



## Evaluation unit

Operating Manual 56925-8BA2

Rev. No.: 00, 07/2019 Embedded software version as of vers. 1.6.0 (CPU) and 1.6.0 (MU)



#### **BERTHOLD TECHNOLOGIES GmbH & Co. KG**

Calmbacher Str. 22 75323 Bad Wildbad, Germany www.berthold.com

Telephone +49 7081 177-0 Fax +49 7081 177-100 industry@berthold.com

## Table of Contents

1	About this Operating Manual	. 7
1.1	Applicable Documents	. 7
1.2	Some Prior Remarks	. 7
1.3	Storage Place	.7
1.4	larget Group	. /
1.5	Validity of the Operating Manual	ŏ. ە
1.0	Structure of the Operating Manual	. 0 0
1.7	Representation	ט. צ
1.0	Warning notes	. 0 9
1.9.1	Symbols Used in the Operating Manual	.9
1.9.2	Symbols Used on the Device	10
1.10	Conformity	11
2	Safety	13
2.1	Dangers and safety measures	13
2.2	Proper Use	13
2.3	Qualification of the Personnel	14
2.4	Operator's Obligations	15
2	System Description	17
31	Overview	17
3.7	Measuring Principle	18
3.3	System Components	18
3.3.1	Front/rear view EVU	19
3.3.2	Type plate	22
3.3.3	Software	22
3.4	Storage	22
		~~
4	Installation	23
4.1	General Instructions	23
4.2	Installation in the 19" subrack	23 71
4.3	Installation with clamp blocks	24
4.5.1		25
5	Electric Installation	27
5.1	General Instructions	27
5.1.1	Circuit Breaker	28
5.1.2	Cables and Lines	28
5.1.3	Cable Glands and Blanking Elements	29
5.1.4	Protective earth and equipotential bonding	29
5.1.5	EIA-485 (RS-485) Network	29
5.2	Wiring diagram	30
5.3 E 4	Electrical connection in the 19 subrack	3 I 22
5.4	Switching current output	22
6	Operation of the Software	35
6.1	System start	25
6.2	······································	35
6.2	EVU standard display	35 36
6.2 6.3	EVU standard display Navigation	35 36 37
6.2 6.3 6.3.1	EVU standard display Navigation Diagram display	35 36 37 37
6.2 6.3 6.3.1 6.4	EVU standard display Navigation Diagram display Status messages	35 36 37 37 37
6.2 6.3 6.3.1 6.4 6.5	EVU standard display Navigation Diagram display Status messages Event Reports	35 36 37 37 37 38
6.2 6.3 6.3.1 6.4 6.5 6.6	EVU standard display Navigation Diagram display Status messages Event Reports Input field	35 36 37 37 37 38 39

7.1	Menu Identification	42
7.1.1	Location	43
7.1.2	Device Information	44
7.2	Menu Access	46
7.3	Menu Setup	48
7.3.1	System (Date / Time, Interfaces, Units, Network, Reset, Repair Det. Software)	48
7.3.2	Sensors (only master EVU)	60
7.3.3	Interphase Densities (only master EVU)	62
7.3.4	Densities Overview (only master EVU)	62
7.3.5	Outputs	63
7.4	Menu Backup/Restore	71
7.4.1	Backup	71
7.4.2	Restore	73
8	Main Menu Diagnostics	75
8.1	Transmitter Temperature	75
8.2	Transmitter Events	
8.2.1	Traspmitter Event Log	76
822	Transm Event Overview	78
83	Change Log	79
84	Data Log (only master EVII)	80
8.5	Export service data (only master EVU)	82
9	Troubleshooting	83
91	Frror Search	05
9.7	Error Codes of the Evaluation Unit	84
9.2	System	07 8/
9.2.1	Application	04
9.2.2	Detector	00
9.2.5	BS /158 Interface	00 87
9.2.4	Process Connection	87
51215		
10	Maintenance and Repair	89
10.1	Replacing of Fuses	90
10.2	Cleaning	92
10.3	Data backup	93
11	Decommissioning	95
11.1	Decommissioning 19" subrack	95
11.2	Disposal of Measurement System	96
12	Appendix	97
12.1	Commissioning Guide	97
12.1.1	Underlying data	98
12.1.2	Implementation on site	100
12.2	Setup Protocol LB 478	103
12.3	Setup Protocol LB480 MPLM	105



## About this Operating Manual

## 1.1 Applicable Documents

This manual contains the following document:

• Technical Information Id. No. 56925TIM (see appendix)

## **1.2** Some Prior Remarks

The product is handed over to you by the manufacturer BERTHOLD TECHNOLOGIES GmbH & Co. KG (designated as Berthold in the following) in a complete and functionally reliable condition.

This operating manual illustrates how to:

- set up/install the product
- make electrical connections
- · perform measurements
- apply software settings
- carry out maintenance on the product
- fix errors
- · disassemble the product
- · dispose of the product.

Read these instructions thoroughly and completely before working with the product. We have tried to compile all information for safe and proper operation for you.

However, should questions arise which are not answered in this operating manual, please refer to Berthold.

## 1.3 Storage Place

This operating manual as well as all product-related documentation relevant to the respective application must be accessible at all times near the device.

## 1.4 Target Group

This operating manual is directed at qualified specialist personnel who are familiar with handling electrical and electronic assemblies as well as with communication and measuring techniques.

Specialist personnel refers to those who can assess the work assigned to them and recognise possible dangers through their specialist training, knowledge and experience as well as knowledge of the relevant regulations.



## 1.5 Validity of the Operating Manual

The operating manual is valid from the delivery of the Berthold product to the user until its disposal. Version and release date of this operating manual can be found in the bottom of each page. Modification services are not performed by the manufacturer Berthold.

The manufacturer reserves the right to make changes to this operating manual at any time without stating reasons.

#### NOTICE

The current revision of this operating manual replaces all previous versions.

## 1.6 Structure of the Operating Manual

This operating manual has been divided into chapters. The series of chapters should help you to familiarise yourself quickly and properly with the operation of the product.

## 1.7 Copyright

This operating manual contains copyright-protected information. None of the chapters may be copied or reproduced in any other form without prior authorisation from the manufacturer.

## 1.8 Representation

Identifier	Meaning	Example
Quotation mark	Field in the soft- wareuser interface	"Calibrate"
Vertical line	Path specification	Settings   Selection
Pointed brackets	Keys and buttons	<update></update>
Round brackets	Image reference	Connect the plug (fig. 1, item 1)

In the software description, the term "clicking" is used if a process is to be activated. This also refers to the pressing of a button or an area on the touch display if a mouse is not used for control.

## 1.9 Warning notes

Signal Word

Source:

Consequence: Prevention:

Warning notes are designed as follows:

Source and consequence Explanation, if required U Prevention In case of emergency

- Warning symbols: (warning triangle) draws attention to the hazard.
- Signal word: Indicates the severity of danger.
  - Specifies the type or source of danger.
    - Describes the consequences of non-compliance.
      - Specifies how the hazard can be avoided.
- In case of emergency: Specifies which actions are required in the event of the occurrence of risk.

## 1.9.1 Symbols Used in the Operating Manual

In this manual, warning instructions before instructions for action refer to risks of injury or damage to property. The hazard-prevention measures described must be observed.

#### 

Indicates an **imminent**, major hazard, which will certainly result in serious injuries or even death if the hazard is not avoided.

#### 

Indicates a **potential** hazard, which can result in serious injuries or even death if the hazard is not avoided.

## 

Refers to a **potentially dangerous** situation, which can result in medium or minor physical injuries or damages to property, if it is not avoided.

## NOTICE

If this information is not observed, deterioration in the operation and/or property damage may occur.

## IMPORTANT

Sections marked with this symbol point out important information on the product or on handling the product.

#### Tip



Provides tips on application and other useful information.



## 1.9.2 Symbols Used on the Device

#### Read the operating manual



#### Electrostatic discharge

Please note the handling instructions. Electrostatically endangered components. Please observe the instructions in this operating manual.

#### Protective earth connection

(느)

At this position, connect the protective earth conductor (PE).

#### Equipotential bonding connection

At this position, connect the equipotential bonding conductor.

#### Direct voltage

The device is operated with direct voltage and may only be connected with a direct voltage source.

#### Alternating voltage

The device is operated with alternating voltage and may only be connected with an alternating voltage source.

#### No domestic waste



The electric product must not be disposed of in domestic waste.



## 1.10 Conformity

The company Berthold hereby declares in its sole responsibility that the design of this product, which is brought to the market by Berthold, complies with relevant EU directives stated in the original declaration of conformity.

This statement shall become void in the case of changes not authorised by Berthold or improper use.

For the original declaration of conformity, please refer to Declaration of Conformity in the document "Technical Information".





## 2 Safety

## 2.1 Dangers and safety measures

- Read these instructions thoroughly and completely before working with the product.
- Store the instructions where they are accessible for all users at all times.

## 2.2 Proper Use

The evaluation unit DuoXpert LB 478 MPLM (EVU) measures the level together with compatible detectors and an appropriate radiation source and may only be used for this purpose.

#### The following constitutes proper use:

- Adhering strictly to the instructions and operation sequences and not undertaking any different, unauthorised practices which could endanger your safety and the operational reliability of the EVU!
- · Observing the given safety instructions!
- · Carrying out the prescribed maintenance measures or having them carried out for you!
- Only use accessories and spare parts from Berthold.



#### Improper use to be prevented:

- Failing to observe the specified safety instructions and instructions for the operation, maintenance and disposal in the operating manual.
- · Any non-compliance with the present operating manual for the supplied products.
- Applying conditions and requirements which do not conform to those stated in the technical documents, data sheets, operation manuals and assembly instructions and other specific guidelines of the manufacturer.
- Use of the product if parts of it are damaged or corroded. This also applies for seals and used cables.
- Restructuring or changing the system components.
- The product is not suitable for use in potentially explosive areas and may therefore not be operated in such areas. The product is not explosion-proof.
- · Operation ...
  - o in a state where live parts are accessible.
  - in a wall housing with inadequately sealed glands and / or insufficiently tightened or damaged cable glands.
- · Operation without the safety precautions provided by the manufacturer.
- · Manipulation or avoidance of existing safety equipment.

Berthold shall only accept liability for / guarantee the correspondence of the device to its publicised specifications.

If the product is used in a way which is not described in the present operating manual, the device's protection is compromised and the warranty claim becomes invalid.



The device is not approved according to IEC 61508 "Functional safety of safety-related electric/electronic/programmable electronic systems".

## 2.3 Qualification of the Personnel

#### NOTICE

A minimum requirement for all work on or with the product would be employees with general knowledge who are instructed by an expert or authorised person.

At different parts in this operating manual, reference is made to groups of people with certain qualifications who can be entrusted with different tasks during installation, operation and maintenance.

These three groups of people are:

- Employees with General Knowledge
- Experts
- Authorised Persons.



#### Employees with General Knowledge

#### NOTICE

i

Employees with general knowledge must always be guided by an expert at the very least. When dealing with radioactive substances, a radiation safety officer must also be consulted.

Employees with general knowledge are e.g. technicians or welders, who can undertake different tasks during the transportation, assembly and installation of the product under the guidance of an authorised person. This can also refer to construction site personnel. The persons in question must have experience in handling the product.

#### **Experts**

Experts are persons who have sufficient knowledge in the required area due to their specialist training and who are familiar with the relevant national health and safety regulations, accident prevention regulations, guidelines and recognised technical rules.

Expert personnel must be capable of safely assessing the results of their work and they must be familiar with the content of this operating manual.

#### **Authorised Persons**

Authorised persons are those who are either designated for the corresponding task due to legal regulations or those who have been authorised by Berthold for particular tasks. When dealing with radioactive materials, a radiation safety officer must also be consulted.

## 2.4 Operator's Obligations

The operator of the product must regularly train his personnel in the following topics:

- Observation and use of the operating manual and the legal provisions.
- · Intended operation of the product.
- Observation of the plant security instructions and the operating instructions of the operator.
- Regular monitoring/maintenance of the product.





# **3** System Description

## 3.1 Overview

Multiphase Level Measurement (MPLM) systems can be used to determine the interface position between liquid phases or the height of different product layers.

Berthold's EmulsionSENS is used to determine multiphase layers and consists of detectors and evaluation units.

A complete measuring system consists of the following components, which are, according to customer's order, individually assembled:

- Evaluation unit DuoXpert LB 478 MPLM (master and extension module)
- EmulsionSENS Control Cabinet
- Source
- · Shields
- Point detector / rod detector(s) LB 480

These instructions concern the operation of the evaluation unit DuoXpert LB 478 MPLM (Fig. 1, item 7 and item 8). The operation of other system components is part of the independent instructions of the respective system components. Depending on the measuring task and the composition of the measuring product and containers, different arrangements and system components are required.



Fig. 1 Example measurement arrangement

BERTHOLD

## 3.2 Measuring Principle

The Multiphase Level Measurement system EmulsionSENS consists of radiometric density devices LB 480 at different elevations through the measurement span (see Figure 1) and evaluation units LB 478 MPLM. Such an arrangement allows accurate measurement of the density at each of these positions and provides a density profile of the fluids as they separate inside the vessel.

The density measurement is based on the through-beam method and uses the physical laws of absorption of radiation by matter. The resulting measurement effect is the ratio I/I0 between the unattenuated radiation I0 and the radiation I weakened by the measured medium.

The density values of the individual detectors LB 480 are sent to the evaluation unit LB 478 via RS485. By means of an algorithm implemented in the evaluation unit LB 478, interface levels can be calculated from the measured density values. This allows the operator to see where the top of a layer is located as well as the bottom of a layer.



Further information on the functional principle of the detectors can be found in the operating manual of the detector.

## 3.3 System Components



Fig. 2 System components

## 3.3.1 Front/rear view EVU

#### Front view EVU

The following control elements are found on the front of the EVU:

- · LEDs for status display of individual operating states
- · 3.5" Touch display
- USB port.



Fig. 3 Front view of the EVU

#### **Operation Display / Touch Screen**

The EVU is operated via the touch screen. Alternatively, the device can be connected to a mouse at the USB port. The device can also be operated by remote control (see chap. 7.3.1).

#### NOTICE

#### Damage to the touch screen

Pointed or sharp objects can damage the plastic surface of the touchscreen.

 Operate the touch screen only with your fingers or with a touch pen or connect a mouse. The mouse pointer automatically becomes visible when a mouse is inserted into the USB port.



i

## Status displays of the master EVU

The LEDs (Fig. 3, items 2-4) below the touch display show the current operating status of the master EVU.

Display LED	Description
RUN WARNING ERROR	<b>RUN</b> This LED lights up green if the device is in operation and fault-free. The current measurement is carried out.
	<b>RUN (flashing)</b> The RUN LED flashes green while the measurement is (held) in the STOP state by user actions (e.g. stop func- tion, simulation mode, plateau recording).
O RUN O WARNING O ERROR	WARNING This LED lights up yellow when a system event of the type "Outside of specification", "Mainte- nance required" or "Function check" is present. All system events are described in chapter 9.2.
O RUN WARNING ERROR	<b>ERROR</b> This LED lights up red if a system event of the type "Fail- ure" is present. The current measurement is retained. Check the Device Setup. All system events are described in chapter 9.2.
- O-RUN - O-WARNING - O-ERROR	<b>RUN / WARNING / ERROR flashing</b> All three LEDS flash during the system test which is per- formed as part of the start up process.



#### Rear view EVU

The following connections are located on the back of the EVU:

- · Master/slave connector, 4-pin
- · RJ45 socket for Ethernet
- · 32-pin plug connector



Fig. 4 Rear view EVU



## 3.3.2Type plate



Fig. 5 Type plate

## 3.3.3 Software

The EVU is delivered with pre-installed software. The revision status (version) of the software can be seen on the screen display when starting up the EVU or in the menu "Device information" (Chapt. 7.1.2).

This operating manual describes the software version 1.6.0 (Control Unit / CU) and 1.6.0 (Measurement Unit / MU).

## 3.4 Storage

Keep devices in a dry (no condensation), dark (no direct sunlight), clean and lockable room. Stay within the temperature range for storage (see document "Technical Information").



## **4** Installation

## 4.1 General Instructions

The applicable national regulations of the country of use have to be observed! Repair and maintenance on the devices may only be performed by experts (see chapter 2.3). In case of doubt, the complete device must be returned to Berthold for repair.

#### NOTICE

The Evaluation unit is not explosion protected and is not designed for hazardous environments.

Only mounting accessories approved by Berthold should be used for installation of the devices. The device should only be operated if firmly installed.

## 4.2 Unpacking/Scope of Delivery

The product will be delivered completely configured according to the purchase order. Check your delivery for completeness and damage according to your order. Please report missing, defective or incorrect parts immediately.



## 4.3 Installation in the 19" subrack

The 19" subrack can be equipped differently, depending on requirements. The rear clamp blocks (Fig. 6, item 3) are used for the electrical connection.

#### NOTICE

The 19" subrack may only be installed in a dry environment.

The subrack is installed in a 19" control cabinet or a control panel (switchboard). The 4 side holes that should be provided with fitting screws are used to fasten the subrack.

## NOTICE



i

The EVU is delivered equipped, depending on the order. The installation of the modules is only necessary if:

- u Another measurement channel is to be fitted
- u A defective module is to be replaced



## 4.3.1 Installation with clamp blocks

#### 🛆 DANGER



- u Installation/maintenance may only be carried out if the device has been deenergised.
- u Test of absence of harmful voltages when the front side is open.

In case of an electric shock, carry out first aid measures and immediately call an emergency service.

#### NOTICE

i

The master's EVUs / extension modules must be secured against pulling out by fixing screws (Fig. 6, item 6).

The device must be disconnected from the mains voltage before it is pulled out.



Fig. 6 19" subrack with clamp block

- 1. Set master EVU / extension module (Fig. 6, item 1, item 2) in the guide rails (Fig. 6, item 5).
- **2.** Carefully slide module into the subrack until the plug connector is inserted into the clamp block.
- **3.** Tighten fixing screws (Fig. 6, item 6).
- u The EVU is correctly inserted and can be connected.





# 5 Electric Installation

## General Instructions

#### 



5.1

Danger to life from electric shock!

- u The installation may only be carried out by a qualified electrician.
- u Please adhere to the relevant safety regulations.
- u Open the housing only in a dry environment and for installation, maintenance and servicing.
- During installation and servicing on the hardware as well as during wiring of the detector, the measuring system, connected relay contacts and all inputs and outputs must be de-energised.
- u Connect only devices onto the product that comply with the applicable safety standards.

In case of an electric shock, carry out first aid measures and immediately call an emergency service.

#### NOTICE

Apply the voltage of the specified and marked range only!

#### NOTICE

Ť

The relay of the LB 478 MPLM can only switch low voltages. Please note the specifications in appendix Technical Information.

The voltage parameters of all devices connected to the outputs of the product (e.g. relay circuit, RS-485, current output) must comply with the limit values of the safety directives for electrical measurement, control, regulation and laboratory devices (DIN EN 61010-1) and be equipped with double or reinforced insulation.

These protective measures are necessary to avoid the risk of contact with life-threatening voltages.

Changing the installation without precise knowledge of this operating manual is not permitted.

#### General important points for installation

- u Connect the earth conductor.
- u Ground the housing.
- u Please observe the information signs on the devices.



## 5.1.1 Circuit Breaker

A circuit breaker according to DIN EN 61010-1

- · must be available,
- must be easily accessible for the maintenance personnel and
- is to be included in the company-internal documentation.

The master EVUs / extension modules are not equipped with a separate ON/OFF switch to connect or disconnect the voltage supply. Make sure that the system can be de-energised via the external power supply.

The circuit breaker can be installed as an automatic fuse or switch and has to comply with the requirements according to IEC 947-1 and IEC 947-3. If a fuse is applied, it must not be triggered under a current strength of 4 A per device.

#### IMPORTANT

The circuit breaker must be located near the device and be properly marked as belonging to it.

## 5.1.2 Cables and Lines

- u Lines are to be connected with special care.
- u Connection lines and routing must comply with the applicable regulations.
- u When routing the cables, make sure that the cable insulation cannot be mechanically damaged by sharp edges or movable metal parts.
- u Use the approved Berthold cable or a cable with equivalent specifications for the connection.

For intrinsically safe systems, the detector must be connected to the equipotential bonding of the system. The detector is connected via a 2-core (0.2 ... 2.5 mm<sup>2</sup>) cable with approx. 5 ... 10 mm diameter. A screened cable can be used in systems with extremely strong electrical noise. The screen may only be laid out on one side of the detector. The maximum cable length depends on the cable resistance, which may not exceed a total (there and back) of 40 ohms. For standard cables from Berthold (Id. no. 32024), this results in a cable length of 1000m, from the evaluation unit to the detector. For intrinsically safe systems, the maximum permissible inductance and capacitance of the cable should be considered to the max. 40 ohm.

When routing the connection lines, make sure that

- u no dirt or moisture reaches the connection room,
- u the conductors are not damaged when the cable insulation is removed,
- u the conductor insulation or the sleeve of the wire end ferrules reach into the housing of the terminal unit,
- u blank, conductive segments of the lines (e.g. wires of a litz wire) do not reach outside the terminal unit,
- u the wire end ferrule or the stripped wire have a length of 8 mm so that the wire is held securely in the clamp,
- u the line insulation reaches into the sleeve of the wire end ferrule if these components are used,
- u the admissible minimum bending radius for the respective line cross-section is not exceeded and



- u the cables are laid out in a strain-relieved and friction-free manner.
- u Only use cables whose diameters are approved for the respective cable gland. The cables must comply with the requirements and cross-sections specified in the technical data.
- u The connected cables must be appropriate for temperatures of at least 10°C above the maximum admissible ambient temperature.

## 5.1.3 Cable Glands and Blanking Elements

- u The feeding of cables into the wall housing is only permitted via a cable entry.
- u Cable glands must be suitable for the respective application.
- u All cable glands must be assembled according to manufacturer's instructions and be tightened to the appropriate tightening torque.
- u Cable glands that are not required for installation must be covered with suitable blanking elements.
- u Line cross-sections must comply with the respectively used cables.
- u Cable bushings and blanking elements must comply with the applicable IP protection class and with the requirements for the operational environment.
- u We recommend ordering missing cable glands, blanking elements or adapters from Berthold.

## 5.1.4 **Protective earth and equipotential bonding**

- u The protective earth conductor has to be connected to the terminals marked with "PE".
- u The housing must be connected to local equipotential bonding.

## 5.1.5 **EIA-485 (RS-485) Network**

For integration of EVU units into an EIA-485 (RS-485) network, all participants must be connected one after the other in the configuration Master-Master. Star connection is not permitted.

The first and last station (physical, independent of the master's position) on the network needs a terminating resistor of 121  $\Omega$ .



## 5.2 Wiring diagram

The detectors LB 480 are connected to the evaluation unit via the RS485 terminals. Connect the communication lines according to assignment to clamp blocks (Fig. 7). Observe the circuit diagrams of the EVU in the document "Technical Information".

#### NOTICE

Observe the circuit diagrams of the control unit and the detectors for connecting.



Fig. 7 Communication connection detector-EVU



## 5.3 Electrical connection in the 19" subrack

The electrical connections are made via clamp blocks. These clamp blocks are already installed in the 19" subracks and are also available as an optional accessory.

#### 



#### Danger to life from electric shock!

- u The installation may only be carried out by a qualified electrician.
- u Please adhere to the relevant safety regulations.
- u Installation/maintenance may only be carried out if the device has been deenergised.
- u Only open the device when free of voltage.

In case of an electric shock, carry out first aid measures and immediately call an emergency service.

#### IMPORTANT

In the case of applications with clamp blocks a contact protection must be provided by the customer when voltage is applied. The cable connections of clamp blocks have to be in accordance with IEC 61010-1 (2010).



Fig. 8 Electrical connection in the 19" component rack

- 1. Connect the lines to the clamp blocks according to assignment (see chapt.5.2 and the document "Technical Information"), note in addition the wiring diagrams of the control unit.
- 2. Open the clamping screw (Fig. 8, item 1) and insert the stripped wire (min. 8 mm).
- u The terminal connections are designed for wires with a conductor cross-section from 0.2 mm<sup>2</sup> to 2.5 mm<sup>2</sup>.
- 3. Screw the clamping screws with a tightening torque of 0.4 0.5 Nm.



- 4. Plug the network plug into the RJ45 socket (Fig. 8, item 3) (optional).
- 5. Check the correct connection of the PE conductor (Fig. 8, item 4).



#### NOTICE

Note the specification relating to Protective earth and equipotential bonding in chapter 5.1.4 .

u The connection was made correctly.



## 5.4 Switching current output

Switching between "SOURCE" (active) and "SINK" (passive) is possible using the slide switch on the I/O board. Factory setting EVU is delivered in "SOURCE" mode.

#### ▲ DANGER



Danger to life from electric shock!

- u The switching may only be carried out by a qualified electrician.
- u Please adhere to the relevant safety regulations.
- u Switching may only be carried out if the device has been de-energised.
- u Only open the device when free of voltage.

In case of an electric shock, carry out first aid measures and immediately call an emergency service.



Fig. 9 Switching of the current output

- 1. Loosen the four sunken screws on the front side of the EVU (Fig. 9, item 1).
- 1. Pull out the housing (Fig. 9, item 2) carefully.
- 2. Slide the switch (Fig. 9, item 3) to position a for "active" (SOURCE), to position p for "passive" (SINK).
- **3.** Carefully insert the front panel into the housing. Pay attention to the correct guide rail!
- **4.** Screw the front panel to the housing (Fig. 9, item 2) with the four screws (Fig. 9, item 1).
- u The switching has been carried out correctly.







## **Operation of the Software**

## 6.1 System start



Fig. 10 Start screens with display of the software version

## System start with invalid application software

A different menu structure is present in this mode.

2015-03 11:32:3	3-20 5	
Nc	o valid application software found.	_
c	complete a software update.	Ľ
		>

Fig. 11 Start screen (Invalid application software)

#### IMPORTANT

The communication between the sensor and EVU is limited to 1200 baud. Accordingly, there is a load time for data that are retrieved in the detector.

All set values as well as the calibration data are stored in a non-volatile memory.

The real-time clock for date and time is buffered via a capacitor and continues to run for up to approx. 4 weeks even when the device is switched off. See also "Date and time" on page 49.



## 6.2 EVU standard display

#### **IMPORTANT**

i

Changing the language of the user interface is changed in menu Device Setup | Setup | System | Interfaces | Languages.



Fig. 12 Standard display of the EVU

Level scale	The scale range of the value is displayed.
Date and Time	The date and time displayed in the menu Device Setup   Setup   System   Date/Time was set.
Process value	The current process value with the set unit is displayed.
Selected Interface	The set interface from the menu Device Setup   Setup   Outputs   Analog Outputs
Unit	The set unit from the menu Device Setup   Setup   System   Units is displayed
Button "Menu"	The main menu opens
Button "Diagram"	The view changes to diagram and monotonized densities
Status information	The current system status is displayed.
Location	Name of the measuring point in the menu De- vice Setup   Identification.


# 6.3 Navigation



Fig. 13 Icons for navigation

# 6.3.1 **Diagram display**

Clicking the symbol (Fig. 12, item 6) changes the view to the display monotonized of densities (Fig. 14, item 4; only available for the master unit). Clicking the symbol (Fig. 13, item 2) changes the view to the diagram display. The diagram of the extension modules shows only the time course of the selected interface, whereas the master unit shows the time c ourse of all interfaces. The arrow keys (Fig. 14, item 2) are used to view the colors (Fig. 14, item 1) of the interfaces lines.



Fig. 14 Diagram display of the EVU

# 6.4 Status messages



#### Fig. 15 Status information



# 6.5 Event Reports

Events are displayed in the standard display and in the submenus as a symbol. All events are displayed on the main screen. A specific "D" (for detector) indicates that a detector has an event, the prefix "M" (for measuring unit) indicates that there is an event in the LB 478 transmitter. In the event of a detector fault, the operating manual of the detector must be observed. Only the event with the highest priority will be displayed. Refer to the menus Transmitter Events (chapter 8.2) and Detector Event Log "Detector-Service" (chapter 7.3.2) for additional information.



Fig. 16 Event reports

- 1. Click on the icon (Fig. 16, item 1, item 2) to display detailed information about the event.
- 2. Click the button <Acknowledge> to confirm an event that requires a manual confirmation.
- u The event description indicates the next event or reports no further events.
- 3. Click <Close> to return to the submenu or to the standard display.
- u The icon disappears from the status information.

#### **IMPORTANT**



If you click the button <Close>, the event message is closed, the icon continues to be displayed.



# 6.6 Input field



The input field appears by clicking on the blue display panels.

Description: 1234567890 qwertyuiop	Description:
asdfghjkl Î=zxcvbnm← 123esc ←→ ↓	$\begin{array}{c c} 4 & 5 & 6 \\ \hline 1 & 2 \\ \end{array} \begin{array}{c} 3 & 0 \\ \hline \end{array} \begin{array}{c} \bullet \\ \bullet \\ \bullet \end{array} \begin{array}{c} \bullet \\ \bullet \\ \bullet \end{array} \begin{array}{c} \bullet \\ \bullet \\ \bullet \\ \bullet \end{array} \begin{array}{c} \bullet \\ \bullet \\ \bullet \\ \bullet \end{array} \begin{array}{c} \bullet \\ \bullet \\ \bullet \\ \bullet \end{array} \begin{array}{c} \bullet \\ \bullet \\ \bullet \\ \bullet \\ \bullet \end{array} \begin{array}{c} \bullet \\ \bullet \\ \bullet \\ \bullet \\ \bullet \\ \bullet \end{array} \begin{array}{c} \bullet \\ \bullet \\ \bullet \\ \bullet \\ \bullet \\ \bullet \end{array} \begin{array}{c} \bullet \\ \bullet $
1 2 3 4     5 6     7 8	9 10 11 12
1 Input line	7 Input key
2 Shift key for numbers	8 Delete key
3 Shift to upper case characters	9 Number field
4 Escape key	10 Home (item1) key
5 Keypad	11 End key
6 Navigation buttons	12 Shift key for letters

Fig. 17 Screen keyboard





# 7 Main Menu Device Setup

#### Menu structure master EVU



#### Menu structure extension modules





Fig. 18 Menu "Main Menu", "Device Setup"

# 7.1 Menu Identification

Device Setup | Identification

You can make the following settings and read information in the Identification menu:

- Display and change the location name
- · Display of hardware and software information
- · Display of the device ID
- · Perform software update



Fig. 19 Menu "Identification"



# 7.1.1 Location

Device Setup | Identification | Location

The location of the evaluation unit is displayed in the "Location" menu. The name can only be edited (7.2 Menu Access) in the access level "Standard". The Location is displayed on the EVU standard display (Fig. 12, item 8).

Device location:	G	Location		
< Edit2		LOCATION	Device location:	
	<	I	Edit	

Fig. 20 Device location

- 1. Click <Edit> (Fig. 20, item 2) to open the input field.
- 2. Enter a location name for the evaluation unit.
- **3.** Confirm with the Enter key.
- u The name has been changed.



# 7.1.2 Device Information

Device Setup | Identification | Device Information

Information about hardware and software of the evaluation unit are displayed in the window "Device Information".



Fig. 21 Device Information (example)

#### NOTICE

During an update where the first or second digit of the version changes, it is necessary to reset the EVU to factory settings.

# NOTICE

Settings are deleted!

- u Carry out a backup of the measuring channel settings before resetting and the update of the EVU (10.3 Data backup).
- u The secured settings should then be imported after the successful software update.

#### Tip



The current software versions can be downloaded from the Berthold website (www.berthold.com).



#### Perform CU update

1. Save the current update file of the CU software on a USB storage device.

#### IMPORTANT

In order for the system to detect the update file it must not be located in an index in the USB storage device.

- 2. Connect a USB storage device to the device (Fig. 3, item 5).
- **3.** The USB storage device is recognised by the system after a few seconds and the <CU Update> (Fig. 21, item 1) button can be clicked.
- **4.** Click on the button <CU Update> (Fig. 21, item 1) and select the appropriate file.
- u The update is performed.
- 5. Measurement is interrupted.
- 6. Click the Button <Restart> to reboot the EVU.

#### Perform MU update

- 1. Save the current update file of the MU software on a USB storage device.
- 2. Connect a USB storage device to the device (Fig. 3, item 5).
- 3. The USB storage device is recognised by the system after a few seconds and the <MU Update> (Fig. 21, item 1) button can be clicked.
- Click on the button <MU Update> (Fig. 21, item 2) and select the appropriate file.
- u The update is performed.
- 5. Measurement is interrupted.
- 6. Click the Button <Restart> to reboot the EVU.

#### NOTICE

Berthold recommends calibrating the current outputs whenever a module has been installed/replaced or if a software update has been carried out.



# 7.2 Menu Access

Device Setup | Access

You can set the user rights via the user levels and assign passwords in the submenu Access. After assigning a password the system is protected against unauthorized manipulation of the parameters.

#### IMPORTANT



Make sure that the password is known to you before you select the "Basic" access level and lock the device. The same applies if you set "Automatically log-out".

If you do not know the password, you will not be able to unlock the device! If in doubt, enter a new password with "Change Password".

(	Access	
	Access level	٦
	O Admin	
	Standard	+
	O Basic	Ļ
		1
	Automatic logour	
	1 2	
1	Check box "Automatic logout"	
2	Change <change password=""></change>	
3	User level "Basic"	
4	User level "Standard"	
5	User level "Admin"	

Fig. 22 Access

```
Automatic logout Activating the selection box (Fig. 22 item 1) au-
tomatically resets the access level Standard to
"Basic" when the system changes to the stand-
ard display after the timeout (chapter 7.3.1).
```

The following user levels are available to you:

User Level Basic	Select "Basic" to lock the device against un- wanted manipulation. After the device has been locked, it is still possible to read all data, but changes to the data are no longer possible.		
	If "Basic" is already set, then the device is al- ready in the locked state.		
	To unlock the device, select the access level "Standard".		
User Level Standard	If the device is in the "Standard" access level, all parameters are accessible and can be changed.		

56925-8BA2 Rev.00, 07/2019



If the device is in the "Basic" access level (locked), you can unlock the device with the "Standard" access level. The password will be asked for. You can unlock the device only if you enter the correct password.

User Level Admin

This access level is only intended for the system management by Berthold.

#### NOTICE

Incorrect measurement and calibration parameters can be set through unauthorised inputs. These can possibly lead to production losses and damage in the system.

Protect the measuring system from unauthorised entries with a password and activate the function "Automatic logout".

#### Assign / change password

To set or change a password, select "Standard" (Fig. 22, item 4) and click on <Change password> (Fig. 22, item 2) to open the input field.



Fig. 23 Change Password

- 1. Click on the text field (Fig. 23, item 1.) to open the input field.
- 2. Enter a password (case-sensitive!).
- **3.** Confirm with the Enter key.
- 4. Click <OK> (Fig. 23, item 2) to confirm.
- u The password has been set / changed.



# 7.3 Menu Setup

Device Setup | Setup

<b>D</b>	Setup
	System
	Sensors
	Interface Densities
	Densities Overview
<	Outputs

Fig. 24 Menu "Setup"

# 7.3.1 System (Date / Time, Interfaces, Units, Network, Reset, Repair Det. Software)

Device Setup | Setup | System

3	System
	 Date / Time
	Interfaces
	Units
	Network
	Reset Device
<	Repair Detector Software

Fig. 25 Submenu "System"



#### Set Date and Time

Device Setup | Setup | System | Date / Time

#### **IMPORTANT**



The date and time must always be set correctly so that all records (log files) have the correct metadata.

The correct date is also indispensable for the decay compensation.



Fig. 26 Date / Time, calendar

- 1. Click on the arrow key (Fig. 26, item 5) in order to set the date.
- u The calendar is opened (Fig. 26, item 6).
- 2. Click on the year number (Fig. 26, item 8) in order to enter the year.
- 3. Set the month (Fig. 26, item 7) by clicking on the arrow keys.
- 4. Set the day by clicking on a number in the calendar.
- 5. Change the time by clicking on the arrow keys (Fig. 26, item 4).
- 6. Click on <Apply> (Fig. 26, item 3), to accept the date and time settings.
- u The date and time are set.

#### NOTICE

The real-time clock for date and time is buffered via a capacitor and continues to run for up to approx. 4 weeks even when the device is switched off.

u If the device has been out of operation for more than 4 weeks, error M116 appears. The date and time must then be reset.



#### Interfaces

Device setup | Setup | System | Interfaces

You can adjust the following settings in the submenu "Interfaces" (Fig. 27):

- · Local Display
  - o Brightness / Touch
  - o Input / Output
- Language
- · CE Remote Control

6	Interfaces
	Local Display
	Language
	CE Remote Control
<	

Fig. 27 Menu "Interfaces"

#### Local Display

Device setup | Setup | System | Interfaces | Local Display



Fig. 28 Submenu "Local Display"



#### **Brightness / Timeout**

Device Setup | Setup | System | interfaces | Local Display | Brightness / Timeout



Fig. 29 Brightness / Time out settings

"Time out" refers to the period of time during which the display is not operated. The value "Time out display dimming" cannot be set greater than the value at "Time out display shutdown".

Display dimming	In the field Display dimming, clicking the input fields allows the entering of the brightness (Fig. 29, item 2) in percent, that is set after expiry of the time (Fig. 29, item 1).
Display shutdown	In the field Display shutdown, clicking the input fields allows the entering of the brightness (Fig. 29, item 4) in percent, that is set after expiry of the time (Fig. 29, item 3).
Menu Timeout	Under "Menu Timeout" clicking on the input field (Fig. 29, item 5) changes the time period (seconds) in which the menu view changes to the standard view.



#### Input / Touch

Device Setup | Setup | System | Interfaces | Local Display | Input / Touch



Fig. 30 Input/ Touch

#### Calibrate touch screen

The calibration may only be carried out with direct skin contact. Take gloves or any other protective equipment off your hands.





- 1. Click on <Calibrate touch screen>.
- u The calibration screen opens.
- 2. Press the middle of the displayed cross with your finger.
- u If you take your finger off the cross again, the cross jumps to the top left corner.
- **3.** Repeat the process until the cross is no longer displayed and the calibration is finished.
- **4.** Confirm the calibration by clicking on the white screen to go back to "Input/Touch"
- 5. Execute a restart of the EVU after prompting.
- u The screen calibration was performed.



#### Language

Device Setup | Setup | System | Interfaces | Language

🔂 Language		
System Language ENGLISH GERMAN ITALIAN JAPANESE KOREAN	——1	
1 Selection arrow		

Fig. 32 Language

#### Change system language

- 1. Click on the selection arrow (Fig. 32, item 1) and select a language.
- u A message window "Restart" appears.
- **2.** Confirm with "OK" to restart the device.
- u The device is restarted and the language has been changed.



# **CE Remote Control**

Device Setup | Setup | System | interfaces | Remote Control

By activating (Fig. 33, item 1) the CE Remote Control, the unit can be operated via the network connection. The software of the remote control (RC software) is stored on the device and can be copied to a USB storage device.

CE Remote Control	
Enabled	—1
Get RC Software	2
<ol> <li>Check box remote control "Enabled"</li> <li>Button <copy rc="" software=""></copy></li> </ol>	

Fig. 33 Remote Control

#### Copy RC software

- 1. Connect a USB storage device to the device (Fig. 3, item 5).
- u The USB storage device is recognised by the system after a few seconds and the button <Copy RC software> (Fig. 33, item 2) can be clicked.
- 2. Click on the button <Copy RC software> (Fig. 33, item 2).
- u The software is copied to the USB storage device.

#### Information



The RC software includes the file "LB47xRemoteControl.exe" and runs without installation. Operation of the RC software is described in Chapter "Remote Control Software" (see next but one chapter).



#### Units

Device Setup | Setup | System | Units

Clicking on the individual selection arrow lists the available units for the measuring value. The selected unit is used in the output settings and is shown in the standard display.

Fig. 34 Units



#### Network

Device Setup | Setup | System | Network

In the Network settings, you can make changes to the network settings. The information can only be edited in the access level "Standard" (see chap. 7.2 Menu Access). You can set the network address either manually or using DHCP (automatic assignment). To do this, check the "DHCP active" in the selection field (Fig. 35, item 1).

network		1
DHCP Enabled		— I
IP Address: 192.1	.68.2.118	—2
Subnet Mask: 255.2	255.255.0	—3
Default Gateway: 192.1	.68.0.1	—4
DNS Server: 192.1	.68.0.1	— 5
MAC: 54-F6	-66-6F-FF-01-	— 6
Apply		—7
1 Check box "DHCP active"	5	DNS server
2 IP address	6	MAC address
<ol> <li>Subnet mask</li> <li>Standard gateway</li> </ol>	7	Button <apply></apply>

Fig. 35 Network Settings

#### **IMPORTANT**

The PC and the LB 478 have to be in the same IP subnet.

 In the event of an automatic assignment of the IP address by a DHCP server, you can only look at the given IP address. A modification of the IP address is not possible. On this side, you can also read the MAC address of the device (Fig. 35, item 6).

#### Manual settings

- 1. Click on the text field (Fig. 35, item 2 5) to open the input field.
- 2. Enter the appropriate network addresses.
- **3.** Confirm with the Enter key.
- 4. Click on <Apply> (Fig. 35, item 7) to adopt the network settings.

#### **IMPORTANT**

All settings performed must be confirmed by clicking on <Apply> so that the settings become real.



#### **Remote Control Software**

If the EVU is connected to a network at the RJ45 socket (Fig. 3, item 2), the EVU can be operated via a computer. The software can be loaded onto a USB storage device (see Chapter "CE Remote Control").

#### **IMPORTANT**

In order for the Remote Control to function, the selection check mark in the menu "CE Remote Control" must be set to "Active" (Fig. 33, item 1).



Fig. 36 Establishing connection to the EVU using the RC software

- 1. Click on "LB47xRemoteControl.exe", to start the program.
- u The program starts (Fig. 36, item 1).
- 2. Click on the <File> tab and then on <Connect...> (Fig. 36, item 2), to establish a connection to the EVU.
- A new window "Connect" is opened (Fig. 36, item 3) and the connected transmitters are listed.

#### **IMPORTANT**

The IP address of the EVU must be in the same sub-network (Fig. 35, item 3) as the network adapter of the computer (see previous Chapter "Network").

- **3.** Click on the identifier of the transmitter (Fig. 36, item 4) or enter the IP address of the EVU in the input box (Fig. 36, item 5) (see Fig. 36, item 2).
- 4. Click on <OK>.
- u The connection to the EVU is established.
- u You can enlarge the view in the "Zoom" menu (2x, 3x).



#### **Reset Device (evaluation unit)**

Device Setup | Setup | System | Reset Device

The evaluation unit can be restarted and reset to factory settings in the window "Reset Device".

Û	Reset Device	Factory Reset Warning Al device settings will be reset back to the factory settings and device will be reboted! Are
_	Reboot	3 2 Reboot Warning
<		Device will be rebooted and measurement will be interrupted! Continue? 4 Yes No
1	Button Reboot	
2	Button Factory settings	
3	Window "Warning: Reboot"	
4	Window "Warning: Factory Reset"	

Fig. 37 Reset Device

#### **Restart the device**

**IMPORTANT** 



The measurement is interrupted during a restart!

- 1. To restart the device, click the button <Restart> (Fig. 37, item 1).
- u A window with a warning "Restart" (Fig. 37, item 3) opens.
- 2. Click on <Yes> to confirm.
- u The device is restarted.

#### **Reset device (Factory Reset)**

#### IMPORTANT



When there is a reset to factory settings, all data logs are deleted and all user-defined configuration settings are reset!

- u If error M102 appear, the device possibly must be reset twice.
- 1. To reset the evaluation unit to the factory settings, click the button <Factory settings> (Fig. 37, item 2).
- u A window with the warning "Factory settings" (Fig. 37, item 4) opens.
- 2. Click on <Yes> to confirm.
- u The device is reset to factory settings and restarts.



#### **Repair Detector Software**

Device Setup | Setup | System | Repair Detector Software

In the window "Repair Detector Software" detectors can be repaired and updated.

#### Tip

The current software versions for the detectors can be downloaded from the berthold website (www.berthold.com).





#### ▲ DANGER



#### Danger to life from electric shock!

- u The repair may only be carried out by a qualified electrician.
- u Please adhere to the relevant safety regulations.
- u During servicing on the hardware as well as during wiring of the detector, the measuring system, connected relay contacts and all inputs and outputs must be de-energised.

In case of an electric shock, carry out first aid measures and immediately call an emergency service.

- 1. Disconnect all detectors from the EVU except for the one you want to repair / update.
- 2. Insert a USB flash drive with the application software of the detector into the USB port of the EVU.
- 3. Click the <Update / Repair> button and follow the installation instructions.
- **4.** After the update is complete, restart the detector by disconnecting and connecting its power supply.
- u The update / repair was done.



# 7.3.2 Sensors (only master EVU)

Device Setup | Setup | Sensors

All settings of the sensors are carried out via the master EVU.

ω	Sensors
	Detector Configuration

Fig. 39 Menu "Sensors" (only Master EVU)



#### **Detector configuration**

Device Setup | Setup | Sensors | Detector configuration

In the window "Detector configuration" the detectors for the measuring system are added and configured. By clicking the button <Edit> (Fig. 40, item 2) the device ID, the position and description can be changed.





- 1. Click on < + > to add a new detector.
- 2. Enter the Detector Device ID (Fig. 40, item 8), located at the socket of the detector (Fig. 41, item 2).
- **3.** Enter the position of the detector (Fig. 40, item 7), depending on measurement arrangement (Fig. 41, item 3). This values must be agreed with an authorised person.
- 4. Enter a Description (Fig. 40, item 6) and click < OK >.
- u The detector was added successfully.



Fig. 41 Detector position / Device ID

BERTHOLD

# 7.3.3 Interphase Densities (only master EVU)

An interphase denotes the layer with the respective density. In this menu the interphases are defined and a density is assigned to each interphase. The entered density should be the upper limit value of the density for this interphase. In addition to the individual interfaces, it is possible to add a total level. The gas density should be specified as the value for the total level.



Fig. 42 Interphase settings

# 7.3.4 **Densities Overview (only master EVU)**

Device Setup | Setup | Densities Overview

The control unit reads the detectors' individual local densities in order to build a density profile of the vessel's content. This density profile might not be perfectly monotonous (see Fig. 43, item 1), due to statistics, noise and vortexes inside the vessel. Thus, the measurement algorithm uses mathematical techniques such as a special averaging / regression to monotonize and slightly smooth the local densities (see Fig. 43, item 2), before determining the interface positions between the different phases. As a result, interface positions calculated by LB 478 will never be ambiguous or inconsistent.



Fig. 43 Densities Overview



# 7.3.5 Outputs

Device Setup | Setup | Outputs

You can make the following settings and read information in the window "Outputs":

- Analog Output (AO)
  - AO Mapping
  - $\circ$  AO monitoring
  - $\circ \ \ \, \text{Failure mode}$
  - o Current limits
  - o Calibrate
- · Additional Outputs
  - $\circ$  Configuration
  - AO Mapping

n Outputs		ω	Analog Ouptut (AO)
	Analog Output (AO)		AO Mapping
	Additional Outputs		AO Monitoring
			AO Failure Mode
			AO Limits
<		<	AO Calibration

Fig. 44 Menu "Outputs" (only Master EVU), Menu "Analog Output (AO)"



#### 7.3.5.1 Analog Output (AO)

#### AO Mapping (Master EVU)

Device Setup | Setup | Outputs | Analog Output (AO) | AO Mapping

An Interface can be assigned to the analog output of the master EVU in the window "AO Mapping". The current output signal is between 4 mA and 20 mA. The corresponding level values can be freely assigned. The value which is assigned to a current output signal of 4 mA must be smaller than the value which is assigned to that of 20 mA.



Fig. 45 Analog Output Mapping (Assignment)

#### AO Monitoring

#### Device Setup | Setup | Output | Analog Output (AO) | AO Monitoring

If "AO Monitoring" is activated (Fig. 46, item 1), the current output will be monitored. It is continuously monitored whether the current value flowing in the current loop is correct. In the event of a variation e.g. owing to an error in the hardware, too large a load or a disruption in the loop, an error message is triggered.



Fig. 46 Analog output monitoring



#### **AO Failure Mode**

Device Setup | Setup | Output | Analog Output (AO) | AO Failure Mode

The alarm function is set when an error is detected at the current output in the window "Error mode".



#### Fig. 47 Analog output (AO Failure Mode)

The following behaviour of the current output can be assigned in case of error:

Namur High	22 mA (in error mode).		
Namur Low	2 mA (in error mode).		
Hold Value	Last value before the error.		
User-defined value	The value can be set manually.		

#### NOTICE

If the value "Hold Value" is set, it is recommended that the error relay is connected in order to allow device errors to be transmitted to the control system.



# **AO Limits**

Device Setup | Setup | Output | Analog Output (AO) | AO Limits

By clicking on the input fields (Fig. 48, item 1, item 2), the values [mA] for the lower and upper current limit can be set. In addition to the default value of 3.8 mA, the lower current limit can be set to 0 mA in order to switch the current output from 4 ... 20 mA to 0 ... 20 mA.



Fig. 48 Analogue output (AO Limit)



If the measurement is operated according to Namur, the standard current values of 3.8 or 20.5 mA must be maintained.



#### Calibrate

Device Setup | Setup | Output | Analog Output (AO) | AO Calibration

If there are any discrepancies between the target value and the actual value of the current signal, then the current output may be calibrated again.

#### NOTICE

For calibration of the current output, an ammeter (not included in the scope of delivery) is required, which is connected to the current output.



Berthold recommends calibrating the current outputs whenever a module has been installed/replaced or if a software update has been carried out.



Fig. 49 Analog Output (Calibration)



#### Perform calibration

#### 



#### Danger to life from electric shock!

- u The repair may only be carried out by a qualified electrician.
- u Please adhere to the relevant safety regulations.
- During servicing on the hardware as well as during wiring of the detector, the measuring system, connected relay contacts and all inputs and outputs must be de-energised.

In case of an electric shock, carry out first aid measures and immediately call an emergency service.

- 1. Connect the measuring lines of the current measuring device on the back of the EVU to the analog output.
- 2. Observe the terminal allocation in the document "Technical information".
- 3. Click on the button <Calibration>.
- u The device switches to test mode and a new window (Analog output calibration) opens.
- u The calibration point 4 mA is displayed and the current measuring instrument shows a value.
- **4.** Enter the indicated value on the current measuring instrument in the input box (Fig. 49, item 4).
- 5. Click on the button <Continue>.
- u The calibration point 20 mA is displayed and the current measuring instrument shows a value.
- 6. Enter the indicated value on the current measuring instrument in the input box (Fig. 49, item 5).
- 7. Click on the button <Continue>.
- u A message appears "Calibration successful".
- 8. Click on the button <Continue>.
- u The calibration of the analog output is completed.



#### 7.3.5.2 Additional Outputs (only master EVU)

Û	Additional Outputs
	Configuration
	Out 1
	Out 2
	Out 3
<	



Already configured outputs are displayed, otherwise only the menu "Configuration" is displayed.

#### Configuration

Device Setup | Setup | Outputs | Analogue Output (AO) | Additional Outputs | Configuration

<b>Onfiguration</b>				+		
	Device ID	Description	Status		Description:	
	0016092	Out 1	ok		Description.	
	0016054	Out 2	ok		Device ID:	
	0016074	Out 3	ok			
<	+	Edit -	4		Ok	Cancel
1	Device ID (	extension mo	odule	4	Button < - >	
ว	2 Putton z L		5	Description of analog output		
2			5			
3	Button < E	dıt >		6	Input field Device ID	extension module

In this menu further outputs of extension modules are added.

Fig. 51 Additional Outputs configuration

- **1.** Click on the button < + >.
- 2. Enter a Description (Fig. 51, item 5).
- 3. Add the Device ID of the extension module (Fig. 51, item 6).
- You can find the extension's Device ID in the menu Device Setup | Identification | Device Information of the respective extension module.
- 4. Click < OK >.
- u "OK" is displayed in the status column and the analog output of the extension module was added successfully.



#### AO Mapping (additional outputs)

Device Setup | Setup | Outputs | Analog Output (AO) | Out # | AO Mapping

An Interface can be assigned to an analogue output of the extension module in the window "AO Mapping". The current output signal is between 4 mA and 20 mA. The corresponding level values can be freely assigned. The value which is assigned to a current output signal of 4 mA must be smaller than the value which is assigned to that of 20 mA.



Fig. 52 Analog Output Mapping (Assignment)



#### Example



# 7.4 Menu Backup/Restore

Device Settings | Backup/Restore

You can make a backup copy of the configuration data, and perform a recovery in the submenu "Backup/Restore".

Fig. 54 Menu "Backup/Restore"

# 7.4.1 Backup

Device Settings | Backup/Restore | Backup



Fig. 55 Backup

#### **Perform Backup**

- 1. Connect a USB storage device to the device (Fig. 3, item 5).
- 2. The USB storage device is recognised by the system after a few seconds and the button <Backup> (Fig. 55, item 1) can be clicked.
- 3. Click on the button <Backup> (Fig. 55, item 1).
- u The window "Enter description" appears.



- u The message "Error free" (Fig. 55, item 3) appears in the field "Backup data" for error-free backup files.
- 4. Click the button <Edit>, enter a description, and confirm with the Enter key.
- 5. Click on the button <Save>.
- u The backup files are copied to the USB storage device.
- u The message "Backup successful!" appears after a successful copy process (Fig. 55, item 6).

#### Information



The backup includes an XML file that is created in the folder "Backup\_LB47x". The file name is derived from "Backup", the date and time (Backup\_YYYYMMDD\_ hr-min-sec).


## 7.4.2 Restore

Device settings | Backup/Restore | Restore

G	Restore	Restore Device Settings			
_	28/08/2017 17:58:04 No Errors 7	Device settings will be replaced with those from the selected Backup. Are you sure?			
<	1 2 3 4	Yes No			
1	Button < "previous"				
2	Number of recovery files on the USB stora	ge device			
3	Button > "next"				
4	Button <restore></restore>				
5	Recording date of the backup file				
6	Time of backup file recording				
7	Information about the backup data (error	/ error free)			
8	Confirmation window				

Fig. 56 Restore

#### **Executing restore**

- 1. Connect a USB storage device to the device (Fig. 3, item 5).
- 2. Select the backup file with the buttons (Fig. 56, item 1,3)
- u The date and time of the backup is displayed (Fig. 56, item 5, 6). Only errorfree (Fig. 56, item 7) backup files can be loaded.
- 3. Click on the button "Restore" (Fig. 56, item 4).
- 4. A confirmation message (Fig. 56, item 8) appears.
- 5. Click on < Yes > to confirm, click on < No > to cancel.
- u The restore of data is carried out.









Fig. 57 Menu "Diagnostics"

### 8.1 Transmitter Temperature

Diagnostics | Transmitter Temperature

Temperature values from the evaluation unit (processor) are displayed in the menu item "Transmitter Temperature".



Fig. 58 Transmitter Temperature



## 8.2 Transmitter Events

Diagnostics | Transm. Events

Fig. 59 Menuu "Transm. Events"

#### Information

Events of the respective detector can be seen at Device Setup | Setup | Sensors | [NAME OF DETECTOR] | Detector Service.

### 8.2.1 Trasnmitter Event Log

Diagnostics | Transmitter Events | Transmitter Event Log

The last 25 events of the detector are displayed in the window "Event Log".



Fig. 60 Event log (Transmitter)



#### **Display Event Description**

6	Transm. Event L	_og		Device event information
	Date / Time	No.	Description 🔺	205
	2017-08-28 05:58:01	311	Backup in prog	
	2017-08-28 05:53:45	309	Measurement s	T smperature Upper Limit out of specification
	2017-08-28 05:51:05	304	Temperature U	Dutector temperature at / above upper limit
	2017-08-28 05:50:51	306	Temperature C	
	2017-08-28 05:48:26	320	PV Lower Limit	
	2017-08-28 05:47:21	322	PV Lower Limit 🚽	
			•	
		ar Ever	nt Log	
	1	2		3456 7
1	Button	-		
2	Button <clear ever<="" th=""><th>nt loa</th><th></th><th></th></clear>	nt loa		
2	Lighlighted event	ntiog		
3	Highlighted event			
4	Event no.			
5	Event title			
6	Event description			
7	Button <close></close>			

Fig. 61 display an event log

- 1. Click on a line in the list (Fig. 61, item 3).
- **2.** Click on <?> (Fig. 61, item 1).
- u The event description appears.
- 3. With the button <Close>, close the event description (Fig. 61, item 7).

#### NOTICE

With the button <Clear event log> (Fig. 61, item 2) all events are deleted irrevocable.



### 8.2.2 Transm. Event Overview

Diagnostics | Transm. Events | Transm. Event Overview

All events that can be logged are chronologically presented in tabular form in the window "Event overview". Activate the check box "Non-zero Counter only" (Fig. 62, item 5) in order to display events that have occurred.



Fig. 62 Transmitter Event Overview

- 1. Click on a line in the list (Fig. 62, item 4).
- **2.** Click on <? > (Fig. 62, item 7).
- u The event description appears.
- 3. Close the event description with the button <Close>.
- **4.** Slide the bar of the horizontal scroll bar (Fig. 62, item 6) to the right to see at what times (date, time) the event occurred.
- u The last 5 time points are displayed.



# 8.3 Change Log

Diagnostics | Change Log

You can track changes that were performed on the device in the window "Change log".

D	ate / Time	De	scripti	Value1	
	2017-08-28 05:58:01	Ba	::kup	1.0	
	2017-08-28 05:54:49	Si	n Mode	0	
_	2017-08-28 05:53:45	Si	n Mode	1.0	
	2017-08-28 05:52:11	D	UL e	0	
_	2017-08-28 05:52:11	D1	LL ev	0	
	2017-08-28 05:51:05	D	UL e	2.0	
_	2017-08-28 05:51:05	D	LL ev	2.0	_
4	2017-08-28 05:50:51	_ D1	UL e	1.0	•
	<b>Ⅰ</b>			▶	
	1	:	2		
	_				
٦	Time of the chan	ge			
5	Short info of the	ch	ange		
``	Jalua 1		<u> </u>		
•					

Fig. 63 Change Log (Transmitter)



# 8.4 Data Log (only master EVU)

#### Diagnostics | Data Log

You can set the log interval as well as delete and export the log data in the submenu "Data log".

#### IMPORTANT

The data cannot be viewed on the EVU Display or via Ethernet. The data must be exported to a USB storage device to view it on a PC.



Fig. 64 Data Log

#### Change log interval

- 1. Click on the button <Stop> (Fig. 64, item 2) to stop the data log process.
- 2. Click on the input field "Log interval" (Fig. 64, item 1) and enter the time in seconds.
- **3.** Confirm with the Enter key.
- u The interval was accepted
- 4. Click on the button <Start> (Fig. 64, item 4) to start the data log process.



#### Export log data

- 1. Click on the button <Stop> (Fig. 64, item 2) to stop the data log process.
- 2. Connect a USB storage device to the device (Fig. 3, item 5).
- 3. Click on the button <Export> (Fig. 64, item 5).
- u The export process is started and can take several minutes to complete under certain circumstances.
- u The message window "Export successful!" appears with a successful export.

#### Information

The export includes a zip file that is created in the folder "ExtendedLogExport". The file name is derived from "ExtendedLogExport", the date and time (ExtendedLogExport\_YYYYMMDD\_hr-min-sec).

#### **Exported data structure**



Fig. 65 Data structure



# 8.5 Export service data (only master EVU)

Diagnostics | Export Service Data



Fig. 66 Export Service Data

- 1. Connect a USB storage device to the device (Fig. 3, item 5).
- 2. The USB storage is recognised by the system after a few seconds and the button <Export service data> (Fig. 66, item 1) can be clicked.
- 3. Click on the button <Export service data> (Fig. 66, item 1).
- u The process window is displayed and the export of service data will be carried out (Fig. 66, item 2).
- u After successful export a confirmation message Fig. 66, item 3) is displayed.

#### Information

Four .txt files are copied to the USB drive when exporting:

- ChangeLog
- ErrorHistory
- ErrorSummary
   SystemInfo

# Information



A new folder "Backup\_LB47x" is created and the Backup-File (.xml) is copied when exporting.



# 9 Troubleshooting

# 9.1 Error Search

Problem	Cause	Measure
Master unit: Screen black; LEDs are not illuminated	EVU does not work	<ul> <li>Check power supply and fuses</li> </ul>
Extension module: LEDs are not illuminated	Extension module not clamped properly	<ul> <li>Check cabling, contact so- ckets</li> </ul>
Extension module: is not detected	Incorrect device ID	u Check Device ID in the menu Identification of the extension module
No signal	Detector does not work	u Check the functioning of the detector
	Shield not opened or not opened correctly	u Check lock and ensure it is in OPEN position
Count rate too low	Incorrect focus of the effec- tive radiation on the detector	<ul> <li>Correct and optimise the alignment</li> </ul>
	Objects in the beam path	u Offset irradiation level
	Source at the end of its usa- ble life span	u Replace source
	Defect in detector	u Check detector
The display deviates	Count rate too low (see above)	<ul> <li>Check source age and irra- diation level, replace detector</li> </ul>
	Incorrect Calibration	u Check calibration values
	Terminals / wiring	<ul> <li>Check terminal connec- tion; check terminal assignment</li> </ul>
Detector is not detected (soft- ware)	Damaged line	u Check cable; examine with measurement device.
	Incorrect ID in the configura- tion	<ul> <li>Check ID of the detector (see type plate on the de- tector)</li> </ul>
Touch panel does not respond	Error in operating system	u Restart EVU
Buttons are missed when you click	Incorrect screen calibration	u Calibrate screen again



## 9.2 Error Codes of the Evaluation Unit

In the following tables you can find the EVU and detector error codes which give you exact information on how to fix them. The error codes of the detectors can be found in the operating instructions of the respective detectors.

System events are classified in

- FAILURE (F)
- OUT OF SPECIFICATION (S)
- FUNCTION CHECK (C)
- MAINTENANCE REQUIRED (M)

#### Failure (F)

Severe device error. The current output emits an error current. The error relay gives alarm (contact opens).

#### Out of specification (S)

The detector, one of its components or the process itself, are out of normal specification.

#### Function Check (C)

Indicates that entries are made at the detector or a function check/simulation is being performed.

Code	Message	NAMUR107	Help Text
M101	HW Module	F	Hardware electronics module corrupt. Restart the device. Contact Berthold ser- vice, if this event occurs repeatedly.
M102	Device data-set	F	Failure of the permanent memory. No parameter set found. Factory reset and / or restart the device. Contact Berthold service, if this event occurs repeatedly. The device possibly must be reset twice.
M103	RAM, Flash or CPU	F	Internal hardware failure. Restart the device. Contact Berthold service, if this event occurs repeatedly.
M104	WD Reboot	М	The Watchdog has caused the device to restart. Contact Berthold service, if this event occurs repeatedly. Check, if mas- sive electromagnetic interferences have caused this event.
M105	WD Failure	F	Watchdog malfunction. Contact Berthold service, if this event occurs re- peatedly.
M106	WD Off	М	Watchdog is inactive. Activate Watch- dog

## 9.2.1 **System**



M107	Error in the internal real time clock	М	Malfunction of the real-time clock. Check Date and Time. If the event oc- curs frequently, contact Berthold Service.
M108	CPU temperature sensor	М	The temperature sensor of the device is defective. Contact Berthold Service. The hardware is defective and, if necessary, must be checked and replaced.
M109	Lower temperature limit: Maintenance required	М	The internal temperature of the device is close to the lower threshold value of the permissible operating temperature.
M110	Temp LL OOS	S	The internal temperature of the device is below the lower limit. The correct function of the device can not be guar- anteed. It is recommend to have the device checked by Berthold Service, even if it seems to work normally.
M111	Temp UL maintenace	М	The internal temperature of the device is close to the upper limit.
M112	Temp UL OOS	S	The internal temperature of the device is above the upper limit. The correct function of the device can not be guar- anteed. It is recommend to have the device checked by Berthold Service, even if it seems to work normally.
M113	Power On Reboot	С	The device was restarted due to un- known reasons.
M114	Software Reboot	С	The device was restarted by user input.
M115	Extern RTC malfunction	М	Failure of the external real time clock. Contact Berthold service, if this event occurs repeatedly.
M116	Corrupt Date	М	The date could not be verified at startup. Check date and time and set if necessary.

# 9.2.2 Application

Code	Message	NAMUR107	Help Text
M310	PV calc not possible	F	Process value could not be calculated.
M311	Backup process	С	Backup in process.
M312	Restore process	С	Restore in process.
M399	Internal program error	F	Internal software failure. Restart the de- vice. Contact Berthold service, if this event occurs repeatedly.

# 9.2.3 Detector

Code	Message	NAMUR107	Help Text
D501	Detector not found	S	Connection to at least one detector lost. Check detector configuration and / or detector connection. Contact Berthold service, if this event persists.
D502	Detector comm error	М	Communication to at least one detector interrupted. Check detector connection. Contact Berthold service, if this event oc- curs repeatedly.
D503	Detector failure	S	At least one detector signals "failure". Check detector events.
D504	Detector out of spec	S	At least one detector signals "out of specification". Check detector events.
D505	Detector func check	с	At least one detector signals "function check". Check detector events.
D506	Detector maintenance	М	At least one detector signals "mainte- nance required". Check detector events.
D507	Ex-i Board failure	S	
D599	Internal program err	F	Internal software failure. Restart the device. Contact Berthold service, if this event occurs repeatedly.



# 9.2.4 RS 458 Interface

Code	Message	NAMUR107	Help Text
D601	Detector not found	S	Connection to at least one "additional output" lost. Check configuration and / or connection. Contact Berthold service, if this event persists.
D602	Detector comm error	М	Communication to at least one "addi- tional output" interrupted. Check connection. Contact Berthold service, if this event occurs repeatedly.
D699	Internal program err	F	Internal software failure. Restart the de- vice. Contact Berthold service, if this event occurs repeatedly.

# 9.2.5 **Process Connection**

Code	Message	NAMUR107	Help Text
M701	Current output	F	Deviation of analog output value from feeback value is too high. Calibrate ana- log output. Contact Berthold service, if this event occurs repeatedly.
M702	Current loop open	F	Current output loop open. Check cable connection.
M703	Software update	с	Indicates that a software update is run- ning. No action necessary. The device automatically returns to measuring mode after the software update is fin- ished.
M799	Internal program error	F	Internal software failure. Restart the de- vice. Contact Berthold service, if this event occurs repeatedly.





# **10** Maintenance and Repair

The replacing of fuses, the cleaning and data storage of the EVU are described in the maintenance and repair chapter.

#### IMPORTANT

- u The applicable national regulations of the respective country of use have to be observed!
  - Repair and servicing on the EVU may only be carried out by experts. In case of doubt, the complete EVU is to be sent to Berthold.

#### NOTICE

u Repair on electronic circuits on the circuit boards of a field device may only be carried out in the manufacturer's factory.



- u When working at electronic components, the relevant safety regulations must always be observed. Particularly observe the safety instructions in the chapter 2 Safety.
- u De-energise the detector and potentially connected relay contacts as well as all inputs and outputs.

#### **IMPORTANT**



To achieve optimum measuring accuracy, we recommend recalibrating the measuring system after a repair (not after changing the housing).



# 10.1 Replacing of Fuses

#### \Lambda DANGER



#### Danger to life from electric shock!

Replacing of fuses may only be carried out by a qualified electrician.

- u Please adhere to the relevant safety regulations.
- u Installation/maintenance may only be carried out if the device has been de-energised.
- u Only open the device when free of voltage.

In case of an electric shock, carry out first aid measures and immediately call an emergency service.

#### NOTICE

Ū.

Damage to the device! Short circuit!

The EVU can be damaged if incorrect fuses are used.



- Only use fuses which correspond to the fuses on the circuit board of the module. Fuses:
  - EVU: 250V 1A T (5x20 mm)
  - EVU: 250V TR5 T80mA (Ø 8,5 mm)



Fig. 67 Replacing fuses master EVU

- **1.** De-energise the device.
- 2. Loosen the four fixing screws (Fig. 67, item 1) and remove the EVU from the wall housing or subrack.
- 3. Loosen the four sunken screws on the front side of the EVU (Fig. 67, item 2).
- 4. Pull out the housing (Fig. 67, item 3) carefully.



- 5. Remove the protective covering of the fuse (Fig. 67, item 4)
- 6. Remove the fuse (Fig. 67, item 5, item 7).
- 7. Insert the new fuses and attach the protective covering again.
- 8. Carefully slide the circuit board into the housing.
- 9. Screw the front panel to the housing with the four screws (Fig. 67, item 2).
- **10.** Set module into the guide rails and push it gently until the plug connector of the module is inserted into the socket board.
- 11. Tighten all fixing screws (Fig. 67, item 1).
- u The fuse change was carried out correctly.



# 10.2 Cleaning

The display is designed for maintenance-free operation. Make sure you keep the touch screen clean. Use a cleaning cloth dampened with a cleaning agent to clean the equipment. Only use water with a little liquid soap or a screen cleaning foam.

#### NOTICE



Unintentional reaction!

When cleaning the touchscreen, touching keys can trigger an unintentional reaction in the EVU.

u When cleaning, make sure that no unintentional reactions are triggered.

#### NOTE

#### Damage caused by unauthorized cleaning products!

The display may be damaged if compressed air, steam jet blowers, aggressive solvents or scouring powders are used for cleaning purposes.

u Do not clean the EVU with compressed air or steam jet blowers. Do not use aggressive solvents or scouring powder.



Fig. 68 Cleaning the display

- 1. Shut down the device.
- 2. Spray the cleaning solution onto a cleaning cloth. Do not spray directly onto the display.
- 3. Clean the display.
- 4. When cleaning the display wipe from the screen edge inwards.
- 5. Let the display dry out.
- u The cleaning of the display was performed correctly.





#### 10.3 Data backup

Activate the data log (see chapter 8.4) so that all data are recorded. Perform a log data and service data backup at regular intervals.

#### **Export service data**

Diagnostics | Export Service Data

Four .txt files are copied to the USB drive when exporting service data:

- ChangeLog
- ErrorHistory
- ErrorSummary
- · SystemInfo

Export Service Data	Export Starting export:
Start Export	Starting export ->export successful Starting backup: ->beckup successful Starting XML2CSV conversion: ->MML2CSV conversion successful All done.
1 Button <export data="" service=""></export>	
2 Process window	
3 Confirmation message	

Fig. 69 Export Service Data

- 6. Connect a USB storage device to the device (Fig. 3, item 5).
- 7. The USB storage is recognised by the system after a few seconds and the button <Export service data> (Fig. 66, item 1) can be clicked.
- 8. Click on the button <Export service data> (Fig. 66, item 1).
- u The process window is displayed and the export of service data will be carried out (Fig. 66, item 2).
- u After successful export a confirmation message Fig. 66, item 3) is displayed.

#### Information



A new folder "Backup\_LB474" is created and the Backup-File (.xml) is copied when exporting.





# **11** Decommissioning

#### ▲ DANGER



Danger to life from electric shock!

Decommissioning may only be carried out by qualified electricians.

- u Please adhere to the relevant safety regulations.
- Decommissioning may only be carried out if the device has been de-energised.
- u Only open the device when free of voltage.

In case of an electric shock, carry out first aid measures and immediately call an emergency service.

# 11.1 Decommissioning 19" subrack



Fig. 70 Decommissioning wall housing

- 1. De-energise the device.
- 2. Remove all lines (Fig. 70, item 1) from the clamp blocks.
- 3. Remove the network plug (Fig. 70, item 2).
- **4.** Remove the PE cable (Fig. 70, item 3).
- 5. Remove the connections (Fig. 70, item 4) and pull the subrack from the 19" rack.
- u The decommissioning is completed.

**BERTHOLD** 

# 11.2 Disposal of Measurement System

#### 

#### Toxic!



The product contains electronic components containing toxic substances that are harmful to health.

u Disposal is to be carried out in accordance with the disposal regulations via a disposal expert.

If the device is to be decommissioned, have it disposed of according to legal regulations (e.g. RL 2002/96/EC) by a specialised waste management company.



# **12** Appendix

# 12.1 Commissioning Guide

This guide is intended to support and document the installation and commissioning of Multiphase Level Measurement (MPLM) systems. The document is divided into the following sections:

- · Underlying data
- · Implementation on site
- Setup protocol

#### **IMPORTANT**

This guide does not replace the instructions for installation, electrical installation and setup in the corresponding operating instructions.



Detailed descriptions of the procedures can be found in the respective operating instructions of the detector or the evaluation unit.

u The safety instructions of the corresponding user manuals must be observed!



# 12.1.1 Underlying data

The data required for an installation of the MPLM system are listed here. Some data have already been defined in Berthold's factory or clarified with the customer in advance. All information must be agreed with the customer or the responsible process engineer.

#### Information about the system

Customer	
Measuring point	

#### Detectors

#	Detector-ID	Position in	Process path in
1			
2			
3			
4			
5			
6			
7			
8			
9			
10			
11			

#### Sources

Total number of sources installed in the dip tube: \_\_\_\_\_ Nuclide: \_\_\_\_\_

#	Number of sources per wave core	Length of the wave core	Source numbers (activities)
1			
2			
3			
4			
5			



#### Interfaces

Interface	Density limit value in <sup>1</sup>	Expected position ranges in1
Total Level (Gas Density)		

#### **Evaluation Unit**

Device-ID Master	Location

### **Extension Module**

#	Device-ID Extension	Description

<sup>&</sup>lt;sup>1</sup> The selected units must be consistent with the settings of the evaluation unit!



## 12.1.2 Implementation on site

At this point, the steps to be carried out on site are listed. A detailed description of the installation, electrical installation, etc. can be found in the respective operating instructions. Safety instructions must be observed!

#### **Detector Positioning**

In most cases, the detectors are installed by the customer. Prerequisite: Mounting for the detectors is provided. Fix detectors to the mounting using appropriate clamps. Positioning of the detectors according to specifications (see Table "Detectors" in chapter 12.1.1). During commissioning, only a fine adjustment of the detectors is usually necessary.

#### **Evaluation Unit**

The evaluation unit is usually preconfigured in the factory. Commissioning is documented by means of a commissioning protocol (see 12.2 Setup Protocol LB 478).

The following settings on the Master Control Unit must be checked:

1. Are all detectors incl. position applied? See Table "Detectors" in chapter 12.1.1



Fig. 71 Example Detector Configuration

2. Are all required interfaces created? See Table "Interfaces" in chapter 12.1.1



Fig. 72 Example Interface Densities



**3.** Are all additional outputs created? An analog output must be created for each additional evaluation unit (extension module).



Fig. 73 Example extension module (outputs) configuration

**4.** Scaling of analog outputs according to the specifications for position ranges (see Table "Interfaces" in chapter 12.1.1.). Note: This data must be agreed with the relevant process engineer.



Fig. 74 Example extension module (outputs) configuration

#### **Calibration of detectors**



- u Each detector must be calibrated individually!
- u As part of the calibration procedure, the container must be emptied or filled, time must be allowed for this.

For the detectors of a MPLM system, a two-point calibration is performed, i.e. first the calibration of all detectors is carried out at density A and then at density B. In individual cases, a 1-point calibration may suffice. The calibration of the detectors is described in detail in the corresponding operating instructions, at this point only assistance is given. Calibration is documented individually for each detector in the form of a commissioning protocol (template see 12.2 Setup Protocol LB 478).

- Connect the HART communicator to detector #1 Note: LB480 Evaluation Units are passive, 24 V auxillary voltage in series to 4 - 20 mA output required.
- 2. Set the basic parameters.
- 3. Calibrate first density.
- 4. Repeat Step 2 and 3 for the remaining detectors.
- 5. Calibrate the second density at all detectors and trigger calibrate.



# 12.2 Setup Protocol LB 478

The commissioning is carried out by means of setup protocols for the evaluation units and the detectors. Note: A separate protocol must be completed for each detector!

General data			
Date			
Measuring point			
Number of dectors	LB480		
Number of sources			
lsotope	Cs-137	Co-60	
Container			
Products			

Device configuration		
Model		
Installation variant	Wall housing	Subrack
Power supply	🗌 100-240 V AC	🗌 18-32 V DC
Number of master EVU		
Device ID master EVU		
Number of extension modules		
Device ID's extension modules		
Software Version		



# Setup Protocol (Continued)

Parameters	
Password	
Language	□ DE □ EN □
CE Remote Control	🗌 enabled
Network	DHCP active IP Adress•• Subnet•• Ggateway•• DNS-Server•• MAC Adress
Process Value Range (Master)	min. Value 0 / 4,00 mA max. Value 20,00 mA
Process Value Range (Extension ID)	min. Value 0 / 4,00 mA max. Value 20,00 mA
Process Value Range (Extension ID)	min. Value  0 / 4,00 mA max. Value  20,00 mA
Process Value Range (Extension ID)	min. Value  0 / 4,00 mA max. Value  20,00 mA
Process Value Range (Extension ID)	min. Value 0 / 4,00 mA max. Value 20,00 mA
Process Value Range (Extension ID)	min. Value  0 / 4,00 mA max. Value  20,00 mA
Process Value Range (Extension ID)	min. Value 0 / 4,00 mA max. Value 20,00 mA
AO Assignment Master	Interface:
AO Assignment Extension ID	Interface:



# 12.3 Setup Protocol LB480 MPLM

Measuring point	Date	
lsotope	Detector Position (Height)	
Detector	HV	

Path	Parameters	Unit	Standard	Setup	
Device Config u Meas Parameter u Meas Data					
	Background	cps	50		
Device Config u l	Meas Parameter ပ M	eas Data u Meas Tal	ble		
Point	Density	Rate			
1					
2					
3					
4					
Meas Parameter	J Meas Settings				
	Nuclide	Cs-137 / Co-60	Cs-137		
	Cal Method	1-POINT LIN MULTIPOINT 1- POINT EXP 2- POINT EXP	MULTIPOINT		
	Cal Curve Type	STANDARD / REVERSED	STANDARD		
	Meas. Path				
	Unit Family				
	Unit				
	ReadIn Time	S	60		
Device Config u Setup u Sensor Configuration u Date - Time					
	Date	MM/DD/YYYY			
	Time	hh:mm:ss			

Device Config u Setup u Sensor Configuration u Sensor Settings					
	Detector Code	0 50	0		
	HV Mode	AUTO / MANUAL			
	HV Live	V			
	HV Average	V			
	HV Manual	V			
	HV Default	V			
Device Config u S	ietup u Signal Condi	ition u Signal Parar	neter		
	Time Const	s	20		
	Error Handling	NORMAL / SENSITIVE	NORMAL		
	Signal Unlocked	OFF / ON	OFF		
Device Config u S	ietup u Signal Condi	tion u Reading Rang	je		
	Lower Range Value	SPG	0		
	Upper Range Value	SPG			
	Factor				
	Offset	SPG			
Device Config u	Setup u Signal Con	dition u Signal Dep	pendency		
	Response Mode	DISABLED RAPID SWITCH RAD INTERFERENCE	DISABLED		
	lo Factor		1,5		
	Waiting Time	S	60		
	RI Sigma		10		
	Meas Delay Time	S	4		



Device Config u Setup u I/O Setup u Current Output				
	Current Loop Montitoring	DISABLED ENABLED	ENABLED	
	Loop Alarm Type	HIGH LOW HOLD LAST VALUE VALUE	HIGH	
	Error Current Va- lue	mA	22,00	
	Current Lower Li- mit	mA	3,80	
	Current Upper Li- mit	mA	20,50	



Modifications due to technical advancement reserved.

© BERTHOLD TECHNOLOGIES GmbH & Co. KG 07/2019

Language: English Rev. no.: 00

Printed in Germany

#### BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Str. 22 75323 Bad Wildbad Germany www.Berthold.com




Unité d'évaluation **Duo** XPERT LB 47x Détecteurs **Duo** SERIES LB 4700

Informations sur la sécurité 56925BA59

Rev. No.: 03, 05/2019

# A propos de ce manuel d'utilisation

## 1.8 Avertissement

Les avertissements sont identifiés comme suit :

#### Signalement



Source et conséquence Explication si requise u Prévention En cas de danger

- Symboles d'alerte :
- · Signalement :
- · Source :
- · Conséquence :
- Prévention :
- En cas de danger :
- (triangle d'alerte) attire l'attention sur le risque. Indique la sévérité du danger.
- Précise le type ou la source de danger.
- Décrit les conséquences d'un non respect.
- Précise comment le risque peut être écarté.
- Précise quelles actions sont requises en cas
  - d'occurrence du risque

## 1.8.1 Symboles employés dans le manuel d'utilisation

Dans ce manuel, les avertissements indiqués avant les instructions d'utilisation se réfèrent aux risques de blessures ou de dégâts matériels. Les mesures de prévention de danger décrites doivent être respectées.



Indique un danger majeur imminent, qui entraînera certainement des blessures sérieuses ou la mort s'il n'est pas évité.

#### **AVERTISSEMENT**

Indique un danger potentiel qui peut entraîner des blessures sérieuses ou la mort s'il n'est pas évité.

#### PRUDENCE



Se réfère à une situation potentiellement dangereuse qui peut entraîner des blessures physiques mineures ou graves, ou des dégâts matériels si elle n'est pas évitée.



## RECOMMANDATION

1

Si cette information n'est pas appliquée, un dysfonctionnement et/ou un dégât matériel peuvent apparaître.

#### **IMPORTANT**

i

Les sections identifiées avec ce symbole signalent des informations importantes du produit ou de son fonctionnement.

Tip



Fournit des conseils sur l'application ou d'autres informations utiles.



## 1.8.2 Symboles utilisés sur l'appareil



#### Lire le manuel d'utilisation

Veuillez suivre les instructions dans ce manuel d'utilisation.

#### Décharge électrostatique

Veuillez noter les instructions de manipulation. Composants sensibles aux décharges électrostatiques. Veuillez suivre les instructions de ce manuel d'utilisation.

Connexion de mise à la terre



Raccorder le conducteur de mise à la terre à cet endroit.

#### **Raccordement equipotentiel**

Raccorder le conducteur d'équipotentialité à cet endroit

#### **Tension continue**

L'appareil fonctionne en tension continue et ne doit être raccordé qu'à une source de tension continue.

#### **Tension alternative**

L'appareil fonctionne en tension alternative et ne doit être raccordé qu'à une source de tension alternative.

#### Déchet non domestique



Cet appareil électrique ne doit pas être éliminé avec les déchets domestiques



## 1.9 Conformité

La société BERTHOLD déclare par la présente, sous son entière responsabilité, que la conception de ce produit mis sur le marché par BERTHOLD est conforme aux directives EU indiquées dans la déclaration de conformité originale.

Cette disposition devient nulle en cas de modifications non autorisées par Berthold ou dans le cas d'une utilisation impropre.

Pour la déclaration de conformité originale, se «Technical information».



# 2 Sécurité

## 2.1 Dangers et mesures de sécurité

- · Lire ces instructions entièrement et avec attention avant d'utiliser l'appareil.
- Stocker ces instructions dans un endroit accessible à tous les utilisateurs en permanence.

### 2.2 Utilisation appropriée

#### Ce qui suit constitue une utilisation appropriée :

- Se conformer strictement aux instructions et séquences d'utilisation mentionnées. Ne pas procéder à des pratiques différentes non autorisées qui pourraient engager votre sécurité et la fiabilité fonctionnelle de l'EVU !
- Suivre les instructions de sécurité mentionnées !
- Effectuer les opérations de maintenance prescrites ou les faire réaliser pour vous !
- Utiliser uniquement les accessoires et pièces de rechange BERTHOLD.



#### Utilisation inappropriée à éviter:

- Ne pas suivre les instructions de sécurité et les instructions pour l'utilisation, la maintenance et la mise au déchet indiquées dans le manuel.
- Un non respect quelconque avec le présent manuel d'utilisation pour le produit délivré.
- Appliquer des dispositions et conditions non conformes à celles mentionnées dans les documents techniques, feuilles de spécifications, manuels d'utilisation et instructions de montage, ou tout autre document spécifique du constructeur.
- Utiliser l'appareil si des éléments sont endommagés ou corrodés. Ceci s'applique aussi aux joints et aux câbles.
- · Modification ou changement des éléments du système.
- L'appareil ne doit pas être installé en atmosphère explosive et de ce fait, ne peut pas être utilisé dans une telle atmosphère. Il n'est pas antidéflagrant.
- Utilisation...
  - o où les éléments sous tension sont accessibles.
  - dans un boîtier mural avec presse-étoupes insuffisamment étanches et/ou non adaptés pour le passage des câbles.
- Utilisation sans les précautions de sécurité recommandées par le constructeur.
- Manoeuvre inappropriée ou oubli des équipements de sécurité présents.

BERTHOLD assume la responsabilité de la garantie seulement dans le cadre de ses spécifications publiées.

Si le produit est utilisé dans des conditions autres que celles décrites dans le présent manuel, la sécurité du produit est compromise et la garantie devient nulle.

#### RECOMMANDATION



L'appareil n'est pas conforme à IEC 61508 « Sureté de fonctionnement des systèmes comportant des composants électriques, électroniques ou électroniques programmables »



## 2.3 Qualification du personnel

#### RECOMMANDATION



Le minimum requis pour intervenir sur nos appareils ou pour les utiliser est un personnel avec des connaissances générales complétées par une formation d'un expert ou d'une personne autorisée.

A plusieurs endroits dans ce manuel d'utilisation, il est fait références à des groupes de personnes avec des qualifications particulières et à qui différentes tâches peuvent être confiées pendant l'installation, l'utilisation et la maintenance.

Les trois groupes de personnes sont :

- Employés avec des connaissances générales
- Experts
- Personnes autorisées

#### Employés avec connaissances générales

#### RECOMMANDATION

1

Les employés avec des connaissances générales doivent être guidés par un expert pour le moins. Lors de la mise en œuvre de matières radioactives, la personne compétente en radioprotection doit être consultée.

Les employés avec connaissances générales sont, par exemple, des techniciens, des soudeurs, qui vont assurer différentes tâches lors du transport, de l'assemblage et de l'installation de l'appareil sous l'encadrement d'une personne autorisée. Il peut s'agir aussi de personnel de montage du site. Les personnes concernées doivent posséder une expérience dans la manipulation du produit.

#### Experts

Les experts sont des personnes avec des compétences suffisantes dans le domaine requis, dues à leur formation spécialisée et qui sont familiers avec les lois nationales relatives à la santé et la sécurité, les règlements concernant la prévention des accidents, et les usages techniques applicables.

Le personnel expert doit être capable de déterminer et d'évaluer le résultat de ses tâches et doit être très familier avec le contenu de ce manuel d'utilisation.

#### Personnes autorisées

Les personnes autorisées sont celles désignées pour les tâches correspondantes dans le cadre de dispositions réglementaires, ou celles dûment autorisées par BERTHOLD pour des tâches particulières. Lors de la mise en œuvre de matières radioactives, la personne compétente en radioprotection doit être consultée



## 2.4 Les obligations de l'opérateur

L'opérateur de ces appareils doit régulièrement former son personnel sur les sujets suivants :

- Connaissance et utilisation du manuel d'utilisation et des clauses légales.
- Utilisation prévue de l'appareil.
- Respect des instructions de sécurité du site et des conditions d'utilisation de l'opérateur.
- Gestion régulière de la maintenance du produit.



Sous réserve de modifications dans le cadre du progrès technique.

 $^{\mbox{\scriptsize C}}$  BERTHOLD TECHNOLOGIES GmbH & Co. KG 05/2019

Sprache: Français Rev.-Nr.: 03

Printed in Germany

BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Str. 22 75323 Bad Wildbad Germany www.Berthold.com





## **EmulsionSENS**

Technical Information Technische Information

56925TI1M Rev. No.: 00, 07/2019



Evaluation unit Auswerteeinheit



## Installation variants 19" subrack

Einbauvarianten 19" Baugruppenträger



Item Pos.	Compontens Komponenten	Connection Anschluss
1	4x EVU (1 Master, 3 extension modules)	4 Terminal blocks 4 Klemmenblöcke
2	3x EVU, 1x blanking panel (1 Master, 2 extension modules)	3 Terminal blocks 3 Klemmenblöcke
3	2x EVU, 2x blanking panel (1 Master, 1 extension module)	2 Terminal blocks 2 Klemmenblöcke
4	1x Master EVU, 3x blanking panel	1 Terminal block 1 Klemmenblock

## Assignment terminal block master EVU

Belegung Klemmenblock Master AWE

		Ċ			A		
Signal	Pin		8	(		Pin	Signal
not assigned nicht belegt	C - 2		$\oslash$		$\oslash$	A - 2	not assigned nicht belegt
not assigned nicht belegt	C - 4		$\oslash$		$\oslash$	A - 4	not assigned nicht belegt
not assigned nicht belegt	C - 6		$\oslash$		$\oslash$	A - 6	not assigned nicht belegt
not assigned nicht belegt	C - 8		$\oslash$		$\oslash$	A - 8	not assigned nicht belegt
RELAY RELAIS 3 COM	C - 10		$\oslash$		$\oslash$	A - 10	RELAY RELAIS 3 NO
RELAY RELAIS 2 COM	C - 12		$\oslash$		$\oslash$	A - 12	RELAY RELAIS 2 NO
RELAY RELAIS 1 NC	C - 14		$\oslash$		$\oslash$	A - 14	RELAY RELAIS 2 NC
RELAY RELAIS 1 COM	C - 16		$\oslash$		$\oslash$	A - 16	RELAY RELAIS 1 NO
DIGITAL IN 1	C - 18		$\oslash$		$\oslash$	A - 18	DIGITAL IN 1 GND
DIGITAL IN 2	C - 20		$\oslash$		$\oslash$	A - 20	+ 24 V
CURRENT IN +	C - 22		$\oslash$		$\oslash$	A - 22	CURRENT IN -
RS 485 B (Detector LB 480)	C - 24		$\oslash$		$\oslash$	A - 24	RS 485 A (Detector LB 480)
CURRENT OUT –	C - 26		$\oslash$		$\oslash$	A - 26	CURRENT OUT +
not assigned nicht belegt	C - 28		$\oslash$		$\oslash$	A - 28	not assigned nicht belegt
Main Netz N, DC 24 V –	C - 30		$\oslash$		$\oslash$	A - 30	Main Netz L1, 24 V DC +
Protective conductor PE Schutzleiter PE	C - 32		$\oslash$		$\bigcirc$	A - 32	Protective conductor PE Schutzleiter PE



BERTHOLD

## 19" subrack

19" Baugruppenträger



Technical Data Technische Daten	
Dimensions Abmessungen	3HE/84TE/5T, 482x132x172mm (WxHxD)
Max. Assembly Max. Bestückung	- 4 Master
Weight (with circuit board, without modules) Gewicht (mit Anschlussplatine, ohne Module)	1.4 kg
Weight terminal block Gewicht Klemmenblock	220 g
Operational temperature Betriebstemperatur	-20°C +50°C, not condensing nicht kondensierend
Storage temperature Lagerungstemperatur	-30°C +60°C
Degree of protection Schutzklasse	IP20

## EVU AWE



Technical Data Technische Daten	
Dimensions Abmessungen	117/128/172mm (WxHxD)
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	-20°C +85°C
Degree of protection Schutzgrad	IP20
$\boldsymbol{\mathcal{C}}$	



Connections	<ul> <li>- USB port for the connection to the USB storage medium</li> <li>- Master/slave connection (4-pin) and plug</li> <li>- RJ45 connection for Ethernet (on back wall)</li> <li>- 32-pin plug connector according to DIN 19465 Series C</li> </ul>
Anschlüsse	- USB-Port zum Anschluss von USB-Speichermedium - Master/Slave Buchse (4-polig) und Stecker - RJ45-Buchse für Ethernet (an Rückwand) - 32 polige Stiftleiste nach DIN 19465 Baureihe C
Display	<ul> <li>graphical LCD display</li> <li>320 x 240 points, 262,000 colours</li> <li>Dimmable LED background lighting</li> <li>Touch screen</li> <li>graphisches LCD-Display</li> <li>320 x 240 Punkte, 262.000 Farben</li> <li>Dimmbare LED Hintergrundbeleuchtung</li> <li>Touchscreen</li> </ul>
Computer core	<ul> <li>Processor: Dual Core DSP/ARM Controller</li> <li>clock frequency: 300 MHz internal (20 MHz external quartz)</li> <li>ROM: 512 KByte</li> <li>RAM: 64 MByte ext. SDRAM, 128 KByte int. shared RAM</li> <li>FLASH: 8 MByte external serial</li> </ul>
Rechnerkern	- 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

Power Supply Stromversorgung	
Voltage Spannung	100-240 V AC 50/60 Hz (wide range input) +/- 10% 21-32 V DC (24 V DC power input)
Power consumption Leistungsaufnahme	22 VA, 15 W
Fuses Sicherungen	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	<ul> <li>4-20mA internally switched from power source to sink current (according to NAMUR recommendation NE 006 and NE 043).</li> <li>Continuous short circuit proof and isolated (500 V).</li> <li>Internal resistance about 105 ohms max.</li> <li>Burden when operating as a power source: 850 ohm.</li> <li>Internal monitoring of the loop current and additional error signalling by hardware on detection of a fault condition.</li> </ul>			
Stromausgang	4-20mA (nach Namur-Empfehlung NE 006 und NE 043) intern von Stromquelle auf Stromsenke umschaltbar. 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 Stromeingang	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 ohm max. input voltage: 24 V DC 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 Impuls-eingang	Frequency 0-100 kHz, Umax = 28 V, right angle signal form, low <1,5 V; high 4 – 28 V. Switchable to current input Frequenz 0-100 kHz, Umax = 28 V, Rechteck-Signalform, Low <1,5V; High 4 – 28 V. Umschaltbar auf Stromeingang			
Digital outputs Digitale Ausgänge	3 relays, Umax = 33 V ACeff, 46 V DC; Imax = 1 A functions: Relay 1: SPDT for error signalling Relay 2: SPDT assignable by software Relay 3: SPST assignable by software			
	3 Relais, Umax = 33 V ACeff, 46V DC; Imax = 1 A Funktionen: Relais 1: SPDT zur Fehlersignalisierung Relais 2: SPDT über Software zuweisbar Relais 3: SPST über Software zuweisbar			
Digital inputs Digitale Eingänge	2 x together electrically isolated (500 V) Switch between DigIn and GND, Uoutmax approx. 24 V Function configurable via software 2 x gemeinsam potentialgetrennt (500 V), Schalter zwischen DigIn und GND, Uoutmax ca. 24 V Funktion über Software konfigurierbar			
External supply Externe ersorgung	Output voltage:24 V DCOutput current:max. 150 mAAusgangsspannung:24 V DCAusgangsstrom:max. 150 mA			
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) für Master/Master Kommunikation und Prüf-und Testzwecke. Nicht potentialgetrennt von Hauptelektronik und USB-Anschluss potentialgetrennt von restlichen I/Os (500 V)			



USB port	1 x USB 2.0 Type A (Host) via front plate to the connection of an ext. mouse, keyboard or storage medium Uout = 5 V, loutmax = 0.5 A 1 x USB 2.0 Typ A (Host) über Frontplatte zum Anschluss einer ext. Maus, Tastatur oder Speichermedium
	Uout = 5 V, $Ioutmax = 0,5 A$
Ethernet	RJ45 connection via back wall, 10 Mbit, DHCP supported, max. 3 m
	RJ45-Buchse über Rückwand, 10 Mbit, DHCP unterstützt, max. 3 m



## Number Key LB 47x

Nummernschlüssel LB 47x



\* used by othe hardware



## Declaration of Conformity Konformitätserklärung



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: radiomatric evaluation system DuoXpert

Type / model: LB 47x

	directiv	/e		applied star	ndards
LVD	2014/35	5/EU		EN 61010-1	2010
RoHS	2011/65	5/EG			
EMC	2014/30	D/EU		EN 61326-1 EN 61000-4- EN 61000-4- EN 61000-4- EN 61000-4- EN 61000-4- EN 61000-4- EN 61000-4-	2013 2 3 4 5 6 6 11
				Namur NE21	2012
This d	leclaration is iss	sued by the manufacturer			
BERTH Calmb	HOLD TECHNOL bacher Str. 22,	.OGIES GmbH & Co. KG D-75323 Bad Wildbad, Germ	nany		
releas Dr. Jü Head Bad W	rgen Briggman of R&D /ildbad, 1 <sup>st</sup> of S	n eptember, 2015			
Pers	Reç sönlich haftende Gese Reç Deutsche	jistergericht / Court of Registration Ilschafterin / Fully liable Associates jistergericht / Court of Registration Geschäftsführung / Management USL-Id-Nr. / VAT Reg. No. e Steuernummer / German Tax No. WEEE-Reg. No.	Stuttgart H BERTHOLD Stuttgart H Horst Knau DE8130505 49038/080 DE9946869	RA 330991 TECHNOLOGIES Verwa RB 331520 ff, Dr. Dirk Mörmann 11 38 0	iltungs-GmbH
PF-CW bank	75323 Bad Wildbad 75119 Pforzheim 75105 Pforzheim	Konto/Account No. 8 045 003 (BLZ 666 5 Konto/Account No. 957 004 (BLZ 666 9 Konto/Account No. 6 511 120 (BLZ 666 8	00 85) SWIF 00 00) SWIF 00 13) SWIF	T-BIC PZHSDE66 F-BIC VBPFDE66 F-BIC DRES DEFF 666	IBAN: DE37 6665 0085 0008 0450 0 IBAN: DE85 6669 0000 0000 9570 0 IBAN: DE05 6668 0013 0651 1120 0



Sparkas

Volksbar Commer



#### 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.

LB 47x

Produktbezeichnung:

#### *radiometrisches Auswertesystem DuoXpert*

Typenbezeichnung / Modell:

angewendete Normen und Richtlinie (Fundstelle) 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 Briggmahn Leiter Entwicklung Bad Wildbad, den 1. September 2015

Registergericht / Court of Registration Persönlich haftende Gesellschafterin / Fully liable Associates Registergericht / Court of Registration Stuttoart HRA 330991 Stuttgart HRA 330991 BERTHOLD TECHNOLOGIES Verwaltungs-GmbH Stuttgart HRB 331520 Horst Knauff, Dr. Dirk Mörmann DE813050511 Registergericht / Court or Kegistrauon Geschäftsführung / Management USL-Id-Nr. / VAT Reg. No. Deutsche Steuernummer / German Tax No. WEEE-Reg. No. 49038/08038 DE99468690 Konto/Account No. 8 045 003 (BLZ 666 500 85) Konto/Account No. 957 004 (BLZ 666 900 00) Konto/Account No. 6 511 120 (BLZ 666 800 13) Sparkasse PF-CW 75323 Bad Wildbad SWIFT-BIC PZHSDE66 IBAN: DE37 6665 0085 0008 0450 03 IBAN: DE85 6669 0000 0000 9570 04 IBAN: DE05 6668 0013 0651 1120 00 Volksbank 75119 Pforzheim SWIFT-BIC VBPEDE66 SWIFT-BIC DRES DEFF 666 Commerzbank 75105 Pforzheim



## Certificates Zertifikate

## NRTL certification US/CAN wall-mounted housing

NTRL Zertifikat US/CAN Wandgehäuse

Nen	nko C	Certificate of Compliance
Nemko-C	CCL, Inc.	
Certificate:	NA201610530	Date Issued: January 20, 2016
Project:	257087-7.1	
Issued to:	Berthold Technologies GmbE Calmbacher Straße 22 75323 Bad Wildbad Germany	I & Co. KG
The produc tl	ts listed below have been certifie he specifications listed and are el	d as being compliant with all applicable requirements of ligible to bear the following certification mark
	cNe	<b>mko</b> <sub>us</sub>
Issued by:	Potent Keller	Robert Keller, Senior Engineer/Safety Supervisor
Authorized	by: The	Thomas Jackson, Certification Manager
PRODUCTS	<u>s</u>	
	MENT. CONTROL. OR LABOR.	ATORY EQUIPMENT – Certified to US and Canada
MEASUREN Standards	,	
MEASUREN Standards Product: Pro Model: Wall different soft Ratings: Wa 2M: 44VA 10	ocess measurement unit I-mounted LB 47x, 1M/3S; Wall- ware versions for the master and ill-mounted LB 47x, 1M/3S: 40V, 00-240V, 50/60Hz, Class I	mounted LB 47x, 2M (x can be 0 to 8 and describes slave modules not affecting safety). A 100-240V, 50/60Hz, Class I; Wall-mounted LB 47x,
MEASUREN Standards Product: Pro Model: Wall different soft Ratings: Wa 2M: 44VA 10	ocess measurement unit I-mounted LB 47x, 1M/3S; Wall- ware versions for the master and ill-mounted LB 47x, 1M/3S: 40V, 00-240V, 50/60Hz, Class I	mounted LB 47x, 2M (x can be 0 to 8 and describes slave modules not affecting safety). A 100-240V, 50/60Hz, Class I; Wall-mounted LB 47x,
MEASUREN Standards Product: Pro Model: Wall different soft Ratings: Wa 2M: 44VA 19	ocess measurement unit I-mounted LB 47x, 1M/3S; Wall- ware versions for the master and ill-mounted LB 47x, 1M/3S: 40V. 00-240V, 50/60Hz, Class I	mounted LB 47x, 2M (x can be 0 to 8 and describes slave modules not affecting safety). A 100-240V, 50/60Hz, Class I; Wall-mounted LB 47x,
MEASUREN Standards Product: Pro Model: Wall different soft Ratings: Wa 2M: 44VA 10	ocess measurement unit I-mounted LB 47x, 1M/3S; Wall- ware versions for the master and ill-mounted LB 47x, 1M/3S: 40V. 00-240V, 50/60Hz, Class I	mounted LB 47x, 2M (x can be 0 to 8 and describes slave modules not affecting safety). A 100-240V, 50/60Hz, Class I; Wall-mounted LB 47x,
MEASUREN Standards Product: Pro Model: Wall different soft Ratings: Wa 2M: 44VA 10 The confifcation system	n, az described in ISO/IEC Guide 67 (Conformity Assessme	mounted LB 47x, 2M (x can be 0 to 8 and describes slave modules not affecting safety). A 100-240V, 50/60Hz, Class I; Wall-mounted LB 47x, nu-Fundamentals of Product Confiftcation), most classly resembles System 3



# NRTL certification US/CAN wall-mounted housing (continued) NTRL Zertifikat US/CAN Wandgehäuse (Fortsetzung)

APPLICA	BLE REQUIREMENTS	
UL Std.	No. 61010-1 2nd Edition	<ul> <li>Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: Gener Requirements</li> </ul>
CAN/CSA	-C22.2 No. 61010-1-04 Secon	d Edition - Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: Gener Requirements
This certificate	is issued on condition that the holder compli- pursuant to the terms	es and will continue to comply with the requirements of the above mentioned specifications a and conditions specified in the Certification Agreement.
_		

# NRTL certification US/CAN wall-mounted housing (continued) NTRL Zertifikat US/CAN Wandgehäuse (Fortsetzung)

	Supplement	to Certificate of Compliance
Certificate:	NA201610530	Project: 257087-7.1
Nemko-CCL	grants a license to the appl and that the mark shall on	icant to apply the Certification Mark to the certified products ily be affixed at the following factory locations
	F	actory Information
Factory Nam	ie	Location
Berthold Tech	hnologies GmbH & Co. KG	Calmbacher Straße 22 75323 Bad Wildbad Germany
	The products listed, in are eligible to be marked	cluding the latest revision described below, in accordance with the referenced Certificate.
	Produ	ct Certification History
Project	Date	Description
257087-7.1	January 20, 2010	Wall-mounted LB 47x, 2M (x can be 0 to 8 and describes different software versions for the master and slave modules no 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 Suppleme	ent forms an integral part of the	Certificate of Compliance
		(compared Freedom take of Brachast Contification) must cleach surroubler Output 2
The certification system	n, as described in ISO/IEC Guide 67 (Conformity 2	Esessment – Franklimentals of Frontici Certylcauony, most closely resembles system 5

# NRTL certification US/CAN DuoXpert LB 47x NTRL Zertifikat US/CAN DuoXpert LB 47x

Nen	nko	Certificate of Compliance
Nemko-C	CL, Inc.	
Certificate:	NA201510498	Date Issued: September 17, 2015
Project:	235982-14.1	
Issued to:	Berthold Technologies G Calmbacher Straße 22 75323 Bad Wildbad Germany	mbH & Co. KG
The produc	ts listed below have been ce he specifications listed and	rtified as being compliant with all applicable requirements of are eligible to bear the following certification mark
	c N	lemko
Issued by:	Robert Keller	Robert Keller, Senior Engineer/Safety Supervisor
Authorized	by: The	Thomas Jackson, Certification Manager
PRODUCTS MEASUREN Standards	<u>š</u> MENT, CONTROL, OR LA	BORATORY EQUIPMENT – Certified to US and Canada
Product: Pro Model: Duol	ocess measurement unit for t Xpert LB47x-02-M; DuoXp sions for the master and slav 47x-02-M: 100-240V AC 22	ouilding-in ert LB47x-02-S (x can be 0 to 8 and describes different e modules not affecting safety) 2VA 50/60Hz; LB47x-02-S: 100-240V AC 6VA 50/60Hz
software vers Ratings: LB		
software vers Ratings: LB <u>APPLICAB</u>	LE REQUIREMENTS	
software vers Ratings: LB <u>APPLICAB</u> UL Std.	LE REQUIREMENTS No. 61010-1 3rd Edition	<ul> <li>Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements</li> </ul>
software vers Ratings: LB <u>APPLICAB</u> UL Std. CAN/CSA-C	LE REQUIREMENTS No. 61010-1 3rd Edition 22.2 No. 61010-1-12 Third	<ul> <li>Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements</li> <li>Edition – Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements</li> </ul>
software vers Ratings: LB <u>APPLICAB</u> UL Std. CAN/CSA-C This certificate is i	LE REQUIREMENTS No. 61010-1 3rd Edition 22.2 No. 61010-1-12 Third ssued on condition that the holder complia pursuant to the terms	<ul> <li>Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements</li> <li>Edition – Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements</li> <li>and will continue to comply with the requirements of the above mentioned specifications and and conditions specified in the Certification Agreement.</li> </ul>
software vers Ratings: LB <u>APPLICAB</u> UL Std. CAN/CSA-C This certificate is i	LE REQUIREMENTS No. 61010-1 3rd Edition 222.2 No. 61010-1-12 Third ssued on condition that the holder complie pursuant to the terms	<ul> <li>Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements</li> <li>Edition – Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 1: General Requirements</li> <li>and will continue to comply with the requirements of the above mentioned specifications and and conditions specified in the Certification Agreement.</li> <li>Assessment – Fundamentals of Product Certification), most closely resembles System 3</li> </ul>

# NRTL certification US/CAN DuoXpert LB 47x (continued) NTRL Zertifikat US/CAN DuoXpert LB 47x (Fortsetzung)

	Supplement	to Certificate of Compliance
Certificate:	NA201510498	Project: 235982-14.1
Nemko-CCL	grants a license to the appl and that the mark shall on	icant to apply the Certification Mark to the certified products nly be affixed at the following factory locations
	F	actory Information
Factory Nam	e	Location
Berthold Tech	mologies GmbH & Co. KG	Calmbacher Straβe 22 75323 Bad Wildbad Germany
	The products listed, in are eligible to be marked	cluding the latest revision described below, in accordance with the referenced Certificate.
	Produ	ct 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 differen software versions for the master and slave modules not affecti safety) Ratings: LB47x-02-M: 100-240V AC 22VA 50/60Hz; LB472 02-S: 100-240V AC 6VA 50/60Hz
This Suppleme	nt forms an integral part of the	Certificate of Compliance
	as described in ISO/IEC Guide 67 /Confermity .	(corement - Fundamentals of Product Cartification) most closely recembles System 3
The certification system.	the description of 1969 and of 1969 mary 2	a second a second construction of a second construction of a second



## Parts overview

Übersicht Zubehör

ID. No. <i>Id. Nr.</i>	Description Beschreibung	
68867	LB 478-01-M MPLMTransmitter (Master, 24 VDC) LB 478-01-M MPLM-Messgerät (Master, 24 VDC)	
68866	LB 478-02-M MPLM Transmitter (Master, 100240 VAC) LB 478-02-M MPLM-Messgerät (Master, 100240 VAC)	
68869	LB 478-01-E Transmitter (Extension, 24 VDC) LB 478-02-E Messgerät (Erweiterung, 100240 VAC)	
68868	LB 47x-02-E Transmitter (Extension, 100240 VAC) LB 47x-02-E -Messgerät (Erweiterung, 100240 VAC)	
56925-8BA1	Operating manual DuoSeries LB 478 MPLM, German Betriebsanleitung DuoSeries LB 478 MPLM, Deutsch	
56925-8BA2	Operating manual DuoSeries LB 478 MPLM, English Betriebsanleitung DuoSeries LB 478 MPLM, Englisch	
64607	19" rack, 84 HP / 3 RU for use with terminal blocks 19"-Baugruppenträger für den Einsatz mit Klemmblöcken	
59477	Terminal block for LB 47x, Master Klemmenblock für LB 47x, Master	
37526	Front Cover Plate 21 HP / 3 RU Blindplatte 21TE / 3 HE	

Modifications due to technical advancement reserved.

Änderungen im Zuge technischer Weiterentwicklung vorbehalten.