



Concentration Meters MicroPolar (++) LB 566

User's Guide Hardware Manual 41986BA2

Rev. Nr.: 08, 07/2023

The units supplied should not be repaired by anyone other than Berthold Technologies Service engineers or technicians by Berthold Technologies.

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

The complete user's guide consists of the hardware manual and the software manual.

The hardware manual comprises:

- mechanical components
- > installation
- electrical installation
- technical data
- electrical and mechanical drawings

The **software manual** comprises:

- > operation of the evaluation unit
- parameter description
- basic setup
- ➤ calibration
- error messages

The present manual is the hardware description.

Subject to change without prior notice.

BERTHOLD TECHNOLOGIES GmbH & Co. KG	
Calmbacher Str. 22 · D-753	23 Bad Wildbad
Switchboard:	Service:
Phone +49 7081 177 0	Phone +49 7081 177 111
Fax +49 7081 177 100	Fax +49 7081 177 339
industry@Berthold.com www.Berthold.com	Service@Berthold.com

Table of Contents

	Page
Chapter 1. Safety Summary	7
1.1 Symbols and Warnings1.2 General Information1.3 General Safety Instructions	7 8 9
Chapter 2. General Information	11
2.1 Use and Function2.2 Frequency License2.3 Intended Use2.4 Explanation of Terms	11 12 14 15
Chapter 3. System Description	17
 3.1 Principle of Measurement 3.2 Calculation of Measured Values 3.3 Temperature Compensation 3.4 Throughput Calculation and Output 3.5 Mechanical Components 3.5.1 Evaluation Units 3.5.2 FlowCell 3.5.3 Container Probe 3.5.4 High-frequency Cable 3.6 Assembly on a Pipeline 3.7 Assembly on a Container 	17 18 19 20 21 23 25 28 32 34 35
Chapter 4. Getting Started	36
 4.1 Transport 4.2 Commissioning the FlowCell 4.2.1 Installing the FlowCell 4.2.2 Installing the Evaluation Unit 4.2.3 Connecting the HF Cable 4.3 Commissioning the Container Probe 4.3.1 Installing the Container Probe 4.3.2 Setting Up the Evaluation Unit 4.3.3 Connecting the HF Cable 4.5 Connecting the Evaluation Unit 4.5.1 Pin Configuration of the Connector Strip 4.5.2 Digital Outputs, Relay 	36 37 40 41 43 43 45 45 45 47 48 50
Chapter 5. Service instructions	53
5.1 General Information 5.2 Wear Parts 5.3 Instrument Cleaning 5.4 Battery 5.5 Fuse Replacement	53 54 56 57 58

Chapter 6. Tec	chnical Data	60
6.1	Technical Data Evaluation Unit	60
6.2	Technical Data Sensors	63
6.3	Technical Data HF Cable	69
6.4	Format of Serial Data Output RS232 and RS485	/1
Chapter 7. Cer	tificates	73
7.1 7.2	EC Declaration of Conformity Frequency License	73 76
Chapter 8. Tec	chnical Drawings	85
8.1 8.2 8.3 8.4 8.4 8.4 8.5	Dimensions Drawings Evaluation Unit Wall Housing Electrical Wiring Diagram Dimensional Drawings FlowCell 8.3.1 Type LB 5660-102-00X FlowCell DN 50 Flange, FDA 8.3.2 Type LB 5660-202-00X FlowCell DN 65 Flange, FDA 8.3.3 Type LB 5660-302-00X FlowCell DN 80 Flange, FDA 8.3.4 Type LB 5660-402-00X FlowCell DN 100 Flange, FDA 8.3.5 Type LB 5660-602-00X FlowCell DN 125 Flange, FDA 8.3.6 Type LB 5660-602-00X FlowCell DN 150 Flange, FDA 8.3.7 Type LB 5660-112-00X FlowCell DN 50 G-BS/M 8.3.8 Type LB 5660-112-00X FlowCell DN 50 G-BS/M 8.3.9 Type LB 5660-112-00X FlowCell DN 50 G-BS/M 8.3.10 Type LB 5660-12-00X FlowCell DN 100 G-BS/M 8.3.11 Type LB 5660-12-00X FlowCell DN 100 G-BS/M 8.3.12 Type LB 5660-12-00X FlowCell DN 100 G-BS/M 8.3.13 Type LB 5660-132-00X DN 50 8.3.14 Type LB 5660-132-00X DN 50 8.3.14 Type LB 5660-322-00X DN 50 8.3.15 Type LB 5660-32-00X DN 50 8.3.16 Type LB 5660-432-00X DN 100 8.3.17 Type LB 5660-432-00X DN 150 Dimensional Drawings Container Probes 8.4.1 Type LB 5650-01 8.4.2 Type LB 5650-02 8.4.3 Type LB 5650-03 8.4.4 Type LB 5650-03 8.4.4 Type LB 5650-03 8.4.4 Type LB 5650-03 8.4.7 Installation Situation in Pipelines Dimensional Drawings Container Flush Probes 8.5.1 Type LB 5651-01 8.5.2 Type LB 5651-01 8.5.2 Type LB 5651-02 8.5.3 Type LB 5651-03 8.5.4 Type LB 5651-04 8.5.5 Type LB 5651-05 8.5.6 Installation Situation in Pipelines Assembly sheets for LB 5650 (Container probe) Accembly cheets for LB 5651 (Container probe)	85 87 88 88 89 90 91 92 93 94 95 96 97 98 90 101 102 103 104 105 106 107 108 109 110 111 112 113 114 115 116 117 118 119 121



Chapter 1. Safety Summary

1.1 Symbols and Warnings

In this user manual, the term Berthold Technologies stands for the company Berthold Technologies GmbH & Co.KG.

To rule out bodily injury and property damage, please keep in mind the warning and safety instructions provided in this user manual. They are identified by the following sings: DANGER, WARNING, CAUTION or NOTICE.

A DANGER	Indicates imminent danger. If it cannot be avoided, death or most severe personal injuries may be the consequence.
	Indicates a possibly dangerous situation. The consequences may be death or most severe personal injuries.
	Indicates a possibly harmful situation The consequences may be minor or medium personal injuries.
NOTICE	Indicates a situation that may cause material damage if the instructions are not followed.
	IMPORTANT Paragraphs with this symbol provide important information on the product and how to handle it.



Contains user tips and other useful information.

Other symbols used in this documentation:



1.2 General Information

The most important safety measures a summarized in this user manual. They supplement the corresponding regulations which *must* be studied by the personnel in charge.

Please pay attention to:

- > the national safety and accident prevention regulations
- > the national assembly and installation directions
- > the generally recognized engineering rules
- the information on transport, assembly, operation, service, maintenance
- the safety instructions and information in these operating instructions
- > the enclosed technical drawings and wiring diagrams
- the characteristic values, limit values and the information on operating and ambient conditions on the type labels and in the data sheets
- the signs on the devices
- > the country-specific licensing schemes

1.3 General Safety Instructions



	Electrical hazards
	Disconnect power to ensure that contact with live part is avoided during installation and when servicing.
	Disconnect the power supply before opening the instrument. Work on open and live instruments is prohibited.
NOTICE	Caution! Potential hazards, material damage! Device type:
	LB 566-12 MicroPolar ++ (ID no. 51833-02) When connecting the 24 V DC power supply, the + and – poles must be connected correctly. There is no reverse polarity pro- tection!
NOTICE	Spare fuses must match the rating specified by the device man- ufacturer. Short-circuiting or manipulation is not permitted.
	IMPORTANT The LB 566 and all additional equipment must be connected to mains via grounded connection.
	I IMPORTANT
	The concentration meter LB 566 may be installed, serviced and repaired only by qualified specialists.
Qualified persons	Qualified specialists are persons who through professional train- ing have acquired sufficient skills in the respective field and who are familiar with the relevant national industrial safety regula- tions, accident prevention directions, guidelines and accepted en- gineering standards. They must be able to safely assess the re- sults of their work and they must be familiar with the contents of these user manual.

Chapter 2. General Information

2.1 Use and Function

	The MicroPolar LB 566 has been designed as a concentration measuring system and may be used only for this purpose. If the devices are used in a manner that are not described in this user manual, the protection of the devices is compromised and the warranty will be lost.
	Berthold Technologies is liable and guarantees only that the devices comply with its published specifications. The LB 566 may only be installed in an undamaged, dry and clean condition. Alterations and modifications to the system components are not permitted.
	The LB 566 is not qualified as a "safety-related measurement".
Conformity to standards	The standards and guidelines the LB 566 complies with are item- ized in these device instructions in <i>chapter 2.2 Frequency License</i> and <i>chapter 7.1 EC Declaration of Conformity</i> .
Pressure equipment directive	The FlowCell and the container probe has been classified as pres- sure equipment acc. to art. 4 sect. 3 of guideline 2014/68/EU. Before use, check whether the case of use corresponds to this classification. In particular, the media compatibility of the com- ponents in contact with the fluid must be checked.
Protection type	The protection type of the LB 566 to IEC 60529 is max. IP 65.
	The following use is inappropriate and has to be prevented:
Warning against misuse	The use under conditions other than the terms and conditions stated by the manufacturer in their technical documentation, data sheets, operating and installation manuals and other specifications.
	The use after repair by persons not authorized by Berthold Technologies.
	The use in a damaged or corroded state.
	 Operation with open or inadequately closed cover.
	 Operating with insufficiently tightened adapters and cable glands.
	 Operation without the manufacturer's recommended safety precautions.
	 Manipulating or bypassing existing safety facilities.
Authorized persons	Authorized persons are those who, by law, are permitted to per- form the respective activity, or who have been approved by Berthold Technologies for certain activities.

2.2 Frequency License

This device complies with Part 15 of the FCC Rules and with Industry Canada licence-exempt RSS standard(s).

Operation is subject to the following two conditions:

- (1) this device may not cause harmful interference, and
- (2) this device must accept any interference received, including interference that may cause undesired operation.

Le présent appareil est conforme aux CNR d'Industrie Canada applicables aux appareils radio exempts de licence.

L'exploitation est autorisée aux deux conditions suivantes:

- (1) l'appareil ne doit pas produire de brouillage, et
- (2) l'appareil doit accepter tout brouillage radioélectrique subi, même si le brouillage est susceptible d'en compromettre le fonctionnement.

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

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

License labels

This equipment has been tested and found to comply with the limits for a Class B digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference in a residential installation. This equipment generates, uses and can radiate radio frequency energy and, if not installed and used in accordance with the instructions, may cause harmful interference to radio communications. However, there is no guarantee that interference will not occur in a particular installation. If this equipment does cause harmful interference to radio or television reception, which can be determined by turning the equipment off and on, the user is encouraged to try to correct the interference by one or more of the following measures:

- Reorient or relocate the receiving antenna.
- Increase the separation between the equipment and receiver.

- Connect the equipment into an outlet on a circuit different from that to which the receiver is connected.
- Consult the dealer or an experienced radio/TV technician for help.

Changes or modifications made to this equipment not expressly approved by BERTHOLD TECHNOLOGIES may void the FCC authorization to operate this equipment.

The MicroPolar and MicroPolar ++ comply with the R&TTE Directive 1999/5/EC and thus meet all the requirements for this type of high-frequency devices. As a mark of conformity in accordance with the CE mark, the devices bear the no. 0682 of the certification body. The certificate can be found in chapter 7.3 Frequency License.

€ € 0682

IMPORTANT

The frequency approvals and conformities are only applicable in combination with the container probes and the FlowCell.

To comply with ETSI (European Telecommunications Standards Institute), the LB 566 ++ in combination with horn and spiral antennae may only be operated with a 10dB attenuator pad in the sending path (MTX).



IMPORTANT

The LB 566 has been manufactured in compliance with the safety requirements for microwave devices. It will be the user's responsibility to adhere to any special legal provisions regarding the use of microwaves.



IMPORTANT

Any change in the frequency or otherwise tampering with the microwave device will lead to a loss of the frequency license and may result in criminal consequences.

The microwave modules do not include any replaceable components and must not be opened.

2.3 Intended Use

The LB 566 can be used to determine the water or moisture content and the concentration of virtually any material. The following sensors and evaluation versions are available:

- 1. The container probes have been designed for installation into containers and pipelines with a nominal width of \geq 200 mm. The probe is mounted so that both measuring rods (transmitter and receiver) are immersed into the product being measured.
- 2. The The flow cell is a tubular probe that is either installed into the existing pipeline system inline or into a bypass.

The evaluation unit is available in two versions: The standard model MicroPolar and the high dynamic version MicroPolar ++. The Micro Polar ++ requires a microwave signal attenuation of at least 40 dB. The MicroPolar must be used for lower microwave attenuation.

During operation, the concentration meters MicroPolar and MicroPolar ++ emit electromagnetic radiation in the frequency range between 2.4 GHz and 2.5 GHz (range limitations depending on local regulations in your country). Microwaves are not dangerous to human beings and the environment (power radiation < 10 mW). The microwaves are emitted directed from the microwave window; the product is not altered by the microwaves.

To ensure proper function of the measuring system, please pay attention to the following:



- The material to be measured may be electrically conductive only to a limited degree.
 - The product must not contain any gas bubbles or gas bubbles have to be compressed with adequate pressure when carrying out measurements in pipelines.
 - The ion concentration, for example, salt content must be nearly constant.

2.4 Explanation of Terms

Attenuation	Weakening of microwave signals, microwaves measuring effect
EVU	Evaluation Unit
Factory setting	In the factory setting all parameters have been set to default values. In most cases this considerably facilitates the calibration of the instrument. Despite factory setting, calibration al-ways has to be performed.
FlowCell	Tubular probe for simple integration into the existing pipeline system.
Flush probe	Container probe with flushing device.
HF cable	High frequency cable
Microwaves	Designation for electromagnetic waves in a specific frequency range.
Phase	Phase or phase shift, microwave measuring effect.
Quad cable	Combination of four HF cables of equal length in a corrugated tube.
Softkeys	Software associated keys.
ТС	Temperature compensation



Chapter 3. System Description

3.1 Principle of Measurement

As the microwaves pass through the product, their propagation velocity is slowed down (= phase shift) and their intensity is attenuated (= attenuation). Figure 3-1 illustrates the principle of measurement: Compared to a reference signal, the propagation velocity of microwaves passing through the product is slowed down (phase shift) and their intensity (attenuation) is reduced.



Figure 3-1: Schematic diagram: Change in microwave by product

The prerequisite is that the product being measured shows dielectric properties. Generally, water is a very distinct dielectric fluid. The water or dry mass concentration can therefore be determined by measuring the phase shift and/or attenuation.

The concentration to be detected in the product is therefore in good approximation linearly dependent on the phase shift and the attenuation. For this reason, we can measure the concentration or the dry matter content of the product using a linear calibration (see *chapter 3.2 Calculation of Measured Values*).

3.2 Calculation of Measured Values

The microwave parameters phase and attenuation are calibrated according to an automatic plausibility analysis.

During calibration, the phase and/or attenuation of a concentration value (or density value) are assigned through sampling. The calibration is done automatically and the sampling process is supported by the evaluation unit.

Which of the parameters (phase, attenuation or both) will be used for the calibration depends on the size and interference of the measuring effect. For example, the attenuation is significantly more sensitive to electrolytic conductivity (salt content).

In many cases, the pure phase measurement is recommended and the measured value is calculated as follows:

Measured value = $A \cdot Phase + C$ Eq. 3-1

where:

Measured valueConcentration / Moisture / Dry matterA, CCoefficients of the respective calibration
function

The LB 566 allows you to calibrate, display and output two concentrations: Con1 and Con2. You have to enter the calibration coefficients separately for concentration 1 and 2. For more information please refer to the Software Manual.

Limitations

- Weakly bound water can be detected depending on the strength of the binding. Thus, the measuring effect may be dependent on the grain size distribution and the chemical properties of the product being measured, provided this changes the binding of water to the solid matter.
- Walls made of plastic, rubber or insulation materials with fairly low dielectricity hardly affect the measurement and are calibrated at a constant level.
- Ice and crystal water cannot be measured because the water molecules cannot rotate freely (ice and crystal water are dry).

Conductive materials such as graphite or coke cannot be transmitted by microwaves. Metal walls can also not be transmitted by microwaves.

Compensation In addition to the water content, the product temperature, product density and a varying material load (varying microwave irradiation path) may have an influence on the phase and attenuation. This influence has to be compensated for during calibration.

3.3 Temperature Compensation

Temperature compensation (TC) is required for fluctuating product temperature. It is generally advisable to provide a temperature compensation, i.e. a temperature signal (0/4...20 mA or PT100) to be connected to the evaluation unit and, optionally, to enable the compensation in the evaluation unit. The evaluation unit is designed so that the required TC's can be calculated automatically. The variation in temperature where TC becomes absolutely essential is dependent on the product and on the water content. In the first approximation, \pm 2° C should be set as fluctuation limit.

For example, if the product temperature is measured via the PT100 input, then Eq. 3-1 is expanded as follows:

Measured value = $A \cdot Phase + D \cdot T_{meas} + C$ Eq. 3-2

where:

Measured value Concentration / Moisture / Dry matter A, D, C Coefficients of the calibration function T_{meas} Product temperature

How to work with the temperature compensation is described in detail in the Software Manual.

3.4 Throughput Calculation and Output

For pipeline applications, the LB 566 allows you to calculate the throughput (mass flow) and to output the result via a current output.

The calculation is based on the microwave measured value; if this value correlates with the product density, one can calculate the throughput, if some additional information is available. The additional information needed is: internal pipe diameter/cross section and the product speed. The product speed has to be fed via current input.

For details, please see the *Software Manual, chapter 4.2.12 Mass Flow.*

3.5 Mechanical Components

The measuring system comprises an evaluation unit, a probe/antenna pair and one set of special high-frequency cables (short HF-cable). The evaluation unit is available in two versions: the standard model MicroPolar LB 566 and the high dynamics version MicroPolar ++ LB 566 (see Figure 3-4 and Figure 3-5).



Figure 3-4: Evaluation unit MicroPolar LB 566



Figure 3-5: Evaluation unit MicroPolar ++ *LB 566* The probes/antenna pairs are available in different versions, as pipeline and container probe with or without flushing device (see Figure 3-6 to Figure 3-7).





Figure 3-7: FlowCell LB 5660-102-00x nominal width 50 mm with V flange



The FlowCell is available in the nominal pipe sizes 50 -150 mm (50, 65, 80, 100, 125 and 150). The following connections types are available:

- Hygiene milk pipe screw connection DIN 11853-1
- V flange EN 1092-1/11
- Welding pipe

3.5.1 Evaluation Units

The evaluation units comprise the evaluation computer and the microwave unit. The microwaves are generated, received and analyzed in the microwave unit. Signal processing and communication take place in the evaluation computer. For simple operation, the measuring system includes a display, 4 softkeys and an alphanumeric keypad. Different functions are assigned to the softkeys on the display.

Differences between MicroPolar ++ und MicroPolar

The MicroPolar ++ evaluation unit has an additional HF amplifier module in comparison to the standard model; therefore, it also has a larger wall housing (dimensions see *chapter 6.1 Technical Data Evaluation Unit*). Otherwise, the evaluation units differ only in their application.

MicroPolar ++ The high dynamics version MicroPolar ++ permits higher product attenuations. Larger measuring paths can be irradiated, i.e. FlowCells with larger nominal diameters can be used. Which type of evaluation unit is used depends on the product attenuation. MicroPolar is used up to an attenuation of 50 dB; MicroPolar ++ is used for higher attenuations. The MicroPolar ++ generally requires an attenuation of 40 dB. If this attenuation is not reached, the software displays an error message.

An RS232 interface is included on the bottom side of the instrument.





BERTHOLD

Interpretation of LEDs

Five LEDs on the instrument front panel indicate the current device status.

- Run
- Error
- Signal 1

Figure 3-10: LEDs on the front panel of the evaluation unit

- Signal 2
- Comm

LED	Function
Run	<u>On:</u> Device in measurement mode <u>Flashes + ERROR LED off:</u> Device in warning state, on hold, paused or low load state. A display message with error code indicates the cause (see <i>Software Manual, chapter 11. Error</i> <i>Lists and Device States).</i>
Error	<u>On:</u> Device in error state. A display message with error code indicates the cause (see <i>Software Manual, chapter 11. Error</i> <i>Lists and Device States).</i> Canceled after reset or if error has been elimi- nated.
Signal 1	Display depending on the selected function of relay 1, possible functions: Error, no product, alarm min., alarm max., measurement stopped, low load
Signal 2	Display depending on the selected function of relay 2, possible functions: Error, no product, alarm min., alarm max., measurement stopped, low load
Comm	Communication active, e.g. via RS232 and RS485

Terminal block

The electrical connections of the LB 566 are located on a connector strip in the wall cabinet. The terminal block can be accessed from the front by opening the cover of the housing. There, you also find the power cut-off switch and the fuses. The highfrequency terminals are located on the outside of the housing. All other elements, especially the voltage-carrying elements (on the motherboard) are provided with a protective cover.

3.5.2 FlowCell

The FlowCell is available in the nominal sizes of 50 to 150 mm (see fig. 3-11). As connection, the versions V flange EN 1092-1/11 or Hygiene milk pipe screw connection DIN 11853-1 are available. For technical data please see *chapter 6.2 Technical Data Sensors*.



Figure 3-11: FlowCell versions

> A: with V flange EN 1092-1/11 B: with Hygiene milk pipe screw connection DIN 11853-1

C: with welding pipe

The FlowCell consists of a sturdy stainless steel body. The microwave transmitter and receiver are firmly welded to the outside of the pipe. The entire product pipe is PTFE-coated and thus meets the special requirements for use in food for use in food.

There are no objects that extend into the pipe (e.g. a measuring sensor). Depending on the version, the FlowCell can be mounted with the V flange, milk pipe screw connection or welding pipe to the piping. For the versions with V flange, ASA adapter flanges are available as accessory.

The FlowCell has two HF ports to feed in and output the microwave signals. The input and output can be allocated as needed (M-Tx, M-Rx). The microwave signals transmit the product over the entire pipeline cross-section.

For all versions, the following accessories are available:

- 1. Pipe-mounted PT100 or Inline PT100
- 2. Conductivity measuring device
- 3. Sampling valve (combination with 1. and 2. possible)

Overview accessories (see also chapter 6.2):



Conductivity measuring device



Inline PT100

Sampling valve





Combination Conductivity measuring device with Sampling valve

Combination Inline PT100 with Sampling valve



3.5.3 Container Probe

The container probe is available in a version with and without flushing device (see Figure 3-12). The technical data are listed in *chapter 6.2 Technical Data Sensors*.





A: High-frequency connections B: Process connection, flanges of different sizes

Container probe type LB 5650 and type LB 5651

The container probe has been specially designed for concentration measurements in containers. Both measuring rods are immersed into the product. Microwaves are emitted from one end of the rod and received by the other end of the rod; they are emitted only towards the opposite end of the rod. This directional characteristic of the probe minimizes the interfering influence of metal parts in the vicinity of the probe and allows installation if only limited space is available. For example, the concentration of sugar strike can be measured continuously to find the suitable inoculation time.

The plastic caps of the measuring rods meet the special requirements for application in foodstuffs.

Two different probe types are available:

- The standard type is the container probe without flushing device.
- The flushing probe is used in processes where incrustations are likely to occur, for example, due to increased depositions. The flushing device prevents deposition on the microwave exit windows.



For the case depicted in Figure 3-13, the flow direction of the product being measured must be perpendicular to the drawing plane. This ensures a representative product between the measuring rods, provided the product is mixed thoroughly.



Figure 3-13: Container probe LB 5650

PT100

The container probe LB 5650 is the only one equipped with a PT100 connection and can be connected to the evaluation unit via a 4-core cable. The PT100 wiring diagram is described in *chapter* 4.6.1 Pin Configuration of the Connector Strip. To minimize the danger of incrustation in the immediate vicinity of the measuring rods, the container flush probe is not provided with a PT100.



Warning: Possible material damage! Do not open the cover screws on the front of the container probes, see Figure 3-14.



Figure 3-14: Front view container probe NOTICE

Container probe type LB 5651 with flushing device

The flushing probe LB 5661 has been designed for processes in which depositions, for example, due to incrustations are likely to occur on the probe.

The flushing probe has two flushing channels which keep the measuring rods free from incrustations; this ensures that the microwaves come into direct contact with the product being measured. All probe parts coming into contact with the product meet the special requirements for application in foodstuffs. Figure 3-15 shows the probe design.

Warning: Disturbed communication signal! Possible material damage!

When installing the probe with flushing, make sure that the HF connector is always dry.



Figure 3-15: Flushing probe LB 5651

2 x 3/8" Flushing connections, internal thread

nectoris, internal triead



The flushing slit width is the same for both probe rods and is depicted in Figure 3-16.



Figure 3-16: Rod head with flushing pipe

3.5.4 High-frequency Cable

High-frequency cables (HF cable) are used to transmit microwave signals.

HF cables change their conductivity (for microwaves) with temperature and would therefore produce measurement errors with varying ambient temperature. This error is compensated for by enabling the automatic cable compensation. The influences of the ambient temperature on the signal cable are compensated for by means of the reference cable. To this end, the sum of the reference cables has to match the length of the sum of the signal cables.

Two different HF cable types of different lengths are available:

Version 1: The so-called HF-cable quad: It consists of four single HF-cables of equal length, whose ends are terminated by one HF-connector (N-type). Available cable lengths: 2, 4, 6 and 10 m (see Figure 3-17).

The reference line of the HF cable on the probe side is shortcircuited using an N-connector (see Figure 3-18).









Version 2: It consists of a single HF-cable whose ends are terminated by an HF-connector (N-type). Available lengths: 2, 2.5, 3, 3.5 and 4 m (see Figure 3-19).



3.6 Assembly on a Pipeline

The evaluation unit is installed close to the FlowCell to keep the HF cable between evaluation unit and probe as short as possible. The shorter the cable connection, the better the stability of the measurement. The standard length is 2 m and the maximum length of the HF cable is 10 m. The HF cable must be at least 2 m long.

The FlowCell is integrated into the existing pipeline system or in a bypass. The orientation of the FlowCell can either be vertical or horizontal. To avoid possible sedimentary deposits, vertical installation in a riser is preferred (see Figure 3-20).

The installation should preferably be close to a sample sampling point to ensure representative sampling for calibration.

A representative temperature signal (current signal or PT100) should be connected to the evaluation unit for product temperature compensation.



Figure 3-20: Typical measurement arrangement on a pipeline

3.7 Assembly on a Container

The evaluation unit is installed in the immediate vicinity of the container probe to keep the HF-cable between evaluation unit and probe fairly short. The shorter the cable connection, the better the stability of the measurement. The standard length is 2 m and the maximum length of the HF cable is 10 m. The HF cable must be at least 2 m long.

The installation should preferably be close to a sampling point to ensure representative sampling for calibration. A representative temperature signal (current signal or PT100) should be connected to the evaluation unit for a possibly required product temperature compensation.

Our example in Figure 3-21 shows the measurement configuration on a process container. The probe is fixed to the container wall using a flange coupling so that both measuring rods are immersed into the product.



Figure 3-21: Example: Pipeline measurement configuration on a process container

Chapter 4. Getting Started

4.1 Transport

NOTICE	Warning: Possible material damage! System parts may get damaged during transportation
	Transport all components in their original packaging. Protect parts against shocks.
	In particular, the measuring rods of the container probes must be protected against mechanical shock!
	After unpacking, make sure all parts listed on the packing list have been delivered and show no sign of damage; if necessary, clean these parts.
	If you detect any damage, please notify the forwarder and the manufacturer immediately.
	The weight of the system components may exceed 25 kg, de- pending on the version. We recommend, therefore, that you

wear safety boots.
4.2 Commissioning the FlowCell

4.2.1 Installing the FlowCell

Da	anger of injury from leaking operating medium.
Th	e FlowCell must be mounted correctly.
In mi	particular, the correct tightening torque for the flange screws ust be observed.
Re	quirements for the hygienic installation of the FlowCell:
۶	The mounting position must guarantee self-draining proper- ties.
۶	The device has been developed for Cleaning in Place (CIP) applications and must not be dismantled for cleaning.
۶	Do not use cleaning equipment which are scraping or abrasive to avoid damage of product contact surfaces.
۶	Do not use aggressive cleaning agents or chemical which can affect the product contact surface.
۶	The customer is obligated to clean the FlowCell appropiately before it comes into contact with food.
	To meet the requirements for EHEDG certification, the sensor must be connected with process connections in accordance with the EHEDG position paper entitled "Easy Cleanable Pipe Couplings and Process Connections" (www.ehedg.org).
No	te the following points when installing the FlowCell:
•	The FlowCell is integrated at a suitable location in the pipeline system. Keep in mind that material sampling directly behind the FlowCell should be possible for calibration.
•	The FlowCell should be installed in a vertical riser, if possible. It must be ensured that material deposits cannot form on the pipe walls and no bubble formation occurs in the product.
	When aligned horizontally, the drainability of the pipeline has to be respected. To ensure this, the FlowCell must be mounted acc. to fig. 4-2.
\checkmark	There should be a straight pipe section of at least 300 mm and equal nominal width before and after the FlowCell to en- sure a fairly homogeneous flow profile and to rule out possibly occurring microwave reflections in the pipeline. Also, not fit- tings must be installed in these pipe sections.

- No gas inclusions should be present in the product. If gas bubbles cannot be ruled out, a pressure of at least 4 bar is required in the pipeline to minimize the influence of gas bubbles. Please observe the max. permissible operating pressure, see *chapter 6.2 Technical Data Sensors*.
- The high-frequency cables should preferably be connected to the FlowCell from below to prevent flowing water from getting to the connection sockets.
- The signal and reference cable should as far as possible follow the same path to make sure both cables are exposed to the same temperature and should not come into contact with hot pipelines. We recommend installing the HF cable through a single protection tube. If you are working with the HF cable quad, this function is taken over in good approximation by the corrugated tube.





4.2.2 Installing the Evaluation Unit

Note the following points when installing the evaluation unit:

- Position the evaluation unit depending on the length of the HF cable in the vicinity of the microwave probe.
- The evaluation unit has to be protected against vibrations. In some cases, it is advisable to set up the evaluation unit on a stand separated from the pipeline system.
- For instrument installation you should foresee a cutoff device to allow easy and quick disconnection of the device from the power supply.
- Provide an automatic separating device (line circuit breaker) that disconnects the unit from power within 0.03 seconds in case of failure. The separating device must be matched to the cable cross-section of the supply line, but at least it must be designed for 1 A continuous current.
- When installing the evaluation unit on a crystallizer, use a distance rail to minimize the thermal radiation and conduction. See Figure 4-3:
- For outdoor applications, the evaluation unit must be protected from direct sunlight and rain, for example by a suitable shelter.



Figure 4-3: View from above: Installation of the evaluation unit on a hot container wall

4.2.3 Connecting the HF Cable

The FlowCell is connected with the evaluation unit via the HF cables. Two different options are available:

Version 1:	1 x HF cable quad and 1 x N-connector
Version 2:	2 x HF cables (as signal cable)
	1 x HF cable (as reference cable)

The decision for a certain cable version is taken by Berthold Technologies in the planning stage. It is subject to the application and the desired distance between evaluation unit and FlowCell.

Prerequisite for a correct measurement is the correct installation of cables! Please keep in mind:

ттр Make sure the cable does not get into contact with hot pipes over the entire length (corrugated tube and single cable section after splitting), e.g. direct contact with the device wall (not insulated). This will ensure that all individual cables are exposed to the same ambient conditions and that the compensation of the cable drift works properly. Never bend the HF cable! The bending radius should not be less NOTICE than 100 mm. Fix the HF cable with cable binders or other suitable means, so that the cable cannot slip anymore! During assembly, ensure that the end of the corrugated tube is NOTICE bent down on the probe side. By doing so, the ingress of fluids into the corrugated tube is prevented. Version 1 The HF cable quad and the HF connectors on the evaluation unit are labeled. Connect the FlowCell to the evaluation unit as shown in Figure 4-4; only cables with identical labeling can be combined. The two connections on the FlowCell are not labeled, the assignment of both cable connectors M-Tx and M-Rx is arbitrary. Connect the cable connectors R-Tx and R-Mx with the Nconnector (short-circuited). Version 2 Connect the FlowCell to the evaluation unit as shown in Figure 4-5; make sure that the reference cable (ring line) is connected to R-Tx and R-Mx. The reference cable must be as long as the sum of both signal cables.





Hand tighten all screwed connections of the HF cable (2 Nm = 0.2 kg/m)! Before tightening, carefully screw on the cable by hand. **Caution! Threaded joint jams easily.**

Check occasionally if the screwed connection is still properly tightened. If the installation is exposed to vibrations, the screwed connection may come loose and this may result in inaccurate measurements or corrosion of the connections.

As long as the cables are not connected, the coaxial sockets have to be covered with plastic caps and the cable connectors have to be protected against moisture and dirt.

4.3 Commissioning the Container Probe

4.3.1 Installing the Container Probe

Danger of injury from leaking operating medium.
The Container probe must be mounted correctly.
in particular, the correct tightening torque for the flange screws must be observed.
or installation of the probes, please keep in mind:
Select the installation site such that good mixing and a ho- mogeneous product are ensured and no bubbles are present in the probe. A sample tap should be available in the direct vicinity to allow representative sampling.
The probe must be flanged to the container so that the mate- rial to be measured flows between the two measuring rods. That means the fork (both measuring rods) must be installed perpendicular to the material flow.
The distance between the measuring rod tips and any metal- ized walls (heating element, stirrer, container wall) should be at least 60 mm .

The following mounting hole dimensions in the fitting flange are required for installation of the probe:

Flange	Minimum mounting hole size \varnothing (mm)
DN 65 / PN 6	102 ± 0.3
others	104 ± 0.5

- For further installation dimensions, such as the required installation depth, please refer to *chapter 8.6* and *8.7 Assembly Sheets*.
- Use the appropriate flat gasket (standard accessory) to compensate for minor surface tolerances in the fitting flange.

Installation on process containers

Figure 3-26 shows the position of the container probe on the container. This position is also valid for the container flush probe. Follow the instructions in *chapter 3.9 Assembly on a Container.*

The assembly sheets in *chapters 8.6 and 8.7* include all the information required for installation.

Dismounting in the event of a stuck flange

Depending on the process, the probe flange may stick to the process flange. In such case, during the dismounting process, remove the two eyebolts first. Then insert two longer screws (M8) into their bore holes up to the process flange. If you continue to rotate the screws, you will push the probe away from the process flange.

Installation in pipelines

Container probes can be installed in pipelines with a nominal width \geq 200 mm using an adapter flange. Note the position and orientation of the container probe, see the technical drawings in *chapter 8.4.7 and 8.5.6 Installation Situation in Pipelines.*

Connecting the flushing pipes

The container probe with flushing device has two flushing connections with G3/8'' female thread (DIN ISO 228-1). The flushing connections should be sealed only at this thread. Sealing the probe cover, for example with silicone, is not permitted.

If the flushing pipe is not connected, the container probe must be closed with a blind plug.

Flushing parameters (only for flushing probe)

The degree of deposition or incrustation is essential for the flush settings, i.e. flush interval and flush duration. The flushing parameters must be matched to the product and the process.

The following product and process independent flushing parameters must be strictly adhered to:

Flush solution	Water, condensate
Temperature of	Maximum 120°C
flush solution	
Pressure	≥3 bar, max. 8 bar
Fittings	2 x G3/8" female thread (DIN ISO 228-1)
Supply pipe	≥ 1/2 inch

In general: The flushing connections can be flushed simultaneously or sequentially. The flushing parameters are valid for each flushing connection.



The flush duration has to take into account possible inertias of the system, e.g. valve openings. The flush supply pipes must be well insulated to prevent that the flush solution is initially colder.

Amounts of water With a 5 bar flushing pressure, the amount of water per flushing connector is approximately 0.8 liters per second.

4.3.2 Setting Up the Evaluation Unit

Set up the evaluation unit as described in *chapter 4.2.2*.

4.3.3 Connecting the HF Cable

Connect the container probe to the evaluation unit via the HF cable quad.

Prerequisite for a correct measurement is the correct installation of cables! Please keep in mind:



Make sure the cable does not get into contact with hot pipes over the entire length (corrugated tube and single cable section after splitting), e.g. direct contact with the device wall (not insulated). This will ensure that all individual cables are exposed to the same ambient conditions and that the compensation of the cable drift works properly.



Never bend the HF cable! The bending radius should not be less than 100 mm. Fix the HF cable with cable binders or other suitable means, so that the cable cannot slip anymore!

The HF cables and the HF connections on the evaluation unit and on the probe are labeled. Connect the probe to the evaluation unit as shown in Figure 4-6; only cables with identical labeling can be combined.



Evaluation unit, underside



Hand tighten all screwed connections of the HF cable (2 Nm = 0.2 kg/m)! Before tightening, carefully screw on the cable by hand. **Caution! Threaded joint jams easily.**

Check occasionally if the screwed connection is still properly tightened. If the installation is exposed to vibrations, the screwed connection may come loose and this may result in inaccurate measurements or corrosion of the connections.

As long as the cables are not connected, the coaxial sockets have to be covered immediately with plastic caps and the cable connectors have to be protected by suitable provisions against moisture and dirt.

4.5 Connecting the Evaluation Unit

	Electrical hazards:
	Disconnect power to rule out any contact with live parts during installation and servicing. Turn off power before opening the instrument. NEVER work on open and live instruments.
NOTICE	Caution! Potential hazards, material damage! Device type: LB 566-12 MicroPolar ++ (ID no. 51833-02)
	When connecting the 24 V DC power supply, the + and – poles must be connected correctly. There is no reverse polarity pro-tection!
	The line cross-section the for power supply must be at least 1.0 mm^2 .
	Connect all desired input and output signals to the terminal strip as shown on the following pages. Use the M feed-through

- strip as shown on the following pages. Use the M feed-through to keep the degree of protection.
- Check if the voltage indicated on the type plate matches your local supply voltage.
- Connect the deenergized power cable to the terminals 3(L1), 2(N) and 1(PE).
- Verify that the test switch (power interruption) is in position "ON" (see Figure 5-1).
- > Close the instrument housing and turn on the power supply.

4.5.1 Pin Configuration of the Connector Strip

The connector strip of the evaluation unit includes the following terminals:



Figure 4-7: LB 566 wiring diagram

Power supply: Terminals 3 (L1, +), 2 (N, -) and 1 (PE,)

For MicroPolar, depending on device type, see name plate on the housing outer wall.

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

For MicroPolar ++, depending on device type, see name plate on the housing outer wall.

- 1.) 100...240 V AC, 45...65 Hz
- 2.) 24 V DC: 18...36 V, no reverse polarity protection

Current input no. 1 (terminals 20+ and 8-), insulated Current input no. 2 (terminals 22+ and 10-), not insulated

Input as 0/4 - 20 mA signal. For example, for temperature compensation or reference signal recording.

Current output no. 1 (terminals 27+ and 15-), insulated

Output as 4 - 20 mA signal. Output options: Concentrations (1 / 2), current input signals (1 / 2), PT100 signal, mass flow

Current output no. 2 (terminals 19+ and 7-), insulated Output as 0/4 - 20 mA signal. Output options same as for current input no. 1.

PT100 (terminals 23+ and 11-)

Connection for temperature measurement.

When commissioning the container probe, connect the 4-core cable of the PT100 to the connector strip of the evaluation (see fig. 4-8). Connect the other end of the cable with the corresponding plug on the container probe (see fig. 4-6).

Figure 4-8: PT100 connection container probe



() Terminal no.



In the delivered state, the PT100 cable has a length of 10 meters. We recommend shortening the cable to the required length.

Combine the two core pairs in accordance with figure 4-8.

After connection of the PT100 cable, the included ferrite core must be wound around the cable. The applicable position is within the housing, as close as possible to the cable bushing.

Figure 4-9: Ferrite core



Digital input 1: DI1 (terminals 24+ and 12-)

Only for potential-free contacts! Configuration options:

- No function
- Measurement: Start (closed) and stop (open)

Digital input 2: DI2 (terminals 25+ and 13-)

Only for potential-free contacts! Configuration options:

- No function
- > Average value: hold (closed) and continue averaging (open)
- Product selection: product 1 (open) and product 2 (closed)

Digital input 3: DI3 (terminals 26+ and 14-)

Only for potential-free contacts! Configuration options:

- No function
- Start sampling, open: no action, closed: unique measurement starts
- Product selection

Relay 1: (Terminals 4, 5 and 6) and Relay 2: (Terminals 16, 17 and 18)

Changeover contacts (SPDT), insulated, configuration option:

- No function
- Error message
- Stop measurement
- Limit value min. and max.
- No product
- Below load limit

RS485 interface (terminals 21 (RS1) and 9 (RS2)) and RS232 interface (on instrument underside)

Serial data interface for output of the live data (all readings for every sweep (measuring cycle), the protocol and data logs. Data format: Data transfer rate 38400 baud, 8 data bits, 1 stop bit, no parity, no handshake

4.5.2 Digital Outputs, Relay

The status of the measurement is output via two relays:

- > Error
- Alarm (alarm min. and max.)
- No product

On the Plausibility menu, you can enter a min. attenuation for pause detection (e.g. for process pause, no product present); if this value is not reached, "No product" is signaled via a relay and the current output drops to 0 or 4 mA.

- Measurement stopped
- Below load

The respective switching state is also signaled via LEDs on the front panel (LEDs signal 1 and 2).

Relay no.	Error, alarm, no product, measurement stopped, Below load, currentless condition	Normal
1	4 0	4 0 5 0 com 6 0
2	16 O	16 O 17 O com 18 O

The relays with changeover contacts can either be operated as make contact, terminals 4 & 5 (open at error, alarm ...) or as break contact, terminals 5 & 6 (closed at error, alarm ...).



Chapter 5. Service instructions

5.1 General Information

A malfunction of the measuring system is not always due to a defect in the instrument. Often the error is caused by incorrect operation, improper installation or irregularities in the product being measured. If a malfunction occurs, anyway, the measuring system helps you to identify and eliminate errors by displaying error messages on the display, indicating operator errors and defects of the electronics.

Defective modules of the evaluation unit cannot usually be repaired but must be replaced. The microwave module is firmly bolted to a screening hood and must not be opened.

For device disposal, please contact the Berthold Service and apply for a recycling passport.

5.2 Wear Parts

The evaluation unit does not include any parts that are subject to wear or components that require special maintenance.

The plastic caps of the measuring rods of the container probes and the PEEK Microwave windows of the FlowCell may be subject to abrasion over the course of time. A low to medium abrasive influences the measurement only very little and can be compensated for by a calibration. Therefore, check the wear parts in intervals of about 2 years. In case of heavy wear, the plastic caps of the measuring rods of the container probe and the PEEK Microwave windows of the FlowCell can be replaced.

Replacing the plastic caps of the measuring rods

- **1** Unscrew the two plastic caps from the measuring rods and remove the four sealing rings (see figure 5-1, yellow arrow).
- 2 Clean the four grooves of the sealing rings with a lint-free (if necessary wet) cloth.
- **3** For each measuring rod, insert two new sealing rings into the grooves.
- 4 Screw on the two new plastic caps.





Figure 5-1: PT100 connection container probe

Caps set for container probe	
IdNr. 66049-S	2 pieces of PEEK plastic caps with 4 sealing rings

Replacing the Microwave windows of the FlowCell

- 1 Open the fixing clamp (see fig. 5-2, item 1).
- **2** Remove the antenna (see fig. 5-2, item 2), the microwave window (see fig. 5-2, item 3) and the sealing O- rings.
- 3 Attach the new microwave window, the new sealing O-rings and the antenna to the FlowCell with the fixing clamp according to fig. 5-3. HNBR o-ring must be used instead of EPDM oring once a fat/grease content of 8 % is exceeded.



Figure 5-2: Antenna and Microwave window







5.3 Instrument Cleaning

Clean all system components exclusively with a damp cloth with no chemical cleaning agent. Parts coming into contact with the product (during regular operation) can be cleaned with warm water, taking into account the temperature limits, see *chapter 6.2 Technical Data Sensors*.

5.4 Battery

If the measuring system LB 566 is a long time without power supply (power failure or disconnected from the mains supply), the system clock is powered by the lithium battery on the motherboard.

If the battery voltage is no longer sufficient, the error message CODE 14 "Battery voltage" appears after a restart of the evaluation unit. After acknowledging the error message, the unit continues to work properly; however, the date and time should be checked and corrected, if necessary. Measurement data that are output via a serial interface can be fatally damaged by incorrect date and time information. We recommend changing the batteries immediately.

The service life of the battery, even under continuous stress, is approximately 8 years. Replacement of batteries must be carried out in a device disconnected from mains.

Battery type: 3 Volt lithium cell (button cell), type CR2032

5.5 Fuse Replacement

The mains fuses of the LB 566 are located in the wall housing. Replace the fuses only if the instrument is disconnected from mains.

Use only fuses with the correct rating, see chapter 6.1



Spare fuses must match the rating specified by the device manufacturer. Short-circuiting or manipulation is not permitted.



Figure 5-4: View with open housing wall MicroPolar

Chapter 6. Technical Data

General specifications	
Method	Microwave transmission measurement
Operating fre- quency	2.4 - 2.5 GHz (ISM band), depending on local regulations
Transmission power	MicroPolar: < 0.1 mW (< - 10 dB) MicroPolar ++: < 10 mW (< 10 dBm) All coaxial line power
Application	Concentration / moisture measurement in containers and pipelines

6.1 Technical Data Evaluation Unit

Evaluation unit		
Housing	Wall housing made of stainless steel, see dimensional drawing in <i>chapter 8</i>	
	MicroPolar: HxWxD: 300 x 323 x 140 mm	
	MicroPolar++: HxWxD: 400 x 338 x 170 mm	
Protection type	IP 65	
Weight	MicroPolar: approx. 6.5 kg	
	MicroPolar ++: approx. 8.0 kg	
Ambient conditions during operation	-20 +60 °C (253 K333 K), no condensation	
	Relative humidity: max. 80 %	
	Altitude: max. 2000 m	
Ambient conditions during storage	-20 +70 °C (253 K343 K), no condensation	
	Relative humidity: max. 80 %	
Achievable accuracy	\leq 0.2 weight % (standard deviation) depending on product and sensor	
Display	Dot matrix LC display, 114 mm x 64 mm, 240 x 128 pixels, with back-lighting, automatic contrast setting	
Keyboard	Freely accessible foil keypad, light-stable and weatherproof: alphanumeric key- board and 4 softkeys (software-assigned buttons)	

Power supply	 For Micro Polar, depending on device type: 1.) 100 240 V AC, 45 65 Hz 2.) 24 V DC: 1836 V; 24 V AC: -20%, +5%, 40 440 Hz For MicroPolar++, depending on device type: 1.) 100240 V AC, 4565 Hz 2.) 24 V DC: 1836 V, no reverse polarity protection
Power consumption	For Micro Polar: max. 30 VA (AC/DC), depending on configuration For MicroPolar ++: max. (48/60) VA (AC/DC), depending on configuration
Fuses	For Micro Polar: 2 x 2.0 A / slow-blow For MicroPolar ++: 2 x 2.0 A / slow-blow at 100 240 V AC or 2 x 6.3 A / slow-blow at 24 V DC
Battery type	3 V Lithium button cell, type CR2032
Measured value	e.g. concentration, dry matter content
Inputs and outputs	
Cable cross-section	min. 1.0 mm ² (mains supply)
Cable feed-through	2 x M20x1.5 for cable 514 mm (depending on application) 4 x M16x1.5 for cable 58 mm (depending on application)
Sensor connection	Inputs and outputs for signal and reference channel, 50 Ω N-socket
HF cable	Cable lengths: 2, 4, 6 and 10 m; 50 Ω ; both sides with 4 N connectors
Current input	2 x current input $0/4 \dots 20$ mA, ohmic resistance 50 Ω , 1x insulated, 1x instrument ground e.g. for temperature compensation
Current output	Current output 1: 420 mA, ohmic resistance max. 800 Ω , insulated Current output 2: 0/420 mA, ohmic resistance max. 800 Ω , insulated e.g. for measured value or temperature out-
PT100 connection	Measuring range: -50 +200 °C (223 473 K); measurement tolerance: < 0.4 °C

Digital input	3 x digital inputs (DI13), for floating con- nectors (do not connect to a power supply).
	<u>Configuration options:</u> DI1: none, measurement start/stop DI2: none, measurement hold, product selec- tion DI3: none, sampling, product selection
	 <u>Function description:</u> Measurement (Start/Stop), <u>open:</u> Measurement stopped, <u>closed:</u> Measurement started and/or measurement running
	 Hold measurement, <u>open:</u> measurement running, <u>closed:</u> measurement stopped, i.e. average values and current output are held
	3. Product selection via a DI: <u>open:</u> Product 1 (P1), <u>closed:</u> P2
	Product selection via two DI's: <u>DI2 & DI3 open</u> : P1 <u>DI2 closed & DI3 open</u> : P2 <u>DI2 open & DI3 closed</u> : P3 <u>DI2 & DI3 closed</u> : P4
	single measurement starts
Relay outputs	2 x relays (SPDT), insulated <u>Configuration options:</u> - Collective failure message - Stop measurement - Limit value (min. and max.) - No product - Low load
	Load capacity: AC: max. 400VA DC: max. 90W AC/DC: max. 250V, max. 2A, non-inductive ≥ 150V: voltage must be grounded
	The cable used at the relay output must corre- spond to a mains cable.
	Restrictions at 24 V AC/DC (DC: 1836 V; AC: 24 V +5 %, -20 %) mains supply, if the ground conductor is not connected to terminal <u>1 (PE):</u> AC: max. 50 V
	DC: max. 70 V
Serial interfaces	RS232 on the bottom side RS485 via terminal strip Data format: 38400 Bd, no handshake, 8 data bits 1 stop bit no parity

6.2 Technical Data Sensors

FlowCell	
Application	Microwave FlowCell with various nominal diame- ters and flanges for concentration measurement on pipelines
Material	Inline housing made of stainless steel 1.4404 pol- ished (AISI 316L)
	Mikrowave windows made of PEEK
	Product touching sealing made of EPDM
Process coupling	 Three versions: 1. Hygiene milk pipe screw connection DIN 11853-1 EHEDG certified 2. Flange according to EN 1092-1/11 (V flange) FDA-approved materials 3. Welding pipe Optional adapter for the V flange version with ASA flange
Process pressure	up to 16 bar (relative)
Temperature range	Product temperature: 10130 °C (283403 K), temporarily up to 140 Ambient temperature: -2060 °C (253333 K) Storage temperature: 1080 °C (283353 K)
Connections	2 x HF connections: N female, 50 Ω for HF cable with max. 10 m length
Versions	Nominal pipe widths from 50 150 mm
Dimensions	See dimensional drawings in chapter 8.

Overview FlowCells with V flange

Designation	ID no.	Nominal width [mm]	Flange	Pressure [bar]
LB 5660-102-00x	66744-001	50	DN 50 / PN 16	
LB 5660-202-00x	66744-002	65	DN 65 / PN 16	
LB 5660-302-00x	66744-003	80	DN 80 / PN 16	
LB 5660-402-00x	66744-004	100	DN 100 / PN 16	
LB 5660-502-00x	66744-005	125	DN 125 / PN 16	
LB 5660-602-00x	66744-006	150	DN 150 / PN 16	
LB 5660-402-200 FlowCell Vfl. FDA Immersion cap	66744-031	100	DN 100	16
LB 5660-502-200 FlowCell Vfl. FDA Immersion cap	66744-032	125	DN 125	
LB 5660-602-200 FlowCell Vfl. FDA Immersion cap	66744-033	150	DN 150	

Overview FlowCells with Hygiene milk pipe screw connection

Designation	ID no.	Nominal width [mm]	Pressure [bar]
LB 5660-112-00x	66744-013	50	
LB 5660-212-00x	66744-014	65	
LB 5660-312-00x	66744-015	80	
LB 5660-412-00x	66744-016	100	
LB 5660-512-00x	66744-017	125	
LB 5660-612-00x	66744-018	150	
LB 5660-412-200 G-BS-M Immersion cap (hygienic)	66744-034	100	16
LB 5660-512-200 G-BS-M Immersion cap (hygienic)	66744-035	125	
LB 5660-612-200 G-BS-M Immersion cap (hygienic)	66744-036	150	

	-		
Designation	ID no.	Nominal width [mm]	Pressure [bar]
LB 5660-132-00X	66744-025	50	
LB 5660-232-00X	66744-026	65	
LB 5660-332-00X	66744-027	80	
LB 5660-432-00X	66744-028	100	
LB 5660-532-00X	66744-029	125	
LB 5660-632-00X	66744-030	150	
LB 5660-432-200 Immersion cap (hy-	66744-037	100	16
		4.25	
LB 5660-532-200	66744-038	125	
Immersion cap (hy- gienic)			
LB 5660-532-200	66744-039	150	
Immersion cap (hy- gienic)			

Overview FlowCells with welding pipe

Overwiew Microwave windows set

The microwave windows of the FlowCell are available in a reinforced glass fibres design for applications with a high grade of abrasion. This design is not approved for the food sector.

ID no.	Description
66624-S	2 pieces of PEEK Microwave windows with 6 seal- ing O-rings
66625-S	2 pieces of PEEK Microwave windows GF30 glass fibre reinforced with 6 sealing O-rings
75514-S	2 pieces of PEEK EHEDG Microwave windows with 6 sealing O-rings

NOTICE

Probes with PEEK EHEDG Microwave windows can only be used from a nominal width of DN100.

Overview ASA flange adapter

Designation	ID no.	
ASA flange adapter set for Flow Cell 50	62324	
ASA flange adapter set for Flow Cell 65	62319	
ASA flange adapter set for Flow Cell 80	62328	
ASA flange adapter set for Flow Cell 100	62331	
ASA flange adapter set for Flow Cell 150	62335	
The kit consists of two adapters, screws and two seals.		

Overview Inline housing, FDA

for temperature or conductivity sensors or sampling valve

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

Overview surface temperature sensor DN 50

Self-adhesive PT100 temperature sensor with fixing material Connection cable 10 m, 4-wire (loose ends) Temperature range: -50 - +200 °C

Designation	IdNr.
Self-adhesive temperature sensor for DN 50	66655
Self-adhesive temperature sensor for DN 65	66656
Self-adhesive temperature sensor for DN 80	66657
Self-adhesive temperature sensor for DN 100	66658
Self-adhesive temperature sensor for DN 125	66659
Self-adhesive temperature sensor for DN 150	66660

Overview sensors

Designation	IdNr.
Conductivity sensor hygienic, Clamp-flange	66693
Inductive conductivity measuring device for liquid media in hygienic applications Measurement range: 0-999 mS/cm Process connection: Clamp-flange Process pressure: Max. 16 bar Power supply: 18-36 V DC, max. 190 mA Output: 4-20 mA	
Temperature sensor EHEDG, Clamp-flange	66694
PT100 temperature sensor for hygienic applica- tions Measurement range: -50 - +250 °C Length 20 mm, diameter 4 mm Process connection: Clamp-flange With connection cable 10 m (loose ends)	

Overview Sampling valve and accessories

Designation	IdNr.
Sampling valve aseptic, Clamp-flange	66738
Aseptic Inline Sampling valve Stainless steel 1.4404 (AISI 316L), bellows PTFE Discharge port S-DN 10 without flushing connection Process connection: Clamp-flange	
Clamp-blind flange	66737
Clamp-blind flange for Inline housing DN 50- 150 Stainless steel 1.4306 (AISI 304L)	
Clamp coupling	66736
1 piece Clamp coupling for Inline housing DN 50-150 Stainless steel 1.4306 (AISI 304L)	

Container probes with and without flushing device for concentration measurement in process containers and pipelines with nominal width ≥ 200 mm.		
Plastic caps, stainless steel PT100 connection cable: Silicon / Teflon		
Flange according to DIN EN 1092 type 05 DN65 / PN6, DN 80, 100, 150 / PN16; ASA flange 2.5" / 150 PSI (More on request)		
Up to 16 bar, depending on model		
Product temperature: 10120 °C (283393 K) Ambient temperature: -2060 °C (253333 K) Storage temperature: 1080 °C (283353 K)		
4 x HF connections: N female, 50 Ω for HF cable with max. 10 m length		
See dimensional drawings in chapter 8.		
washer		
Klingersil C-4400		
3 mm		
Caps set for container probe		
2 pieces of PEEK plastic caps with 4 sealing rings		

Overview container probes and sealing washers

Designation	ID no.	Flange	Pres- sure [bar]	ID no. Seals
LB 5650-01	65464-01	DN 65 / PN6	6	32175
LB 5650-02	65464-02	DN 80 / PN16	16	33717
LB 5650-03	65464-03	DN 100 / PN16	16	46661
LB 5650-04	65464-04	DN 150 / PN16	16	46664
LB 5650-05	65464-05	ASA 2.5" / 150 PSI	16	46665
LB 5650-09	65464-09	ASA 3" / 150 PSI	16	50659
LB 5651-01	65937-01	DN 65 / PN6	6	32175
LB 5651-02	65937-02	DN 80 / PN16	16	33717
LB 5651-03	65937-03	DN 100 / PN16	16	46661
LB 5651-04	65937-04	DN 150 / PN16	16	46664
LB 5651-05	65937-05	ASA 2.5" / 150 PSI	16	46665

6.3 Technical Data HF Cable

HF cable Quad		
Material	Corrugated tube: Polyamide (PA6) Cable sheath: Polyethylene (PE)	
Protection type	IP 66	
Temperature	In operation: -30 +70 °C When installing: -20 +70 °C	

Cable length [m]	ID no.
2	43431
4	43432
6	43433
8	43434
10	43435

HF-Kabel Quad, hygienic			
Material	Corrugated tube: Polyamide (PA6) Cable sheath: Polyethylene (PE)		
Protection type	IP 66		
Temperature	In operation: -30 +70 °C When installing: -20 +70 °C		

Cable length [m]	ID no.
2	67048
4	67049

HF cable Quad (solid cable)			
Material	Cable sheath: Polyethylene (PE)		
Protection type	IP 68 when unscrewed		
Temperature	In operation: -40 +85 °C When installing: -40 +85 °C		
Attenuation coefficient	about 0.3 dB/m		



Cable length [m]	ID no.
2.0	11476
2.5	11477
3.0	11478
3.5	11479
4.0	11480

6.4 Format of Serial Data Output RS232 and RS485

Headline

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

Following lines

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

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

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

13 14 15 16 17 18

Column no.	Description	Format	
1	Date and time	DD.MM.YY·HH:MM:SS	
2	State	4 digits, HEX	
3	Status: Information about the quality of the last measurement	0 : Measurement OK < 0 : Error	
4	Product synchronization	5: not active 1: still asynchronous 0: all values synchronous -1: Error -2: Time too short for syn. -3: Speed outside range	
5	Product number	X (1 to 4)	
6	Attenuation [dB]	X.XX	
7	Phase [°/GHz]	X.XX	
8	Dispersion of the phase regression	X.XX	
9	Device temperature [temperature unit]	X.X	
10	Current input 1 [unit of current input]	X.X	
11	Current input 2 [unit of current input]	X.X	
12	PT100 temperature [temperature unit]	X.X	
	[] with selection of the unit g/cm ³		
13	Concentration 1 live	X.XX [X.XXXX]	
14	Concentration 1 averaged	X.XX [X.XXXX]	
15	Concentration 2 live	X.XX [X.XXXX]	
16	Concentration 2 averaged	X.XX [X.XXXX]	
17	Mass flow for concentration 1	X.XXX	
18	Mass flow for concentration 2	X.XXX	

Special characters

" \rightarrow " Tabulation"¶" Carriage return + Line feed

"·" Blank character
Chapter 7. Certificates

7.1 EC Declaration of Conformity



BERTHOLD TECHNOLOGIES GmbH & Co.KG

Caimbacher Str. 22 75323 Bad Wildbad, Germany

Phone +49 7081 177-0 Fax +49 7081 177-100 mfo@BertboldTech.com www.BertholdTech.com

EC – Declaration of Conformity

We herewith confirm that the constraction of the following indicated products / systems / units is brought into circulation to comply with the relevant EC regulations.

This declaration is declared void should alterations or unintended use take place without our authorisation,

Title: Concentration-Measuring System Micro-Polar 1

Type: LB 566-XX

Relevant EC regulations:

89/336/EWG (electromagnetic compatibility) reviewed: 91/263/EWG, 92/31/EWG, 93/68/EWG, 93/97/EWG

73/23/EWG (low voltage guidelines) reviewed: 93/68/EWG

The following norms were considered for the assessment of the pruducts:

EN 55011:1998 + A1:1999 + A2:2002 EN 61010-1:2002-08 EN 61006-6-2:2001 EN 61000-4-2:1995 + A1:1998 + A2:2001 EN 61000-4-3:2002 + A1:2002 EN 61000-4-4:1995 + A1:2001 + A2:2001 EN 61000-4-5:1995 + A1:2001 EN 61000-4-6:1996 + A1:2001

This declaration is issued by the manufacturer:

BERTHOLD TECHNOLOGIES GmbH & Co. KG P.O. Box 100163 D-75312 Bad Wildbad / Germany

by

Dr. J. Briggmann Development Manager Process Control

Bad Wildbad, 04.05.2004

H. Tetzlaff specialitat CE Certification Illerer T. Adaff

Registergesicht / Court of Registration Calw HRA 992 Persönlich baltende Gesellschafterin / Fully kable Associates BERTHOLD TECHNOLOGIES Verwaltungs-GrebH Registergeschi / Court of Registration Calw HRB 1520 Geschaftsführung / Managemens Hans 1. Oberhofer (Vers./CED), Dr. Wilfried Reuter Beinatsvorstzenkler / Chalman of the Boind Dr. Frtz Berthold USL-Johr, VAT Reg. No DEB33050511 Deutsche Steuermernmer / German 74k Ng Persähler Bank, 75105 Pforzheim Konto/account no.6 511 120 (BL2 666 800 13)/ SWIFT-BIC DRES DE FF 665 18AN: DE05 6668 9013 0651 1120 00 SperkassePf-CW 75323 8ad Wilfbad Konto/account no.6 351 05 (SL2 666 500 13)/ SWIFT-BIC DZH50E66 18AN: DE05 6005 10/03 0008 045005





BERTHOLD TECHNOLOGIES GmbH & Co.KG

Calmbacher Str. 22 75323 Bad Wildbad, Germany

Phone +49 7081 177-0 Fax +49 7081 177-100 mfo@Berthold.com www.Berthold.com

EC-Declaration of Conformity

We herewith conform that the construction of the following indicated products / systems / units is brought into circulation to comply with the relevant EC regulations.

This declaration is declared vold should alteration or unintended use take place without our authorisation.

Description:

Concentration-Measuring Systems Micro-Polar ++

Туре:	LB 566-XX			
	EC-Regulation	and Reviews	considered Norms	
EMC	2004/108/EG		EN 55011	1998
				+A1:1999
	1.0			+A2:2002
			EN 61326-1	2006-05
			EN 61000-4-2	1995
	· .		에는 물고 가장 가지 않는 것이 같이 많이 봐.	+A1:1998
	· .			+A2:2001
:			EN 61000-4-3	2006-12
	· .		EN 61000-4-4	2004
			EN 61000-4-5	1995
				+A1:2001
	i an thi a		EN 61000-4-6	1996
	부모가 같아 가지?			+A1:2001
		•	EN 61000-4-11	1994-08
	14 - 14 - 14 - 14 - 14 - 14 - 14 - 14 -	·		+A1:2001-02
			Namur NE21	¹¹ 2004
Low Voltage Directive	a 73/23/EWG	93/68/EWG	EN 61010 Part 1	2002-08
R&TTE	1999/5/EC		ETSI EN 300 440-1	2007-08
			ETSI EN 300 440-2	2007-08

This declaration is issued by the manufacturer BERTHOLD TECHNOLOGIES GmbH & Co. KG Calmbacher Str. 22 75323 Bad Wildbad, Germany

updated by Δ

Dr. Wilfried Reuter **Technical** Director Bad Wildbad, 06.06.2008

 Registergericht / Court of Registration
 Stuttgart H8A 330091

 Personitch haltende Gesellschaftenin / Fully liable Associates
 BERTHOLD TECHROLOGIES Verwaßungs-GmbH

 Registergericht / Court of Registration
 Stuttgart H8A 330091

 Bersonitch haltende Gesellschaftenin / Fully liable Associates
 BERTHOLD TECHROLOGIES Verwaßungs-GmbH

 Registergericht / Court of Registration
 Stuttgart H8A 330091

 Geschaftsfahrung / Management
 Hans J. Oberholer, Dr. Wiltried Reuter

 Beizetsvorsitzender / Charman of the Boar
 Dis Fritz Berthuld

 USt-1d-Nr, / VAT Reg. No.
 DE813050511

 Deutsche Steuernummer / Germän Tax No.
 1509468690

 WEE-Reg. No.
 DE99468690

 Drewiner Bank
 75105 Pforbleim
 Kontry/Account No. 6 511 120 (BLZ 666 800 13) /SWIFT-BIC DRES DEFF 666 TBAN: DE05 6668 0013 3651 1123 30

 Spärkasse 9F-CW 75323 Bad Wildbad Kontry/Account No. 60 450 03 (BLZ 566 500 85) /SWIFT-BIC PZHSDE66
 TBAN: DE37 6665 0085 0008 0450 03



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

7.2 Frequency License

CETECOM ICT Services GmbH

EC Identification number 0682

authorized by the German Gos crument



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

CERTIFICATE EXPERT OPINION

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

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

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



The scope of this evaluation relates to the submitted documents only γ . The exclusions only valid in can anchor with the following principly of annexes.

1

Number of annexes:

Suarbrücken, 24.06.2004 Place, Date of Issue

Signed by Ernst Hussinger Notified Body



CETECOM ICT Services Gmb11. Unterturkheimer Straße 6-10. D-66117 Starbrucken, Germany http://www.cetecom.de

÷



CETECOM ICT Services GmbH



CERTIFICATE OF CONFORMITY

Registration-No.:	E814059R-CC	Number of annexes:
Certificate Bolder:	Berthold Technologies GmbH & Co KG Calmbacher Strasse 22	
	D-75323 Bad Wildbad	
Product Designation.	LB 465xx, LB 466xx, LB 565xx, LB 56	6 6- 33
Product Description:	Short Range Devices (humidity sensor)	
Product Manufacturer:	Berthold Technologies GmbH & Co KG Ceimbacher Strasse 22	
	D-75323 Bad Wildbad	

Specifications and fest reports:

S	pecification	Test re	port no. & date	Name of test laboratory	Notes
E	N 55011:1998+A1:1999 (c	(lass A) 2003-73	31-1182-REN dated Sept. 9, 2003	ELMAC GmbH	conform
E	N 61000-6-2:2001				
E	N 300 440-1 V1.3.1 (2001	-09) 2-3389-	01-01/03 dated May 14, 2004	CETECOM ICT	conform
E	N 300 440-2 V1.1.1 (2001	-09)			L

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

CETECOM ICT Services is authorized to act no Notified Body in accordance with the R&TTE Directive 1999/5/EC of 09. March 1999

Saarbräcken, 24.06.2004 Place, Date offissie



Signed by Ernst Hüssinger Notified Body

CETECOM ICT Services (imbH, Unionid), heimer Stieße 6-10, U-66117 Saarbrücken, Germany

CETECOM ICT Services GmbH



CERTIFICATE OF CONFORMITY

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

Specifications and test reports:

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

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

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

	<i>///</i> ·	AND NOTIFICO B
Saarbrücken, 2008-12-05	CAH &	19. De
Place, Date of Issue	Signed by Ernst Hussinger Notified Body	OCHE BENAMOTE

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

CETECOM ICT Services GmbH

EC Identification Number 0682

authorized by the German Government



BNetzA-bS-02/51-52

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

EXPERT OPINION

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

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

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

The scope of this evaluation relates to the submitted documents only. The certificate is only valid in conjunction with the following number of apne:

1

Number of annexes:



<€ 0682

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

Signed by Ernst Hussinger Notified Body

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



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



Certificate Holder:	BERTHOLD TECHNOLOGIES GMBH & CO. KG CALMBACHER STR. 22 75323 BAD WILDBAD
	Germany BNetzA-CAB-03/22-51
ISED Certification Number:	4777A-IC01X01 authorized by the German Bovernment to act as CAB
CTC Registration Number:	1865 EU Canada of 1st November 1965
OATS Facility ID Number:	3462C-1
OATS Facility:	CTC advanced GmbH Untertuerkheimer Str. 6 -10 66117 Saarbruecken Germany Phone: +49 681 598-0 Fax: +49 681 598-8775 Email: info@ctcadvanced.com
Product Description:	Concentration, Dry Mass, Water, Density measuring system

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

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

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

> CTC advanced GmbH CTC advanced GmbH CTC advanced GmbH Ch-Gerald Schmidt, o-CTC advanced GmbH, ou=SCT-141216 meschemidtgetcadvan ced.com, c=DE 2017.09.08 10:20:11 +02/00'

Saarbrücken

BERTHOLD



CTC I advanced

Certificate Holder:	BERTHOLD TECHNOLOGIES GMBH & CO. KG CALMBACHER STR. 22 75323 BAD WILDBAD	۲	Bundesnetzegentur
	Germany		BNetzA-CAB-03/22-51
ISED Certification Number:	4777A-IC01X12		authorized by the German Government to act as CAB
CTC Registration Number:	1866		EU Canada of 1st November 1998.
OATS Facility ID Number:	3462C-1		
OATS Facility:	CTC advanced GmbH		
Western Constant .	Untertuerkheimer Str. 6 -10		
	66117 Saarbruecken		
	Germany		
	Phone: +49 681 598-0		
	FIGHE. +40.001.000-0		
	Fax. info@ctcadvanced.com		
	Entall, info@cleadvanced.com		
Product Description:	Concentration, Dry Mass, Water, Density measuring	syste	m
		a	
The equipment for which this certi- sale or sold unless the equipment La certification du matériel signifie Les demandes de licences nécess de délivrance d'ISDE et dépender présent certificat est délivré à la procédures d'ISDE. Le matériel à distribué loué mis en verte ou	ficate is issued shall not be manufactured, imported, dist complies with the applicable technical specifications and p seulement que le matériel a satisfait aux exigences de la saires pour l'utilisation du matériel certifié sont traitées en o nt des conditions radio ambiantes, du service et de l'empl condition que le titulaire satisfasse et continue de satisf à l'égard duquel le présent certificat est délivré ne doit p vendu à moins d'être confirme aux procédures et aux	norme onséd aceme aire au oas êtr	I, leased, offered for ires issued by ISED. e indiquée ci dessus. juénce par le bureau ent d'exploitation. Le ux exigences et aux e fabriqué, importé, isations techniques
applicables publiées par ISDE.	venda o mono o cap comonica das procedares es das	opeon	ioniono teorinqueo
I have by attact that the subject	imment was tested and found in semalish south the should	n netr	d encoification
J'atteste par la présente que le ma	ipment was tested and found in compliance with the above itériel a fait l'objet d'essai et jugé conforme à la spécificatio	e-noted	specification.
	and the second second second second second		Street Street Street
			Arrive The Tay
			OTIFIC
	and the second secon		N CERTIFICATIO
	مستعمل معمول فليحو والمراجع والمراجع فالمتحر والمتحر والمتحر والمتحر والمتحر والمتحر والمتحر والمتحر والمتحر والمتحر	1	N.C.
		¥ FO	
	11 - ANN		98 ID NO DE0001
		1 11:	

CTC advanced GmbH

Certification Body

0

Q.

cn=Gerald Schmidt, o=CTC advanced GmbH, ou=SCT-141216,

email=gerald.schmidt@ctcadva nced.com, c=DE 2017.09.08 09:10:01 +02'00

Saarbrücken



GRANT OF EQUIPMENT AUTHORIZATION

тсв

Certification

Issued Under the Authority of the Federal Communications Commission

By:

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

Date of Grant: 09/08/2017

Application Dated: 09/08/2017

BERTHOLD

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

Attention: Dirk Moermann , Dr.

NOT TRANSFERABLE

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



TCB

GRANT OF EQUIPMENT AUTHORIZATION

Certification Issued Under the Authority of the Federal Communications Commission By:

NOT TRANSFERABLE

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

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

TCB

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

Attention: Dirk Moermann , Dr.

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

R9ZFCC01X12 Berthold Technologies Equipment Class: Part 15 Field Disturbance Sensor Concentration, Dry Mass, Water, Density measuring system



Grant Notes



Chapter 8. Technical Drawings



8.1 Dimensions Drawings Evaluation Unit Wall Housing





BERTHOLD

8.2 Electrical Wiring Diagram



8.3 Dimensional Drawings FlowCell

1.00 5 5660-102-00X FlowCell 50 VFL. FDA 10 С Σ 66744 ççz.o) ۳S 6 EN GIOHTREE R 164 Ø 165 Ø 125 Ø18 († ë Aussi kante 齨 1 K20140 anzen ohne Toleranz-sich DIN ISO 2768-mK 4E201800013 5 Dateiname Nodell: 66744-001-NI Dateiname Zeichnung: 66744-001-NI Disse Zeichnung darf ohne schriftliche Zustimmung weder Kopiert noch driften Personen Disse Zeichnung darf ohne schriftlich genütztich genützt werden. Copyrights reserved

8.3.1 Type LB 5660-102-00X FlowCell DN 50 Flange, FDA





8.3.2 Type LB 5660-202-00X FlowCell DN 65 Flange, FDA



8.3.3 Type LB 5660-302-00X FlowCell DN 80 Flange, FDA



8.3.4 Type LB 5660-402-00X FlowCell DN 100 Flange, FDA





8.3.5 Type LB 5660-502-00X FlowCell DN 125 Flange, FDA



8.3.6 Type LB 5660-602-00X FlowCell DN 150 Flange, FDA









8.3.8 Type LB 5660-212-00X FlowCell DN 65 G-BS/M



ş 5 (Tewich!) LB 5660-312-00X FlowCell DN 80 G-BS/M 9 С 66744 M15 705 .00 ta/breuu) Ma fishab 節 (a. 195 🕕 BERTHOLD Υ -AUSSER kanten 囱 Pfen en-, Winkelmasse, Form- u. Vieranzen ohne Toleranz-e nach DIN ISO 2768-mK agefol ы Dateiname Nodell: 66744-015-N15 Dateiname Zeichnung: 66744-015-N15 Discs žeichnung dam fahr schriftlich Zustimmung weder kopieri nach draft desersed Discs zeich anderweitig midzichlich genuizt werden. Copyrights reserved

8.3.9 Type LB 5660-312-00X FlowCell DN 80 G-BS/M





8.3.10 Type LB 5660-412-00X FlowCell DN 100 G-BS/M

8.3.11 Type LB 5660-512-00X FlowCell DN 125 G-BS/M







8.3.12 Type LB 5660-612-00X FlowCell DN 150 G-BS/M



8.3.13 Type LB 5660-132-00X DN 50



8.3.14 Type LB 5660-232-00X DN 65



BERTHOLD





8.3.16 Type LB 5660-432-00X DN 100





8.3.17 Type LB 5660-532-00X DN 125



8.3.18 Type LB 5660-632-00X DN 150





8.4 Dimensional Drawings Container Probes

8.4.1 Type LB 5650-01



8.4.2 Type LB 5650-02





8.4.3 Type LB 5650-03


8.4.4 Type LB 5650-04





8.4.5 Type LB 5650-05



8.4.6 Type LB 5650-09









8.5 Dimensional Drawings Container Flush Probes

8.5.1 Type LB 5651-01





8.5.2 Type LB 5651-02



8.5.3 Type LB 5651-03





8.5.4 Type LB 5651-04



8.5.5 Type LB 5651-05





8.5.6 Installation Situation in Pipelines



8.6Assembly sheets for LB 5650 (Container probe)





8.7 Assembly sheets for LB 5651 (Container flush probe)





Index

A

Accuracy · 60 Adapter flange · 44 ASA flange adapter · 66 Assembly on a container · 35 Assembly on a pipeline · 34

B

Battery · 57

С

Cable installation · 41, 45 Calculation of measured values · 18 CE mark · 13 Compatible with foodstuffs · 28 Compensation · 19 Components · 21 Conductive materials · 19 Connector strip · 48 Container probe · 28 Container probe installation · 43

D

Data format RS232 · 71 Data transfer rate · 50 Device disposal · 53 Digital outputs · 50 Distance rail · 40 Distance to metal · 43

E

Evaluation unit \cdot 23 Evaluation unit installation, flow cell \cdot 40 EVU \cdot 15

F

Factory setting · 15 Fitting flange · 43 Flat gasket · 43 Flow cell · 25 Flush probe · 15, 30 Flushing parameters · 44 Format RS232 · 71 *Frequency license* · 12, 76 Fuses · 58

G

Gas inclusions · 14, 38

Η

HF cable · 32 HF-cable quad · 32

Ι

Incrustation · 30 Installation depth · 43 Installation situation in pipelines · 112, 118

L

LED's \cdot 24, 50 Limitations \cdot 18

М

Mains fuses · 58

0

Overview container probes · 68 Overview flow cells · 64, 65 Overview sealing washers · 68

Р

Pipeline pressure · 38 Power radiation · 14 Power supply · 47

R

Recycling passport · 53 Relay · 50 RS232 interface · 50

S

Safety summary · 7 Salt content · 14 Service instructions · 53 Setting up the evaluation unit, container probe · 45 Symbols · 7 Synchronization · 20

T

technical data \cdot Technical data HF cable \cdot Technical data sensors \cdot Technical drawings \cdot Temperature compensation \cdot Throughput calculation \cdot Transmission power \cdot Transport \cdot

V

Vertical riser \cdot 37

W

Warning · 7 Wear parts · 54 Wiring diagram · 87

Notes



 	 	 			 	 	 	 	 		 	 		 	 		_	
 	 	 		 -	 	 	 	 		_		 		 	 	_		
 	 	 	 	 	 	 	 	 		_		 	_	 	 	_	_	
 	 	 	 	 	 	 	 	 	 			 		 	 	_		



Notes



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





Concentration Meters MicroPolar (++) LB 566

User's Guide Software Manual 41986BA2

Rev. Nr.: 07, 03/2018



The units supplied should not be repaired by anyone other than Berthold Service engineers or technicians by Berthold.

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

The complete user's guide consists of the hardware manual and the software manual.

The hardware manual comprises the

- component description
- > assembly instructions
- electrical installation description
- technical data
- ➤ certificates
- dimensional drawings

The **software manual** comprises the description of the

- ➤ operation
- software functions
- ➤ calibration
- > error messages

The present manual is the software description.

Subject to changes without prior notice.

& Co. KG									
Calmbacher Str. 22 · D-75323 Bad Wildbad									
Service:									
Phone +49 7081 177 111									
Fax +79 7081 177 339									
Service@Berthold.com									

Table of Contents

	Page
Chapter 1. Safety Summary	7
Chapter 2. Communication with MicroPolar	9
Chapter 3. Getting Started Guide	11
Chapter 4. Software Functions	13
4.1 Information on the Menu Structure	13
4.2 Menu Structure	14
4.2.1 Start Menu	16
4.2.2 Diagnostic	16
4.2.3 Setup	18
4.2.4 Access Level	19
4.2.5 Language	19
4.2.0 Configuration 4.2.7 General Data	20
4.2.8 Measurement	21
4.2.9 Plausibility	22
4.2.10 Phase Measurement	23
4.2.11 Offline mode	23
4.2.12 Massflow	24
4.2.13 Calibration	25
4.2.14 System Adjust	25
4.2.15 Calibrate Concentration 4.2.16 Sampling	20
4.2.17 Sample Data (expanded)	27
4.2.18 Advanced	28
4.2.19 Calibration	29
4.2.20 Inputs / Outputs	30
4.2.21 Current Output	31
4.2.22 Current Out 1	31
4.2.23 Current Out 2	32
4.2.24 Current Input	32
4.2.25 Current Input 1 4.2.26 Current Input 2	22
4.2.27 PT100	34
4.2.28 Digital Output	35
4.2.29 Digital Input	35
4.2.30 Service	37
4.3 Trend Display	39
Chapter 5. Configuration	41
5.1 Configuration Setup	41
5.1.1 General Data	41
5.1.2 Measurement	42
5.1.3 Plausibility	42
5.1.4 MICrowave	43
5.1.5 UTILS 5.1.6 Markor	43
5.2 Start Calibration Coefficients	44
5.2 Start Cambration Coencients	45 46
Sis compare nausibility	40

Chapter 6	. Calibration	47
	6.1 System Adjust	47
	6.1.1 Verifying the Reference Values	49
	6.2 Sampling	49
	6.2.1 Entering the Lab Values	51
		52
Chapter 7	'. Start-up of MicroPolar ++	53
Chapter 8	8. Calibration and Advanced	55
	8.1 Plausibility Configuring Phi/Att Ratio	55
	8.1.1 Phi/att Ratio	55
	8.1.2 Process Recording 8.2 Adjusting the Calibration	58 58
	8.3 Output of the Start-up Protocol	60
	8.4 Calibration	61
	8.4.1 Calibration with Two Concentrations	64
	8.4.2 Calibration with Split Value	68
	8.5 Typical Calibration Coefficients/Start Values	70
Chapter 9	. Password	71
	9.1 Password Forgotten	71
Chapter 1	0. Error Lists and Device States	73
	10.1 Power failures	73
	10.2 Temperature errors	73
	10.3 Hardware failures	74
	10.4 Sensor errors	75
	10.5 Analog input range errors	75 76
	10.7 Auxiliary measurement errors	70
	10.8 Analog output range errors	70
	10.9 Watchdog error	77
	10.10 System errors	77
	10.11 Density errors	77
	10.12 Input Error	78
	10.13 Device States	/9
Chapter 1	1. Start-up Protocol	81
	11.1 Example of Start-up Protocol	84
	11.2 Sampling	89



Chapter 1. Safety Summary

Please observe all safety instructions in the *Hardware Manual*, especially those in *chapter 1 Safety Summary*.

NOTICE

Parameter settings

Never change the installation and the parameter settings without a full knowledge of these operating instructions, as well as a full knowledge of the behavior of the connected controller and the possible influence on the operating process to be controlled.

Chapter 2. Communication with MicroPolar

The communication with MicroPolar and MicroPolar ++ is carried out via 4 softkey buttons. The function of the individual buttons changes relative to the position in the menu. Values and texts are entered via an alphanumeric keyboard. The instrument status is indicated by 5 LEDs.



Click on the help button ? in the display footer to view useful information.



Chapter 3. Getting Started Guide

To get started, please carry out the steps described below one after the other.

Please read *chapter 8 Start-up of the MicroPolar* ++ before you take the high dynamics version into operation.

Chapter 4 Software Functions describes all software functions and also serves as a reference guide.

1. Step

Configure the analog inputs as needed: Current inputs 1, 2 and PT100. *See chapter 4.2.22 Input / Output.*



All analog inputs and outputs have already been set in the factory. Therefore, no adjustment work is required during commissioning.

2. Step

Review and edit the software parameters of the application. Some parameters have already been set in the factory. Carry out the steps described in *chapter 5. Configuration*.

3. Step

Carry out the calibration with sampling, *chapter 6. Calibration Flow Cell / Container Probe* or *chapter 7. Calibration Conveyor Belt/Chute*

Temperature compensation is required only if the phase/attenuation will be clearly influenced by the product temperature; this is dependent on the product and water content.

4. Step

Configure the current outputs, digital in- and outputs as needed.

Chapter 4. Software Functions

4.1 Information on the Menu Structure

The menu structure on the following pages provides an overview of all functions of the LB 566. Using the **page numbers** indicated you can look up the function of the depicted window.

You have to enter a password to change from the **Read only** level to **User Mode**. The **Service level** is not accessible due to licensing regulations.

NOTICE

The analysis unit and the software contain functions which are not relevant for the measuring system LB 566 (++) and their measuring application. It concerns the following functions: - Load compensation - Synchronizing the measuring devices

In the operation manual, these functions and the corresponding menus are not described.



4.2 Menu Structure













4.2.1 Start Menu

1 -	LB 566	07.05	5 - 13:25
Live Diag Setu Acce	Display nostic ip ess Level		
Lang	guage		
RUN			

Live Display

Shows the live display.

Diagnostic:

This menu item contains the submenu items data logger, error log, device information and print setup.

Setup:

All necessary inputs for operation of the measuring system can be entered here.

Access level:

Select the access level. Areas protected by passwords can be unlocked.

Language:

Select the dialog language.

4.2.2 Diagnostic



Datalog:

Datalog records the data corresponding to the contents of the serial data output RS232 and RS485 (see *Hardware Manual, chapter 6.6*).

All measured data of a measurement (sweep) are averaged over the averaging time (see below) and stored. This time is dependent on the selected log time. The contents of the datalog can be displayed on the live display, see *chapter 4.3 Trend Display*. Output as a text file is also possible by using RS232 and RS485, or the Memory Tool (optional accessory).

•	Log type	Disable
		single
		continuous
		stop at error
•	Log time	Logging period
		15 minutes to 3 days
•	Restart log	Clears the datalog and starts
		with the above setting
•	Averaging time	Obtained from log time
•	Print log	Printout of tables, output via
		RS232 and RS485, format see chapter 6.6 Hardware Manual
		•

Change datalog settings:

If you change the log type from any to "single", the datalog will be cleared and you start again with the current setting. If you change all other log types and log times, the datalog will not be cleared and you continue with the new settings.

Behavior with stopped measurement:

If the measurement is stopped for some time during the datalog, then the measurement pause will be interpreted as log time in the log type "single". For all other log types, the measurement pause will be added to the log time.

Error log:

Shows the logged error The last 20 error messages will be stored with date and time.

Info:

- Tag
 - Device type : LB 566
- Supplier : Berthold Technologies

:...

- Manufacturer : Berthold Technologies
- Device no. : ...
- Production no. : ...-...
- Software Ver. : V...
- SW rev. date : ...

Print Setup:

Printout of the start-up protocol via RS232 and RS485. Format, contents and example see *chapter 11. Start-up Protocol.*



4.2.3 Setup



Configuration:

Setup of

- General data
- Measurement-specific data
- Plausibility data
- Microwave data
- Units
- Marker

Calibration:

- System Adjust
- Calibrate Conc
- Advanced

Input / Output:

- Current output
- Current input
- PT100
- Digital output
- Digital input

Service:

- Factory Setting
- General Reset
- Memory Tool (operation of the memory tool, optional accessory)
- Data Output (via RS232 and RS485, data contents can be selected)

Product:

Product selection (1-4); if you select another product, the product-specific data will be loaded: outputs, inputs and calibration.

When you select the products 2 to 4 for the first time, all settings and contents (e.g. system calibration, sampling table, datalog and calibration) of the current product will be copied to the new product.

Change password:

The password for the User Mode access level can be changed here.

For more information see *chapter 9. Password*.
4.2.4 Access Level



Read only:

In this mode, the measuring system can be protected against unauthorized access. You can exit this level only by entering a password. The measuring system cannot be started and stopped. You can go to Diagnostic and to Access Level only in the main menu.

User Mode:

- The user mode is the default mode and provides access to all user-relevant parameters.
- On the *Read only* level you have to enter a password.
- The password can be changed.

Service:

• This level is reserved for the service personnel.

4.2.5 Language



Language:

• Select the dialog language

4.2.6 Configuration



General Data:

• Enter date, time and tag

Measurement:

- Meas. Mode (batch/continuous)
- Start Mode (keyboard/external)
- Averaging (number of measured values used for averaging)
- Reset averaging (yes/no)
- Current output at stop

For more information please see *chapter 4.2.8 Measurement*

Plausibility:

- The process limits define the permissible range within which the actual concentration must be.
- The phase measurement is subject to a plausibility analysis, which can be set here.
- Enable and define the pause function

For more information please see chapter 4.2.9 Plausibility

Microwave:

Cable (enter the reference and signal cable length) For example, for 4 m HF cable quad, you have to enter 8 m for both lengths.

Mass flow:

If the density is measured in a pipeline application (unit of concentration = g/m^3 selected), the mass flow (throughput) can be output via current output 1 / 2 (in tons per hour).

Details on the mass flow see *chapter 4.2.12 Mass Flow*.

Units:

Depending on the configuration, different units can be selected for concentrations, current inputs and temperature. For the concentration (1 and 2) you can select: none, specific, %, %TS, °BX, g/L, g/cm³, °Be

For current input 1 you can select: none, specific, °C, °F, g/cm³, kg, t/h

For current input 2 you can select: none, specific, °C, °F, cm, m/s

For the PT100 input you can select: none, °C, °F

Markers:

Enter a value and a name (up to 5 characters) for the marker here. The presentation takes place in the live display and refers to the bar chart. To disable the marker, select a marker value outside of the chart limits or the current output limits.

4.2.7 General Data



Date:

• Enter the current date

Time:

• Enter the current time

Tag:

• Enter the name of the measuring point. The tag (max. 8 characters) is displayed in the header on the display.

4.2.8 Measurement



Meas. Mode:

Select continuous or batch. In Batch mode, an average value is calculated between start and stop. In Continuous mode, a moving averaging is calculated depending on the adjusted averaging number.

Start Mode:

The measurement device can be started or stopped via external terminals (digital input) or via keyboard.

Averaging:

Enter the number of averaging processes. This number indicates over how many measurements the concentration value is to be averaged (moving average). This is true only for the measuring mode Continuous.

Reset Averaging:

Reset averaging (yes/no) This refers to Batch and Continuous.

Current output at stop:

Select "0/4 mA" or "Hold". The selection defines how the current outputs behave with stopped measurement. This is true only for the measuring mode Continuous.



4.2.9 Plausibility



Process Limits:

Enter a permissible measuring range exceeding. If the concentration exceeds the range, the concentration average is put on hold and an error message is displayed (error state). The process limits are independent of the current output limits.

Phase Measure:

The phase is subject to a plausibility analysis. For more information please see *chapter 4.2.10 Phase Measurement*.

Offline mode:

Can be enabled or disabled. Switching variable is the attenuation; if the entered min. attenuation is not reached, the evaluation unit switches to the Offline mode:

- Current output drops to the lower current output limit (0/4 mA)
- Message on display
- RUN LED is flashing

For more information on the offline function see *chapter* 4.2.11 Offline Mode.

4.2.10 Phase Measurement



Sigma max.:

Here you set the maximum sigma of the regression Phase vs. Frequency. During normal measurement operation, sigma lies between 0 and 100. Default: Sigma = 100. With Sigma = 0 the plausibility is turned off.

Phi/att ratio:

The correlation between Phase and Attenuation is another plausibility criterion. Enter a fixed ratio value Phi/att. If you do not know it, you have to record it once, see description below "Auto set". The exact procedure is described in *chapter 8.1.*

Auto set:

The automatic measurement Phi/att ratio is turned on and off here. For more information about the function, *see* chapter *8.1.*

4.2.11 Offline mode

Description:

The offline mode is a software feature with allows detection of a certain product state (e.g. no product present). If the measured attenuation falls below a defined minimum attenuation, the evaluation unit switches to the offline mode. The device status for this mode is described in *chapter 10.4 Device States*.



Enabled:

The Offline mode is turned on and off.

Attenuation min:

Enter the minimum attenuation; if this value is not reached, the measurement switches to the offline mode and will quit this mode again as soon as the value is exceeded.



4.2.12 Massflow

The MASSFLOW menu appears only if two prerequisites have been fulfilled:

- 1. The density unit g/cm^3 is selected for the concentration.
- 2. Current input 2 is enabled and m/s has been selected as unit for current input 2.

For this case, the mass flow (throughput) in tons per hour can be displayed and output via the current output, based on the density reading with indication of the pipe cross-section or the internal pipe diameter.



Massflow calculation:

Here the calculation is enabled or disabled.

Cross-sectional area:

Enter the cross-sectional area of the pipe.

Pipe inside diameter:

Enter the internal pipe diameter.

Comment: Only one entry is required: either the cross-sectional area or the internal pipe diameter, the other parameter will be calculated automatically.

IMPORTANT

The massflow calculation is performed only at varying product speed which must be entered via current input 2.

4.2.13 Calibration



System Adjust:

System calibration is started here. For details see *chapter 4.2.14 System Adjust*.

Calibrate Conc:

Opens the calibration menu of concentration 1.

Calibrate Conc2:

Opens the calibration menu of concentration 2. The second concentration is displayed only if a second concentration is selected under menu | ADCANCED | PROCESS TYPE |.

Advanced:

Here you set the tare values, the number of sweeps when recording samples, the process type and the split value. For more details see *chapter 4.2.18 Advanced*.

4.2.14 System Adjust



Adjust:

System adjustment is started. Phase and attenuation are set to zero, and thus, for example, all cable parameters are considered. This adjustment also forms the reference for the measurement.

The system adjustment (= reference measurement) must be carried out once.

Ref. values:

Upon completion of the reference measurement, the reference values for phase, attenuation, slope and Sigma can be output.

Chart Phi:

Shows the phase versus the frequency.

Chart Atten.:

Shows the attenuation versus the frequency.

A system adjustment will not delete the datalog (see *chapter* 4.2.2 *Diagnostic*).

4.2.15 Calibrate Concentration



Sampling:

Shows all measured samples and entered lab values.

Calibration:

Here

- you select the calibration parameters, the temperature and loading compensation
- the calibration coefficients are calculated automatically
- the calibration coefficients are displayed

For more information see *chapter 4.2.19 Calibration*.

Tuning:

Subsequent correction of the reading is possible by entering a factor and an offset.

Calculation is carried out according to the following formula:

Eq. 4-1:

Corrected display = Display * Factor + Offset

Result:

Presentation of calibration curve, display of correlation and coefficients.

4.2.16 Sampling

1 1/1 Sample # 1 07.05 –13:25		
Next sample		
Active	Yes	
Measured value	65.50%	
Lab value	0.00 %	
Advanced		

The header includes the following information (from left to right):

- Product no.
- Current table position / Total number of entries
- Sample no. of current table position
- Date and time of sampling

Up to 30 sample entries are possible. The sample can be assigned to the lab value either via the sample no. or through data/time. The sample no. is assigned on a continuous basis. If a sample is deleted, the sample no. will not be assigned a second time. Up to 999 sample numbers are available. Only if all numbers have been assigned, you may assign a number for the second time; you will be alerted accordingly by a message on the display.

Next sample:

Continue with the next sample.

Active:

You can choose if this sample should be taken into account in the calibration.

Measured value:

Display of the measured values, calculated with the actual coefficient.

Lab value:

Entry position for the lab value.

Advanced:

Switches to the next data page.

Delete:

Briefly push the softkey to delete the indicated sample entry. Push this key for a longer time to delete all sample entries.

4.2.17 Sample Data (expanded)



Current In 1:

Editable display of the first compensation input.

Current In 2:

Editable display of the second compensation input.

PT100:

Editable display of the PT100 input.

Phi(fm):

Not editable display of the measured phase.

Attenuation:

Not editable display of the measured attenuation.

4.2.18 Advanced

1 - Advanced	07.05 – 13:25
Tare values	
Num. Cal. Sw	eeps 20
Process type	Split Conc
Split Value	75.00 %

Tare values:

Option to enter tare values for phase and attenuation. The tare values are added to the phase and/or the attenuation prior to calibration. The calculation is carried out as follows:

Eq. 4-2 and 4-3

Phase = Phase_{meas} - Phi Tare

Attenuation = Attenuation_{meas} - Phi Tare

Number of Calibration Sweeps:

Freely adjustable number of sweeps over which a calibration point (in the course of automatic sample measurement) will be averaged.

Process Type:

Select the operation mode:

- one concentration [1 measuring range]
- two concentrations [2 measuring ranges]
- split concentration [1 measuring range with switching point (split value) for coefficient switchover].

Split Value:

Setting of the switching point on a value basis.

4.2.19 Calibration

Calibration is performed using the following formula:

Eq. 4-4

Measured value = $A \cdot Phase + B \cdot Attenuation$ + C + D · PT100 + E · Input1 + F · Input2 + G · Load

where:

Meas. value	Concentration / Moisture / Dry mass / Density
A	Phase coefficient
В	Attenuation coefficient
С	Offset
D	Compensation coefficient for PT100 input
E	Compensation coefficient for current input 1
F	Compensation coefficient for current input 2
G	Compensation coefficient for loading

The coefficients can be entered manually or calculated automatically from the entries of the sample table.

Start Calibr.

Starts the calibration using the parameters set and the coefficients are calculated automatically from the entries of the sample table.

Cal. Base

Selection of microwave signals, which are taken into account for the calibration. The following parameters can be set:

- Phase
- Attenuation
- Phase and attenuation

Default: Phase





Compensation Input

Here you can select the analog inputs (PT100, current input 1 and 2) required for compensation. Depending on the enabled analog inputs, the following options can be selected:
None

- In1
- In1 + In2
- In1 + PT100
- In1 + In2 + PT100
- In2
- In2 + PT100
- PT100

Coefficients:

Here all coefficients can be entered directly, e.g. start coefficient.

The automatically calculated coefficients are also stored here. Coefficients that are not used are set to zero.

4.2.20 Inputs / Outputs



Current Output:

Both outputs can be adjusted, assigned and set up on the selected level.

Current Input:

Activation level of current input, calibration and display of the live current signal.

PT100:

Here you can enable and adjust a connected PT100. Display of the actual temperature signal.

Digital Output:

Allocation of relays 1 and 2 and test function.

Digital Input:

Status control and assignment of the digital inputs.

4.2.21 Current Output



IMPORTANT

If a measurement is running, enabling a current input which is not used or not adjusted may cause an error.

4.2.22 Current Out 1

1 - Current Out 1 07.05 –13:25		
Assignment	None	
4 mĀ	60.00	
20 mA	95.00	
Test/Adjust		
Error current		
Current at load underc.		

Assignment:

The following signals can be assigned to the current output:

- None
- Concentration
- Concentration 2 (if active)
- Current input 1 or 2 (if active)
- PT100 (if active)

4 mA:

Display value assigned to the 4 mA value.

20 mA:

Display value assigned to the 20 mA value.



Current output 1 only 4 - 20mA possible

If the current output limit is exceeded, the measurement switches to the warning state, *see chapter* 10.4 *Device States.*

Test/Adjust:

Current test, calibration and display of live current.



The measurement should be stopped for test function.

THOLE

To check the current loop and possibly connected remote displays, you can set a current between 4 and 20 mA via the test function. If you quit the test function, the system automatically switches back to the live current.

Error current:

If the measurement switches to the fault state, a fault current is output via the current output; this can be set here. • 22 mA

- 22 mA
 3.5mA
- Bold
- Value (selectable)

4.2.23 Current Out 2



All functions same as current output 1

Current output 2 can either be set to 0/4 or to 20 mA.

Range:

Change the current output

- 0 20mA
- 4 20mA

4.2.24 Current input



Current In 1:

When selected, change to the activation and calibration menu.

Current In 2:

As described above.

4.2.25 Current Input 1

1 - Current Input 1 07.05 –13:2	5	
Status		
Range		
0/4 mA		
20 mA		
Adjust		
Live current		

Status:

Select yes/no to enable or disable the current input.

Range:

Change the current output

- 0 20mA
 - 4 20mA

0/4 mA:

Display value assigned to 0/4 mA value.

20 mA:

Display value assigned to the 20 mA value.

Adjust:

Follow the instructions on the display.

Live current:

Display of the live current signal.

4.2.26 Current Input 2

Settings correspond to current input 1.



4.2.27 PT100



Enabled:

If a PT100 is connected, the input has to be enabled first.



If a measurement is running, enabling a PT100 input which is not used or not adjusted may cause an error.

PT100 Adjust:

You need a 100 Ohm and a 138.5 Ohm resistance. Follow the instructions on the display.

PT100 Live:

Display of the live temperature.

Set and enabled same as input 1.

4.2.28 Digital Output



The meter has two relays. Relay 1 is linked with LED signal 1 and relay 2 with LED signal 2.

Relay 1:

Different functions can be assigned to relay 1:

- None
- Error
- Hold
- No product
- Alarm min
- Alarm max
- Current at load undercut

Function	Description	
None	Relay and LED function disabled	
Error	In case of error, relay and LED will be set.	
Hold	If Hold function is enabled, relay and LED will be set.	
Offline	If the evaluation unit is in the Offline mode, this will be signaled via the relay and LED.	
Alarm min.	The relay switches if the value falls below the limit value to be set.	
Alarm max.	The relay switches if the value exceeds the limit value to be set.	

Relay 2:

Same assignments possible as above.

Test:

The switching status of the relays can be set here and checked at the respective terminals.

4.2.29 Digital Input



The meter has 3 digital inputs to which different functions can be assigned.

Status:

- Shows the status of the input circuit
- open/closed

DI 1 Function

- The following functions can be assigned to DI 1:
- None
- Start (external start)

THOLD

DI 2 Function

- The following functions can be assigned to DI 2:
- None
- Hold (averaging is stopped)
- Product (external product selection)

DI 3 Function

- Assignments for DI 3:
- None
- Sample (external control of sampling)
- Product (external product selection)

For external start function, the start function has to be set to *External* in the *Measurement* menu window.

Hold means that averaging is stopped, but the measurement continues to run.

Sample means that sampling is started by closing the contact.

Product means that another product is selected by closing the contact (product 1 to 4).



If you select a product for the first time (product 2 to 4), all settings and contents of the current product will be copied to the new product, including:

- Configuration data
- System adjust
- Calibration data (including sampling table)
- Input/Output definitions

To switch over all 4 products, DI 3 also has to be set to product. Please take the terminal configuration from the table below.

Terminals	DI 2	DI 3
	13 / 25	14 / 26
Product 1	open	open
Product 2	closed	open
Product 3	open	closed
Product 4	closed	closed

4.2.30 Service



Factory setting and General reset:

See table on the next side.

Memory Tool:

Refers to the communication with the external memory tool (optional accessory). Data transfer takes place via the 9-pole SubD-connector on the bottom of the instrument.

- Save parameters: All instrument parameters for all products will be saved to the memory tool.
- Load parameters: All instrument parameters stored on the memory tool will be loaded onto the evaluation unit. All operating parameters in the evaluation unit will be deleted.
- Save datalog: The datalog will be saved to the memory tool.
- Save log: The start-up log will be saved to the memory tool.

NOTICE

The concentration average value is put on hold during communication with the memory tool. Thus, the measured value via current output is also put on hold!

Data printout:

All measured values are output for each measurement via the serial data interfaces RS232 and RS485. The output can be set as follows:

- None (disabled)
- Row (data transfer, see *Hardware Manual, chapter 6.4*)
 - Table (microwave data for each frequency point)
- Row and table

"Row" is defaulted.



	Factory setting	General reset
Language selection	unchanged	unchanged
Access level	unchanged	default: User mode
Measurement	stopped	stopped
Password	unchanged	default: PASS1
Product selection	unchanged	all products deleted
Error log	not deleted	deleted
Data log	not deleted, default settings	deleted, default set- tings
System Adjust	not deleted	deleted
Cable length	unchanged	default
Sampling	not deleted	deleted
TAG label	default	default
All parameters on menu: Measurement Plausibility Marker Units	default	default
Calibration coefficients	default	default
All settings for the analog and digital inputs and out- puts	default	default
Adjustment of the analog in- puts and outputs	unchanged	unchanged
Comment:	affects only the cur- rent product	affects all products (P1 to P4)

*Default: Default values, see chapter 11.1 Example Start-up Protocol

4.3 Trend Display



Push the **ZOOM** button to enlarge the measurement value which is surrounded by a frame.



By pushing the **ZOOM** button for a longer time, the enlarged measurement value will be displayed as trend over the entire display.

The trend display corresponds to the contents of the datalog. The datalog has to be enabled for the trend display.



As long as the trend builds up, the measured value and/or the current output are put on hold!



Chapter 5. Configuration

Before doing any calibration work, you have to enable and configure the required analog inputs and check and, if necessary, correct the configuration parameters.

If the required inputs are not enabled, some menus are not displayed and a proper configuration and calibration is not possible under certain circumstances. The current outputs, digital outputs can be enabled and configured after the calibration.

The measuring system includes two separate floating current outputs.

5.1 Configuration Setup



Starting from the main menu, go to the display shown on the left via | SETUP |

> CONFIGURATION

5.1.1 General Data



➢ GENERAL DATA



Date

ESC

07.05.2015

?

Example:

Select the respective entry, edit and store it.

DATE

Push **DEL** to delete the entry and then enter the new date. Push **V** to confirm and store the changed date.

TIP The colon for the time input (e.g. 13:25) is invoked by pushing the button [.].

BERTHOLD

5.1.2 Measurement

DEL



You have to check the settings on this display and adapt them to the measurement conditions.

Averaging over 20 measurements is a good choice as a rule.

5.1.3 Plausibility



The **process limits** need to be adjusted. Allow exceeding of the measuring by \pm 5% absolute. Example: The measuring range is 65-95% TS. Enter 60-100% TS as process limit.

The process limits are independent of the current output limits.

For setup of the **phase measurement**, see *chapter 5.3 Configuring Plausibility*.

The **Offline mode** is deactivated by default. Take this function into operation, when necessary, only when the measuring system was put into operation without errors. Refer to the description in *chapter 4.2.11 Offline Mode*.

5.1.4 Microwave



CABLES



If the factory-set cable lengths do not match the actual geometry conditions, you have to correct the values.

Example: For a 4 m long HF quad cable, enter 8 m for the reference and signal cable length. The input value corresponds to twice the quad cable length.

5.1.5 Units



Set the units to the desired dimension.

> UNITS



The units of the concentrations (conc 1 and 2) and those of the enabled analog inputs can be selected.

CONC / CONC 2



Different units can be set for both concentrations.

≻ %







➢ CURRENT IN 1

≻ °C

The temperature input can be set to °C, °F, none or specific.

5.1.6 Marker



You can set a marker comprising max. 5 characters which identify the value set in the live display.

> MARKERS

5.2 Start Calibration Coefficients

1 Calibrate Conc 07.05	-13:25
Sampling	
Calibration	
Result	
1 Calibration 07.0	5 12.25
1 - Calibration 07.0	5 –13:25
1 - Calibration 07.03 Start Calibr.	5–13:25 PHI
1 - Calibration 07.03 Start Calibr. Cal. Base Loading comp.	5–13:25 PHI

Coefficients

Starting from the main menu, go to the display shown on the left via: ESC | SETUP | CALIBRATION | CALIBRATE CONC |

CALIBRATION

➢ COEFFICIENTS

For the default setting, the concentration is calculated as follows:

Measured value = $A \cdot Phase + C$ Eq. 5-1

where: A, C: Calibration coefficients



Check the coefficients A and C and correct them, if necessary, as follows:

C = average measuring range value (concentration value) A = 0

All coefficients that are not needed are automatically set to zero.

Note: With these calibration coefficients the concentration average value and thus the current output is put on hold during startup.

5.3 Configure Plausibility



Starting from the main menu, go to the display shown on the left via:

| SETUP | CONFIGURATION | PLAUSIBILITY | PHASE MEASURE |

The display on the left shows the default settings, which may have to be entered.

During normal measurement operation, **Sigma** lies between 0 to 100. Therefore, Sigma_{max} = 100 is a good choice for most applications. However, if you constantly get sigma values up to 300 during the measurement, you may increase the limit up to 500. Higher sigma values than that usually indicate a fault, such as continuous air bubbles which have to be eliminated.

For most applications, the **Phi/att ratio** is not known; therefore, it has to be determined once. For this, you need the Auto setting, which otherwise must always be set to OFF. The exact procedure is described in *chapter 9.1*.



Chapter 6. Calibration

Note: The measuring system must have reached normal operating temperature (approx. 45 min. warm-up time)).

The flow cell must be completely filled with product or the container probe is completely surrounded by product.

The measuring system has to operate at a normal throughput and/or material flow and the usual material under actual operating conditions.

Prerequisite: Chapter 5. Configuration have been completed.

6.1 System Adjust



MicroPolar (++) LB 566





6.1.1 Verifying the Reference Values



Starting from the main menu, go to the display shown on the left via | SETUP | CALIBRATION | SYSTEM ADJUST |

➢ REFERENCE VALUES

Limits of important parameters for the reference measurement:

1 - Rei	f. Values	s 07.05	- 13:25
Phi(Atte Slop Sign	fm) nuation ne120.55 na	125.00 2 5 °/GHz	0 °/GHz 22.5 dB 0.00

Sigma: <300 (reliable microwave irradiation) Attenuation: <45 dB with standard evaluation unit MicroPolar <65 dB with high dynamic evaluation unit MicroPolar ++ (Cables connected properly, product irradiatable)

6.2 Sampling





For temperature compensation, the product temperature must be entered via one of the analog inputs and in addition the corresponding input has to be enabled. If not, the product temperature is not stored in the device during sampling.

If the measuring system is not yet in the measurement mode, start the measurement now.

Push **RUN** to start the measuring system.

Push \checkmark to confirm the safety prompt and the device switches to the measurement mode.





The display to the left appears if you push RUN.

Note: Push the **SAMPLE** button to start measurement of the raw data. At the same time, the laboratory sample has to be taken and marked. The analysis may be performed later, provided the product is not changed by this.



Sampling is in process.....

Push the X button to stop the sampling process any time..



If the sampling process has been completed without any problem, push the \checkmark button to save the sample in the table and the measurement continues.

The process previously described must be repeated for each additional sample.

The moisture/concentration of the samples should be distributed over the entire measuring range. For additional temperature compensation, the temperature of the samples should be distributed over the entire temperature range.

The minimum number of samples required is dependent on the selected calibration modes. If the sample size is too low, an error message is displayed after you have attempted to run a calibration.

About six samples suffice for a rough calculation of the calibration coefficients, provided the concentration differs by at least 5%. At least 15 samples are required for fine calibration and temperature compensation.

6.2.1 Entering the Lab Values



Chapter 6. Calibration



Sampling Calibration Tuning Result	l
1 1/1 Sample # 1 07.05 – 13:25 Next sample Active Yes Measured value 65.50% Laborwert 0.00 % Advanced ■ DEL ▲ ■	

> SAMPLING

LAB VALUE



Delete default value with $\overline{\text{DEL}}$ and enter new value and confirm with $\sqrt{}$

1 1/1 Sample # 1 07.05 –13:25		
Next sample		
Active	Yes	
Measured value	65.50%	
Lab value	72.40 %	
Advanced		
	\checkmark	

➢ NEXT SAMPLE

and repeat the step described above with the next sample.

After you have entered the last sample by pushing the der dutton you get back to the Calibration menu. (Short push – one page, longer push of the button – you get back to the Calibration menu immediately)

6.3 Calibration

Proceed as described in *chapter 8.4*.



Chapter 7. Start-up of MicroPolar ++

System adjustment and calibration are carried out in just the same way for MicroPolar ++, as they are for the standard system MicroPolar. However, please keep in mind that the ++ unit requires a minimum attenuation of 40 dB over the entire concentration range and during system adjustment. When falling below, the measurement is not precise anymore.

The total attenuation is calculated as follows:

 $dB_{total} = dB_{adjust} + dB_{live} + 0.4 x signal cable length$ **Eq. 7-1**

where:	
dB _{total} :	Total attenuation
dBadjustment:	Attenuation during system calibration
dBlive:	Current attenuation in the measure-
ment mode	
Signal cable length:	e.g. 4 m HF-cable quad results in 8 m signal cable length (to and from)

The evaluation unit monitors the entire attenuation automatically and reports a falling below by an error message (error no. 55).

Remedy when falling short of the attenuation:

If the required overall attenuation is not reached, you have the option to install a 10 dB fixed attenuator into the transmitting branch (see Fig. 7-1). The standard model MicroPolar should be used if the attenuation is clearly below the required value.



MicroPolar (++) LB 566
Chapter 8. Calibration and Advanced

8.1 Plausibility | Configuring Phi/Att Ratio



Starting from the main menu, go to the display shown on the left via | SETUP | CONFIGURATION | PLAUSIBILITY | PHASE MEASURE |

Check or correct corresponding to the settings shown on the left.

> PHI/ATT RATIO

If you know the ratio value, enter it here; otherwise go to *chapter 8.1.1 Phi/att Ratio*.



Delete default value with $\overline{\text{DEL}}$ and enter new value and confirm with \checkmark

8.1.1 Phi/att Ratio

If you do not know the ratio value, proceed as follows:

- 1. Perform a process recording, see *chapter 8.1.2 Process Recording*. A prerequisite is that the process covers the entire concentration range during the recording.
- If the process recording is currently not possible or if the concentration range is small anyway (< ± 5%), then enter Phi/att = 1 as start value.

Subsequent adjustment of the process recording is possible.



8.1.2 Process Recording

Prerequisite:

The steps described in chapter

5. Configuration

6.1 System Calibration

have been completed.

The process recording is used to determine the ratio of phase and attenuation (Phi/att), a parameter of the plausibility analysis for correct determination of the phase.

If you already know the ratio from other measurements, you may enter it directly on the PLAUSIBILITY menu (see *chapter* 4.2.10 *Phase Measurement*); in this case, that process need not be recorded.

IMPORTANT

The measurement takes place automatically; you only have to start and stop it again. During measurement, please keep in mind:

- Do not stop the measurement
- Do not change the concentration erratically (max. 1 %).
- Cover the entire measuring range, if possible



1 | Phase Measure | 07.05 – 13:25 Sigma max. 100.00 Phi/att ratio 1.00 Auto set OFF Starting from the main menu, go to the display shown on the left via:

| SETUP | CONFIGURATION | PLAUSIBILITY | PHASE MEASURE

PHASE MEASUREMENT

AUTO SET

B<u>E</u>RT<u>H</u>QL<u></u>D



> ON

1 | Phase Measure | 07.05 – 13:25

Pause recording:

You have the option to pause recording by turning the recording off. The recording is paused and starts again only after it is turned on again.

Start new recording:

Prerequisite: Recording is turned off. Stop and start the measurement before you start a new recording. The result of older recordings will be deleted.

Stop recording:

After recording, simply turn off the automatic recording; the recorded Phi/att is stored automatically.

Carry out sampling while the process recording is running. Do not forget to enable the process recording again as described above!



8.2 Adjusting the Calibration

A correction factor and an offset may be entered later to obtain a subsequent adjustment of the calibration (fine tuning).

Below please find an example for an offset adjustment.

The display to the left appears if you push RUN.

The display reading is now compared with the analysis value of the lab sample. The difference has to be entered as offset with the correct algebraic sign.

Calculation:

Analysis value – Display = Offset

Push **ESC** to return to the main menu.

Eq. 8-1

 1 | - | LB 566 | 07.05 – 13:25

 Live Display

 Diagnostic

 Setup

 Access Level
 User Mode

 Language
 English

1 | - | Live Display | 07.05 – 13:25

ESC SAMPLE ▲▼ ZOOM

Conc. av.

Concentration av.

65.50 %

Conc. act. 64.35%

> SETUP





CALIBRATION

CALIBRATE CONC





> TUNING

1 - Calibration	07.05 –13:25
Factor	1.00000
Offset	0.000

> OFFSET

Calculation formulas see *chapter 4.2.16 Calibrate Concentration*.



Enter the calculated offset value, confirm with \checkmark button and push the Home button \checkmark four times to return to the main menu.

1 - LB 566 0	7.05 – 13:25
Live Display	
Diagnostic	
Setup	
Access Level	User Mode
Language	English
STOP 🔺	

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

 Concentration av.

 75.50 %

 Conc. av.

 Conc. av.

 Conc. act.

 64.35%

 ESC

 SAMPLE

Select

LIVE DISPLAY

to get back to the display.

The reading value should now correspond to the actual value.



8.3 Output of the Start-up Protocol



Starting from the main menu, go to the display on the left by selecting | DIAGNOSTIC |

➢ PRINT SETUP



Push the \checkmark button to print the setup via RS232 and RS485. Push X to go back one page without printout.

The start-up protocol includes all parameters, system adjustment data, calibration data and entries in the sample table.

For further information on the start-up protocol see *chapter 11*. Start-up Protocol

8.4 Calibration

There are several setup options for calibration; for details see *chapter 4.2.20 Calibration*. Furthermore, the default settings are displayed, which, as a rule, are the best choice.

Prerequisite: The steps described in chapter 5. Configuration 6.1 System Calibration 6.2 Sampling have been completed.

1 Calib	orate Co	nc 07.0	5 –13:25
Sa	npling		
Ca	libration		
Tui	ning		
Re	sult		

-13:25
PHI
None

1 - Calibration 07.05 –13:25 Cal. Base
Phase
Attenuation
Both

Starting from the main menu, go to the display shown on the left via: | SETUP | CALIBRATION | CALIBRATE CONC |

> CALIBRATION

➢ CAL. BASE

PHASE (Phase measurement)

Standard for all applications: Phase



Set the desired compensation:

1 - C	alibratio	n 07.0	5 – 13:25
Sta Ca Lo	art Calib I. Base ading co	r. omp.	PHI
Cc	mp. inpl	ut	None
Co	efficient	s	
	\land		

For compensation such as **temperature compensation** proceed as follows:

COMPENSATION

Here you can select the analog inputs (PT100, current input 1 and 2) required for compensation (e.g. temperature compensation). You can select:

- None
- In1
- In1 + In2
- In1 + PT100
- In1 + In2 + PT100
- In2
- In2 + PT100
- PT100

Select "None" if no compensation is required.

The automatic calculation of the calibration coefficients starts as soon as you have set the parameters for the compensation.

START CALIBRATION





Push the \checkmark button to start the calibration; push X to go back one page without calibration.

OK accepts the calibration and changes to the next display.

When calculating the new coefficient set, the Factor will be reset to 1 and the Offset to 0.







The graph on the left shows the measured value versus the lab value.



Output of the correlation between measured value and lab value.

OK



As soon as you confirm this prompt, the calibration display appears again; from there you get back to the main menu by pushing \triangle four times.



8.4.1 Calibration with Two Concentrations

1 | - | Calibration | 07.05 – 13:25 System Adjust Calibrate Conc Advanced







1 - Calibrate Conc1 07.05 –13:25
Sampling
Calibration
Tuning
Result



left via

cess type as described below.

| SETUP | CALIBRATION |

PROCESS TYPE

> 2 CONC

Push the \checkmark button to accept the selected process type and push the $\triangle \triangleleft$ once to go to the display depicted below.

Calibration for two concentrations starts by changing the pro-

Starting from the main menu, go to the display shown on the

> CALIBRATE CONC (corresponding to concentration 1)

> SAMPLING



There is only one sample table for both calibrations.

The lab values have to be entered for all samples used for calibration of concentration 1. All other samples have to be disabled (Active: Yes/No).

1 | 1/4 | Sample # 1 | 07.05 – 13:25 Next sample Active Yes Measured value 65.50% Lab value 0.00 % Advanced ■ DEL

LAB VALUE



Delete default value with $\ensuremath{\overline{\text{DEL}}}$ and enter new value and confirm with $\ensuremath{\overline{\sqrt{}}}$

1 1/4 Sample # 1	07.05 – 13:25
Next sample	
Active	Yes
Measured value	e 65.50 %
Lab value	60.40 %
Advanced	
	√

- NEXT SAMPLE
- Continue with next sample

1 2/4 Sample # 2 07.05 –13:25		
Next sample		
Active	Yes	
Measured value	74.35 %	
Lab value	67.80 %	
Advanced		
	\checkmark	

\triangleright	ACTI	VE
-	/	

Disable sample

1 2/4 Sample # 2 07.05 – 13:25 Active
No
Yes
ESC ? DEL √

> NO



1 2/4 Sample # 2 07.05 –13:25				
Next sample				
Active	No			
Measured value	74.35 %			
Lab value	67.80 %			
Advanced				
	\checkmark			

Make sure that all samples have been processed and only those samples are active which are relevant for this calibration.

Push d to get to the Calibration page.



1 - Calibration 07.05	- 13:25
Start Calibr.	
Cal. Base	PHI
Loading comp.	
Comp. input	None
Coefficients	

1 | - | Calibration | 07.05 –13:25 Calibrate Now?



➢ START CALIBRATION

Push the \checkmark button to start the calibration; push \times to go back one page without calibration.

OK	accepts	the	calibration	and	changes	to	the	next	display.
----	---------	-----	-------------	-----	---------	----	-----	------	----------





Push $\Delta \triangleleft$ twice to return two pages.

Repeat the steps as described above for concentration 2; all samples have to be enabled again in the sample table. Now you have to disable all samples which are not used for concentration 2.



CALIBRATE CONC 2

1 - Calibrate Conc2 07.05 –13:25		
Sampling		
Calibration		
Tuning		
Result		

> SAMPLING



8.4.2 Calibration with Split Value

 1 | - | Calibration | 07.05 – 13:25

 System Adjust

 Calibrate Conc

 Advanced

 I | - | Advanced | 07.05 – 13:25







With this type of calibration, two characteristic curves (concentrations) are combined in one measuring range; their point of intersection defines the split value.

Conc 1 for the lower and conc 2 for the upper measuring range can be output only together via current output.

> ADVANCED

➢ PROCESS TYPE

> SPLIT CONC

Push the \checkmark button to accept the selected process type and push the $\triangle \triangleleft$ button once to go to the display depicted below. The displayed split value has been set by the manufacturer, but has to be adapted to the respective application.

The sample measurement should be selected such that the last sample of the lower concentration is fairly close to the first sample measurement of the upper concentration. Ideally, the last sample of the initial concentration is the first sample of the final concentration.



Sample measurements are carried out continuously over the entire measuring range with the display depicted to the left. See *chapter 6.3 Sampling*

After completion of sampling, the individual samples will be enabled or disabled during input of the laboratory values, relative to the set split values. All samples smaller or equal to the split value will be assigned to the lower concentration range and all samples above to the upper concentration range.

The assignment of the samples is carried out automatically, for example, by setting the split value or by entering the lab values have been entered (e.g. after re-sampling). The assignment depends on the split value and the lab value.

IIMPORTANT

The split value entry allows you to enable samples that have been disabled earlier through automatic assignment! In these cases, disabled samples should better be deleted or disabled again after a split value entry!

1 | - | Advanced | 07.05 – 13:25 Tare values Num. Cal. Sweeps 20 Process type Split Conc Split Value 75.00 %

1 - Advanced 07.05 – 13:25	
Split Value	
75.00 %	
ESC ? DEL √	

The split value to be set must correspond to the point of intersection of both calibration curves. This will be corrected automatically after the calibration (within certain limits).

SPLIT VALUE

Enter the split value and confirm with \checkmark .

Push Δ to get to the Calibration page.





CALIBRATE CONC

> CALIBRATION

The lower concentration is now calibrated. Then select CONC2 and repeat the calibration process. Go back to the main menu and start the measurement.

8.5 Typical Calibration Coefficients/Start Values

C: Concentration value at system calibration

For applications with container probe or flow cell nominal width DN 50 and without temperature compensation

- A = 0.19 to determine the concentration of dry matter
- A = + 0.19 to determine the concentration of moisture

THOLD

Chapter 9. Password

The measuring system can be protected against unauthorized access by passwords.

The access levels are as follows.

Read only

The measuring system cannot be started and stopped. You can only switch from the live display to Diagnostic and to Access Level.

User mode

The user mode is the default mode and provides access to all user-relevant parameters.

Service

The service level is reserved to service personnel.

You have to enter a password to change from **Read only** to **User Mode**.

At the time of delivery, this password is

PASS1

The password can be changed in the menu \mid SETUP \mid CHANGE PASSWORD \mid .

9.1 Password Forgotten

The device is in the "Read only" mode and the user has forgotten the password. Please proceed as follows to carry out a "Reset" of the user level:

Turn off device.

Turn on device; as soon as all 5 LEDs light up when powering up, press 0 (zero) and keep it depressed for 8 seconds.

Device powers up in the "User Mode". You can now enter a new password.

IMPORTANT

Check your process before turning off the device. The current outputs drop to 0 mA.

Chapter 10. Error Lists and Device States

The LEDs indicate the device status. Once the errors have been corrected, the measurement returns to the state before the error occurred. An acknowledgment is not required.

10.1 Power failures

Code	Error	<i>Probable cause / correction</i>
10	24V power fail	<i>Please contact the Berthold Technologies Service.</i>
11	9V power fail	<i>Please contact the Berthold Technologies Service.</i>
12	5V power fail	<i>Please contact the Berthold Technologies Service.</i>
13	3V power fail	<i>Please contact the Berthold Technologies Service.</i>
14	Battery fail	Battery power is low, re- place immediately, see Hardware Manual, chap- ter 5.4

10.2 Temperature errors

Code	Error	<i>Probable cause / correction</i>
20	Attention: Ambient temperature too high!	<i>Check operating tempera- ture of the evaluation unit, permissible range: -20 to 50° C or 45° C</i>
21	RF temperature out of range	<i>Check operating tempera- ture of the evaluation unit, permissible range: -20 to 50° C or 45° C</i>



10.3 Hardware failures

Code	Error	<i>Probable cause / correction</i>
30	Program memory corrupted	<i>Please contact the Berthold Technologies Service.</i>
31	Data memory corrupted	<i>Please contact the Berthold Technologies Service.</i>
32	Parameter memory corrupted	<i>Compatibility check after software download: A general reset must be carried out.</i>
33	I2C-bus communica- tion	<i>Please contact the Berthold Technologies Service.</i>
34	DAC update failure	<i>Please contact the Berthold Technologies Service.</i>
35	LCD contrast	<i>Please contact the Berthold Technologies Service.</i>
36	LCD controller	<i>Please contact the Berthold Technologies Service.</i>
37	Keypad error	<i>Please contact the Berthold Technologies Service.</i>
38	RF communication er- ror	<i>Please contact the Berthold Technologies Service.</i>
39	RF hardware failure	Faulty cable connection between the motherboard and HF unit. Check con- nector on the mother- board. Caution! First, disconnect the evaluation unit from the power supply!
40	I/O communication er- ror	<i>Please contact the Berthold Technologies Service.</i>
41	I/O module error	<i>Please contact the Berthold Technologies Service.</i>
42	RF Board startup error	<i>Please contact the Berthold Technologies Service.</i>

10.4 Sensor errors

Code	Error	<i>Probable cause / correction</i>
50	Phase variance too high	<i>The measured phase ex- ceeds the allowed Sigma limit.</i>
53	No product	<i>The evaluation unit is in the Offline state (no product present).</i>
54	No system adjustment available	The system calibration has not yet been carried out.
55	Insertion loss under- run	See Software Manual, chapter 8

10.5 Analog input range errors

Code	Error	<i>Probable cause / correction</i>
60	Current input 1 out of range	The enabled current input has not yet been cali- brated or is not occupied.
61	Current input 2 out of range	The enabled current input has not yet been cali- brated or is not occupied.
62	Pt100 temperature out of range	The enabled PT100 input has not yet been cali- brated or is not yet occu- pied.

10.6 Measurement range errors

Code	Error	<i>Probable cause / correction</i>
70	Concentration out of range	The concentration is out- side the process limits.
71	Concentration 2 out of range	Concentration 2 is outside the process limits.
72	Loading value 1 below the limit	Below the minimum load for concentration 1
73	Loading value 2 below the limit	Below the minimum load for concentration 1
74	Loading comp. disa- bled. Cur. Input upper & lower value invalid	Current input is outside the range.
75	Sync. time too long	<i>Review settings for syn- chronization, see chapter 7.3 in Hardware Manual.</i>
76	Synchronization: speed out of range	<i>Review settings for syn- chronization, see chapter 7.3 in Hardware Manual.</i>
77	Waiting for sync. value	<i>The measurement has not yet been synchronized, please wait.</i>

10.7 Auxiliary measurement errors

Code	Error	<i>Probable cause / correction</i>
78	Mass flow 1 calc. disa- bled. Product speed in- valid	<i>Please contact the Berthold Technologies Service.</i>
79	Mass flow 1 calc. disa- bled. Density invalid	<i>Please contact the Berthold Technologies Service.</i>
80	Mass flow 2 calc. disa- bled. Product speed in- valid	<i>Please contact the Berthold Technologies Service.</i>
81	Mass flow 2 calc. disa- bled. Density invalid	<i>Please contact the Berthold Technologies Service.</i>

10.8 Analog output range errors

Code	Error	Probable cause / correction
90	Current output 1 out of range	The concentration calcu- lated on the basis of the current is outside the cur- rent range
91	Current output 2 out of range	The concentration calcu- lated on the basis of the current is outside the cur- rent range

10.9 Watchdog error

Code	Error	<i>Probable cause / correction</i>
92	Watchdog	<i>Please contact the Berthold Technologies</i>
		Service.

10.10 System errors

Code	Error	<i>Probable cause / correction</i>		
120	<i>No time / date setting</i>	<i>Please enter the date and the time.</i>		

10.11 Density errors

Code	Error	<i>Probable cause / correction</i>
150	Density calc.: Radio- metric MPUA out of range	<i>Check the measured value of the radiometric MPUA</i>
151	Density calc.: Height signal out of range	<i>Check the measured value of the layer thickness sensor</i>



10.12 Input Error

Error	Probable Cause
Value too	Input value is too large
large	
Value too	Input value is too small
small	
Table is	Sampling has been selected without
empty	previous sample measurement
Chart data	The measuring system has determined
faulty	faulty chart data during calibration.
No chart data	The calculated chart data have been de-
available	leted or calibration has not been com-
	pleted.
Sampling full	You have tried to measure more than 30
	samples.

10.13 Device States

Error state:

This state occurs also in error codes 50 to 56, 60 to 62 and 70 to 71 (see table above). The evaluation units behave as follows:

<u>LLD3</u> .	Ron hashes, Error on, signal I and 2
	depending on the configuration.
Current outputs:	Fault current, as selected
Display:	Error message with error code

Warning state:

This state occurs also in error codes 14, 21, 90 and 91 (see table above). The evaluation units behave as follows: <u>LEDs</u>: RUN flashes, ERROR off, signal 1 and 2

no connection.Current outputs:liveDisplay:Error message with error code

Hold state:

Measurement stopped via digital input. The evaluation units behave as follows:

The averaged concentration value is frozen. The measurement continues, however, so that a measurement error can cause the fault condition also from the hold state.

<u>LEDs</u> :	RUN flashes, ERROR off, signal 1 and 2
	depending on the configuration.
Current outputs:	frozen
Display:	No display message

Offline state:

The evaluation unit is in the Offline state, no product is present. The evaluation units behave as follows: LEDs: RUN flashes, ERROR off, signal 1 and 2

<u></u> .	
	depending on the configuration.
Current outputs:	lower current output value 0/4 mA
Display:	Error message with code no. 053 spec-
	ification

Chapter 11. Start-up Protocol

The log can be output via RS232 and RS485. The printout takes place on the menu \mid DIAGNOSTIC \mid PRINT SETUP \mid .

The serial interfaces RS232 and RS485 have the following port settings:

Data transfer rate 38400 Bd, 8 data bits, no parity, 1 stop bit

The log is saved to a TXT file using a terminal program. To view the log (e.g. in $Excel^{(R)}$), the following data format must be considered.

Separator:TabulatorDecimal separator:.Thousand separator,

The following code list helps you to interpret the start-up protocol, see example in *chapter 12.1*.

Parameter	Code	Information
	no.	
Log type		Log type:
	0	Disabled
	1	Single
	2	Continuous
	3	Stop on error
Log time		Log time:
-	0	15 mins
	1	1 hour
	2	4 hours
	3	8 hours
	4	1 day
	5	3 days
Measuring		Meas. Mode:
mode	0	Continuous
	1	Batch
Start mode		Start mode (Start/Stop):
	0	Keyboard
	1	Extern

Parameter	Code	Information	
	no.		
Calibration		Exp. cal. input selection:	
input	0	None	
selection	1	Input 1	
	2	Input 1 + Input 2	
	3	Input 1 + PT100	
	4	Input 1 + Input 2 + PT100	
	5	Input 2	
	6	Input 2 + PT100	
	7	PT100	
Calibration		Calibration base:	
variable	0	Phase	
	1	Attenuation	
	2	Phase and attenuation	
Massflow		Throughput calculation:	
calculation	0	Off	
mode	1	On	
Measure		Process type:	
configuration	0	1 concentration	
-	1	2 concentrations	
	2	Split concentration	
AO Assign		Assignment of current output:	
Code	0	None	
	1	Concentration	
	2	Concentration 2	
	3	Current input 1	
	4	Current input 2	
	5	PT100	
AO Alarm se-		Error current for current output:	
lect code	0	22 mA	
	1	3.5 mA	
	2	Hold	
	3	Value	
Range		Measuring range for current out-	
selection	0	put:	
	1	0 20 mA	
		4 20 mA	
AI Range		Measuring range for current input:	
selection	0	0 20 mA	
	1	4 20 mA	

T E C H N O L O G	E	S	

Parameter	Code no.	Information
AI Ena- bled[2]		State current in 2
DO Function		Function of digital outputs:
	0	None
	1	Error
	2	Hold
	3	No product
	4	Alarm min.
	5	Alarm max.
DO Assign-		Digital output: the min./max.
ment		alarm is assigned to the following:
	0	Concentration
	1	Concentration 2
	2	Current input 1
	3	Current input 2
	4	PT100
DI Function		Function of digital inputs:
selection	0	None
	1	Start/Stop
	2	Hold
	3	Sampling
	4	Product selection
Printout		Form of data printout:
mode	0	None
	1	Line
	2	Table
	3	Line + Table
Access level		Access level:
	0	Read only
	1	User mode
	2	Service
Language		Language selection
	0	English
	1	German
	2	French

11.1 Example of Start-up Protocol

Menu:	Start of Setup:	Start-up Prote	ocol		Interpretation:							
					(* Only relevant for service)							
Due du et	Entry.	Des duet4	Prod-	Prod.	Prod.							
Product		Pioducti	uciz	3	4	Les tracies ende list						
Data log	Log type .	0				Log type. see code list						
	Log lime.	2				Log time. see code list						
	NTC temperature :	45.2 °C										
		45.3 C				*						
	max. NTC temperature :	46.7 C				*						
	9v power supply :	8.94 V										
Info	Tag :	-				Tag						
		LB 500				Device type						
	Unique device ID number :	761										
	Serial number :	4294967000										
	Final assembly number :	000-000										
	Software version :	2.00										
	Software release date :	02.02.2016				Software revision date						
	Actual date :	10.02.2016				Date of logging						
Measure-	Actual time :	12:15				Time of logging						
ment	Measuring mode :	0				Measuring mode: see code list						
	Start mode :	0				Start mode: see code list						
	Filter damning value[2] ·	20				Current out-						
	Filter damping value[3]	20				Averaging number when sampling						
	Reset average :	EALSE				Reset Averaging: Yes/No						
Plausibil-	Treset average .	TALOL										
ity	Lower limit :	0.00				Min. process limit:						
	Upper limit :	100.00				Max. process limit						
	Raw data average value	5				*						
	Max. phase sigma :	100				Sigma max.						
	Correlation Phi/Att :	1:00				Ratio phase/attenuation						
	Auto-set mode :	FALSE				Auto set: On/Off						
	Offline Mode detection :	FALSE				Offline function: On/Off						
Miero	Minimum attenuation :	-15 dB				Limit for offline mode						
wave	Ref. cable length :	8.00 m				Reference cable length						
	Meas. cable length :	8.00 m				Signal cable length						
	Wave band selection :	1				*						
	Start frequency :	2				*						
	Frequency step :	1				*						
	Nbr of freq. points :	5				*						
Marker	Marker name :	Mark1				Marker name for concentration						
	Marker value :	75.00 %				Marker value for concentration						
	Marker name[2] :	Mark2				Marker name for concentration 2						
	Marker value[2] :	75.00 %				Marker value for concentration 2						
System adjust	Nbr of sweeps for reference	1				*						
EV/II type	HE amplifier mode -	i				*						
- vo type	Minimal insertion loss :	40.00 dB				*						



Calibrate	Calibration input selection :	0	Exp. Cal. input selection: see code list
CONC	Calibration mode	0	*
	Calibration variable :	0	Calibration basis: see code list
	Phase coefficients :	0	Phase coefficient A
	Attenuation coefficients :	0	Attenuation Coefficient B
	Constant coefficient :	0	Constant C
	d coefficient	0	Comp-coefficient for PT100 input
	e coefficient	0	Comp-coefficient for current input 1
	f coefficient	0	Comp-coefficient for current input 2
	g coefficient	0	Comp-coefficient for loading
	Adjust factor :	1	Factor
	Adjust offset :	0	Offset
	Massflow calculation mode:	0	Throughput calculation: see code list
	Loading comp. Selection:	0	*
	Loading comp. Lower limit:	0	*
Synchroni- zation	Synchronizer mode:	0	*
	Current input1 distance to uWave	0	*
	Current input1 distance to uWave	0	*
Calibrate	Calibration input selection :	0	Exp. Cal. input selection: see code list
2	Lower limit :		
	Upper limit :		
	Calibration mode :	0	*
	Calibration variable :	0	Calibration basis: see code list
	Phase coefficients :	0	Phase coefficient A
	Attenuation coefficients :	0	Attenuation Coefficient B
	Constant coefficient :	0	Constant C
	d coefficient	0	Comp-coefficient for PT100 input
	e coefficient	0	Comp-coefficient for current input 1
	f coefficient	0	Comp-coefficient for current input 2
	g coefficient	0	Comp-coefficient for loading
	Adjust factor :	1	Factor
	Adjust offset :	0	Offset
	Massflow calculation mode:	0	Throughput calculation: see code list
	Loading comp. Selection:	0	*
	Loading comp. Lower limit:	0	*
Advanced	Tare Phase (°/GHz) :	0.00 °/GHz	
	Tare Attenuation (dB):	0.00 dB	
	Measure configuration :	0	Process type: see code list
	Range split value :	75	Split value
Current	AO Assign code :	0	Assignment: see code list
output 1	AO Upper range value :	95	Upper value
	AO Lower range value :	60	Lower limit
	AO Alarm select code :	2	Error current: see code list
	AO Error current value :	22.00 mA	Error current value
Current	AO Assign code[2]	0	Assignment: see code list
output 2	AO Upper range value[2]	95	Upper value
	AO Lower range value[2]	60	Lower limit
	Range selection[2]	1	Range
	AO Alarm select code[2]	2	Fror current: see code list
	AO Error current value[2]	- 22 00 m∆	Error current value

Chapter 11. Start-up Protocol



Current			Enabled:
output 1	AI Enabled :	0	Disabled: 0 1
	AI Range selection :	1	Range: see code list
	AI Upper range value :	100	Upper value
	AI Lower range value :	0	Lower limit
	Analog input filter constant :	5	*
Current		0	Enabled:
output 2		0	
	AI Range selection[2] :	1	Range: see code list
	AI Upper range value[2] :	100	Upper value
	AI Lower range value[2] :	0	Lower limit
	Analog input filter constant :	5	*
PT100 in-	AL Enchlod [2] -	0	Enabled:
P	AI Ellabled[5] .	0	
Relay 1	DO Function :	1	Function: see code list
	DO Assignment :	0	Assignment: see code list
	DO Threshold :	0.00%	*
	DO Hysteresis :	5.00%	*
Relay 2	DO Function[2] :	2	Function: see code list
	DO Assignment[2] :	0	Assignment: see code list
	DO Threshold[2] :	0.00%	*
	DO Hysteresis[2] :	5.00%	*
Digital In- put	DI Function selection :	0	Function digital input 1: see code list
	DI Function selection[2] :	0	Function digital input 2: see code list
	DI Function selection[3] :	0	Function digital input 3: see code list
	Printout mode :	1	Data output: see code list
	Access level :	2	Access level: see code list
	Language :	1	Language: see code list
	End of Setup		End

Start of Reference Data		Syste	m	adjustme	nt data:									
Product 1:														
Mean Atten.:		46.8509 dB												
Phase at fm:		42.6285 deg/GHz												
Phase slope:		380,984 deg/GHz												
Phase sigma:		0.24575	0.24575											
Frequency[GHz]	ency[GHz]		Transformed Phase[Deg]	Atten.[dB]										
	2.42	35.64	35.64	21.98										
	2.43	361.81	361.81	21.95										
	2.44	689.04	689.04 689.04											
	2.45	1014.44	1014.44	22.36										
	2.46	1339.01	1339.01	22.37										
Start of Sample Data:							Samp	lin	ig:					
Product 1: Sample Data for	r Concent	ration 1:												
Sample:	Active:	Con.(%):	Lab.(%):	AIN1:	AIN2:		Temp. (°C):		Phi. (°/GHz):	Att.(dB):				
1 17.08 - 12:37	TRUE	85	40	C		0		0	-0.35	0.0)2			
2 17.08 - 12:37	TRUE	80	35	C		0		0	30.33	5.0	08			
3 17.08 - 12:45	TRUE	70	25	C		0		0	59.02	18.98				
Correlation factor between														
lab and meas values:		1												
End of Sample Data														
Do not use following data!														



11.2 Sampling

No.	Ac- tive	Measured value	Lab value	Current in 1	Current In 2	PT100	Phi(fm):	Attenua- tion	
1									
2									
3									
4									
5									
6									
7									
8									
9									
10									
11									
12									
13									
14									
15									
16									
17									
18									
19									
20									



Notes



Notes



		 											_				
	 _	 					 					 	 	 _			
	 _	 					 					 _		 _			
		 											 _				
	_	 							_				 _				
													_				_
	_	 					 					 		 _			
		 							_				 _				
		 							_				 _				
	 	 			 		 					 		 _			
		 							_				 _	_			
		 			 		 		_								
																\neg	
																\uparrow	
																\neg	
																\neg	
																\neg	
																\neg	
																\neg	