

  
 Dr. Jürgen Briggmann

Head of R&D  
 Bad Wildbad, 1<sup>st</sup> of September, 2015

Registergericht / Court of Registration  
 Persönlich haftende Gesellschafterin / Fully liable Associates  
 Registergericht / Court of Registration  
 Geschäftsführung / Management  
 USt.-Id.-Nr. / VAT Reg. No.  
 Deutsche Steuernummer / German Tax No.  
 WEEE-Reg. Nr.

Stuttgart HRA 330981  
 BERTHOLD TECHNOLOGIES Verwaltungs-GmbH  
 Stuttgart HRG 331520  
 Horst Knauß, Dr. Dirk Mörmann  
 DE813050511  
 45038/06038  
 DE95468690

SparKasse  -CW	75323 Bad Wildbad	Konto/Account No. 8 043 003 (BLZ 566 500 88)	SWIFT-BIC: PZHSDE33	IBAN: DE37 0660 3085 0008 0450 03
Volksbank	75119 Pforzheim	Konto/Account No. 557 004 (BLZ 666 800 00)	SWIFT-BIC: VBP03333	IBAN: DE85 6666 3000 0000 9520 04
Gemeinschaftsbank	75105 Pforzheim	Konto/Account No. 8 511 120 (BLZ 566 880 12)	SWIFT-BIC: ORESDE33	IBAN: DE05 6668 3013 0651 1120 00

## 8. NRTL Zertifikat US/CAN Wandgehäuse

		<h3>Certificate of Compliance</h3>	
<b>Nemko-CCL, Inc.</b>			
<b>Certificate:</b>	NA201610530	<b>Date Issued:</b>	January 20, 2016
<b>Project:</b>	257087-7.1		
<b>Issued to:</b>	Berthold Technologies GmbH & Co. KG Calmbacher Straße 22 75323 Bad Wildbad Germany		
<i>The products listed below have been certified as being compliant with all applicable requirements of the specifications listed and are eligible to bear the following certification mark</i>			
			
<b>Issued by:</b>		Robert Keller, Senior Engineer/Safety Supervisor	
<b>Authorized by:</b>		Thomas Jackson, Certification Manager	
<b><u>PRODUCTS</u></b>			
MEASUREMENT, CONTROL, OR LABORATORY EQUIPMENT – Certified to US and Canada Standards			
<b>Product:</b> Process measurement unit <b>Model:</b> Wall-mounted LB 47x, 1M/3S; Wall-mounted LB 47x, 2M (x can be 0 to 8 and describes different software versions for the master and slave modules not affecting safety). <b>Ratings:</b> Wall-mounted LB 47x, 1M/3S: 40VA 100-240V, 50/60Hz, Class I; Wall-mounted LB 47x, 2M: 44VA 100-240V, 50/60Hz, Class I			
<small>The certification system, as described in ISO/IEC Guide 67 (Conformity Assessment – Fundamentals of Product Certification), most closely resembles System 3</small>			
<small>Nemko-CCL, Inc. 1940 West Alexander Street Salt Lake City, Utah 84119-2039 Tel (801) 972-6146 Fax (801) 972-8432</small>			
<small>NFCC-002 Issue 2 May 2014</small>			<small>Page 1 of 3</small>

**APPLICABLE REQUIREMENTS**

UL Std. No. 61010-1 2nd Edition - Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements

CAN/CSA-C22.2 No. 61010-1-04 Second Edition - Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements

This certificate is issued on condition that the holder complies and will continue to comply with the requirements of the above mentioned specifications and pursuant to the terms and conditions specified in the Certification Agreement.

*The certification system, as described in ISO/IEC Guide 67 (Conformity Assessment – Fundamentals of Product Certification), most closely resembles System 3*

Nemko-CCL, Inc. 1940 West Alexander Street Salt Lake City, Utah 84119-2039 Tel (801) 972-6146 Fax (801) 972-8432



NFCC-002 Issue 2 May 2014

Page 2 of 3

### *Supplement to Certificate of Compliance*

**Certificate:** NA201610530

**Project:** 257087-7.1

*Nemko-CCL grants a license to the applicant to apply the Certification Mark to the certified products and that the mark shall only be affixed at the following factory locations*

#### **Factory Information**

Factory Name	Location
Berthold Technologies GmbH & Co. KG	Calmbacher Straße 22 75323 Bad Wildbad Germany

*The products listed, including the latest revision described below, are eligible to be marked in accordance with the referenced Certificate.*

#### **Product Certification History**

Project	Date	Description
257087-7.1	January 20, 2016	<b>Original Certification:</b> <b>Model:</b> Wall-mounted LB 47x, 1M/3S; Wall-mounted LB 47x, 2M (x can be 0 to 8 and describes different software versions for the master and slave modules not affecting safety). <b>Ratings:</b> Wall-mounted LB 47x, 1M/3S: 40VA 100-240V, 50/60Hz, Class I; Wall-mounted LB 47x, 2M: 44VA 100-240V, 50/60Hz, Class I

This Supplement forms an integral part of the Certificate of Compliance

The certification system, as described in ISO/IEC Guide 67 (Conformity Assessment – Fundamentals of Product Certification), most closely resembles System 3

Nemko-CCL, Inc. 1940 West Alexander Street Salt Lake City, Utah 84119-2039 Tel (801) 972-6146 Fax (801) 972-8432

NFCC-002 Issue 2 May 2014



Page 3 of 3

## 9. NTRL Zertifikat US/CAN DuoXpert LB 47x

		<h3>Certificate of Compliance</h3>	
<b>Nemko-CCL, Inc.</b>			
<b>Certificate:</b>	NA201510498	<b>Date Issued:</b>	September 17, 2015
<b>Project:</b>	235982-14.1		
<b>Issued to:</b>	Berthold Technologies GmbH & Co. KG Calmbacher Straße 22 75323 Bad Wildbad Germany		
<i>The products listed below have been certified as being compliant with all applicable requirements of the specifications listed and are eligible to bear the following certification mark</i>			
			
<b>Issued by:</b>		<b>Robert Keller, Senior Engineer/Safety Supervisor</b>	
<b>Authorized by:</b>		<b>Thomas Jackson, Certification Manager</b>	
<b><u>PRODUCTS</u></b>			
MEASUREMENT, CONTROL, OR LABORATORY EQUIPMENT – Certified to US and Canada Standards			
<b>Product:</b> Process measurement unit for building-in			
<b>Model:</b> DuoXpert LB47x-02-M; DuoXpert LB47x-02-S (x can be 0 to 8 and describes different software versions for the master and slave modules not affecting safety)			
<b>Ratings:</b> LB47x-02-M: 100-240V AC 22VA 50/60Hz; LB47x-02-S: 100-240V AC 6VA 50/60Hz			
<b><u>APPLICABLE REQUIREMENTS</u></b>			
UL Std.	No. 61010-1 3rd Edition	- Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements	
	CAN/CSA-C22.2 No. 61010-1-12 Third Edition	- Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements	
This certificate is issued on condition that the holder complies and will continue to comply with the requirements of the above mentioned specifications and pursuant to the terms and conditions specified in the Certification Agreement.			
<small>The certification system, as described in ISO/IEC Guide 67 (Conformity Assessment – Fundamentals of Product Certification), most closely resembles System 3</small>			
<small>Nemko-CCL, Inc. 1940 West Alexander Street Salt Lake City, Utah 84119-2039 Tel (801) 972-6146 Fax (801) 972-8432</small>			
<small>NFCC-002 Issue 2 May 2014</small>			 <small>Page 1 of 2</small>

### *Supplement to Certificate of Compliance*

**Certificate:** NA201510498

**Project:** 235982-14.1

*Nemko-CCL grants a license to the applicant to apply the Certification Mark to the certified products and that the mark shall only be affixed at the following factory locations*

#### Factory Information

Factory Name	Location
Berthold Technologies GmbH & Co. KG	Calmbacher Straße 22 75323 Bad Wildbad Germany

*The products listed, including the latest revision described below, are eligible to be marked in accordance with the referenced Certificate.*

#### Product Certification History

Project	Date	Description
235982-14.1	September 17, 2015	<b>Original Certification:</b> Model: DuoXpert LB47x-02-M; DuoXpert LB47x-02-S (x can be 0 to 8 and describes different software versions for the master and slave modules not affecting safety) <b>Ratings:</b> LB47x-02-M: 100-240V AC 22VA 50/60Hz; LB47x-02-S: 100-240V AC 6VA 50/60Hz

This Supplement forms an integral part of the Certificate of Compliance

The certification system, as described in ISO/IEC Guide 67 (Conformity Assessment – Fundamentals of Product Certification), most closely resembles System 3

Nemko-CCL, Inc. 1940 West Alexander Street Salt Lake City, Utah 84119-2039 Tel (801) 972-6146 Fax (801) 972-8432

NFCC-002 Issue 2 May 2014



Page 2 of 2

## 10. Zubehör

Mat.-Nr.	Beschreibung
64653	LB 473-01-M0 Grenzstandmessgerät (Master, 24 VDC)
64652	LB 473-02-M0 Grenzstandmessgerät (Master, 100...240 VAC)
72063	LB 473-21-M0-0 Grenzstandmessgerät Ex-i (Master, 24 VDC, Stromquelle)
72064	LB 473-22-M0-0 Grenzstandmessgerät Ex-i (Master, 100...240 VAC, Stromquelle)
72370	LB 473-21-M0-1 Grenzstandmessgerät Ex-i (Master, 24 VDC, Stromsenke)
72369	LB 473-22-M0-1 Grenzstandmessgerät Ex-i (Master, 100...240 VAC, Stromsenke)
56925-3BA1	Betriebsanleitung DuoSeries LB 473 Grenzstand, Deutsch
56925-3BA2	Betriebsanleitung DuoSeries LB 473 Grenzstand, Englisch
56925BA16	Sicherheitshandbuch / Explosionsschutzhandbuch LB 47x Ex-i (ATEX / IECEx), Deutsch
56925BA26	Sicherheitshandbuch / Explosionsschutzhandbuch LB 47x Ex-i (ATEX / IECEx), Englisch
63783	Wandgehäuse für 2x Master LB 47x (24 VDC)
63784	Wandgehäuse für 2x LB 47x Master (110...240 VAC)
64402	Wandgehäuse für 2x LB 47x Master (Klemmblöcke)
72812	Wandgehäuse für 2x LB 47x Ex-i Master (mit Ex-i Klemmblöcke)
59484	19"-Baugruppenträger für LB 47x, 4x Master
64607	19"-Baugruppenträger, 84 TE / 3 HE für den Einsatz mit Klemmblöcken
72051	19"-Baugruppenträger Ex-i, 84 TE / 3 HE mit 4x Zugfederklemmblöcken
37526	Blindplatte 21TE / 3 HE (Master)