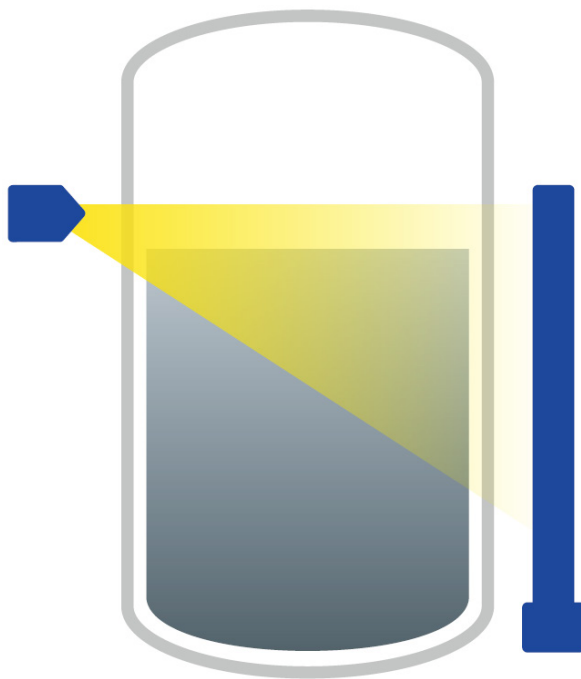




Level

Füllstand

Technical Information

Technische Information

56925TI1L

Rev. No.: 04, 05/2021

1. Information on 2-Wire Technology

The DuoSeries/DuoXPERT measuring system consists of a scintillation detector and a sophisticated evaluation unit (DuoXPERT) for display and operation.

The evaluation unit is a state-of-the-art control unit with robust 3.5" TFT touch panel, powerful Dual Core CPU and diverse operator interfaces. Advanced self-diagnostics and monitoring features ensure a safe function of the system. Furthermore, the data logging functionality allows operators to analyze their processes in depth, e.g. develop trends, track process changes etc.

Sophisticated Measuring System in 2-Wire Technology

- Unique: Radiometric system with intrinsically safe power supply (Full Ex-i) for detectors
- Real 2-wire technology, only 2 wires in the field
- Advanced self-diagnostics and monitoring features
- Easy to use touch screen panel for local display and operation
- Integrated gas density compensation feature
- Direct replacement of predecessor model LB 440
- Interfaces with all 2-wire detectors LB 44xx, LB 54xx and LB 47xx
- Optional Radiation Interference Discrimination (LB 470RID)

1. Informationen zur 2-Leiter Technologie

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

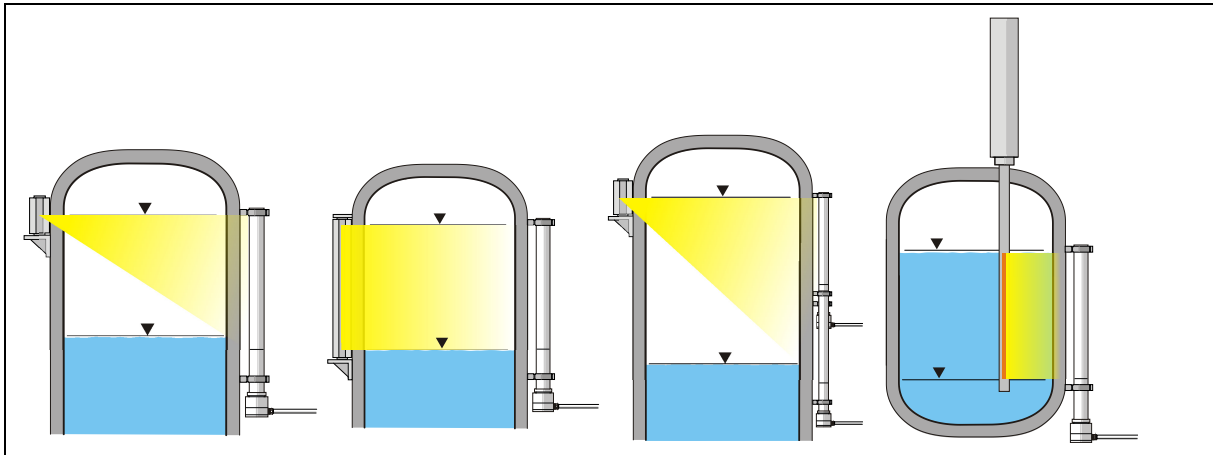
Die moderne Auswerteeinheit verfügt über ein 3,5" Touch Panel, eine starke Dual Core CPU und verschiedenen Bedien-Optionen. Erweiterte Funktionen zur Selbstdiagnose und Überwachung sorgen zudem für höchste funktionale Sicherheit der Messung im Betrieb. Darüber hinaus können die Betreiber die Daten-Log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

Hochentwickeltes Messsystem in 2-Leiter Technologie

- Einzigartig: Radiometrische Messung mit eigensicherer Spannungsversorgung (Voll Ex-i) für den Detektor
- Echte 2-Leiter Technik, nur 2 Adern im Feld
- Verbesserte Diagnosefunktionen und Selbstüberwachung
- Einfache, intuitive Bedienung über Touchscreen
- Integriertes Feature zur Kompensation von Gas-Phasen Schwankungen
- Volle Kompatibilität zum Vorgängermodell LB 440
- Kompatibel zu alle 2-Leiter Detektoren LB 44xx, LB 54xx und LB 47xx
- Optionale Störstrahlungserkennung (LB 470RID)

1.1. Measurement Arrangements with Rod Detector

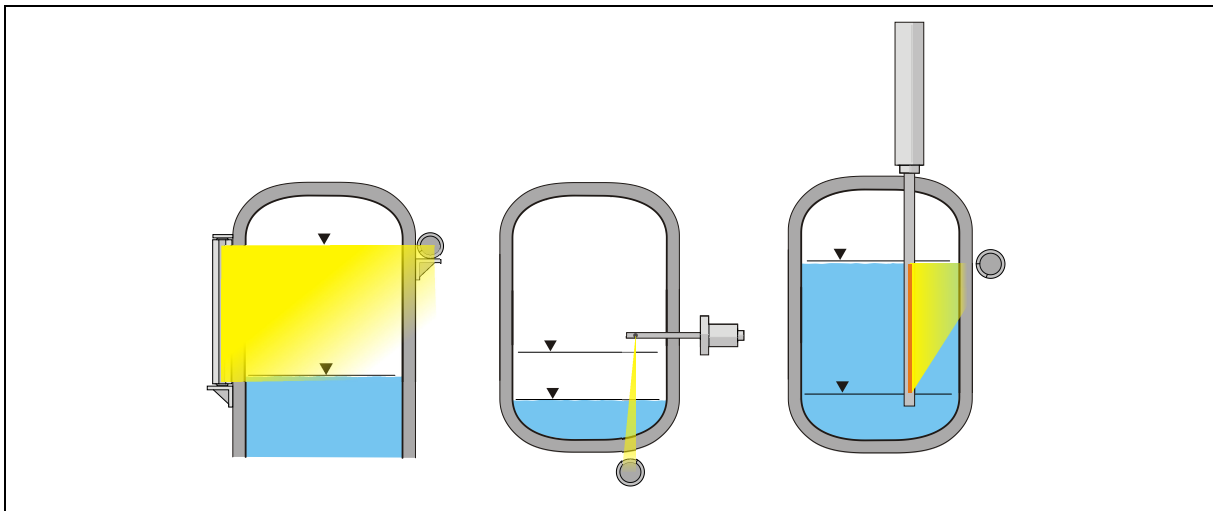
1.1. *Messanordnungen mit Stabdetektor*



TI-Abb. 1 Beispielhafte Messanordnungen mit Stabdetektor
Exemplary measurement arrangements with rod detector

1.2. Measurement Arrangements with Point Detector

1.2. *Messanordnungen mit Punktdetektor*



TI-Abb. 2 Exemplary measurement arrangements with point detector
Beispielhafte Messanordnungen mit Punktdetektor

2. Evaluation Unit

The modules can be installed either in wall housings or 19" subracks. It can be equipped differently, depending on requirements. The rear clamp blocks or terminal panels 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.
- ▶ Please note the applicable documents for versions of the evaluation unit for intrinsically safe operation, see operating manual chapter 1.1.

NOTICE



Note the compatibility!

- ▶ Detector of the type LB 44xx and LB 54xx can capture measurement data only with master EVU.

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 oder Anschlussplatinen 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 .

HINWEIS

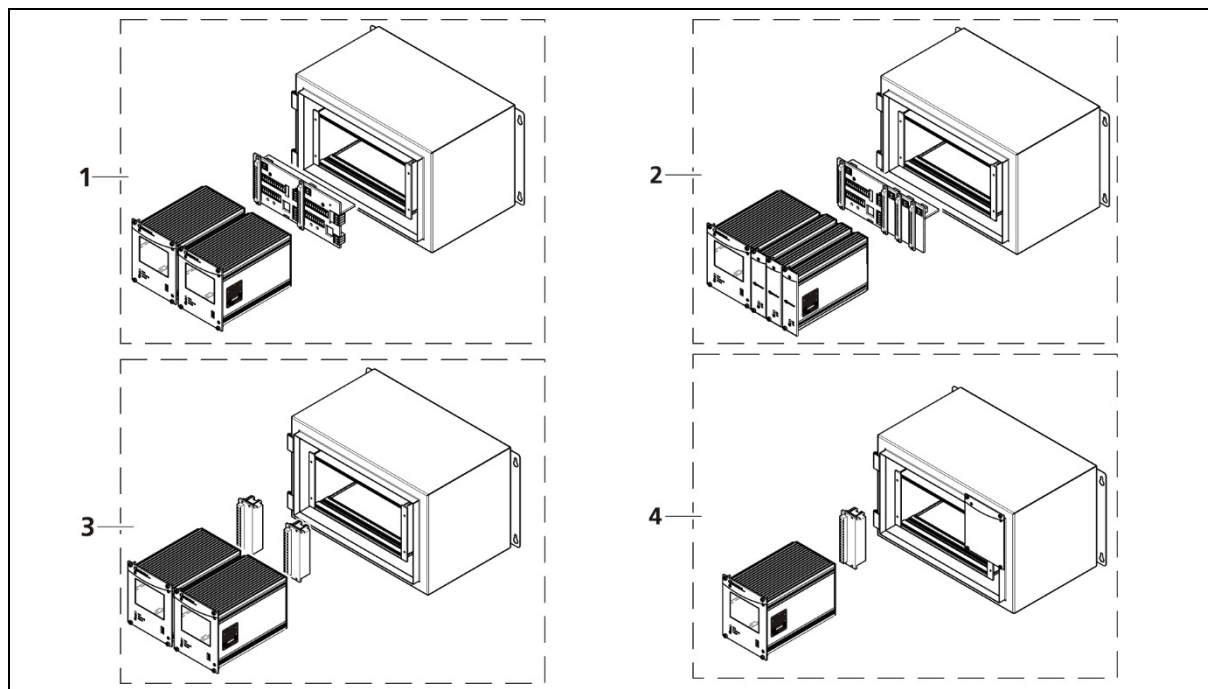


Kompatibilität beachten!

- ▶ Messdaten der Detektoren vom Typ LB 44xx und LB 54xx können nur mit einem Master-Modul erfasst werden.

2.1. Installation Variants Wall Housing

2.1. Installationsvarianten Wandgehäuse



TI-Abb. 3 Installation variants wall housing
Installationsvarianten Wandgehäuse

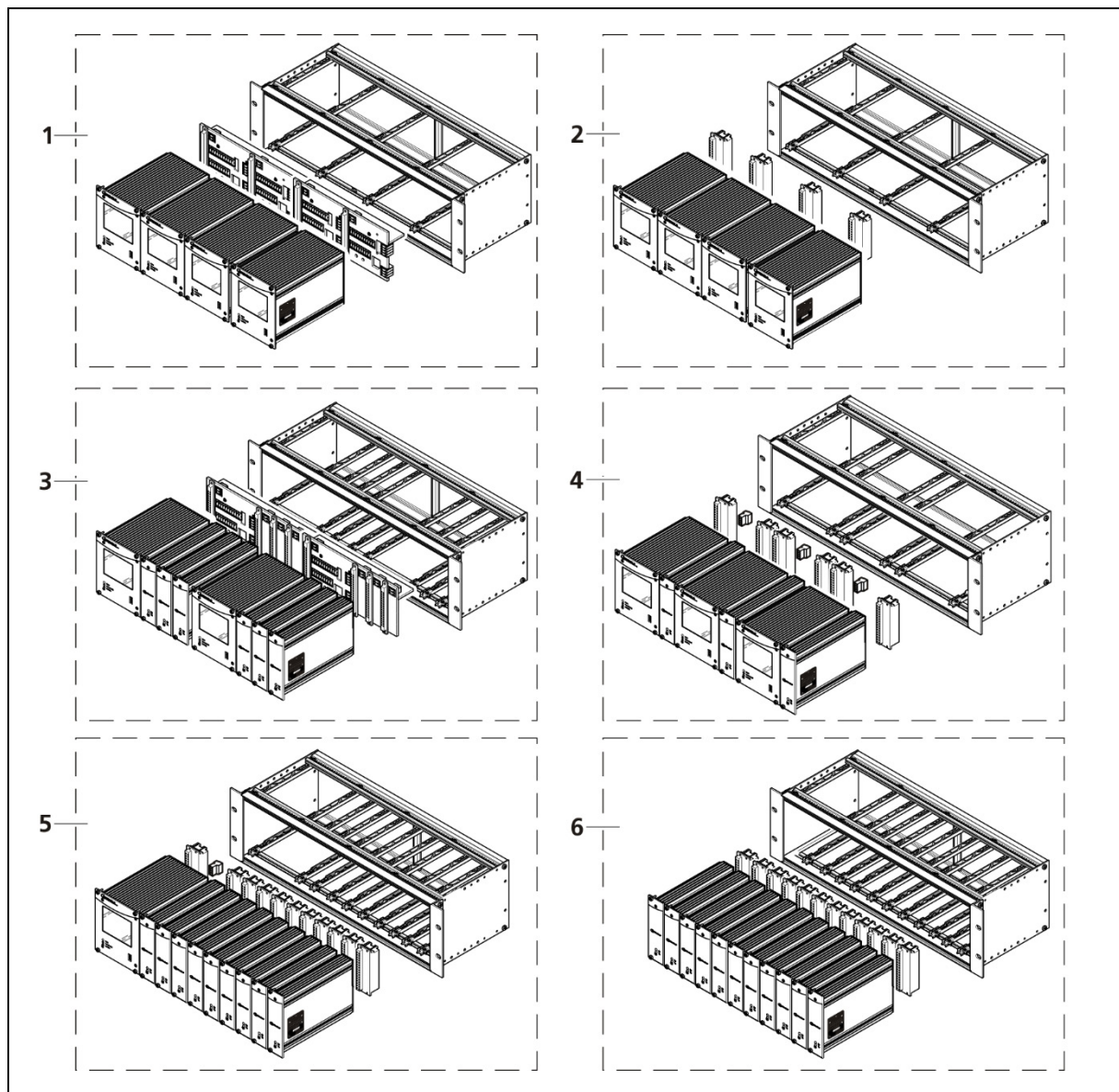
Item Pos.	Componsens Komponenten	Connection Anschluss
1	2 Master	1 Terminal panel master/master ¹ 1 Anschlussplatine Master/Master ¹
2	1 Master, 3 Slaves	1 Terminal panel master/slave ¹ 1 Anschlussplatine Master/Slave ¹
3	2 Master	2 Terminal blocks 2 Klemmenblöcke
4	1 Master, 0 – 3 Slaves	1 Terminal block for master, 0 – 3 Terminal block for slave module 1 Klemmenblock für Master, 0 – 3 Klemmenblöcke für Slave Modul

¹ NRTL certification US/CAN

¹ NRTL Zertifikat US/CAN

2.2. Installation Variants 19" Subrack

2.2. Einbauvarianten 19" Baugruppenträger



TI-Abb. 4 Installation variants 19" subrack
Installationsvarianten 19" Baugruppenträger

Item Pos.	Componsens Komponenten	Connection Anschluss
1	4 Master	2 Terminal panel master 2 Anschlussplatine Master
2	4 Master ²	4 Terminal blocks 4 Klemmenblöcke
3	2x (1 Master, 3 Slaves)	2 Terminal panel master/slave 2 Anschlussplatten Master/Slave
4	4x (1 Master, 1 Slave) ²	6 Terminal blocks; master/slave plugs 6 Klemmenblöcke; Master/Slave Stecker

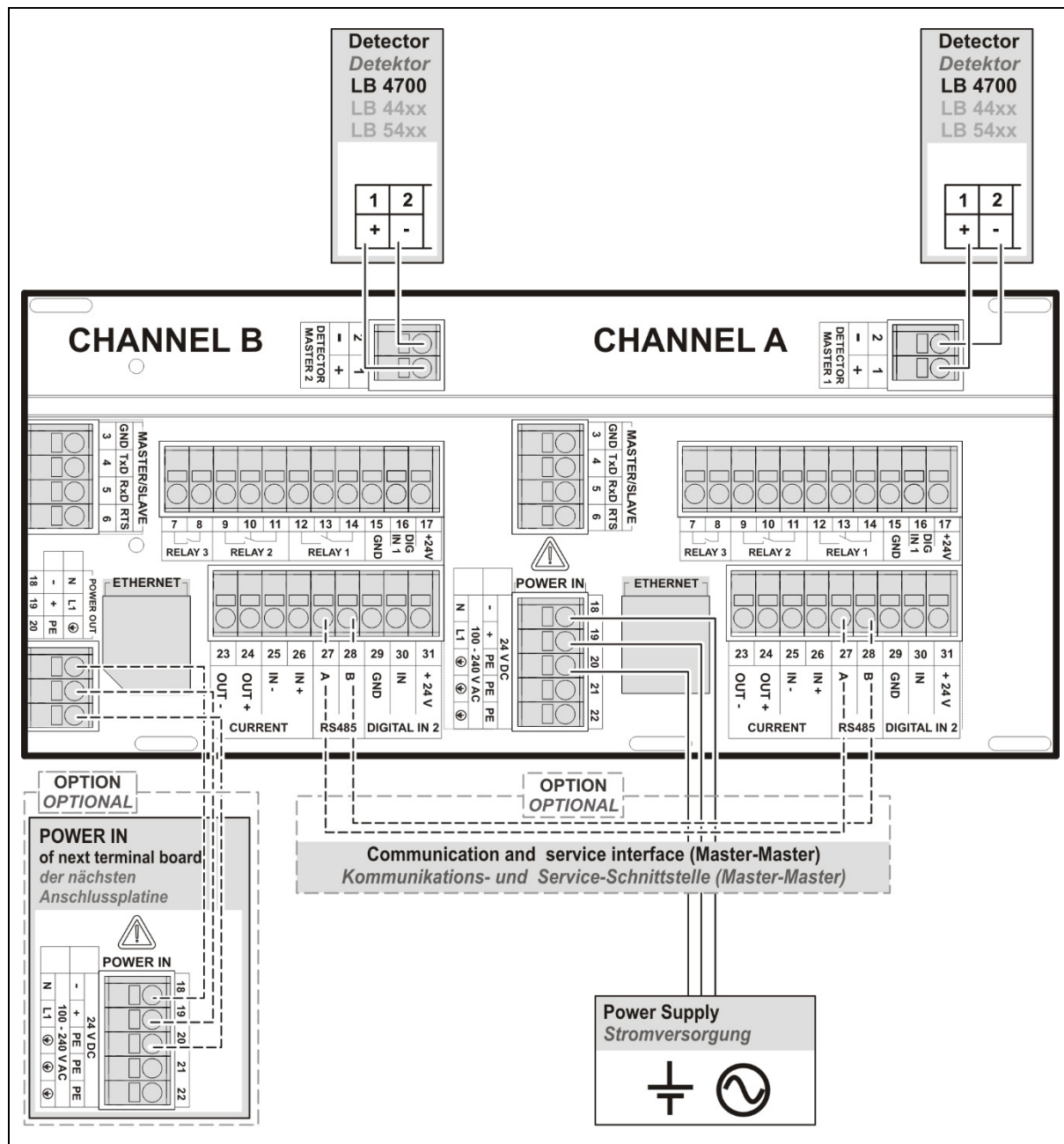
5	1 Master, 9 Slaves ²	10 Terminal blocks; master/slave plug <i>10 Klemmenblöcke; Master/Slave Stecker</i>
6	11 Slaves ²	11 Terminal blocks <i>11 Klemmenblöcke</i>

² Application example. The modules can be arranged arbitrarily with terminal blocks.

² Anwendungsbeispiele. Mit Klemmenblöcken können Module frei zusammengestellt werden.

2.3. Connection Diagram Terminal Board Master/Master

2.3. Anschlussplan Anschlussplatine Master/Master



TI-Abb. 5 Connection Diagram Terminal Board Master/Master
Anschlussplan Anschlussplatine Master/Master

IMPORTANT



In a 19" subrack for 4 masters (Mat. No. 59484), there is another one with Channel C / D next to the connector board for Channel A / B. The channel assignment of Channel C / D is identical to that of Channel A / B.

LB 44xx and LB 54xx detectors can only be operated with master evaluation units, which can be used as slave units in cascaded systems.

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.

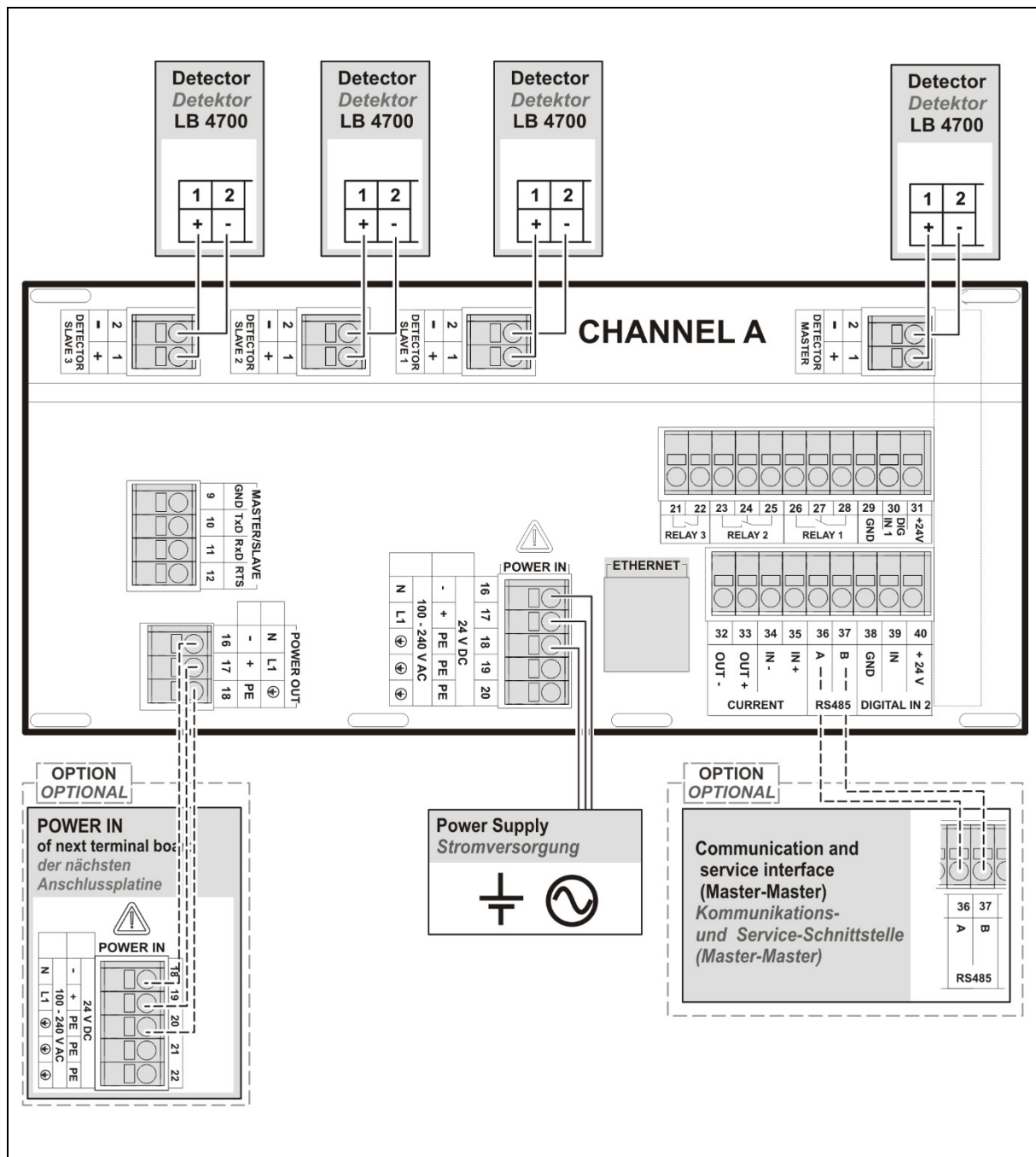
Detektoren vom Typ LB 44xx und LB 54xx können nur mit Master-Auswerteeinheiten betrieben werden, welche bei kaskadierten Systemen als Slaves eingesetzt werden können.

Terminals terminal board master/master
Klemmen Anschlussplatine Master/Master

#	Connection <i>Anschluss</i>	Function <i>Funktion</i>
1	DETECTOR MASTER +	Connection Detector LB 4700 / LB 44xx / LB 54xx <i>Verbindung Detektor LB 4700 / LB 44xx / LB 54xx</i>
2	DETECTOR MASTER -	
3	MASTER/SLAVE GND	Connection of additional slave units <i>Anschluss von weiteren Slave-Einheiten</i>
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	Error DIGITAL OUT <i>Fehler DIGITAL OUT</i>
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 –	Not used for LB 470 / LB 470RID <i>Keine Verwendung bei LB 470 / LB 470RID</i>
26	CURRENT IN +	
27	RS 485 A	Communication and service interface (Master-Master) <i>Kommunikations- und Serviceschnittstelle (Master-Master)</i>
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. Connection Diagram Terminal Board Master/Slave

2.4. Anschlussplan Anschlussplatine Master/Slave



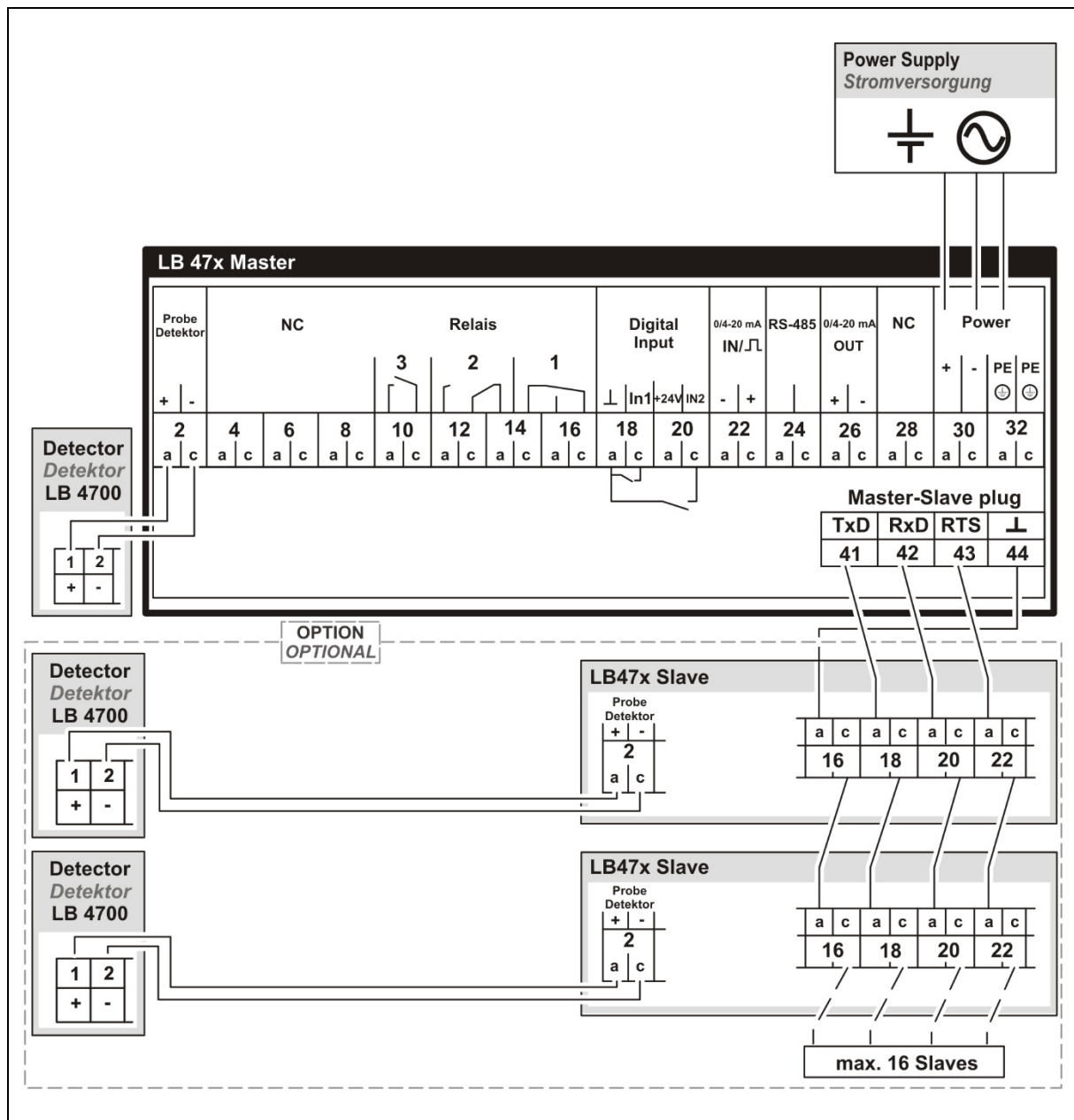
Ti-Abb. 6 Connection Diagram Terminal Board Master/Slave
Anschlussplan Anschlussplatine Master/Slave

Terminals terminal board master/slave
Klemmen Anschlussplatine Master/Slave

#	Connection <i>Anschluss</i>	Function <i>Funktion</i>
1	DETECTOR MASTER +	Connection Detector LB 4700 <i>Verbindung Detektor LB 4700</i>
2	DETECTOR MASTER -	
1	DETECTOR SLAVE 1 +	Connection Detector LB 4700 <i>Verbindung Detektor LB 4700</i>
2	DETECTOR SLAVE 1 -	
1	DETECTOR SLAVE 2 +	
2	DETECTOR SLAVE 2 -	
1	DETECTOR SLAVE 3 +	
2	DETECTOR SLAVE 3 -	
9	MASTER/SLAVE GND	Connection of additional slave units <i>Anschluss von weiteren Slave-Einheiten</i>
10	MASTER/SLAVE TxD	
11	MASTER/SLAVE RxD	
12	MASTER/SLAVE RTS	
16	POWER DC 24 V – / AC N	24 V DC / 100-240 V AC
17	POWER DC 24 V + / AC L1	
18	PE	
19	PE	
20	PE	
21	RELAIS 3 NC	DIGITAL OUT
22	RELAIS 3 COM	
23	RELAIS 2 NC	DIGITAL OUT
24	RELAIS 2 NO	
25	RELAIS 2 COM	
26	RELAIS 1 NC	Error DIGITAL OUT <i>Fehler DIGITAL OUT</i>
27	RELAIS 1 NO	
28	RELAIS 1 COM	
29	DIGITAL IN 1 GND	GND
30	DIGITAL IN 1 IN	Logic Input
31	+ 24 V (GND --> 29)	24 V out (max. 200 mA)
32	CURRENT OUT –	4 mA ... 20 mA
33	CURRENT OUT +	
34	CURRENT IN –	Not used for LB 470 / LB 470RID <i>Keine Verwendung bei LB 470 / LB 470RID</i>
35	CURRENT IN +	
36	RS 485 A	Communication and service interface (Master-Master) <i>Kommunikations- und Service-Schnittstelle (Master-Master)</i>
37	RS 485 B	
38	DIGITAL IN 2 GND	GND
39	DIGITAL IN 2 IN	Logic Input
40	+24 V (GND --> 38)	24 V out (max. 200 mA)

2.5. Assignment Terminal Block Master EVU

2.5. Belegung Klemmenblock Master AWE



TI-Abb. 7 Assignment Terminal Block Master EVU
Belegung Klemmenblock Master AWE

Terminal Block Master EVU Klemmenblock Master AWE

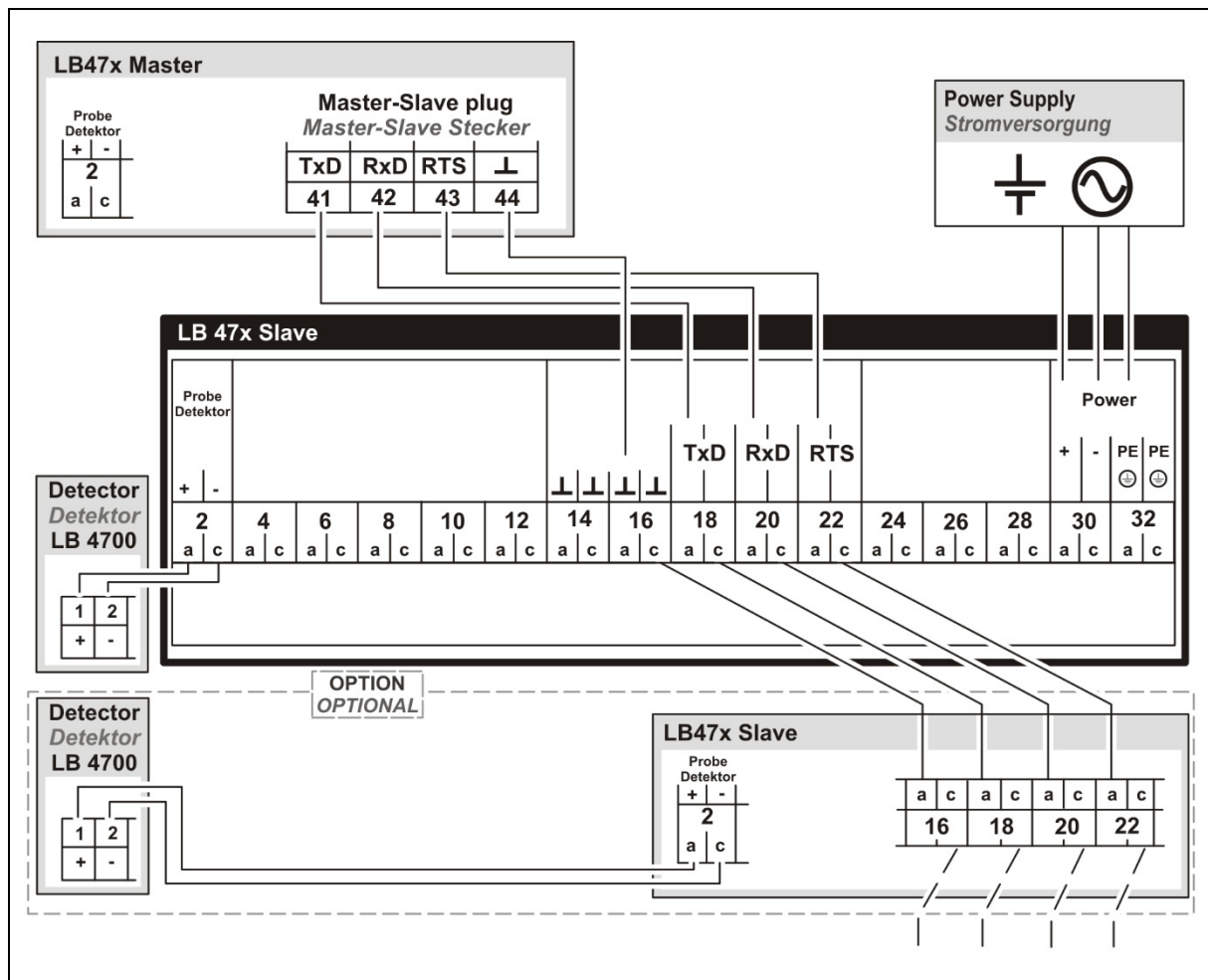
Signal	Pin					Pin	Signal
DETECTOR GND	C - 2					A - 2	DETECTOR +
n.a. *	C - 4					A - 4	n.a. *
n.a. *	C - 6					A - 6	n.a. *
n.a. *	C - 8					A - 8	n.a. *
RELAIS 3 COM	C - 10					A - 10	RELAIS 3 NO
RELAIS 2 COM	C - 12					A - 12	RELAIS 2 NO
RELAIS 1 NC	C - 14					A - 14	RELAIS 2 NC
RELAIS 1 COM	C - 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. *
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

2.6. Assignment Terminal Block Slave

2.6. Belegung Klemmenblock Slave



TI-Abb. 8 Assignment terminal block slave
Belegung Klemmenblock Slave

Terminal Block Slave Klemmenblock Slave

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 ***

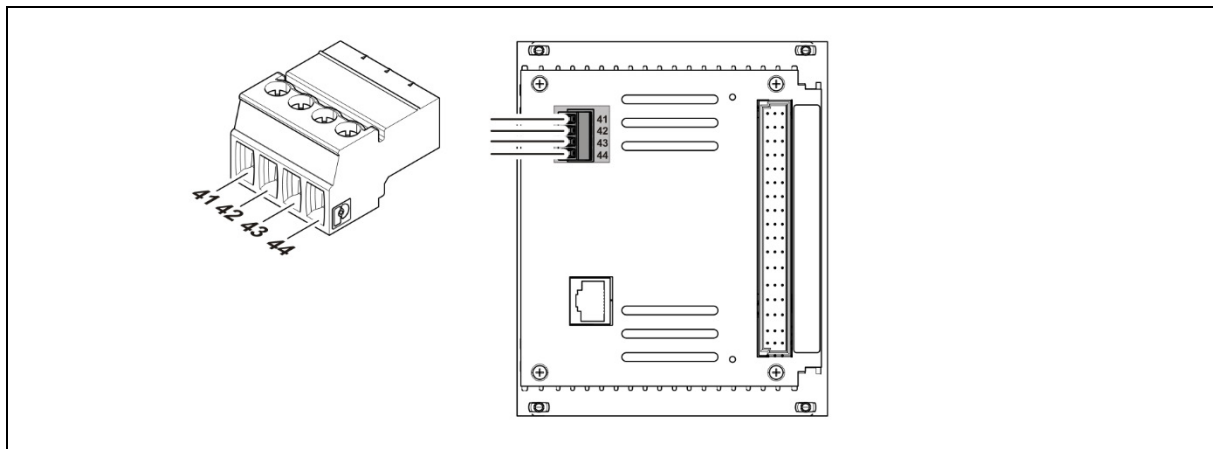
* option
optional

** not assigned
nicht belegt

*** Protective conductor
Schutzleiter

2.7. Assignment Terminals Master/Slave Plug

2.7. Klemmenbelegung Master/Slave Stecker



TI-Abb. 9 Assignment Terminal Master/Slave Plug
Klemmenbelegung Master/Slave Stecker

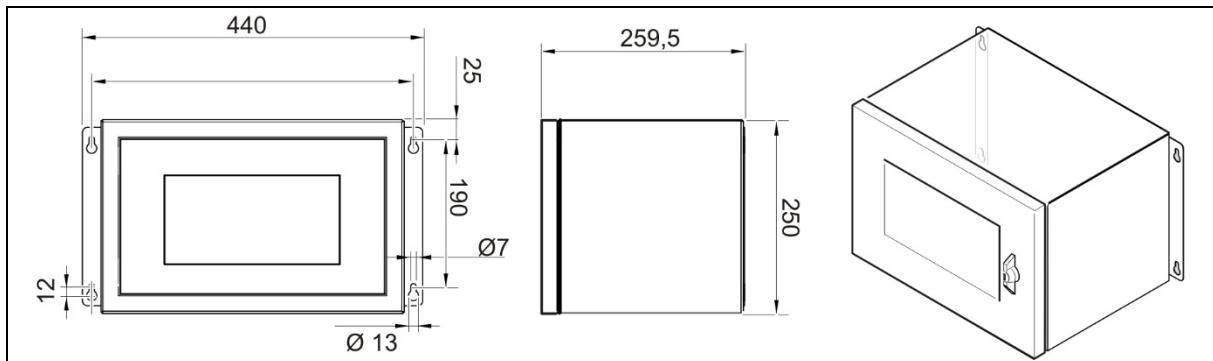
Signal	Pin
TxD	41
RxD	42
RTS	43
GND	44

The master/slave plug is not used by applications with terminal panels. The master/slave plug is contained in the purchase order terminal block (Mat. No. 59477). In the case of existing 19" sub-rack and retrofitting to LB 470, the master-slave plug (Part No. 64608) must be ordered separately.

Der Master/Slave Stecker wird bei Einbauvarianten mit Anschlussplatinen nicht benötigt. Der Master/Slave Stecker ist im Lieferumfang des Klemmenblocks (Mat. Nr. 59477) enthalten. Bei der Nachrüstung eines 19" Baugruppenträgers mit LB 470 Modulen muss der Master-Slave Stecker (Mat. Nr. 64608) gesondert bestellt werden.

3. Wall Housing

3. Wandgehäuse



TI-Abb. 10 Drawing wall housing
Zeichnung Wandgehäuse

Technical Data Technische Daten	
Max. assembly <i>Max. Bestückung</i>	2 Master with terminal board (master/master) ¹ 1 Master, 3 Slave with terminal board (master/slave) ¹ 2 Master with calmp blocks ² 2 Master mit Anschlussplatine (Master/Master) ¹ 1 Master, 3 Slave mit Anschlussplatine (Master/Slave) ¹ 2 Master mit Klemmenblöcken ²
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 ... 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 NRTL certification US/CAN

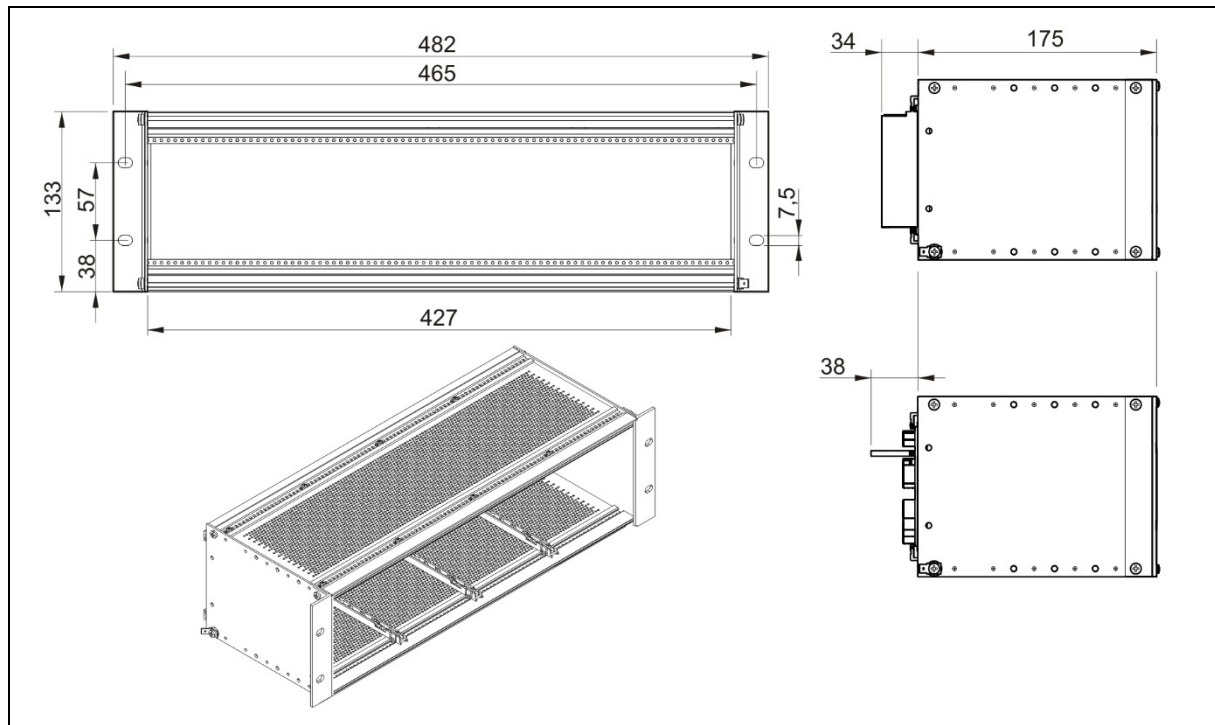
² No certification

¹ NRTL Zertifikat US/CAN

² Kein Zertifikat

4. 19" Subrack

4. 19" Baugruppenträger



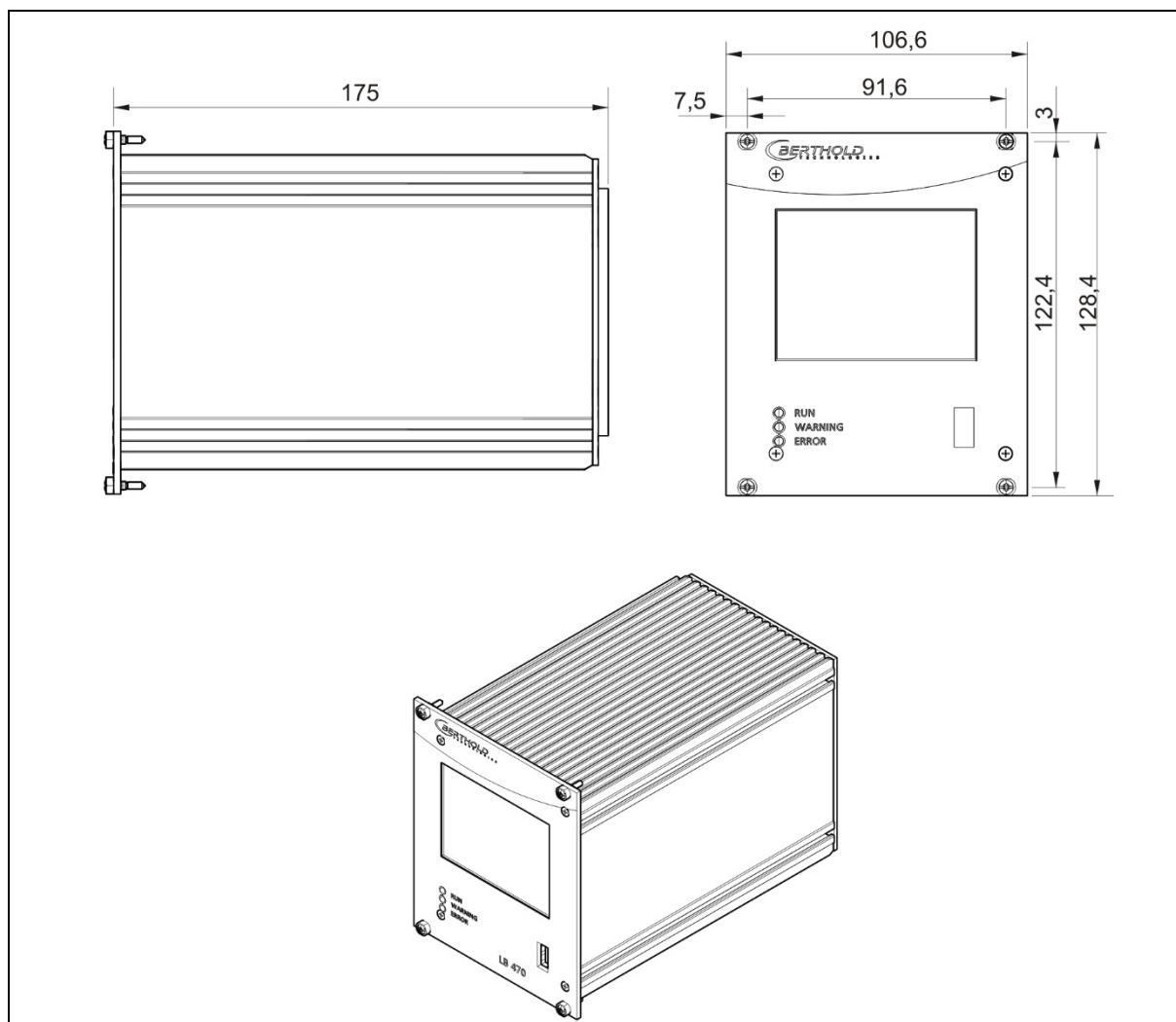
TI-Abb. 11 Drawing 19" subrack
Zeichnung 19" Baugruppenträger

Technical Data Technische Daten

Standard units <i>Normeinheiten</i>	3HE/84TE/5T
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. 12 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 sub-rack. -20°C ... +50°C nicht kondensierend. Direkte Sonneneinstrahlung ist zu vermeiden. Für eine ungehinderte Luftzirkulation ist zu sorgen.
Storage temperature Lagerungstemperatur	-30° C ... +80° C
Degree of protection Schutzart	IP20

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>
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 V AC 50/60 Hz (wide range input) +/- 10% 21 ... 32 V DC (24 V DC 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)

Interfaces

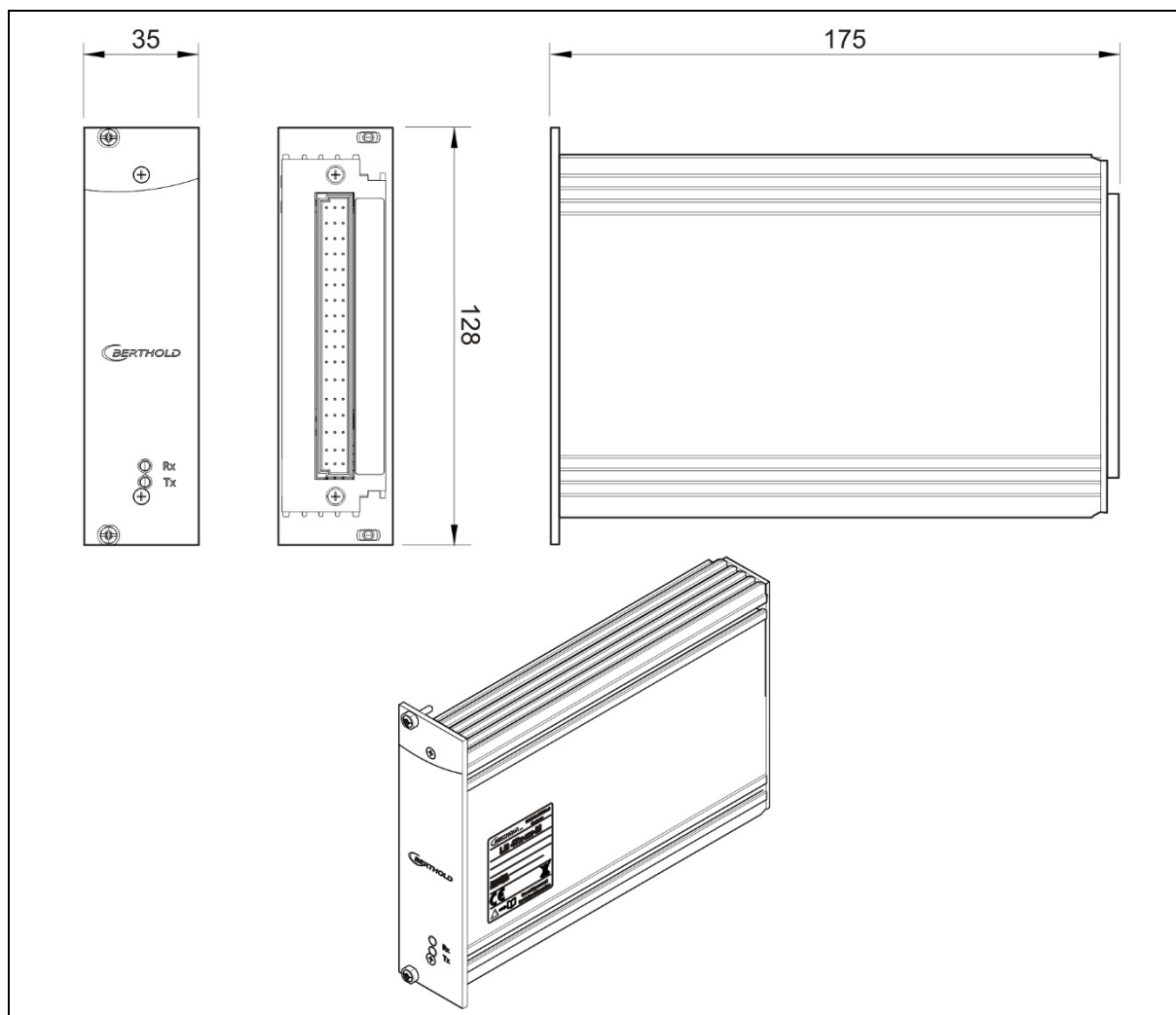
Schnittstellen

Current Output	4-20mA 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 ohms max. Burden when operating as a power source: 850 ohms. Internal monitoring of the loop current and additional error signalling by hardware on detection of a fault condition.
Stromausgang	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. 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.
Current input	4-20 mA (according to NAMUR recommendation NE 006 and NE 043) switchable via software on frequency input, electrically isolated (500 V). Internal resistance approx. 300 ohms max. input voltage: 24 V DC
Stromeingang	4-20 mA (nach Namur-Empfehlung NE 006 und NE 043) per Software umschaltbar auf Frequenzeingang, potentialgetrennt (500 V). Innenwiderstand ca. 300 Ohm max. Eingangsspannung: 24 V DC
Impulse input	Frequency 0 ... 100 kHz, $U_{\max} = 28 \text{ V}$, right angle signal form, low <1,5 V; high 4 - 28 V. Switchable to current input
Impulseingang	Frequenz 0 ... 100 kHz, $U_{\max} = 28 \text{ V}$, Rechteck-Signalform, Low <1,5 V; High 4 - 28 V. Umschaltbar auf Stromeingang
Digital outputs	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
Digitale Ausgänge	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
Digital inputs	2 x together electrically isolated (500 V) Switch between DigIn and GND, U_{outmax} approx. 24 V Function configurable via software
Digitale Eingänge	2 x gemeinsam potentialgetrennt (500 V), Schalter zwischen DigIn und GND, U_{outmax} ca. 24 V Funktion über Software konfigurierbar

External supply	Output voltage: 24 V DC Output current: max. 150 mA
<i>Externe Versorgung</i>	<i>Ausgangsspannung: 24 V DC Ausgangsstrom: max. 150 mA</i>
RS485	For master/master communication, 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>für Master/Master Kommunikation und Prüf-und Testzwecke. Nicht po- tentialgetrennt 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 opera- tion.
<i>Ethernet</i>	<i>RJ45-Buchse über Rückwand, 10 Mbit, DHCP unterstützt, max. 3 m Vorgesehen für Wartungszwecke. Nicht vorgesehen für den Dauerbe- trieb.</i>

6. Slave Module

6. Slave Modul



TI-Abb. 13 Drawing Slave Module
Zeichnung Slave Modul

Mechanical Data

Mechanische 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
Degree of protection <i>Schutzart</i>	IP20

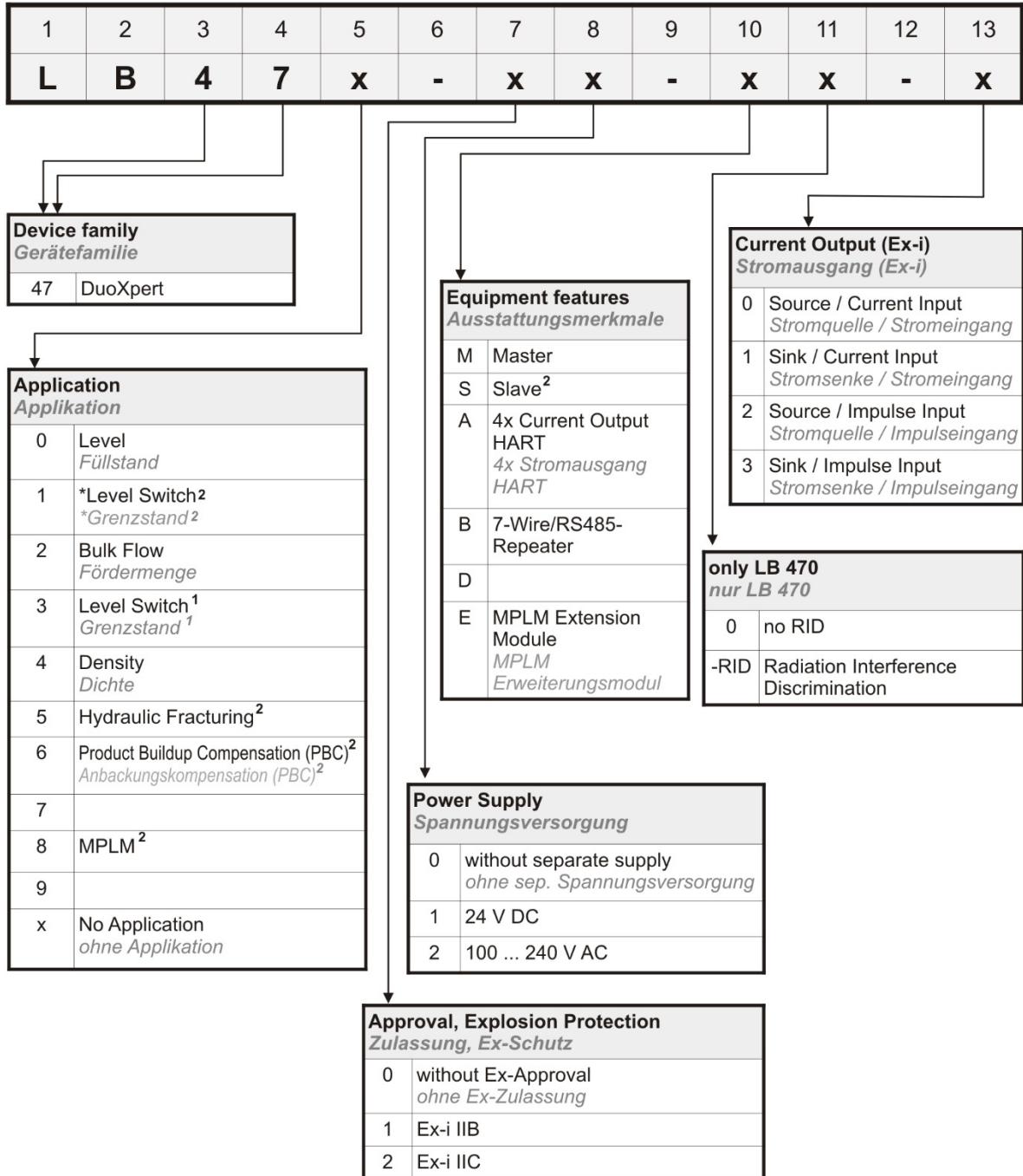
General ambient conditions <i>Allgemeine Umgebungsbedingungen</i>	Overvoltage category: II Pollution Degree: 2 Altitude: up to 2000 m Rel. humidity: 93% or less <i>Überspannungskategorie: II Verschmutzungsgrad: 2 Höhenlage: bis zu 2000 m Rel. Luftfeuchtigkeit: 93% oder weniger</i>
--	--

Electrical Data
Elektrische Daten

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 <i>Intern, 2 x 250 V, 1 A träge, 5x20 mm, 1500 A Abschaltvermögen IEC 60127-2</i>
Connections <i>Anschlüsse</i>	32-pin plug connector <i>32-polige Stiftleiste</i>

7. Number Key LB 47x

7. 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

8. 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-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.
WGL-Reg. No.

Stuttgart HRA 330991
BERTHOLD TECHNOLOGIES Verwaltungs-GmbH
Stuttgart HRB 331520
Herrn/Knaut, Dr. Dirk Mörmann
DE613050511
49038/03038
DE99468690

Sparkasse PF-CW
Volksbank
Commerzbank

75323 Bad Wildbad
75119 Pforzheim
75105 Pforzheim

Konto/Account No. 8 045 033 (BLZ 656 503 85)
Konto/Account No. 957 904 (BLZ 656 900 301)
Konto/Account No. 8 511 128 (BLZ 656 803 13)

SWIFT-BIC PZHS 3333
SWIFT-BIC VPM 2106
SWIFT-BIC DRES 2222

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

8. 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-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
Geschäftsleitung / Management
USt.-Id.-Nr. / VAT Reg. No.
Deutsche Steuernummer / German Tax No.
WEEE-Reg. Nr.

Stuttgart HRA 330981
BERTHOLD TECHNOLOGIES Verwaltungs-GmbH
Stuttgart HRB 331520
Horst Knaut, Dr. Dirk Mörmann
DE813050511
45038/08038
DE95468690

SparKasse R-CW 75323 Bad Wildbad
Volksbank 75119 Pforzheim
Commerzbank 75105 Pforzheim

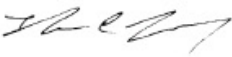
Konto/Account No. 8 043 003 (BLZ 566 500 85)
Konto/Account No. 557 004 (BLZ 666 800 00)
Konto/Account No. 8 511 120 (BLZ 566 800 12)

SWIFT-BIC: PZHSDE33
SWIFT-BIC: VPM0DE33
SWIFT-BIC: ORESDE33

IBAN: DE37 6665 9085 0008 0450 03
IBAN: DE85 6665 9000 0000 9520 04
IBAN: DE95 6668 9013 0651 1120 00

9. NTRL Certification US/CAN wall-mounted housing

9. 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 84119-2039 Tel (801) 972-6146 Fax (801) 972-8432

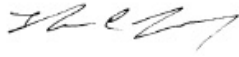
NFCC-002 Issue 2 May 2014



Page 3 of 3

10.NTRL Certifikate US/CAN DuoXpert LB 47x

10.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

11. Parts Overview

11. Übersicht Zubehör

Mat. No. Mat.-Nr.	Description Beschreibung
63284	LB 470-01-M0 Level Transmitter (Master, 24 VDC) LB 470-01-M0 Füllstandsmessgerät (Master, 24 VDC)
63283	LB 470-02-M0 Level Transmitter (Master, 100...240 VAC) LB 470-02-M0 Füllstandsmessgerät (Master, 100...240 VAC)
65092	LB 470-21-M0-0 Level Transmitter Ex-i (Master, 24 VDC, Source) LB 470-21-M0-0 Füllstandsmessgerät Ex-i (Master, 24 VDC, Stromquelle)
65091	LB 470-22-M0-0 Level Transmitter Ex-i (Master, 100...240 VAC, Source) LB 470-22-M0-0 Füllstandsmessgerät Ex-i (Master, 100...240 VAC, Stromquelle)
72364	LB 470-21-M0-0 Level Transmitter Ex-i (Master, 24 VDC, Sink) LB 470-21-M0-0 Füllstandsmessgerät Ex-i (Master, 24 VDC, Stromsenke)
72363	LB 470-22-M0-0 Level Transmitter Ex-i (Master, 100...240 VAC, Sink) LB 470-22-M0-0 Füllstandsmessgerät Ex-i (Master, 100...240 VAC, Stromsenke)
68640	LB 470-01-M-RID Level Transmitter with RID (Master, 24 VDC) LB 470-01-M-RID Füllstandsmessgerät mit RID (Master, 24 VDC)
68639	LB 470-02-M-RID Level Transmitter with RID (Master, 100...240 VAC) LB 470-02-M-RID Füllstandsmessgerät mit RID (Master, 100...240 VAC)
72059	LB 470-21-M-RID-0 Level Transmitter Ex-i with RID (Master, 24 VDC, Source) LB 470-21-M-RID-0 Füllstandsmessgerät Ex-i mit RID (Master, 24 VDC, Stromquelle)
72060	LB 470-22-M-RID-0 Level Transmitter Ex-i with RID (Master, 100...240 VAC, Source) LB 470-22-M-RID-0 Füllstandsmessgerät Ex-i mit RID (Master, 100...240 VAC, Stromquelle)
72366	LB 470-21-M-RID-0 Level Transmitter Ex-i with RID (Master, 24 VDC, Sink) LB 470-21-M-RID-0 Füllstandsmessgerät Ex-i mit RID (Master, 24 VDC, Stromsenke)
72365	LB 470-22-M-RID-0 Level Transmitter Ex-i with RID (Master, 100...240 VAC, Sink) LB 470-22-M-RID-0 Füllstandsmessgerät Ex-i mit RID (Master, 100...240 VAC, Stromsenke)
63286	LB 47x-01-S0 Slave (24 VDC) LB 47x-01-S0 Slave (24 VDC)
63285	LB 47x-02-S0 Slave (100...240 VAC) LB 47x-02-S0 Slave (100...240 VAC)

56925BA1	Operating Manual DuoSeries LB 470 Level, German <i>Betriebsanleitung Füllstand (deutsch)</i>
56925BA2	Operating Manual DuoSeries LB 470 Level, English <i>Betriebsanleitung Füllstand (englisch)</i>
56925-1BA1	Operating Manual DuoSeries LB 470RID Level, German <i>Betriebsanleitung DuoSeries LB 470RID Füllstand, Deutsch</i>
56925-1BA2	Operating Manual DuoSeries LB 470RID Level, English <i>Betriebsanleitung DuoSeries LB 470RID Füllstand, Englisch</i>
56925BA16	Safety Manual / Explosion Protection Manual LB 47x Ex-i (ATEX / IECEx), German <i>Sicherheitshandbuch / Explosionsschutzhandbuch LB 47x Ex-i (ATEX / IECEx), Deutsch</i>
56925BA26	Safety Manual / Explosion Protection Manual LB 47x Ex-i (ATEX / IECEx), English <i>Sicherheitshandbuch / Explosionsschutzhandbuch LB 47x Ex-i (ATEX / IECEx), Englisch</i>
63781	Wall-mounted Housing for LB 47x 1x Master / 3x Slave (24 VDC) <i>Wandgehäuse für LB 47x 1x Master / 3x Slave (24 VDC)</i>
63782	Wall-mounted Housing for LB 47x 1x Master / 3x Slave (110...240 VAC) <i>Wandgehäuse für LB 47x 1x Master / 3x Slave (110...240 VAC)</i>
63783	Wall-mounted Housing for 2x LB 47x Master (24 VDC) <i>Wandgehäuse für 2x Master LB 47x (24 VDC)</i>
63784	Wall-mounted Housing for 2x LB 47x Master (110...240 VAC) <i>Wandgehäuse für 2x LB 47x Master (110...240 VAC)</i>
64402	Wall-mounted Housing for 2x LB 47x Master (terminal blocks) <i>Wandgehäuse für 2x LB 47x Master (Klemmblöcke)</i>
72812	Wall-mounted Housing for 2x LB 47x Ex-i Master (with Ex-i terminal blocks) <i>Wandgehäuse für 2x LB 47x Ex-i Master (mit Ex-i Klemmblöcke)</i>
59493	19" rack for LB 47x, 2x (1x Master & up to 3x Slaves) <i>19"-Baugruppenträger für LB 47x, 2x (je 1x Master & bis zu 3x Slaves)</i>
59484	19" rack for LB 47x, 4x Master <i>19"-Baugruppenträger für LB 47x, 4x Master</i>
59481	19" rack for LB 47x, 3x (1x Master & 1x Slave) <i>19"-Baugruppenträger für LB 47x, 3x (je 1x Master & 1x Slave)</i>
64607	19" rack, 84 HP / 3 RU for use with terminal blocks <i>19"-Baugruppenträger, 84 TE / 3 HE für den Einsatz mit Klemmblöcken</i>
72051	19" rack Ex-i, 84 HP / 3 RU with 4x tension spring terminal blocks <i>19"-Baugruppenträger Ex-i, 84 TE / 3 HE mit 4x Zugfederklemmblöcken</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 for LB 47x Slave (with guide rails) <i>Klemmenblock für LB 47x Slave (mit Führungsschienen)</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>
64608	Master-slave plug and terminal assignment for LB 47x slaves e.g. when changing from LB 44x to LB 47x slaves <i>Master-Slave Stecker und Klemmenbelegung für LB 47x Slaves bei Umrüstung von LB 44x auf LB 47x Slaves</i>