

DuoXpert LB 476 Level+

Technical Information Technische Information



Table of Contents

1.	Evaluation Unit and Extension Modules	3
2.	Connection Diagrams.....	4
2.1.	Assignment Master EVU with LB 4700 Detectors.....	4
2.2.	Assignment Slave Module.....	6
2.3.	Assignment Master EVU with LB 480 Detectors.....	8
2.4.	Assignment 7-Wire/RS485-Repeater Module	10
2.5.	Assignment Current Output/HART module	12
2.6.	Connection Drawing of the Wall Housing	14
2.7.	Assignment Terminals Master/Slave Plug	15
3.	Wall Housing	16
4.	19" Subrack	17
5.	Master EVU	18
6.	Slave Module.....	22
7.	7-Wire/RS485-Repeater Module ¹	24
8.	Current Output/HART Module ¹	26
9.	Number Key LB 47x	28
10.	EG-Declaration of Conformity LB 47x	29
11.	EU-Declaration of Conformity 7-wire/RS-485 repeater	31
12.	EU-Declaration of Conformity 4x current output/HART	32
13.	UK Declaration of Conformity 7-wire/RS-485 repeater	33
14.	UK Declaration of Conformity 4x current output/HART	34
15.	NTRL Certification US/CAN wall-mounted housing.....	35
16.	NTRL Certificate US/CAN DuoXpert LB 47x.....	38
17.	Parts Overview	40

Inhaltsverzeichnis

1.	Auswerteeinheit und Erweiterungsmodule	3
2.	Anschlusspläne	4
2.1.	Belegung Master AWE mit LB 4700 Detektoren	4
2.2.	Belegung Slave Module	6
2.3.	Belegung Master AWE mit LB 480 Detektoren	8
2.4.	Belegung 7-Leiter/RS485-Repeater Modul	10
2.5.	Belegung Stromausgang/HART Modul	12
2.6.	Anschlusszeichnung des Wandgehäuses	14
2.7.	Belegung Master/Slave Stecker	15
3.	Wandgehäuse	16
4.	19" Baugruppenträger	17
5.	Master AWE	18
6.	Slave Modul	22
7.	7-Leiter/RS485-Repeater Modul'	24
8.	Stromausgang/HART Modul'	26
9.	Nummernschlüssel LB 47x	28
10.	EG-Konformitätserklärung LB 47x	30
11.	EU-Konformitätserklärung 7-Leiter/RS-485 Repeater	31
12.	EU-Konformitätserklärung 4x Spannungsausgang/HART	32
13.	UK Konformitätserklärung 7-Leiter/RS-485 Repeater	33
14.	UK Konformitätserklärung 4x Spannungsausgang/HART	34
15.	NRTL Zertifikat US/CAN Wandgehäuse	35
16.	NRTL Zertifikat US/CAN DuoXpert LB 47x	38
17.	Übersicht Zubehör	40

1. Evaluation Unit and Extension Modules

The modules can be installed either in wall housings or 19" subracks. It can be equipped differently, depending on requirements. The rear terminal blocks are used for the electrical connection.

WARNING



Danger to life due to explosion!

- ▶ This version of the evaluation unit is not explosion protected and is not designed for hazardous environments.

1. Auswerteeinheit und Erweiterungsmodule

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

WARNUNG



Lebensgefahr durch Explosion!

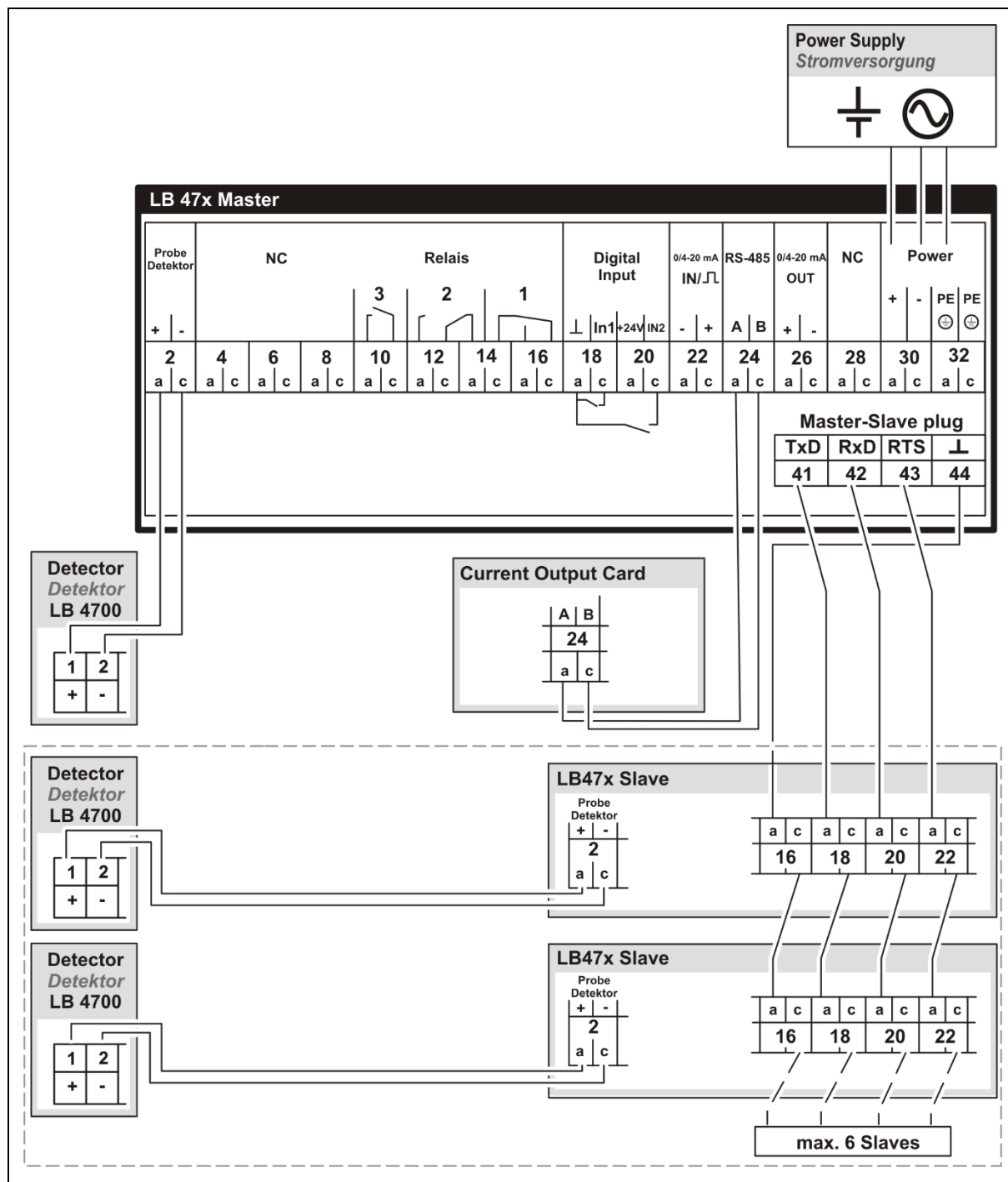
- ▶ Diese Version der Auswerteeinheit ist nicht ex-geschützt ausgeführt und darf nicht in explosionsgefährdete Bereiche verwendet werden.

2. Connection Diagrams

2. Anschlusspläne

2.1. Assignment Master EVU with LB 4700 Detectors

2.1. Belegung Master AWE mit LB 4700 Detektoren



TI-Abb. 1 Assignment master EVU with LB 4700 detectors
Belegung Master AWE mit LB 4700 Detektoren

Terminal Block Master EVU with LB 4700 Detectors

Klemmenblock Master AWE mit LB 4700 Detektoren

Signal	Pin					Pin	Signal
DETECTOR LB 4700 GND	C - 2					A - 2	DETECTOR LB 4700+
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 -
RS485 B (Curr.Output/HART)	C - 24					A - 24	RS485 A (Curr.Output/HART)
CURRENT OUT -	C - 26					A - 26	CURRENT OUT +
n.a. *	C - 28					A - 28	n.a. *
Main <i>Netz</i> N AC, DC 24 V (-)	C - 30					A - 30	Main <i>Netz</i> L1 AC, DC 24 V (+)
PE**	C - 32					A - 32	PE**

* not assigned
nicht belegt

** Protective conductor
Schutzleiter

NOTICE HINWEIS



Since the master EVU and the extension modules (7-Wire/RS485-Repeater module, Current Output/HART module) create a RS485 network, ensure that the last bus participants (first and last) are terminated and thus require a 120 Ω terminating wire resistor between terminal A-24 and C-24.

Da die Master AWE und die Erweiterungsmodule (7-Leiter/RS485-Repeater Modul, Stromausgang/HART Modul) ein RS485-Netzwerk bilden, muss darauf geachtet werden, dass die letzten Busteilnehmer (erster und letzter) terminiert werden und somit einen 120 Ω Abschlussdrahtwiderstand zwischen Klemme A-24 und C-24 benötigen.

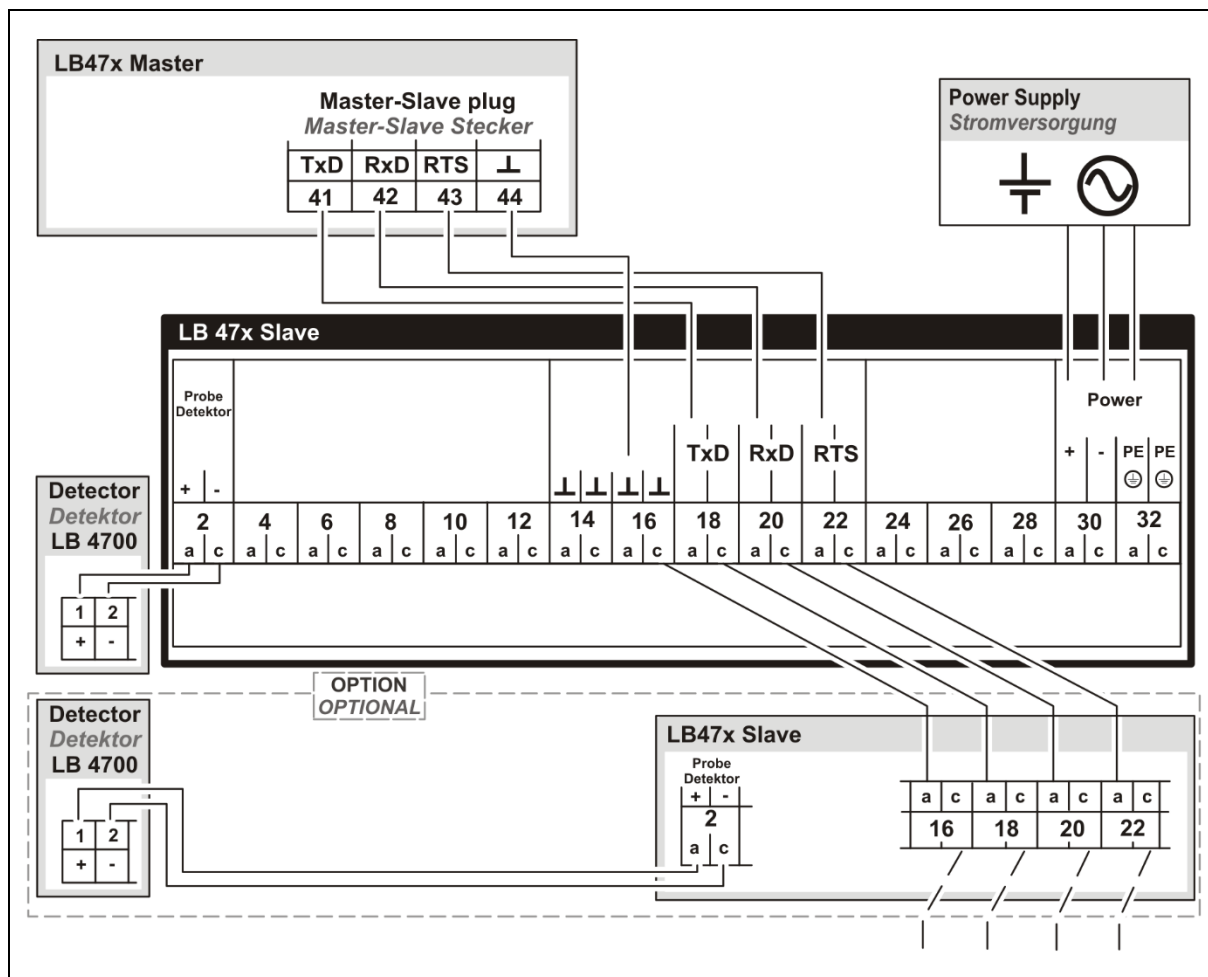
2.2. Assignment Slave Module

2.2. Belegung Slave Module

NOTICE
HINWEIS

Slave modules are only used with LB 4700 detectors.

Slave Module werden nur bei Detektoren des Typs LB 4700 verwendet.



TI-Abb. 2 Assignment slave module

Belegung Slave Modul

Terminal Block Slave Module Klemmenblock Slave Modul

Signal	Pin			Pin	Signal
DETECTOR SLAVE GND	C - 2			A - 2	DETECTOR SLAVE +15 V
n.a**	C - 4			A - 4	n.a**
n.a**	C - 6			A - 6	n.a**
n.a**	C - 8			A - 8	n.a**
n.a**	C - 10			A - 10	n.a**
n.a**	C - 12			A - 12	n.a**
GND	C - 14			A - 14	GND
GND	C - 16			A - 16	GND
TxD to <i>zu</i> SLAVE *	C - 18			A - 18	TxD from <i>von</i> MASTER/SLAVE
RxD to <i>zu</i> SLAVE *	C - 20			A - 20	RxD from <i>von</i> MASTER/SLAVE
RTS to <i>zu</i> SLAVE *	C - 22			A - 22	RTS from <i>von</i> MASTER/SLAVE
n.a**	C - 24			A - 24	n.a**
n.a**	C - 26			A - 26	n.a**
n.a**	C - 28			A - 28	n.a**
Main <i>Netz</i> N AC, DC 24 V (-)	C - 30			A - 30	Main <i>Netz</i> L1 AC, DC 24 V (+)
PE ***	C - 32			A - 32	PE ***

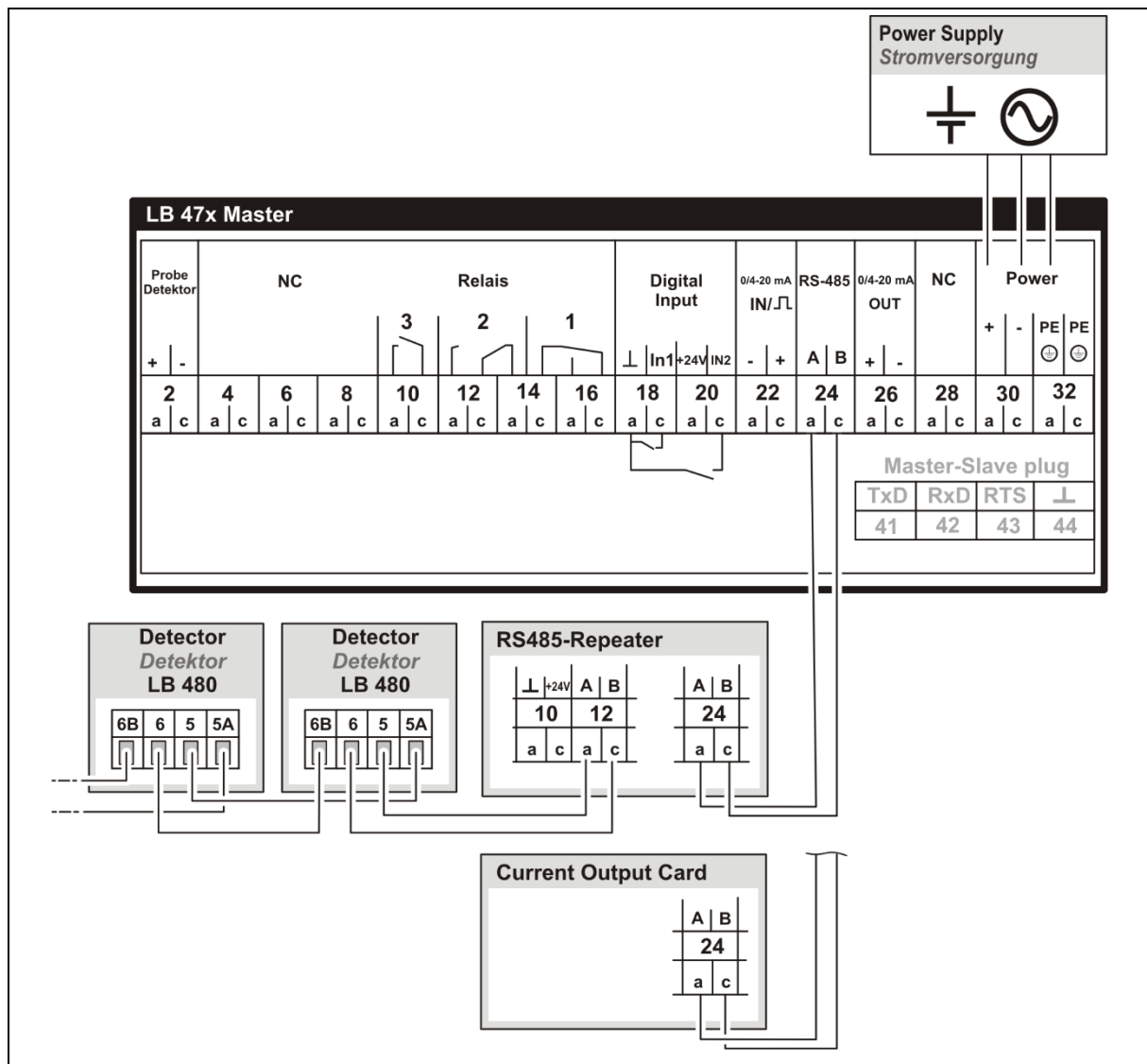
* option
optional

** not assigned
nicht belegt

*** Protective conductor
Schutzleiter

2.3. Assignment Master EVU with LB 480 Detectors

2.3. Belegung Master AWE mit LB 480 Detektoren



TI-Abb. 3 Assignment EVU with LB 480 Detectors

Belegung Master AWE mit LB 480 Detektoren

Terminal Block Master EVU with LB 480

Klemmenblock Master AWE mit LB 480

Signal	Pin					Pin	Signal
n.u. ***	C - 2					A - 2	n.u. ***
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 -
RS485 B ****	C - 24					A - 24	RS485 A ****
CURRENT OUT -	C - 26					A - 26	CURRENT OUT +
n.a. *	C - 28					A - 28	n.a. *
Main <i>Netz</i> N AC, DC 24 V (-)	C - 30					A - 30	Main <i>Netz</i> L1 AC, DC 24 V (+)
PE**	C - 32					A - 32	PE**

* not assigned
nicht belegt

** Protective conductor
Schutzleiter

*** not used for this case
wird für diesen Fall nicht verwendet

**** RS485-Repeater module/Current Output/HART module
RS485-Repeater Module/Stromausgangskarte

NOTICE HINWEIS



Since the master EVU and the extension modules (7-Wire/RS485-Repeater module, Current Output/HART module) create a RS485 network, ensure that the last bus participants (first and last) are terminated and thus require a 120 Ω terminating wire resistor between terminal A-24 and C-24.

Da die Master AWE und die Erweiterungskarten (7-Leiter/RS485-Repeater Modul, Stromausgang/HART Modul) ein RS485-Netzwerk bilden, muss darauf geachtet werden, dass die letzten Busteilnehmer (erster und letzter) terminiert werden und somit einen 120 Ω Abschlussdrahtwiderstand zwischen Klemme A-24 und C-24 benötigen.

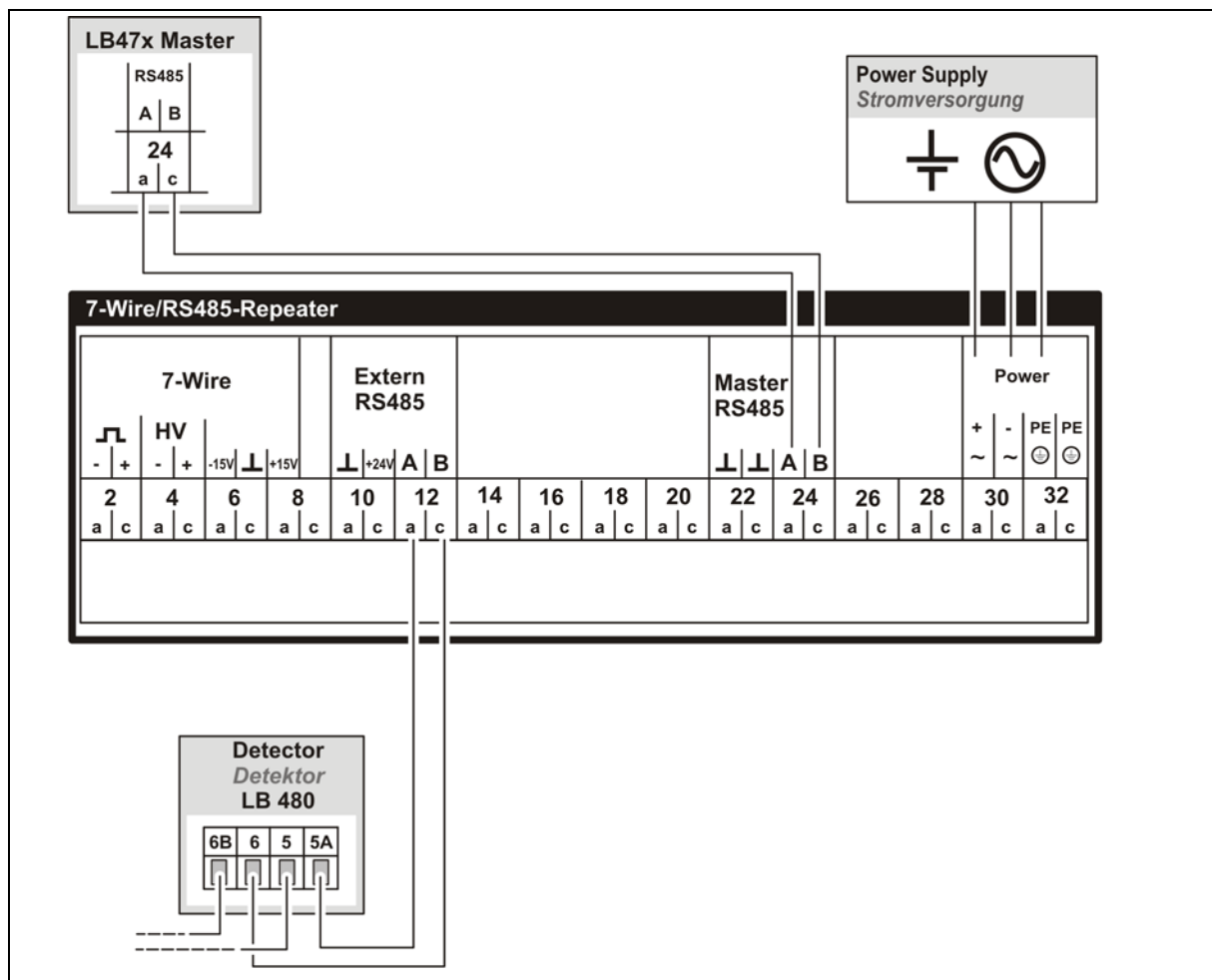
2.4. Assignment 7-Wire/RS485-Repeater Module

2.4. Belegung 7-Leiter/RS485-Repeater Modul

NOTICE
HINWEIS

7-Wire/RS485-Repeater modules are only used with LB 480 detectors.

7-Leiter/RS485-Repeater Module werden nur bei Detektoren des Typs LB 480 verwendet.



TI-Abb. 4 Assignment 7-Wire/RS485-Repeater module
Belegung 7-Leiter/RS485-Repeater Modul

Terminal Block 7-Wire/RS485-Repeater Klemmenblock 7-Leiter/RS485-Repeater

Signal	Pin			Pin	Signal
n.a. *	C - 2			A - 2	n.a. *
n.a. *	C - 4			A - 4	n.a. *
n.a. *	C - 6			A - 6	n.a. *
n.a. *	C - 8			A - 8	n.a. *
+24 V	C - 10			A - 10	GND
RS485 B to Detector LB 480	C - 12			A - 12	RS485 A to Detector LB 480
n.a. *	C - 14			A - 14	n.a. *
n.a. *	C - 16			A - 16	n.a. *
n.a. *	C - 18			A - 18	n.a. *
n.a. *	C - 20			A - 20	n.a. *
GND	C - 22			A - 22	GND
RS485 B to Master	C - 24			A - 24	RS485 A to Master
n.a. *	C - 26			A - 26	n.a. *
n.a. *	C - 28			A - 28	n.a. *
Main <i>Netz</i> N AC, DC 24 V (-)	C - 30			A - 30	Main <i>Netz</i> L1 AC, DC 24 V (+)
PE **	C - 32			A - 32	PE **

* not assigned
nicht belegt

** Protective conductor
Schutzleiter

NOTICE HINWEIS

i Since the master EVU and the extension modules (7-Wire/RS485-Repeater module, Current Output/HART module) create a RS485 network, ensure that the last bus participants (first and last) are terminated and thus require a 120 Ω terminating wire resistor between terminal A-24 and C-24.

Da die Master AWE und die Erweiterungskarten (7-Leiter/RS485-Repeater Modul, Stromausgang/HART Modul) ein RS485-Netzwerk bilden, muss darauf geachtet werden, dass die letzten Busteilnehmer (erster und letzter) terminiert werden und somit einen 120 Ω Abschlussdrahtwiderstand zwischen Klemme A-24 und C-24 benötigen.

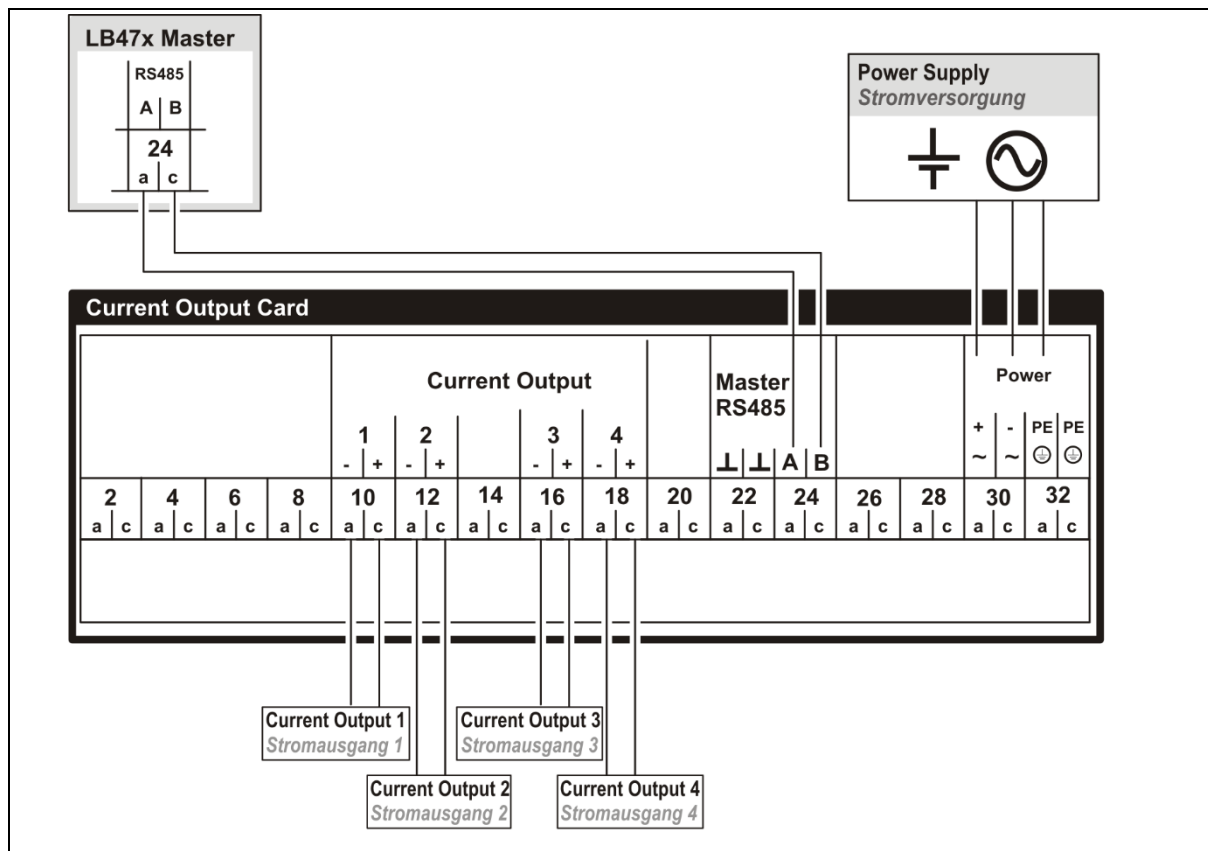
NOTICE HINWEIS

i For the RS485 network between 7-Wire/RS485-Repeater module and LB 480 detector, no termination is necessary as this is already activated internally. It must also be noted that in this case the 7-Wire/RS485-Repeater module is physically placed at the end/beginning of the bus and the LB 480 detector placed at the other end of the bus must have a termination.

Für das RS485-Netzwerk zwischen 7-Leiter/RS485-Repeater Modul und LB 480 Detektor ist kein Abschluss notwendig, da dieser bereits intern aktiviert ist. Außerdem ist darauf zu achten, dass in diesem Falle der 7-Leiter/RS485-Repeater Modul physikalisch am Busende-anfang platziert wird und der am anderen Busende platzierte Detektor LB 480 einen Abschluss haben muss.

2.5. Assignment Current Output/HART module

2.5. Belegung Stromausgang/HART Modul



TI-Abb. 5 Assignment Current Output/HART module
Belegung Stromausgang/HART Modul

Terminal Block Current Output/HART Klemmenblock Stromausgang/HART

Signal	Pin					Pin	Signal
n.a. *	C - 2					A - 2	n.a. *
n.a. *	C - 4					A - 4	n.a. *
n.a. *	C - 6					A - 6	n.a. *
n.a. *	C - 8					A - 8	n.a. *
Current Out 1 +	C - 10					A - 10	Current Out 1 -
Current Out 2 +	C - 12					A - 12	Current Out 2 -
n.a. *	C - 14					A - 14	n.a. *
Current Out 3 +	C - 16					A - 16	Current Out 3 -
Current Out 4 +	C - 18					A - 18	Current Out 4 -
n.a. *	C - 20					A - 20	n.a. *
GND	C - 22					A - 22	GND
RS485 B to Master	C - 24					A - 24	RS485 A to Master
n.a. *	C - 26					A - 26	n.a. *
n.a. *	C - 28					A - 28	n.a. *
Main <i>Netz</i> N AC, DC 24 V (-)	C - 30					A - 30	Main <i>Netz</i> L1 AC, DC 24 V (+)
PE **	C - 32					A - 32	PE **

* not assigned
nicht belegt

** Protective conductor
Schutzleiter

NOTICE HINWEIS

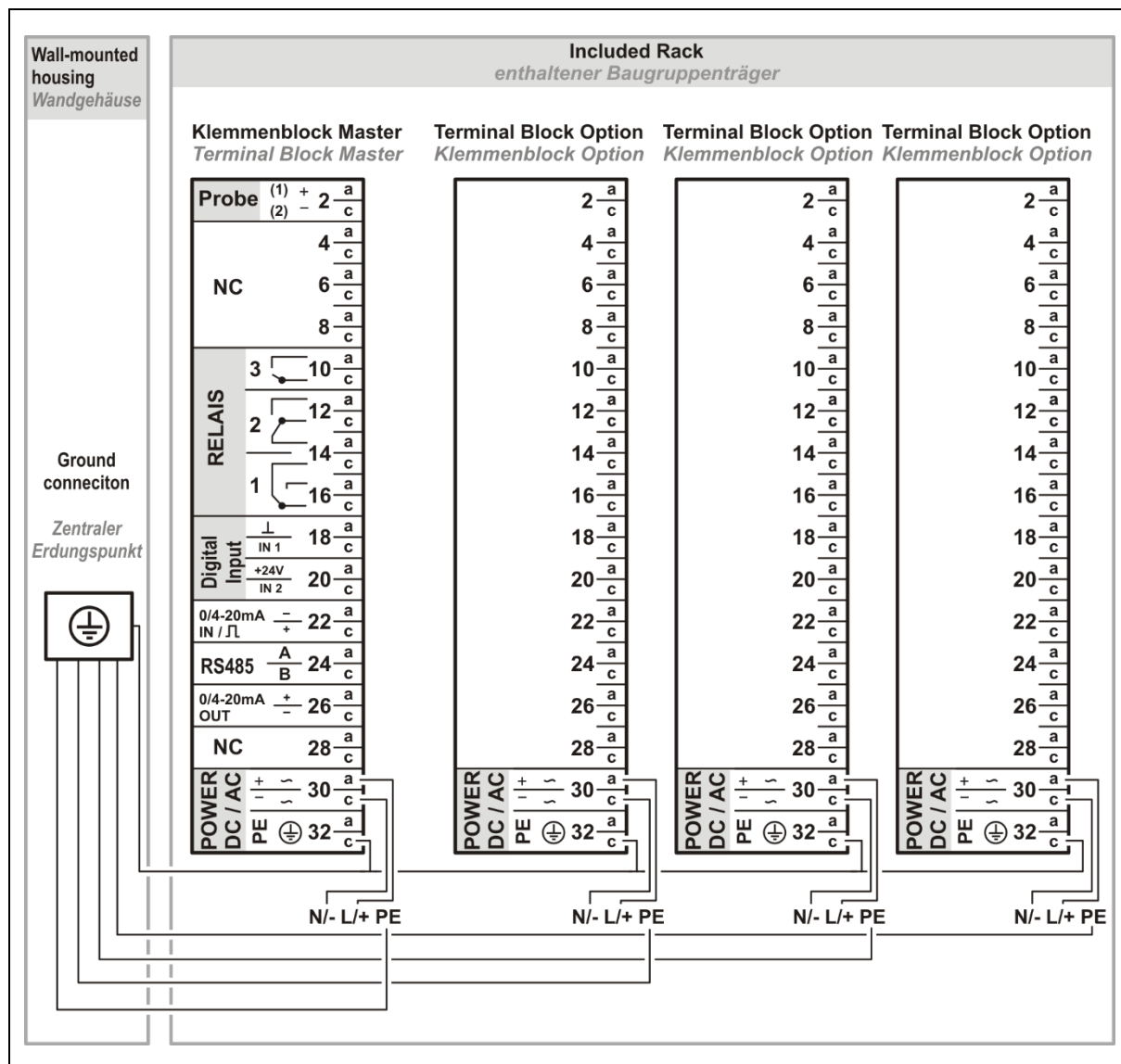


Since the master EVU and the extension modules (7-Wire/RS485-Repeater module, Current Output/HART module) create a RS485 network, ensure that the last bus participants (first and last) are terminated and thus require a 120 Ω terminating wire resistor between terminal A-24 and C-24.

Da die Master AWE und die Erweiterungskarten (7-Leiter/RS485-Repeater Modul, Stromausgang/HART Modul) ein RS485-Netzwerk bilden, muss darauf geachtet werden, dass die letzten Busteilnehmer (erster und letzter) terminiert werden und somit einen 120 Ω Abschlussdrahtwiderstand zwischen Klemme A-24 und C-24 benötigen.

2.6. Assignment of the Wall-mounted Housing for arbitrary Extension Modules

2.6. Belegung des Wandgehäuses für beliebige Erweiterungsmodule



TI-Abb. 6 Assignment drawing of the wall-mounted housing
Belegung des Wandgehäuses

NOTICE HINWEIS



The wall housing includes a sticker as shown in TI-Abb. 6. The terminal assignments of the three possible extension modules are also shown on the sticker.

Das Wandgehäuse enthält einen Aufkleber wie in TI-Abb. 6 dargestellt. Auf dem Aufkleber werden zudem die Klemmenbelegungen der drei möglichen Erweiterungsmodule abgebildet.

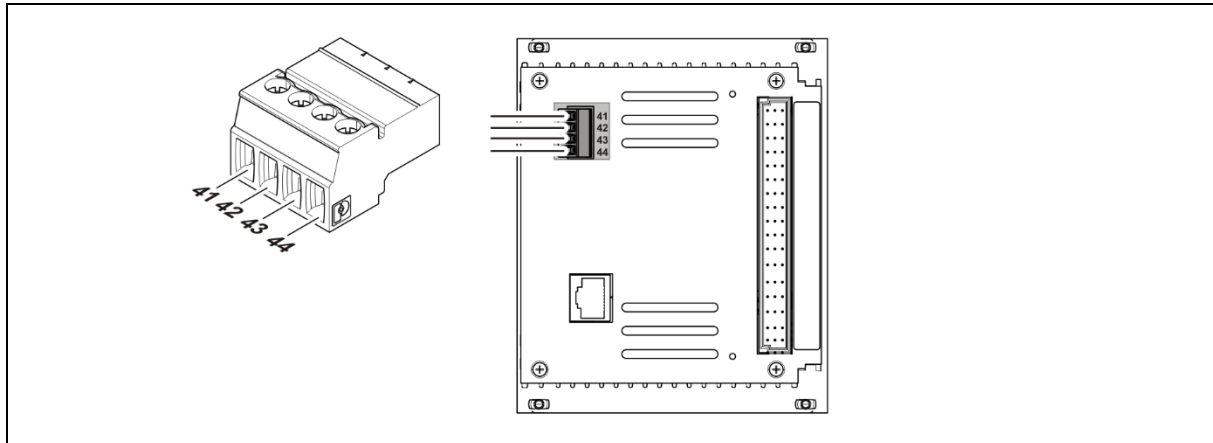
2.7. Assignment Terminals Master/Slave Plug

2.7. Belegung Master/Slave Stecker

NOTICE HINWEIS



The Master/Slave plug is only used in Master/Slave configurations with LB 4700 detectors.
Der Master/Slave Stecker wird nur bei Master/Slave Anordnungen mit LB 4700 Detektoren verwendet.

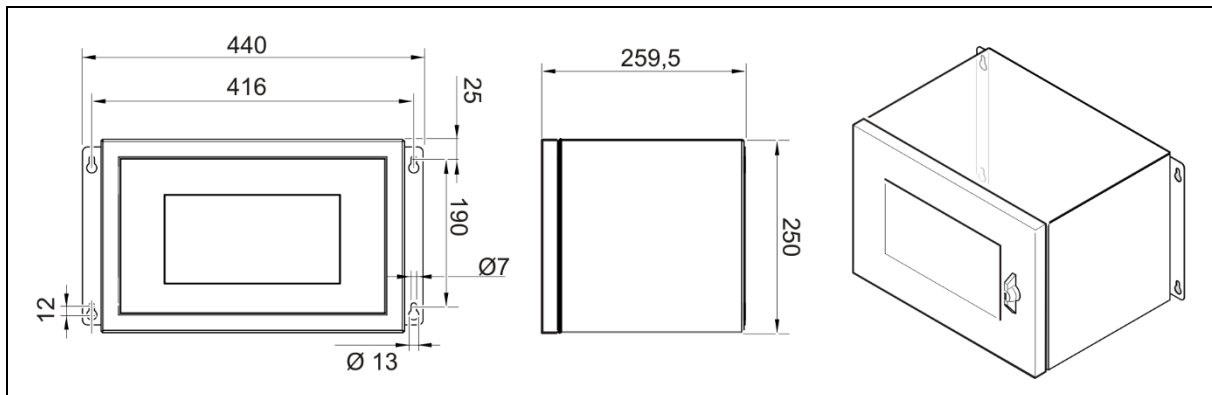


TI-Abb. 7 Assignment Terminal Master/Slave Plug
Klemmenbelegung Master/Slave Stecker

Signal	Pin
TxD	41
RxD	42
RTS	43
GND	44

3. Wall Housing

3. Wandgehäuse



TI-Abb. 8 Drawing wall housing
Zeichnung Wandgehäuse

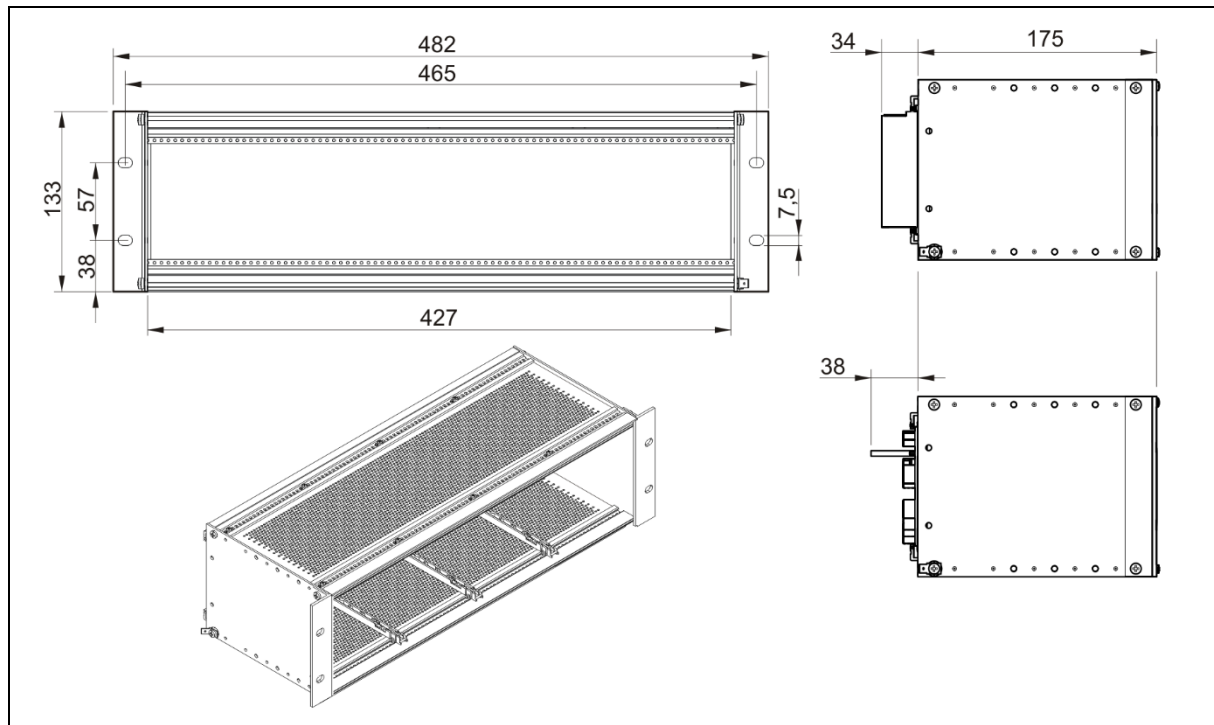
Technical Data Technische Daten	
Max. assembly <i>Max. Bestückung</i>	- 1 Master, 3 arbitrary extension modules with terminal blocks ¹ - 1 Master, 3 beliebige Erweiterungsmodule mit Klemmenblöcken ¹
Weight (with circuit board, without modules) <i>Gewicht (mit Anschlussplatine, ohne Module)</i>	8.8 kg 8,8 kg
Degree of protection <i>Schutzart</i>	IP65
Operational temperature <i>Betriebstemperatur</i>	-20 °C ... +40 °C
Storage temperature <i>Lagertemperatur</i>	-25 °C ... +80 °C
General ambient conditions <i>Allgemeine Umgebungsbedingungen</i>	Overvoltage category: II Pollution Degree: 2 Altitude: up to 2000 m Rel. humidity: 93% or less Überspannungskategorie: II Verschmutzungsgrad: 2 Höhenlage: bis zu 2000 m Rel. Luftfeuchtigkeit: 93% oder weniger
User interface, colours <i>Oberfläche, Farbe</i>	Powder coated, grey Pulverbeschichtet, grau
Cable entry <i>Kabeleinführung</i>	8 x M16, 2 x M32

¹ No NRTL certification

¹ Kein NRTL-Zertifikat

4. 19" Subrack

4. 19" Baugruppenträger



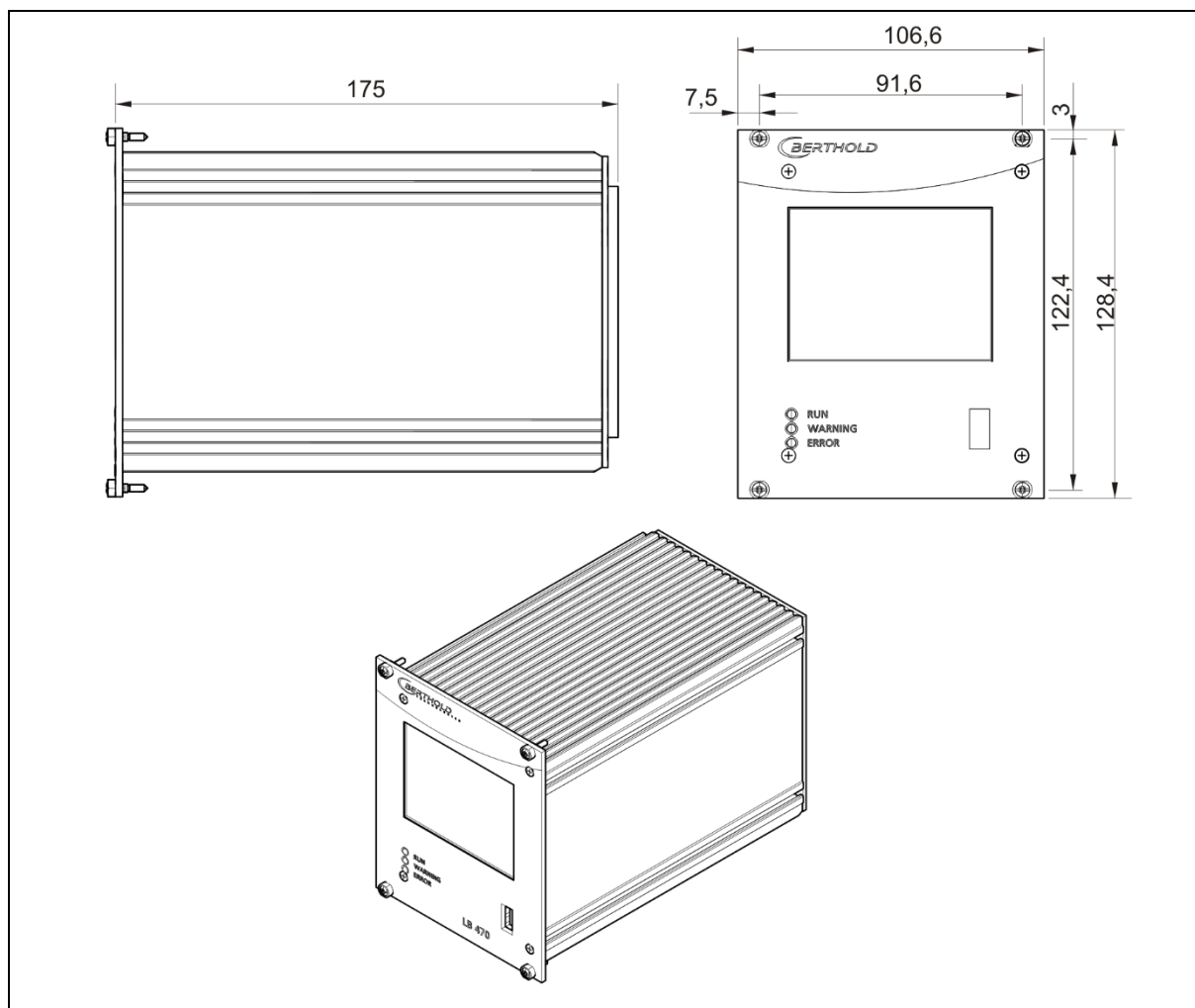
TI-Abb. 9 Drawing 19" subrack
Zeichnung 19" Baugruppenträger

Technical Data Technische Daten

Standard units <i>Normeinheiten</i>	3HE/84TE
Max. Assembly <i>Max. Bestückung</i>	3 Master + 3 Slave / 2 Master + 6 Slave / 4 Master / 1 Master+ 9 Slave / 12 Slave
Weight (with circuit board, without modules) <i>Gewicht (mit Anschlussplatine, ohne Module)</i>	1.4 kg 1,4 kg
Weight terminal block <i>Gewicht Klemmenblock</i>	220 g
Operational temperature <i>Betriebstemperatur</i>	-20 °C ... +50 °C, not condensing -20°C ... +50°C, nicht kondensierend
Storage temperature <i>Lagerungstemperatur</i>	-25°C ... +80°C
Degree of protection <i>Schutzart</i>	IP20

5. Master EVU

5. Master AWE



TI-Abb. 10 Drawing master EVU
Zeichnung Master AWE

Technical Data Technische Daten

Weight Gewicht	1200 g
Operational temperature Betriebstemperatur	-20 °C ... +50 °C not condensing. Avoid direct sunlight. Unobstructed air circulation must be provided to the subrack. -20 °C ... +50 °C nicht kondensierend. Direkte Sonneneinstrahlung ist zu vermeiden. Für eine ungehinderte Luftzirkulation um den Baugruppenträger ist zu sorgen.
Storage temperature Lagerungstemperatur	-30 °C ... +80 °C

General ambient conditions <i>Allgemeine Umgebungsbedingungen</i>	Overvoltage category: II Pollution Degree: 2 Altitude: up to 2000 m Rel. humidity: 93% or less <i>Überspannungskategorie: II Verschmutzungsgrad: 2 Höhenlage: bis zu 2000 m Rel. Luftfeuchtigkeit: 93% oder weniger</i>
Degree of protection <i>Schutzart</i>	IP20 ¹
Connections <i>Anschlüsse</i>	- USB port for the connection to the USB storage medium - Master/slave connection (4-pin) and plug - RJ45 connection for Ethernet (on back wall) - 32-pin plug connector according to DIN 19465 Series C <i>- 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</i>
Display <i>Display</i>	- graphical LCD display - 320 x 240 points, 262,000 colours - Dimmable LED background lighting - Touchscreen <i>- graphisches LCD-Display - 320 x 240 Punkte, 262.000 Farben - Dimmbare LED Hintergrundbeleuchtung - Touchscreen</i>
Computer core <i>Rechnerkern</i>	- Processor: Dual Core DSP/ARM Controller - clock frequency: 300 MHz internal (20 MHz external quartz) - ROM: 512 KByte - RAM: 64 MByte ext. SDRAM, 128 KByte int. shared RAM - FLASH: 8 MByte external serial <i>- Prozessor: Dual Core DSP/ARM Controller - Taktfrequenz: 300 MHz intern (20 MHz externer Quarz) - ROM: 512 KByte - RAM: 64 MByte ext. SDRAM, 128 KByte int. shared RAM - FLASH: 8 MByte extern seriell</i>

Power Supply

Stromversorgung

Voltage <i>Spannung</i>	100...240 VAC 50/60 Hz (wide range input) +/- 10% 21...32 VDC (24 VDC power input)
Power consumption <i>Leistungsaufnahme</i>	22 VA, 15 W
Fuses <i>Sicherungen</i>	Internal, 2 x 250 V, 1A delayed, 5x20 mm, 1500 A breaking capacity IEC 60127-2, 1x 250 V TR5 T80 mA (Ø 8,5 mm)

¹ Designed for indoor use. Outdoor use only possible if installed in wall housing.

¹ Für den Einsatz in Innenräumen geeignet. Außeneinsatz nur bei Installation im Wandgehäuse möglich.

Interfaces

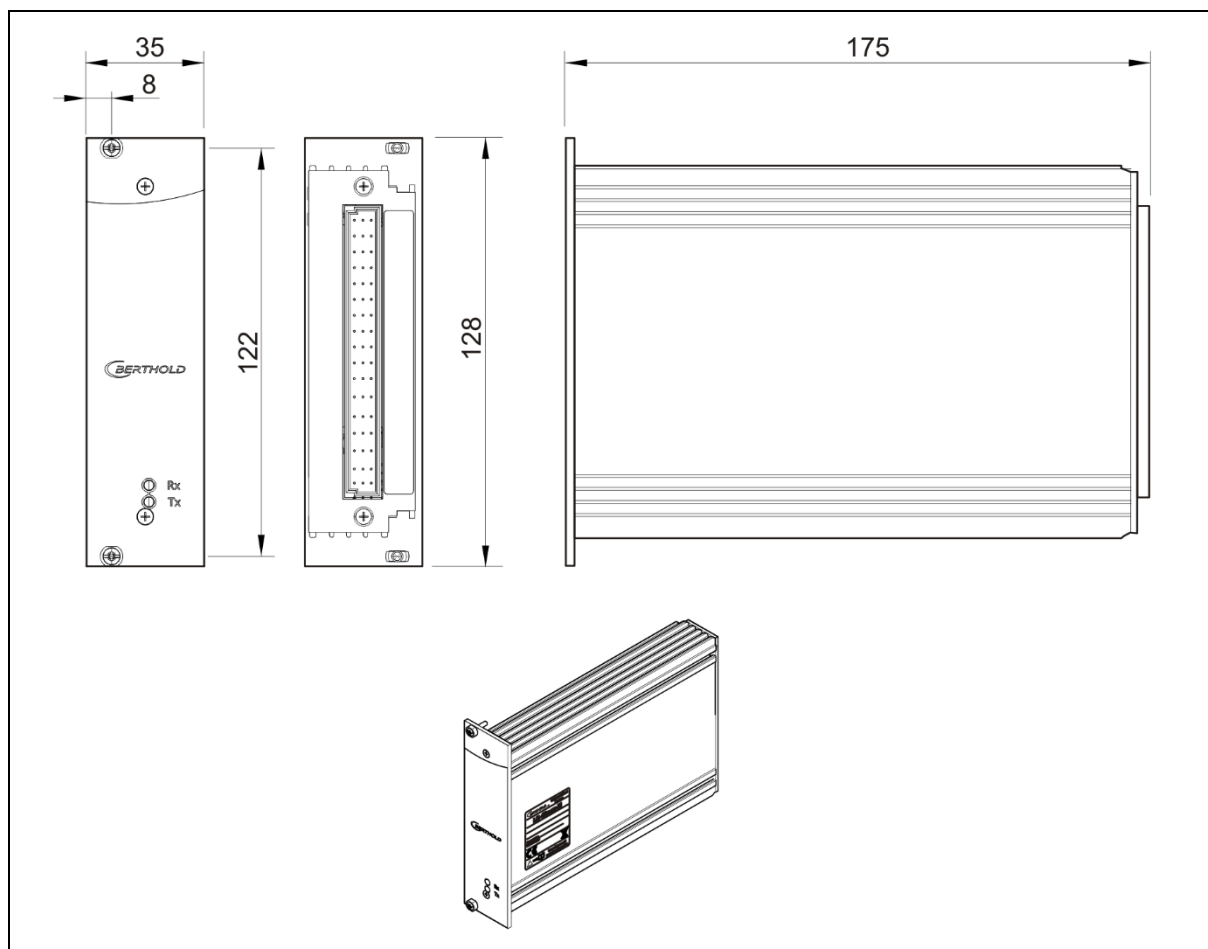
Schnittstellen

Current output <i>Stromausgang</i>	4...20 mA internally switched from power source to sink current (according to NAMUR recommendation NE 006 and NE 043). Dip switch source/sink on the electronic board of the LB 47x. Standard setting is source current. Continuous short circuit proof and galvanically isolated (500 V). Internal resistance about 105 ohm max. Burden when operating as a power source: 850 ohm. Internal monitoring of the loop current and additional error signalling by hardware on detection of a fault condition. <i>4...20 mA (nach Namur-Empfehlung NE 006 und NE 043) intern von Stromquelle auf Stromsenke umschaltbar. Dip-Schalter auf der Elektronik-Platine in der Auswerteeinheit. Standard-Einstellung ist aktiver Stromausgang. 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.</i>
Digital outputs <i>Digitale Ausgänge</i>	3 relays, $U_{\max} = 33 \text{ V AC}_{\text{eff}}, 46 \text{ V DC}; I_{\max} = 1 \text{ A}$ functions: Relay 1: SPDT for error signalling Relay 2: SPDT assignable by software Relay 3: SPST assignable by software <i>3 Relais, $U_{\max} = 33 \text{ V AC}_{\text{eff}}, 46 \text{ V DC}; I_{\max} = 1 \text{ A}$ Funktionen: Relais 1: SPDT zur Fehlersignalisierung Relais 2: SPDT über Software zuweisbar Relais 3: SPST über Software zuweisbar</i>
Digital inputs <i>Digitale Eingänge</i>	2 x together electrically isolated (500 V) Switch between DigIn and GND, U_{outmax} approx. 24 V Function configurable via software <i>2 x gemeinsam potentialgetrennt (500 V), Schalter zwischen DigIn und GND, U_{outmax} ca. 24 V Funktion über Software konfigurierbar</i>
External supply <i>Externe Versorgung</i>	Output voltage: 24 V DC Output current: max. 150 mA <i>Ausgangsspannung: 24 V DC Ausgangsstrom: max. 150 mA</i>

RS485	For communication between Master EVU, 7-Wire/RS485-Repeater module, Current Output/HART module and testing and evaluation purposes. Not isolated from main electronics and USB port, electrically isolated from remaining I/Os (500 V).
<i>RS485</i>	<i>Zur Kommunikation zwischen Master AWE, 7-Leiter/RS485 Repeatermodul und Stromausgang/HART Modul, sowie zu Prüf- und Testzwecken. Nicht potentialgetrennt von Hauptelektronik und USB-Anschluss, potentialgetrennt von restlichen I/Os (500 V).</i>
USB port	1 x USB 2.0 Type A (Host) via front plate to the connection of an ext. mouse, keyboard or storage medium $U_{out} = 5\text{ V}$, $I_{outmax} = 0.5\text{ A}$
<i>USB Anschluss</i>	<i>1 x USB 2.0 Typ A (Host) über Frontplatte zum Anschluss einer ext. Maus, Tastatur oder Speichermedium $U_{out} = 5\text{ V}$, $I_{outmax} = 0,5\text{ A}$</i>
Ethernet	RJ45 connection via back wall, 10 Mbit, DHCP supported, max. 3 m Designed for maintenance purposes. Not designed for long-term operation.
<i>Ethernet</i>	<i>RJ45-Buchse über Rückwand, 10 Mbit, DHCP unterstützt, max. 3 m Vorgesehen für Wartungszwecke. Nicht vorgesehen für den Dauerbetrieb.</i>

6. Slave Module

6. Slave Modul



TI-Abb. 11 Drawing slave module
Zeichnung Slave Modul

Technical Data Technische Daten

Weight <i>Gewicht</i>	600 g
Operational temperature <i>Betriebstemperatur</i>	-20 °C ... +50 °C, not condensing. Avoid direct sunlight. Unobstructed air circulation must be provided to the sub-rack. <i>-20 °C ... +50 °C nicht kondensierend. Direkte Sonneneinstrahlung ist zu vermeiden. Für eine ungehinderte Luftzirkulation um den Baugruppenträger ist zu sorgen.</i>
Storage temperature <i>Lagertemperatur</i>	-30 °C ... +80 °C

General ambient conditions <i>Allgemeine Umgebungsbedingungen</i>	Overvoltage category: II Pollution Degree: 2 Altitude: up to 2000 m Rel. humidity: 93% or less <i>Überspannungskategorie: II Verschmutzungsgrad: 2 Höhenlage: bis zu 2000 m Rel. Luftfeuchtigkeit: 93% oder weniger</i>
Degree of protection <i>Schutzart</i>	IP20 ¹

Power Supply *Stromversorgung*

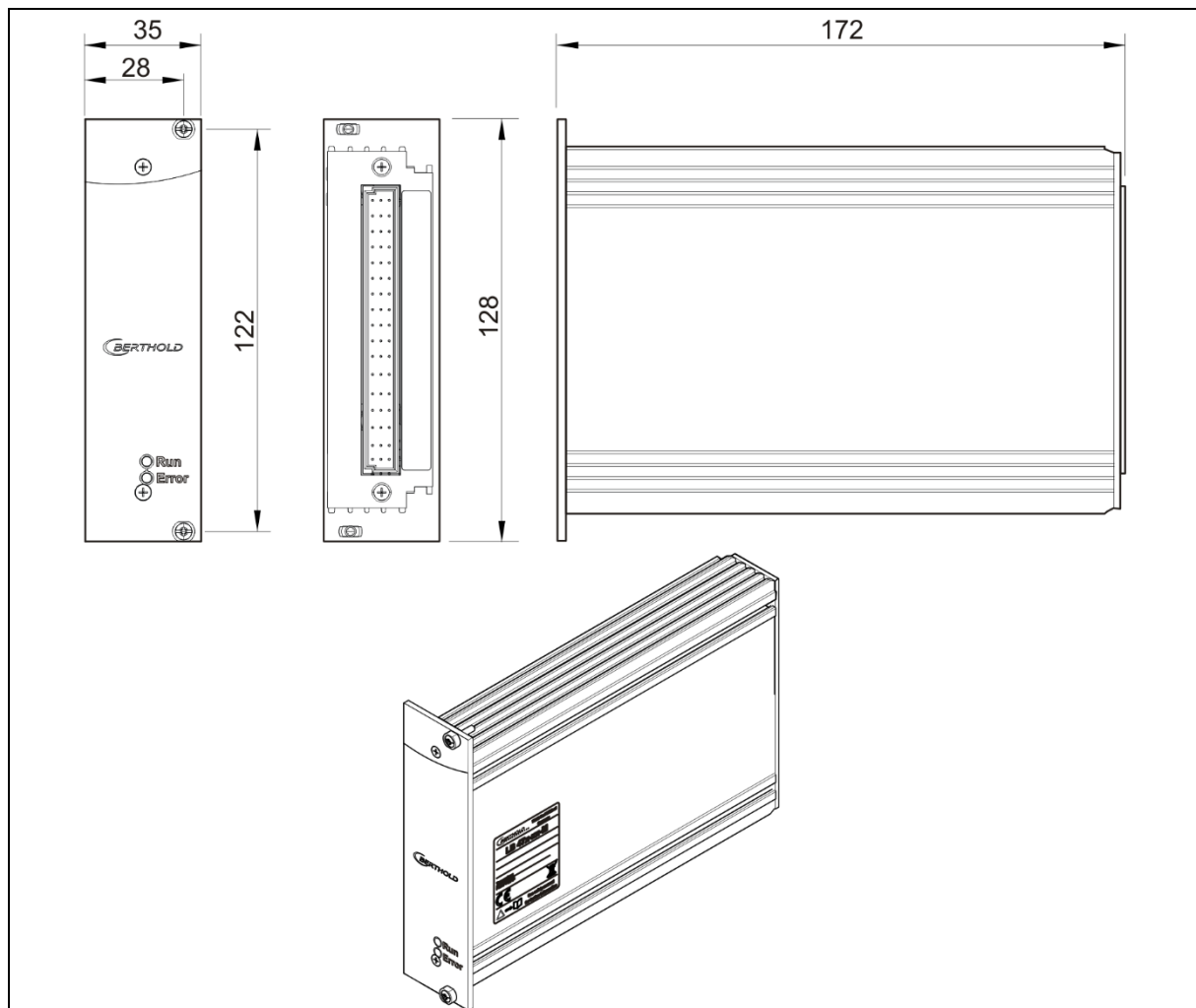
Voltage <i>Spannung</i>	100...240 V AC 50/60 Hz (wide range input) +/- 10% 21...32 VDC (24 VDC power input)
Power consumption <i>Leistungsaufnahme</i>	6 VA, 5 W
Fuses <i>Sicherungen</i>	Internal, 2 x 250 V, 1 A delayed, 5x20 mm, 1500 A breaking capacity IEC 60127-2
Connections <i>Anschlüsse</i>	32-pin plug connector <i>32-polige Stiftleiste</i>

¹ Designed for indoor use. Outdoor use only possible if installed in wall housing.

¹ Für den Einsatz in Innenräumen geeignet. Außeneinsatz nur bei Installation im Wandgehäuse möglich.

7. 7-Wire/RS485-Repeater Module¹

7. 7-Leiter/RS485-Repeater Modul¹



TI-Abb. 12 7-Wire/RS485-Repeater module
7-Leiter/RS485-Repeater Modul

Technical Data

Technische Daten

Weight Gewicht	500 g
Operational temperature Betriebstemperatur	-20 °C ... +50 °C not condensing. Avoid direct sunlight. Unobstructed air circulation must be provided to the subrack. -20 °C ... +50 °C nicht kondensierend. Direkte Sonneneinstrahlung ist zu vermeiden. Für eine ungehinderte Luftzirkulation um den Baugruppenträger ist zu sorgen.
Storage temperature Lagerungstemperatur	-30 °C ... +80 °C

¹ No NRTL certification US/CAN

¹ Kein NRTL-Zertifikat US/CAN

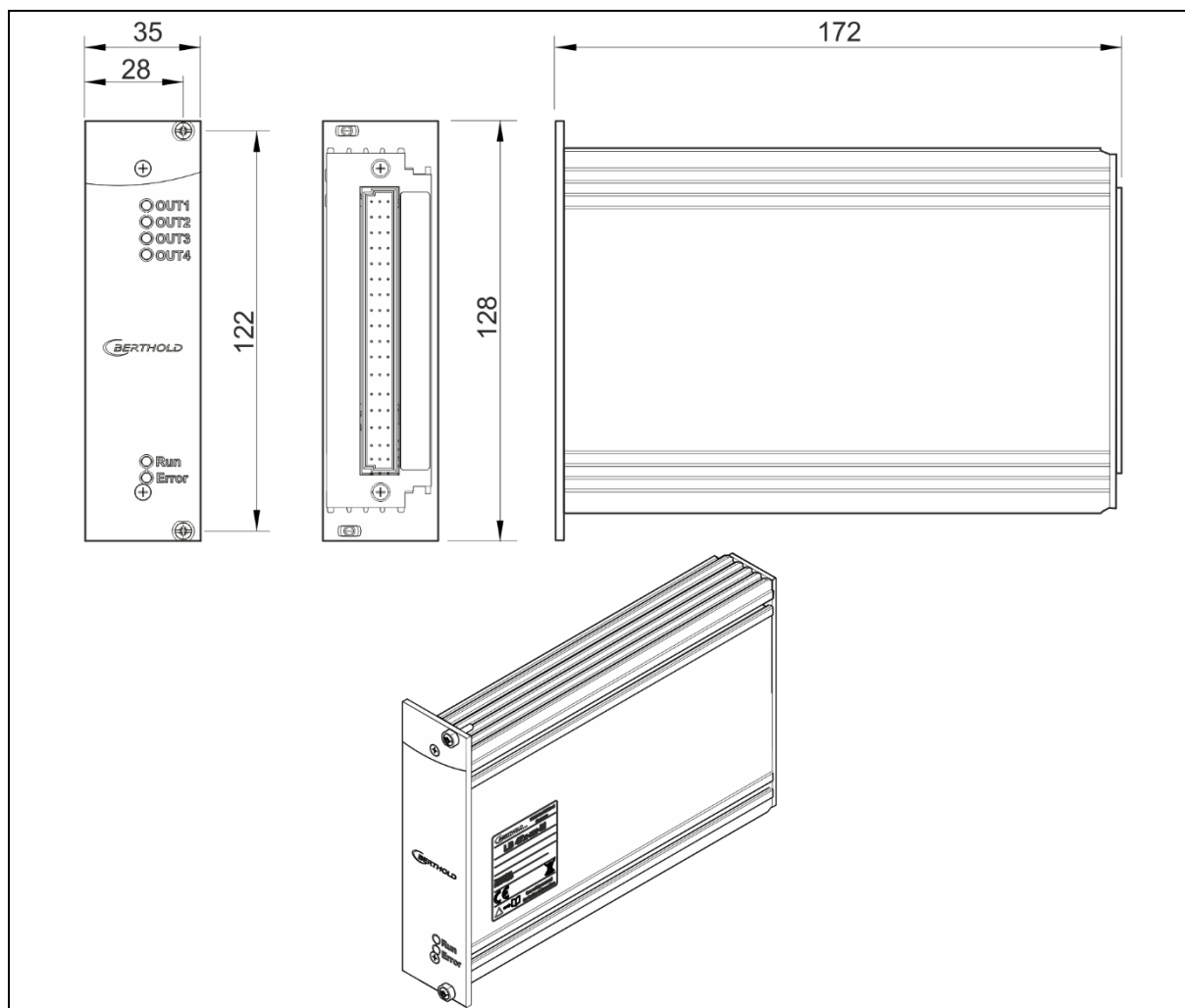
General ambient conditions <i>Allgemeine Umgebungsbedingungen</i>	Overvoltage category: II Pollution Degree: 2 Altitude: up to 2000 m Rel. humidity: 93% or less <i>Überspannungskategorie: II Verschmutzungsgrad: 2 Höhenlage: bis zu 2000 m Rel. Luftfeuchtigkeit: 93% oder weniger</i>
Degree of protection <i>Schutzart</i>	IP20 ²
Power Supply Voltage <i>Versorgungsspannung</i>	100...240 VAC 50/60 Hz (wide range input) +/- 10% 21...32 VDC (24 VDC power input)
Power Consumption <i>Leistungsaufnahme</i>	8 VA, 4 W
Fuses <i>Sicherungen</i>	Internal, 2 x 250 V, 1 A delayed, 5x20 mm, 1500 A breaking capacity IEC 60127-2

² Designed for indoor use. Outdoor use only possible if installed in wall housing.

² Für den Einsatz in Innenräumen geeignet. Außeneinsatz nur bei Installation im Wandgehäuse möglich.

8. Current Output/HART Module¹

8. Stromausgang/HART Modul¹



TI-Abb. 13 Current Output/HART module
Stromausgang/HART Modul

Technical Data Technische Daten

Weight Gewicht	500 g
Operational temperature Betriebstemperatur	-20 °C ... +50 °C not condensing. Avoid direct sunlight. Unobstructed air circulation must be provided to the subrack. -20 °C ... +50 °C nicht kondensierend. Direkte Sonneneinstrahlung ist zu vermeiden. Für eine ungehinderte Luftzirkulation um den Baugruppenträger ist zu sorgen.
Storage temperature Lagerungstemperatur	-30 °C ... +80 °C

¹ No NRTL certification US/CAN

¹ Kein NRTL-Zertifikat US/CAN

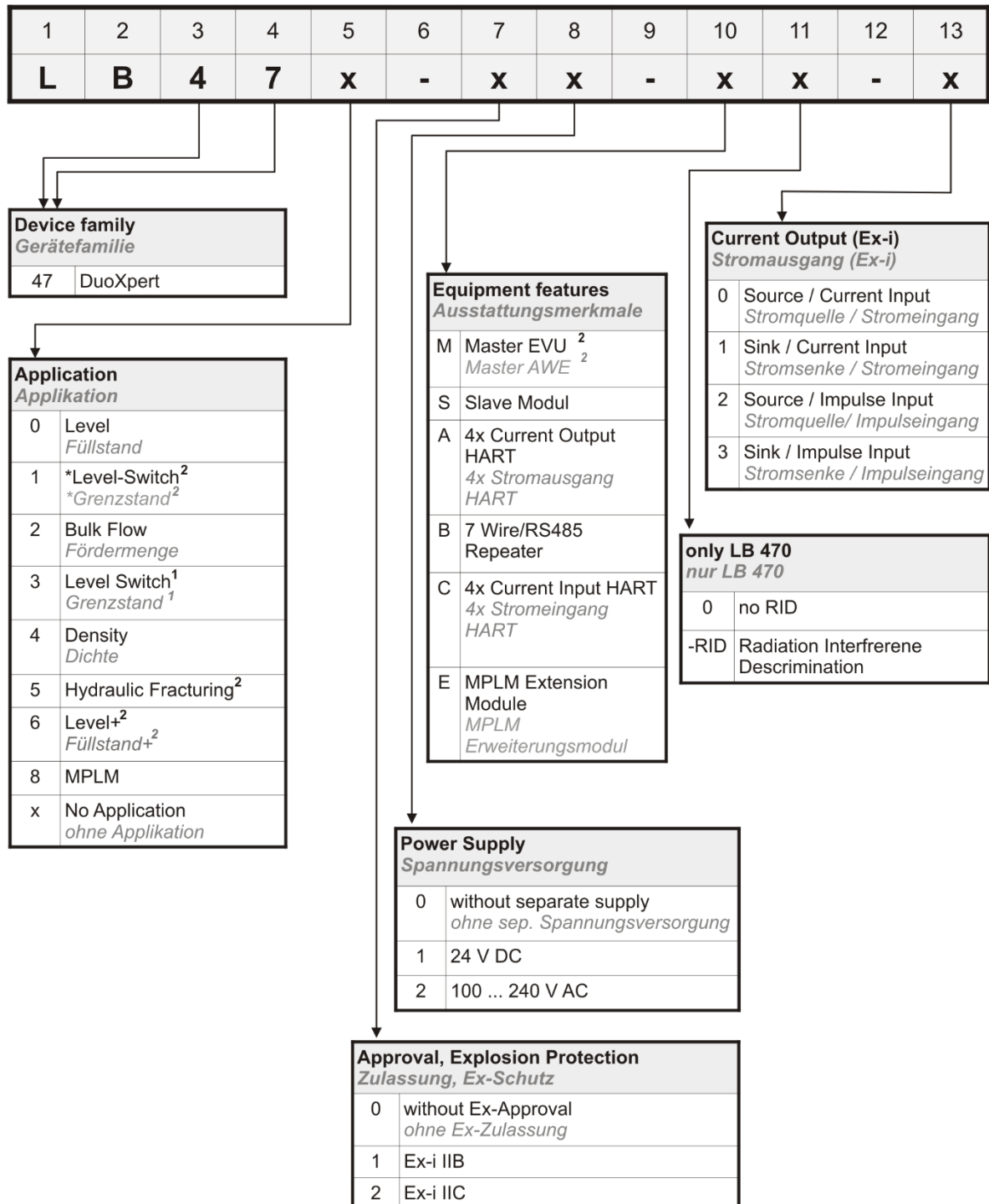
General ambient conditions <i>Allgemeine Umgebungsbedingungen</i>	Overvoltage category: II Pollution Degree: 2 Altitude: up to 2000 m Rel. humidity: 93% or less <i>Überspannungskategorie: II Verschmutzungsgrad: 2 Höhenlage: bis zu 2000 m Rel. Luftfeuchtigkeit: 93% oder weniger</i>
Degree of protection <i>Schutzart</i>	IP20 ²
Power Supply Voltage <i>Versorgungsspannung</i>	100...240 VAC 50/60 Hz (wide range input) +/- 10% 21...32 V DC (24 VDC power input)
Power Consumption <i>Leistungsaufnahme</i>	16 VA, 5 W
Fuses <i>Sicherungen</i>	Internal, 2 x 250 V, 1 A delayed, 5x20 mm, 1500 A breaking capacity IEC 60127-2
Current Outputs <i>Stromausgänge</i>	4x 4...20 mA as power source. Operation as power sink not possible. Continuous short circuit proof and galvanically isolated (500 V). Burden when operating as a power source: 600 ohm. <i>4x 4...20 mA als Stromquelle. Betrieb als Stromsenke nicht möglich. Dauerhaft kurzschlussfest und potentialgetrennt (500 V). Bürde bei Betrieb als Stromquelle: 600 Ohm.</i>

² Designed for indoor use. Outdoor use only possible if installed in wall housing.

² Für den Einsatz in Innenräumen geeignet. Außeneinsatz nur bei Installation im Wandgehäuse möglich.

9. Number Key LB 47x

9. Nummernschlüssel LB 47x



TI-Abb. 14 Number key
Nummernschlüssel

¹ Only available as Ex-i version

² Only available as a standard version

*Other Hardware

¹ Nur verfügbar als Ex-i Version

² Nur verfügbar als Standard-Version

*Andere Hardware

10. EG-Declaration of Conformity 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-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**

	directive	applied standards
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, 1st 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. No.

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

Sparkasse -CW 75323 Bad Wildbad
Volksbank 75119 Pforzheim
Commerzbank 75105 Pforzheim

Konto/Account No. 8 043 003 (BLZ 566 500 00)
Konto/Account No. 557 004 (BLZ 666 800 00)
Konto/Account No. 8 511 120 (BLZ 566 800 12)

SWIFT-BIC: PZHSDE33
SWIFT-BIC: VPM0DE33
SWIFT-BIC: COBSDE33

IBAN: DE37 0660 0000 0000 0450 03
IBAN: DE85 6666 0000 0000 9520 04
IBAN: DE05 6666 0013 0651 1120 00

10. EG-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**

	Richtlinie (Fundstelle)	angewendete Normen und weitere Spezifikationen
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
Geschäftsführung / Management
USt-Id-Nr. / VAT Reg. No.
Deutsche Steuernummer / German Tax No.
WLL-Reg. No.

Stuttgart HRA 330991
BERTHOLD TECHNOLOGIES Verwaltungs-GmbH
Stuttgart HRB 331520
Herrn/Knaut, Dr. Dirk Mörmann
CE613050511
49038/03038
DE59468690

Sparkasse PF-CW
Vollbank
Commerzbank

75323 Bad Wildbad
75119 Pfaffenheim
75105 Pfaffenheim

Konto/Account No. 8 045 003 (RI 7 666 903 85)
Konto/Account No. 957 004 (BLZ 656 900 001)
Konto/Account No. 6 511 128 (BLZ 656 800 13)

SWIFT-BIC PFHS3333
SWIFT-BIC VDPD3333
SWIFT-BIC DRES3333

IBAN: DE37 6565 0085 0038 0450 03
IBAN: DE85 6565 0003 0038 0570 04
IBAN: DE05 6565 0013 0651 1120 00

11. EU-Declaration of Conformity 7-wire/RS-485 repeater

11. EU-Konformitätserklärung 7-Leiter/RS-485 Repeater



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

EU-Konformitätserklärung

EU-Declaration of Conformity

Nr./No.: CE20041-01

Hiermit erklären wir in alleiniger Verantwortung, dass die Bauart des nachfolgend bezeichneten Geräts, in der von uns in den Verkehr gebrachten Ausführung den unten genannten einschlägigen Harmonisierungsvorschriften der EU entspricht.

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

We hereby declare, under our sole responsibility, that the design of the following device placed on the market by us complies with the relevant harmonized rules of the EU.

Unauthorized modifications or unintended use of the product makes this declaration invalid

Hersteller
Manufacturer

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

Produktbezeichnung
Product name

7-Leiter/RS-485 Repeater
7-wire/RS-485 repeater

Typenbezeichnung / Modell
Type / model

LB47x-01-B0
LB47x-02-B0

Richtlinie/Verordnung directive/regulation

angewendete Normen applied standards

*Niederspannungsrichtlinie 2014/35/EU
low voltage directive

EN 61010-1:2010 + A1:2019

EMV / EMC 2014/30/EU

EN IEC 61326-1:2021

Namur NE 021:2017

RoHS 2011/65/EU

* nur die Variante LB47x-02-B0 (100-240 V AC) fällt in den Geltungsbereich der LVD – die Variante LB47x-01-B0 (24 V) nicht.
Only the variant LB47x-02-B0 (100-240 V AC) is within the scope of the LVD – but not the variant LB47x-01-B0 (24 V).

Diese Erklärung wird in Verantwortung für den Hersteller ausgestellt durch
This declaration is issued by the manufacturer released by


Dr. Jürgen Briggmann

Head of R&D
Bad Wildbad, 13.01.2023

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. No.

Bankverbindungen / Bank Details
Sparkasse Pforzheim-Calw
Volksbank Pforzheim
Commerzbank Pforzheim

IBAN BLZ Konto / Account
DE37 6665 0085 0008 0450 03
DE85 6669 0000 0000 9570 04
DE05 6668 0013 0651 1120 00

Stuttgart HRA 330991
Berthold Technologies Verwaltungs-GmbH
Stuttgart HRB 331520
Thomas Bogner
DE813050511
49038/08038
DE99468690

Swift BIC
PZHSDE66
VBPFD66
DRESEFF666

12. EU-Declaration of Conformity 4x Current Output/HART

12. EU-Konformitätserklärung 4x Spannungsausgang/HART



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

EU-Konformitätserklärung

EU-Declaration of Conformity

Nr./No.: CE20040-01

Hiermit erklären wir in alleiniger Verantwortung, dass die Bauart des nachfolgend bezeichneten Geräts, in der von uns in den Verkehr gebrachten Ausführung den unten genannten einschlägigen Harmonisierungs-vorschriften der EU entspricht.

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

We hereby declare, under our sole responsibility, that the design of the following device placed on the market by us complies with the relevant harmonized rules of the EU.

Unauthorized modifications or unintended use of the product makes this declaration invalid

Hersteller
Manufacturer

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

Produktbezeichnung
Product name

4x Spannungsausgang/HART
4x current output/HART

Typenbezeichnung / Modell
Type / model

LB47x-01-A0
LB47x-02-A0

Richtlinie/Verordnung directive/regulation	angewendete Normen applied standards
*Niederspannungsrichtlinie 2014/35/EU low voltage directive	EN 61010-1:2010 + A1:2019
EMV / EMC 2014/30/EU	EN IEC 61326-1:2021 Namur NE 021:2017
RoHS 2011/65/EU	

* nur die Variante LB47x-02-A0 (100-240 V AC) fällt in den Geltungsbereich den LVD – die Variante LB47x-01-A0 (24 V) nicht.
Only the variant LB47x-02-A0 (100-240 V AC) is within the scope of the LVD – but not the variant LB47x-01-A0 (24 V).

Diese Erklärung wird in Verantwortung für den Hersteller ausgestellt durch
This declaration is issued by the manufacturer released by


Dr. Jürgen Briggmann

Head of R&D
Bad Wildbad, 13.01.2023

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. No.

Bankverbindungen / Bank Details
Sparkasse Pforzheim-Calw
Volksbank Pforzheim
Commerzbank Pforzheim

IBAN BLZ Konto / Account
DE37 6665 0085 0008 0450 03
DE85 6669 0000 0000 9570 04
DE05 6668 0013 0651 1120 00

Stuttgart HRA 330991
Berthold Technologies Verwaltungs-GmbH
Stuttgart HRB 331520
Thomas Bogner
DE813050511
49038/08038
DE99468690

Swift BIC
PZHSDE66
VBPFDE66
DRESDEF666

13. UK Declaration of Conformity 7-wire/RS-485 repeater

13. UK Konformitätserklärung 7-Leiter/RS-485 Repeater



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

UK Declaration of Conformity

File No.: UK20041-01

We hereby declare, under our sole responsibility, that the design of the following device placed on the market by us complies with the relevant U.K. legislation for UKCA-marking.

Unauthorized modifications or unintended use of the product make the declaration invalid.

Product name: **7-wire/RS-485 repeater**

Type / model: **LB47x-01-B0**
LB47x-02-B0

Regulation		applied standards
*Electrical Equipment (Safety) Regulations 2016	SI 2016/1101	EN 61010-1:2010+A1:2019
Electromagnetic Compatibility Regulations 2016	SI 2016/1091	EN IEC 61326-1:2021 Namur NE 021:2017
The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012	SI 2012/3032	

* Only the variant LB47x-02-B0 (100-240 V AC) is within the scope of the Electrical Equipment (Safety) Regulations – but not the variant LB47x-01-B0 (24 V).

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, 13th of January 2023

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. No.

Bankverbindungen / Bank Details
Sparkasse Pforzheim-Calw
Volksbank Pforzheim
Commerzbank Pforzheim

IBAN BLZ Konto / AccountSwift
DE37 6565 0085 0008 0450 03
DE85 6669 0000 0000 9570 04
DE05 6668 0013 0651 1120 00

Stuttgart HRA 330991
Berthold Technologies Verwaltungs-GmbH
Stuttgart HRB 331520
Thomas Bogner
DE813050511
49038/08038
DE99468690

BIC
PZHSDE66
VBPFD66
DRESDEFF666

14. UK Declaration of Conformity 4x Current Output/HART

14. UK Konformitätserklärung 4x Spannungsausgang/HART



BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Straße 22
75323 Bad Wildbad, GermanyPhone +49 7081 177-0
Fax +49 7081 177-100
info@Berthold.com
www.Berthold.com

UK Declaration of Conformity

File No.: UK20040-01

We hereby declare, under our sole responsibility, that the design of the following device placed on the market by us complies with the relevant U.K. legislation for UKCA-marking.

Unauthorized modifications or unintended use of the product make the declaration invalid.

Product name: **4x current output/HART**Type / model: **LB47x-01-A0**
LB47x-02-A0

Regulation		applied standards
*Electrical Equipment (Safety) Regulations 2016	SI 2016/1101	EN 61010-1:2010+A1:2019
Electromagnetic Compatibility Regulations 2016	SI 2016/1091	EN IEC 61326-1:2021 Namur NE 021:2017
The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012	SI 2012/3032	

* Only the variant LB47x-02-A0 (100-240 V AC) is within the scope of the Electrical Equipment (Safety) Regulations - but not the variant LB47x-01-A0 (24 V).

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, 13th of January 2023

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. No.

Bankverbindungen / Bank Details
Sparkasse Pforzheim-Calw
Volksbank Pforzheim
Commerzbank Pforzheim

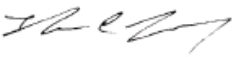
IBAN BLZ Konto / AccountSwift
DE37 6665 0085 0008 0450 03
DE85 6669 0000 0000 9570 04
DE05 6668 0013 0651 1120 00

Stuttgart HRA 330991
Berthold Technologies Verwaltungs-GmbH
Stuttgart HRB 331520
Thomas Bogner
DE813050511
49038/08038
DE99468690

BIC
PZHSDE66
VBPFDE66
DRESDEFF666

15.NTRL Certification US/CAN wall-mounted housing

15.NRTL Zertifikat US/CAN Wandgehäuse

 Nemko-CCL, Inc.	Certificate of Compliance
Certificate: NA201610530	Date Issued: January 20, 2016
Project: 257087-7.1	
Issued to: 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>	
	
Issued by: 	Robert Keller, Senior Engineer/Safety Supervisor
Authorized by: 	Thomas Jackson, Certification Manager
<u>PRODUCTS</u>	
MEASUREMENT, CONTROL, OR LABORATORY EQUIPMENT – Certified to US and Canada Standards	
Product: Process measurement unit Model: 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). Ratings: 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	
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 1 of 3

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 84119-3039 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	Original Certification: Model: 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). Ratings: 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 84119-3039 Tel (801) 972-6146 Fax (801) 972-8432

NFCC-002 Issue 2 May 2014



Page 3 of 3

16.NTRL Certifikate US/CAN DuoXpert LB 47x

16.NTRL Zertifikat US/CAN DuoXpert LB 47x

		Certificate of Compliance	
Nemko-CCL, Inc.			
Certificate: NA201510498		Date Issued: September 17, 2015	
Project: 235982-14.1			
Issued to: 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>			
			
Issued by:		 Robert Keller, Senior Engineer/Safety Supervisor	
Authorized by:		 Thomas Jackson, Certification Manager	
<u>PRODUCTS</u>			
MEASUREMENT, CONTROL, OR LABORATORY EQUIPMENT – Certified to US and Canada Standards			
Product: Process measurement unit for building-in 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) Ratings: LB47x-02-M: 100-240V AC 22VA 50/60Hz; LB47x-02-S: 100-240V AC 6VA 50/60Hz			
<u>APPLICABLE REQUIREMENTS</u>			
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.			
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 1 of 2	

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	Original Certification: 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) Ratings: 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 84119-3039 Tel (801) 972-6146 Fax (801) 972-8432

NFCC-002 Issue 2 May 2014



Page 2 of 2

17. Parts Overview

17. Übersicht Zubehör

Mat. No. Mat.-Nr.	Description Beschreibung
64656	LB 476-02-M0 Level+ Transmitter (100...240 VAC) <i>LB 476-02-M0 Füllstand + Auswerteeinheit (100...240 VAC)</i>
64657	LB 476-01-M0 Level+ Transmitter (24 VDC) <i>LB 476-01-M0 Füllstand + Auswerteeinheit (24 VDC)</i>
73033	LB 47x-02-A0 4x Current Output/HART (100...240 VAC) <i>LB 47x-02-A0 4x Stromausgang/HART (100...240 VAC)</i>
73072	LB 47x-01-A0 4x Current Output/HART (24 VDC) <i>LB 47x-01-A0 4x Stromausgang/HART (24 VDC)</i>
73148	LB47x-02-B0 7-Wire/RS485-Repeater (100...240 VAC) Only required when using LB 480 (1x in total) <i>LB47x-02-B0 7-Leiter/RS485-Repeater (100...240 VAC)</i> <i>Nur bei Verwendung von LB 480 benötigt (insgesamt 1x)</i>
73147	LB47x-01-B0 7-Wire/RS485-Repeater (24 VDC) Only required when using LB 480 (1x in total) <i>LB47x-01-B0 7-Leiter/RS485-Repeater (24 VDC)</i> <i>Nur bei Verwendung von LB 480 benötigt (insgesamt 1x)</i>
63285	LB 47x-02-S0 Slave (100...240 VAC) Only required when using LB 4700 (per detector) <i>LB 47x-02-S0 Slave (100...240 VAC)</i> <i>Nur bei Verwendung von LB 4700 benötigt (pro Detektor)</i>
63286	LB 47x-01-S0 Slave (24 VDC) Only required when using LB 4700 (per detector) <i>LB 47x-01-S0 Slave (24 VDC)</i> <i>Nur bei Verwendung von LB 4700 benötigt (pro Detektor)</i>
64607	19" subrack for use with terminal blocks <i>19"-Baugruppenträger für den Einsatz mit Klemmblöcken</i>
74818	Wall-mounted housing with preinstalled terminal blocks for 1x Master and 3x arbitrary extension module <i>Wandgehäuse mit vorinstallierten Klemmenblöcken für 1x Master und 3x beliebiges Erweiterungsmodul</i>
59477	Terminal block for LB 47x Master and Master/Slave plug <i>Klemmenblock für LB 47x Master und Master/Slave Stecker</i>
59478	Terminal block für LB 47x Slave <i>Klemmenblock für LB 47x Slave</i>
73149	Terminal block for Current Output/HART <i>Klemmenblock für Current Output/HART</i>
73150	Terminal block for 7-Wire/RS485-Repeater <i>Klemmenblock für 7-Leiter/RS485-Repeater</i>

37526	Front Cover Plate 21 HP / 3 RU (Master) <i>Blindplatte 21TE / 3 HE (Master)</i>
59501	Front Cover Plate 7 HP / 3 RU (Slave) <i>Blindplatte 7TE / 3 HE (Slave)</i>
56925-6BA1	Operating Manual LB 476 Level+, German <i>Betriebsanleitung LB 476 Füllstand +, Deutsch</i>
56925-6BA2	Operating Manual LB 476 Level+, English <i>Betriebsanleitung LB 476 Füllstand +, Englisch</i>