



Detectors Concentration Meters
MicroPolar (++)
LB 566

Konzentrations-Messsysteme
MicroPolar (++)
LB 566

Technical Information
Technische Information

41986TI
Rev. No.: 00, 06/2018

Technical Data

Technische Daten

General specifications <i>Allgemeine Spezifikationen</i>	
Method <i>Verfahren</i>	Microwave transmission measurement <i>Mikrowellen-Transmissionsmessung</i>
Operating frequency <i>Arbeitsfrequenz</i>	2.4 - 2.5 GHz (ISM band), depending on local regulations <i>2.4 – 2.5 GHz (ISM-Band), je nach Landesbestimmung</i>
Transmission power <i>Sendeleistung</i>	MicroPolar: < 0.1 mW (< - 10 dBm) MicroPolar ++: < 10 mW (< 10 dBm) All coaxial line power <i>MicroPolar : < 0.1 mW (< - 10 dBm)</i> <i>MicroPolar ++: < 10 mW (< 10 dBm)</i> <i>Alles koaxiale Leitungsleistung</i>
Applications <i>Applikationen</i>	Concentration measurement in containers and pipes <i>Konzentrationsmessung in Behältern und Rohren</i>

Evaluation Unit

Auswertereinheit

Technical Data Evaluation Unit <i>Technische Daten Auswertereinheit</i>	
Housing <i>Gehäuse</i>	Wall housing made of stainless steel, see dimensional drawing MicroPolar: HxWxD: 300 x 323 x 140 mm MicroPolar ++: HxWxD: 400 x 338 x 170 mm <i>Wandgehäuse aus Edelstahl, siehe Maßbild.</i> <i>MicroPolar: HxBxT: 300 x 323 x 140 mm</i> <i>MicroPolar ++: HxBxT: 400 x 338 x 170 mm</i>
Protection type <i>Schutzart</i>	IP 65 <i>IP 65</i>
Weight <i>Gewicht</i>	MicroPolar: about 6.5 kg MicroPolar ++: about 8.0 kg <i>MicroPolar: ca. 6.5 kg</i> <i>MicroPolar ++: ca. 8.0 kg</i>
Ambient conditions during operation <i>Umgebungsbedingungen im Betrieb</i>	-20 ... +60 °C (253 K ...333 K), no condensation Relative humidity: max. 80 % Altitude: max. 2000 m <i>-20 ... +60 °C (253...333 K), keine Betauung</i> <i>Relative Luftfeuchte: max. 80 %</i> <i>Höhenlage: max. 2000 m</i>
Ambient conditions during storage <i>Umgebungsbedingung bei Lagerung</i>	-20 ... +70 °C (253 K ...343 K), no condensation Relative humidity: max. 80 % -20 ... +70 °C (253...343 K), keine Betauung, Relative Luftfeuchte: max. 80 %
Achievable accuracy <i>Erreichbare Genauigkeit</i>	≤ 0.2 weight % (standard deviation) depending on product and sensor ≤ 0.2 Gewichts % (Standardabweichung) je nach Produkt und Sensor

Display <i>Anzeige</i>	Dot matrix LC display, 114 mm x 64 mm, 240 x 128 pixels, with back-lighting, automatic contrast setting <i>Dot-Matrix LC Display, 114 mm x 64 mm, 240 x 128 Pixel, mit Hintergrundbeleuchtung, automatische Kontrasteinstellung</i>
Keyboard <i>Tastatur</i>	Freely accessible foil keypad, light-stable and weatherproof: alphanumeric keyboard and 4 softkeys (software-assigned buttons) <i>Freizugängliche Folientastatur, lichtstabil und witterungsfest: alphanumerische Tastatur und 4 Softkeys (software zugeordneten Tasten)</i>
Power supply <i>Hilfsenergie</i>	For MicroPolar, depending on device type: 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: 1.) 100 ... 240 V AC, 45 ... 65 Hz 2.) 24 V DC: 18 ... 36 V, no reverse polarity protection <i>Für MicroPolar, je nach Geräteausführung:</i> 1.) 100 ... 240 V AC, 45 ... 65 Hz 2.) 24 V DC: 18 ... 36 V; 24 V AC: -20%, +5%, 40 ... 440 Hz <i>Für MicroPolar ++, je nach Geräteausführung:</i> 1.) 100 ... 240 V AC, 45 ... 65 Hz 2.) 24 V DC: 18 ... 36 V, kein Verpolungsschutz
Power consumption <i>Leistungsaufnahme</i>	For MicroPolar: max. 30 VA (AC/DC), depending on configuration For MicroPolar ++: max. (48/60) VA (AC/DC), depending on configuration <i>Für MicroPolar: max. 30 VA (AC/DC), je nach Konfiguration</i> <i>Für MicroPolar ++: max. (48/60) VA (AC/DC), je nach Konfiguration</i>
Fuses <i>Sicherungen</i>	For MicroPolar: 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 <i>Für MicroPolar: 2 x 2.0 A / Tr.</i> <i>Für MicroPolar ++: 2 x 2.0 A / Tr. bei 100 ... 240 V AC oder 2 x 6.3 A / Tr. bei 24 V DC</i>
Battery type <i>Batterietyp</i>	3 V Lithium button cell, type CR2032 <i>3 V Lithium Knopfzelle, Typ CR2032</i>
Measured value <i>Messwert</i>	e.g. concentration, dry content <i>Z.B. Konzentration, Trockengehalt</i>

Inputs and outputs Evaluation Unit *Eingänge und Ausgänge Auswertereinheit*

Cable cross-section <i>Leiterquerschnitt</i>	min. 1.0 mm ² (mains supply) <i>min. 1.0 mm² (Netzanschluss)</i>
Cable feed-through <i>Kabeleinführung</i>	2 x M20x1.5 for cable 5...14 mm (depending on application) 4 x M16x1.5 for cable 5 ...8 mm (depending on application) <i>2 x M20x1.5 für Kabel 5...14 mm (je nach Einsatz)</i> <i>4 x M16x1.5 für Kabel 5...8 mm (je nach Einsatz)</i>
Sensor connection <i>Sensoranschluss</i>	Inputs and outputs for signal and reference channel, 50 Ω N-socket <i>Ein- und Ausgänge für Mess- und Referenzkanal, 50 Ω N-Buchse</i>
HF cable <i>HF-Kabel</i>	Cable lengths: 2, 4, 6 and 10 m; 50 Ω; both sides with 4 N connectors <i>Kabellängen: 2, 4, 6 und 10 m; 50 Ω; beidseitig mit 4 N-Stecker</i>
Current input <i>Stromeingang</i>	2 x current input 0/4 ...20 mA, ohmic resistance 50 Ω, 1x insulated, 1x instrument ground e.g. for temperature compensation <i>2 x Stromeingang 0/4 ...20 mA, Bürde 50 Ω, 1x isoliert, 1x Gerätemasse Z.B. für Temperaturkompensation</i>
Current output <i>Stromausgang</i>	Current output 1: 4...20 mA, ohmic resistance max. 800 Ω, insulated Current output 2: 0/4...20 mA, ohmic resistance max. 800 Ω, insulated e.g. for result or temperature output <i>Stromausgang 1: 4...20 mA, Bürde max. 800 Ω, isoliert</i> <i>Stromausgang 2: 0/4...20 mA, Bürde max. 800 Ω, isoliert Z.B. zur Messwert- oder Temperaturausgabe</i>
PT100 connection <i>PT100 Anschluss</i>	Measuring range: -50 ... +200 °C (223 ... 473 K) measurement tolerance: < 0.4 °C <i>Messbereich: -50 ... +200 °C (223 ... 473 K)</i> <i>Messtoleranz: < 0.4 °C</i>
Digital input <i>Digitaler Eingang</i>	3 x digital inputs (DI1..3), for floating connectors (Do not connect to a power supply). Configuration options: DI1: none, measurement start/stop DI2: none, measurement hold, product selection DI3: none, sampling, product selection Function description: 1. Measurement (Start/Stop), <u>open</u> : Measurement stopped, <u>closed</u> : Measurement started and/or measurement running 2. 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 4. Start sampling: <u>open</u> : no actions, <u>closed</u> : single measurement starts <i>3 x digitale Eingänge (DI1..3), für potential-freie Kontakte (nicht an einer Spannungsversorgung anschließen).</i> <i>Konfigurationsmöglichkeiten:</i> <i>DI1: Keine, Messung Start/Stopp</i> <i>DI2: Keine, Messung halten. Produktauswahl</i>

	DI3: Keine, Probeaufnahme, Produktauswahl Funktionsbeschreibung: 1. Messung (Start/Stopp), <u>offen</u>: Messung gestoppt, <u>geschlossen</u>: Messung gestartet bzw. Messung läuft 2. Messung anhalten, <u>offen</u>: Messung läuft, <u>geschlossen</u>: Messung angehalten, d.h. Mittelwerte und Stromausgang sind eingefroren 3. Produktauswahl über ein DI: <u>offen</u>: Produkt 1 (P1), <u>geschlossen</u>: P2 Produktauswahl über zwei DI's: DI2 & DI3 <u>offen</u>: P1 DI2 <u>geschlossen</u> & DI3 <u>offen</u>: P2 DI2 <u>offen</u> & DI3 <u>geschlossen</u>: P3 DI2 & DI3 <u>geschlossen</u>: P4 4. Probeaufnahme: <u>offen</u>: keine Aktionen, <u>geschlossen</u>: einmalige Aufnahme startet
Relay outputs	2 x relays (SPDT), insulated <u>Configuration options:</u> - Collective failure message - Stop measurement - Limit value (min. and max.) - No product <u>Load capacity:</u> 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 correspond to a mains cable. <u>Restrictions at 24 V AC/DC (DC: 18 ...36 V; AC: 24 V +5 %, -20 %) mains supply, if the ground conductor is not connected to terminal 1 (PE):</u> AC: max. 50 V DC: max. 70 V
Relaisausgänge	2 x Relais (SPDT), isoliert <u>Konfigurationsmöglichkeiten:</u> - Sammelstörmeldung - Messung anhalten - Grenzwert (min. und max.) - kein Produkt <u>Belastbarkeit:</u> AC: max. 400VA DC: max. 90W AC / DC: max. 250V, max. 2A, induktionsfrei ≥ 150V: Spannung muss erdbezogen sein Das an den Relaisausgang verwendete Kabel muss einer Netzanschlussleitung entsprechen. <u>Einschränkungen bei 24 V AC/DC (DC: 18 ...36 V; AC: 24 V +5 %, -20 %) Netzversorgung, wenn der Schutzleiter nicht an die Klemme 1 (PE) angeschlossen ist:</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
Serielle Schnittstellen	RS232 auf der Geräteunterseite, RS485 über Klemmenleiste Datenformat: 38400 Bd, kein Handshake, 8 Datenbits, 1 Stoppbit, keine Parität

Sensors

Sensoren

Technical Data Flow Cells <i>Technische Daten Flow Cell</i>	
Application <i>Anwendung</i>	Microwave FlowCell with various nominal diameters and flanges for concentration measurement on pipelines <i>Mikrowellen-FlowCell verschiedener Nennweiten und Flansche zur Konzentrationsmessung an Rohrleitungen</i>
Material <i>Material</i>	Inline housing made of stainless steel 1.4404 polished (AISI 316L) Microwave windows made of PEEK Product touching sealing made of EPDM <i>Inline-Gehäuse aus Edelstahl 1.4404 poliert (AISI 316L)</i> <i>Mikrowellenfenster aus PEEK</i> <i>Produktberührende Dichtung aus EPDM</i>
Process coupling <i>Prozessankopplung</i>	Two 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 Optional adapter for the V flange version with ASA flange <i>Zwei Varianten:</i> <i>1. Hygiene Milchrohrverschraubung DIN 11853-1</i> <i>EHEDG-Zulassung</i> <i>2. Flansch nach EN 1092-1/11 (V-Flansch)</i> <i>FDA-zugelassene Materialien</i> <i>Optionale Adapter für die V-Flansch Variante mit ASA-Flansch</i>
Process pressure <i>Prozessdruck</i>	up to 16 bar (relative) <i>Bis 16 bar (relativ), je nach Nennweite und Flanschtyp, siehe Tabelle unten</i>
Temperature range <i>Temperaturbereich</i>	Product temperature: 10...130 °C (283...403 K), temporarily up to 140 Ambient temperature: -20...60 °C (253...333 K) Storage temperature: 10...80 °C (283...353 K) <i>Produkttemperatur: 10...130 °C (283...403 K), kurzzeitig bis 140°C (413 K)</i> <i>Umgebungstemperatur: -20...60 °C (253...333 K)</i> <i>Lagertemperatur: 10...80 °C (283...353 K)</i>
Connections <i>Anschlüsse</i>	2 x HF connections: N female, 50 Ω for HF cable with max. 10 m length <i>2 x HF-Anschlüsse: N-Buchse, 50 Ω</i> <i>Für HF-Kabel mit max. 10 m Länge</i>
Versions <i>Varainten</i>	Nominal pipe widths from 50 ... 150 mm <i>Rohrnennweiten von 50 ... 150 mm</i>
Dimensions <i>Maße</i>	See dimensional drawings. <i>Siehe Maßzeichnungen</i>

Overview FlowCells with V flange Übersicht FlowCell mit V-Flansch

Designation Bezeichnung	ID no. Id.-Nr.	Nominal width [mm] Nennweite [mm]	Flange Flansch	Pressure [bar] Druck [bar]
LB 5660-102-00x	66744-001	50	DN 50 / PN 16	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	

Overview FlowCells with Hygiene milk pipe screw connection Übersicht FlowCell mit Milchrohrverschraubung

Designation Bezeichnung	ID no. Id.-Nr.	Nominal width [mm] Nennweite [mm]	Flange Flansch	Pressure [bar] Druck [bar]
LB 5660-112-00x	66744-013	50	LB 5660-112-00x	16
LB 5660-212-00x	66744-014	65	LB 5660-212-00x	
LB 5660-312-00x	66744-015	80	LB 5660-312-00x	
LB 5660-412-00x	66744-016	100	LB 5660-412-00x	
LB 5660-512-00x	66744-017	125	LB 5660-512-00x	
LB 5660-612-00x	66744-018	150	LB 5660-612-00x	

Overview Microwave windows set Übersicht Mikrowellenfenster-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.

Die Mikrowellenfenster der FlowCell sind für Anwendungen mit ho-hem Abrieb auch in einer glasfaserverstärkten Ausführung erhältlich. Diese ist nicht für den Lebensmittelbereich zugelassen.

ID no. Id.-Nr.	Designation Bezeichnung
66624-S	2 pieces of PEEK Microwave windows with 6 seal-ing O-rings 2 Stück PEEK Mikrowellenfenster mit 6 Dichtungs-O-Ringen
66625-S	2 pieces of PEEK Microwave windows GF30 glass fibre reinforced with 6 sealing O-rings 2 Stück PEEK Mikrowellenfenster GF30 glasfaserverstärkt mit 6 Dichtungs-O-Ringen

Overview ASA flange adapter Übersicht ASA-Flansch-Adapter

Designation Bezeichnung	ID no. Id.-Nr.
ASA flange adapter set for Flow Cell 50 <i>ASA-Flansch-Adapter-Set für Flowcell 50</i>	62324
ASA flange adapter set for Flow Cell 65 <i>ASA-Flansch-Adapter-Set für Flowcell 65</i>	62319
ASA flange adapter set for Flow Cell 80 <i>ASA-Flansch-Adapter-Set für Flowcell 80</i>	62319
ASA flange adapter set for Flow Cell 100 <i>ASA-Flansch-Adapter-Set für Flowcell 100</i>	62331
ASA flange adapter set for Flow Cell 150 <i>ASA-Flansch-Adapter-Set für Flowcell 150</i>	62335
The kit consists of two adapters, screws and two seals. <i>Das Set besteht aus zwei Adaptern, Schrauben und zwei Dichtungen.</i>	

Overview Inline housing, FDA Übersicht Inline-Gehäuse, FDA

for temperature or conductivity sensors or sampling valve
 für Temperatur-, Leitfähigkeitssensor oder Probenahmeventil

Designation Bezeichnung	VFL Id.-Nr.	G-BS/M Id.-Nr.
Inline housing for Flowcell 50 <i>Inline-Gehäuse für Flowcell 50</i>	67078	67084
Inline housing for Flowcell 65 <i>Inline-Gehäuse für Flowcell 65</i>	67079	67085
Inline housing for Flowcell 80 <i>Inline-Gehäuse für Flowcell 80</i>	67080	67086
Inline housing for Flowcell 100 <i>Inline-Gehäuse für Flowcell 100</i>	67081	67087
Inline housing for Flowcell 125 <i>Inline-Gehäuse für Flowcell 125</i>	67082	67088
Inline housing for Flowcell 150 <i>Inline-Gehäuse für Flowcell 150</i>	67083	67089

Overview surface temperature sensor Übersicht Oberflächen-Temperaturfühler

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

Selbstklebender PT100 Temperaturfühler mit Befestigungsmaterial, Anschlusskabel 10 m, 4-Leiter (lose Enden) Temperaturbereich: -50 - +200 °C.

Designation Bezeichnung	ID no. Id.-Nr.
Self-adhesive temperature sensor for DN 50 <i>Selbstklebender Temperaturfühler für DN 50</i>	66655
Self-adhesive temperature sensor for DN 65 <i>Selbstklebender Temperaturfühler für DN 65</i>	66656
Self-adhesive temperature sensor for DN 80 <i>Selbstklebender Temperaturfühler für DN 80</i>	66657
Self-adhesive temperature sensor for DN 100 <i>Selbstklebender Temperaturfühler für DN 100</i>	66658
Self-adhesive temperature sensor for DN 125 <i>Selbstklebender Temperaturfühler für DN 125</i>	66659
Self-adhesive temperature sensor for DN 150 <i>Selbstklebender Temperaturfühler für DN 150</i>	66660

Overview sensors Übersicht Sensoren

Designation Bezeichnung	ID no. Id.-Nr.
Conductivity sensor hygienic, Clamp-flange 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 <i>Leitfähigkeitssensor hygienisch, Clamp-Flansch</i> <i>Induktives Leitfähigkeitsmessgerät für flüssige Medien in hygienischen Anwendungen</i> <i>Messbereich: 0-999 mS/cm</i> <i>Prozessanschluss: Clamp-Flansch</i> <i>Betriebsdruck: max. 16 bar</i> <i>Hilfsspannung: 18-36 V DC max. 190 mA</i> <i>Ausgang: 4-20 mA</i>	66693
Temperature sensor EHEDG, Clamp-flange PT100 temperature sensor for hygienic applications Measurement range: -50 - +250 °C Length 20 mm, diameter 4 mm Process connection: Clamp-flange With connection cable 10 m (loose ends)	66694

<i>Temperatursensor EHEDG, Clamp-Flansch</i> <i>PT100 Temperatursensor für hygienische Anwendungen</i> <i>Messbereich: -50 - +250 °C</i> <i>Länge 20 mm, Durchmesser 4 mm</i> <i>Prozessanschluss: Clamp-Flansch</i> <i>Mit Anschlusskabel 10 m (lose Enden)</i>	66694
--	-------

Overview Sampling valve and accessories Übersicht Probenahmeventil und Zubehör

Designation Bezeichnung	ID no. Id.-Nr.
Sampling valve aseptic, Clamp-flange Aseptic Inline Sampling valve Stainless steel 1.4404 (AISI 316L), bellows PTFE Discharge port S-DN 10 without flushing connection Process connection: Clamp-flange <i>Probenahmeventil aseptisch, Clamp-Flansch</i> <i>Aseptisches Inline Probenahmeventil</i> <i>Edelstahl 1.4404 (AISI 316L), Faltenbalg PTFE</i> <i>Auslaufstutzen S-DN 10</i> <i>ohne Spülanschluss</i> <i>Prozessanschluss: Clamp-Flansch</i>	66738 66738
Clamp-blind flange Clamp-blind flange for Inline housing DN 50-150 Stainless steel 1.4306 (AISI 304L) <i>Clamp-Blindflansch</i> <i>Clamp-Blindflansch für Inline-Gehäuse DN 50-150</i> <i>Edelstahl 1.4306 (AISI 304L)</i>	66737 66737
Clamp coupling 1 piece Clamp coupling for Inline housing DN 50-150 Stainless steel 1.4306 (AISI 304L) Clamp-Klammer 1 Stück Clamp-Klammer für Inline-Gehäuse DN 50-150 Edelstahl 1.4306 (AISI 304L)	66736 66736

Container probes
Behältersonden

Application <i>Anwendung</i>	Container probes with and without flushing device for concentration measurement in process containers and pipelines with nominal width ≥ 200 mm. <i>Behältersonden mit und ohne Spülvorrichtung zur Konzentrationsbestimmung in Prozessbehältern und in Rohrleitungen mit NW ≥ 200 mm.</i>
Material <i>Material</i>	Plastic caps, stainless steel PT100 connection cable: Silicon / Teflon <i>Kunststoffkappen, Edelstahl</i> <i>PT100 Anschlusskabel: Silicon / Teflon</i>
Process coupling <i>Prozessankopplung</i>	Flange according to DIN EN 1092 type 05 DN65 / PN6, DN 80, 100, 150 / PN16; ASA flange 2.5" / 150 PSI (More more on request) <i>Flansch nach DIN EN 1092 Typ 05 DN65 / PN6, DN 80, 100, 150 / PN16; ASA Flansch 2.5" / 150 PSI (weitere auf Anfrage)</i>
Process pressure <i>Prozessdruck</i>	Up to 16 bar, depending on model <i>Bis 16 bar, je nach Ausführung</i>
Temperature range <i>Temperaturbereich</i>	Product temperature: 10...120 °C (283...393 K) Ambient temperature: -20...60 °C (253...333 K) Storage temperature: 10...80 °C (283...353 K) <i>Produkttemperatur: 10...120 °C (283...393 K)</i> <i>Umgebungstemperatur: -20...60 °C (253...333 K)</i> <i>Lagertemperatur: 10...80 °C (283...353 K)</i>
Connections <i>Anschlüsse</i>	4 x HF connections: N female, 50 Ω for HF cable with max. 10 m length <i>4 x HF-Anschlüsse: N-Buchse, 50 Ω</i> <i>Für HF-Kabel mit max. 10 m Länge</i>
Dimensions <i>Maße</i>	See dimensional drawings. <i>Siehe Maßzeichnungen.</i>

Accessory sealing washer
Zubehör Dichtungsscheibe

Material <i>Material</i>	Klingersil C-4400
Thickness <i>Dicke</i>	3 mm

Caps set for container probe
Kappenset für Behältersonde

Id.-Nr. 66049-S	2 PEEK Plastic caps with 4 sealing rings <i>2 Stück PEEK Kunststoffkappen mit 4 Dichtungsringen</i>
-----------------	--

Overview container probes and sealing washers Übersicht Behältersonden und Dichtungsscheiben

Designation <i>Bezeichnung</i>	ID no. <i>Id.-Nr.</i>	Flange <i>Flansch</i>	Pressure [bar] <i>Druck [bar]</i>	ID no. seals <i>Id.-Nr. Dichtungen</i>
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

Technical Data HF Cable Technische Daten HF Kabel

HF cable Quad <i>HF-Kabel Quad</i>	
Material	Corrugated tube: Polyamide (PA6) Cable sheath: Polyethylene (PE)
<i>Material</i>	<i>Wellrohr: Polyamid (PA6)</i> <i>Kabelmantel: Polyethylen (PE)</i>
Protection type <i>Schutzart</i>	IP 66 <i>IP 66</i>
Temperature <i>Temperatur</i>	In operation: -30 ... +70 °C When installing: -20 ... +70 °C <i>Bei Betrieb: -30 ... +70 °C</i> <i>Bei Installation: -20 ... +70 °C</i>
Cable length [m] <i>Kabellänge [m]</i>	ID no. <i>Ident.-Nr.</i>
2.0	43431
4.0	43432
6.0	43433
8.0	43434
10.0	43435

HF-Kabel Quad, hygienic <i>HF-Kabel Quad, hygienisch</i>	
Material	Corrugated tube: Polyamide (PA6) Cable sheath: Polyethylene (PE)
<i>Material</i>	<i>Wellrohr: Polyamid (PA6)</i> <i>Kabelmantel: Polyethylen (PE)</i>
Protection type	IP 66

<i>Schutzart</i>	<i>IP 66</i>
Temperature <i>Temperatur</i>	In operation: -30 ... +70 °C When installing: -20 ... +70 °C <i>Bei Betrieb:</i> -30 ... +70 °C <i>Bei Installation:</i> -20 ... +70 °C
Cable length [m] <i>Kabellänge [m]</i>	ID no. <i>Ident.-Nr.</i>
2.0	67048
4.0	67049

HF-Kabel Quad, solid cable
HF-Kabel, Festmantelkabel

Material <i>Material</i>	Cable sheath: Polyethylene (PE) <i>Kabelmantel: Polyethylen (PE)</i>
Protection type <i>Schutzart</i>	IP 68 when unscrewed <i>IP 68 im aufgeschraubten Zustand</i>
Temperature <i>Temperatur</i>	In operation: -40 ... +85 °C When installing: -40 ... +85 °C <i>Bei Betrieb:</i> -40 ... +85 °C <i>Bei Installation:</i> -40 ... +85 °C
Cable length [m] <i>Kabellänge [m]</i>	ID no. <i>Ident.-Nr.</i>
2.0	11476
2.5	11477
3.0	11478
3.5	11479
4.0	11480

Format of Serial Data Output RS232 and RS485

Headline

Date·Time→State→Status→Synchronizer→Product→Att→Phi→R2→Tint→IN1→IN2→PT100→
 C→Cm→C2→C2m→MF1→MF2

Following lines

01.01.2005-00:00:00→0000→0→0→1→0.43→5.30→0.07→0.0→0.0→0.0→0.0→
 1 2 3 4 5 6 7 8 9 10 11 12

75.36→75.00→0.00→0.00→0.000→0.000¶
 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

“→” Tabulation

“¶” Carriage return + Line feed

“.” Spaces

Format der seriellen Datenausgabe RS232 und RS485

Titelzeile

Date·Time→State→Status→Synchronizer→Product→Att→Phi→R2→Tint→IN1→IN2→PT100→
C→Cm→C2→C2m→MF1→MF2

¶

Folgezeilen

01.01.2005·00:00:00→0000→0→0→1→0.43→5.30→0.07→0.0→0.0→0.0→0.0→
1 2 3 4 5 6 7 8 9 10 11 12

75.36→75.00→0.00→0.00→0.000→0.000¶
13 14 15 16 17 18

Spalte Nr.	Beschreibung	Format
1	Datum und Uhrzeit	TT.MM.JJ·HH:MM:SS
2	State	4 –Stellig, HEX
3	Status: Information über die Qualität der letzten Messung	0 : Messung OK < 0 : Fehler
4	Produktsynchronisation	5: nicht aktiv 1: noch unsynchron 0: alle Werte synchron -1: Fehler -2: Zeit zu klein für syn. -3: Geschw. außer Ber.
5	Produktnummer	X (1 bis 4)
6	Dämpfung [dB]	X.XX
7	Phase [°/GHz]	X.XX
8	Streuung der Phasenregression	X.XX
9	Gerätetemperatur [Temperatureinheit]	X.X
10	Stromeingang 1 [Einheit des Stromeingangs]	X.X
11	Stromeingang 2 [Einheit des Stromeingangs]	X.X
12	PT100 Temperatur [Temperatureinheit]	X.X
		[...] mit Auswahl der Einheit g/cm
13	Konzentration 1 live	X.XX [X.XXXX]
14	Konzentration 1 gemittelt	X.XX [X.XXXX]
15	Konzentration 2 live	X.XX [X.XXXX]
16	Konzentration 2 gemittelt	X.XX [X.XXXX]
17	Massenstrom für Konzentration 1	X.XXX
18	Massenstrom für Konzentration 2	X.XXX

Sonderzeichen

“→” Tabulation

“¶” Carriage return + Line feed

“.” Leerzeichen

EC Declaration of Conformity



BERTHOLD TECHNOLOGIES GmbH & Co.KG

Calmbacher Str. 22
75323 Bad Wildbad, Germany

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

EC – Declaration of Conformity

We herewith confirm 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 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/EEG (electromagnetic compatibility)
reviewed: 91/263/EEG, 92/31/EEG, 93/68/EEG, 93/97/EEG

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

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

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
EN 61000-4-11:1994 + A1:2001

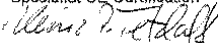
This declaration is issued by the manufacturer:

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

by


Dr. J. Briggmann
Development Manager
Process Control

H. Tetzlaff
Specialist CE Certification



Bad Wildbad, 04.05.2004

Registergericht / Court of Registration
Persönlich haftende Gesellschafterin / Fully liable Associates
Registergericht / Court of Registration
Geschäftsführung / Management
Beiratsvorsitzender / Chairman of the Board
USt.-IdNr. / VAT Reg.No
Deutsche Steuernummer / German Tax No

Calw HRA 991
BERTHOLD TECHNOLOGIES Verwaltungs-GmbH
Calw HRB 1520
Hans J. Oberhofer (Vors./CEO), Dr. Wilfried Reuter
Dr. Fritz Berthold
DEB13050511
49038/08038

Dresdner Bank 75105 Pforzheim Konto/account no.6 511 120 (BLZ 666 800 13)/ SWIFT-BIC DRES DE FF 666 IBAN: DE05 6668 0013 0651 1120 00
Sparkasse PF-CW 75323 Bad Wildbad Konto/account no.80 450 05 (BLZ 606 510 70)/ SWIFT-BIC PZHSDE66 IBAN: DE76 6065 1070 0008 0450 05

EC Declaration of Conformity (continued)



BERTHOLD TECHNOLOGIES GmbH & Co.KG

Calmbacher Str. 22
75323 Bad Wildbad, Germany

Phone +49 7081 177-0
Fax +49 7081 177-100
info@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 void should alteration or unintended use take place without our authorisation.

Description: **Concentration-Measuring Systems
Micro-Polar ++**

Type: **LB 566-XX**

	EC-Regulation	and Reviews	considered Norms	
EMC	2004/108/EG		EN 55011	1998 +A1:1999 +A2:2002
			EN 61326-1	2006-05
			EN 61000-4-2	1995 +A1:1998 +A2:2001
			EN 61000-4-3	2006-12
			EN 61000-4-4	2004
			EN 61000-4-5	1995 +A1:2001
			EN 61000-4-6	1996 +A1:2001
			EN 61000-4-11	1994-08 +A1:2001-02
			Namur NE21	2004
Low Voltage Directive	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

Dr. Wilfried Reuter
Technical Director
Bad Wildbad, 06.06.2008

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

Stuttgart HRA 330991
BERTHOLD TECHNOLOGIES Verwaltungs-GmbH
Stuttgart HRB 331520
Hans J. Oberhofer, Dr. Wilfried Reuter
Dr. Fritz Berthold
DE813050511
49038/08038
DE99468690

Dresdner Bank 75105 Pforzheim Konto/Account No. 6 511 120 (BLZ 666 800 13) /SWIFT-BIC DRES DEFF 666 IBAN: DE05 6668 0013 0651 1120 00
Sparkasse PF-CW 75323 Bad Wildbad Konto/Account No. 80 450 03 (BLZ 666 500 85) /SWIFT-BIC PZHSDE66 IBAN: DE37 6665 0085 0008 0450 03

EG-Konformitätserklärung



BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Str. 22
75323 Bad Wildbad, Germany

Phone: +49 7081 377 0
Fax: +49 7081 377 100
info@bertholdtech.com
www.BertholdTech.com

EG – Konformitätserklärung

Hiermit erklären wir, daß die Bauart des(r) nachfolgend bezeichneten Gerätes / Systems / Anlage in der von uns in den Verkehr gebrachten Ausführung den unten genannten einschlägigen EG-Richtlinien entspricht.

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

Bezeichnung: **Konzentrations-Messsystem Micro-Polar 1**

Typ: **LB 566-XX**

Einschlägige EG-Richtlinien:

89/336/EWG (Elektromagnetische Verträglichkeit)
geändert: 91/263/EWG, 92/31/EWG, 93/68/EWG, 93/97/EWG

73/23/EWG (Niederspannungsrichtlinie)
geändert: 93/68/EWG

Zur Beurteilung des Erzeugnisses wurden folgende Normen herangezogen:

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
EN 61000-4-11:1994 + A1:2001

Diese Erklärung wird verantwortlich für den Hersteller:

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

abgegeben durch

Dr. J. Briggmann
Entwicklungsleiter
Prozessmesstechnik

H. Tetzlaff
Fachkraft CE Zertifizierung

Bad Wildbad, den 04.05.2004

Reg.-Einsicht / Court of Registration
Personen betreffende Gesellschaften / Publicly traded Companies
Registrierung / Court of Registration
Geschäftsführung / Directors
Beratungsvorstand / Chairman of the Board
USA, China / V.A. Reg. No.
Deutsche Steuernummer / German Tax No.

CE-Verfahren
BERTHOLD Technologies Verwaltungs GmbH
Calb. Str. 22
75323 Bad Wildbad
Deutschland
DE 42 00000
DE 42 00000
49036/08038

Dresdner Bank – 75105 Pforzheim – Konto/account no. 53 1 120 (BIC 2666 100) / SWIFT BIC 2666 100
Sparkasse Pforzheim – 75323 Bad Wildbad – Konto/account no. 20 450 05 (BIC 2666 100) / SWIFT BIC 2666 100
IBAN: DE 05 5056 0012 0001 11 70 00
IBAN: DE 20 6005 0070 0003 0450 05

EG-Konformitätserklärung (Fortsetzung)



BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Str. 22
75323 Bad Wildbad, Germany

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

EG-Konformitätserklärung

Hiermit erklären wir, dass die Bauart des(r) nachfolgend bezeichneten Geräte / Systems / Anlage in der von uns in den Verkehr gebrachten Ausführung den unten genannten einschlägigen EG-Richtlinien entspricht.

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

Produktbezeichnung: **Konzentrations-Messsysteme
Micro-Polar ++**

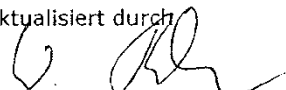
Typenbezeichnung: **LB 566-XX**

Richtlinie und Änderungen		angewendete Normen	
EMV	2004/108/EG	EN 55011	1998
			+A1:1999
			+A2:2002
		EN 61326-1	2006-05
		EN 61000-4-2	1995
			+A1:1998
			+A2:2001
		EN 61000-4-3	2006-12
		EN 61000-4-4	2004
		EN 61000-4-5	1995
			+A1:2001
		EN 61000-4-6	1996
			+A1:2001
Niederspg.	73/23/EWG 93/68/EWG	EN 61010 Teil 1	2002-08
R&TTE	1999/5/EG	ETSI EN 300 440-1	2007-08
		ETSI EN 300 440-2	2007-08

Diese Erklärung wird verantwortlich für den Hersteller

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

aktualisiert durch


Dr. Wilfried Reuter
Technischer Geschäftsführer
Bad Wildbad, den 06.06.2008

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

Stuttgart HRA 330991
BERTHOLD TECHNOLOGIES Verwaltungs-GmbH
Stuttgart HRB 331520
Hans J. Oberhofer, Dr. Wilfried Reuter
Dr. Fritz Berthold
DE91305351:
49038/08038
DE99468690

Dresdner Bank 75105 Pforzheim Konto/Account No. 6 511 120 (BLZ 666 800 13) / SWIFT-BIC DRES DEFF 666 IBAN: DE05 6668 0013 0651 1120 00
Sparkasse PF-CW 75323 Bad Wildbad Konto/Account No. 80 450 03 (BLZ 666 500 85) / SWIFT-BIC PZHSDE66 IBAN: DE37 6665 0065 0006 0450 03

EHEDG Certificate EHEDG Zertifikat

CERTIFICATE OF COMPLIANCE



Valid until end: December, 2018

TUM Certification
hereby declares that the product

microwave flow cell with PEEK windows and EPDM O-ring seals in size DN50 –
DN150

from

Berthold Technologies GmbH & Co. KG, 75323 Bad Wildbad, Germany

has been evaluated for compliance with the
Hygienic Equipment Design Criteria of the EHEDG, by:

TUM (Forschungszentrum für Brau- u. Lebensmittelqualität) at Weihenstephan, Germany
and meets the criteria as demonstrated by:

Evaluation Report No. 596TUM2017

Signed Jürgen Hofmann
Dr. Jürgen Hofmann

Evaluation Officer

Signed Fritz Jacob
Prof. Dr. Fritz Jacob

Head of Department

Date of issue: 13 November 2017

Certificate No. 17/2017



85354 Freising-Weihenstephan, Germany
©EHEDG

Frequency License Frequenzzulassung

CETECOM ICT Services GmbH

EC Identification number 0682

authorized by the German Government



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
Certificate 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 A) 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

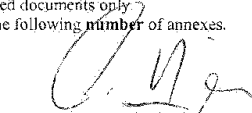
Marking: The product shall be signed with CE and our notified body number as shown right hand.

CE 0682

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

Number of annexes: 1

Saarbrücken, 24.06.2004
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.de>

Frequency License (continued) Frequenzzulassung (Fortsetzung)

CETECOM ICT Services GmbH



CERTIFICATE OF CONFORMITY

Registration-No.: E814059R-CC Number of annexes: ---

Certificate 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
(humidity sensor)

Product Manufacturer: Berthold Technologies GmbH & Co KG
Calmbacher Strasse 22
D-75323 Bad Wildbad

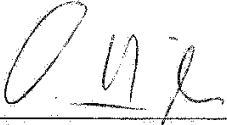
Specifications and test reports:

Specification	Test report no. & date	Name of test laboratory	Notes
EN 55011:1998+A1:1999 (class A) EN 61000-6-2:2001	2003-731-1182-REN dated Sept. 9, 2003	ELMAC GmbH	conform
EN 300 440-1 V1.3.1 (2001-09) EN 300 440-2 V1.1.1 (2001-09)	2-3389-01-01/03 dated May 14, 2004	CETECOM ICT	conform

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

Saarbrücken, 24.06.2004
Place, Date of Issue


Signed by Ernst Hussinger
Notified Body



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

Frequency License (continued) Frequenzzulassung (Fortsetzung)

CETECOM ICT Services GmbH



CERTIFICATE OF CONFORMITY

Registration-No.: E815580V-CC Number of annexes: ---

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
(humidity sensor)

Product Manufacturer: BERTHOLD TECHNOLOGIES GmbH & Co KG
Calmbacher Strasse 22
75323 Bad Wildbad
Germany

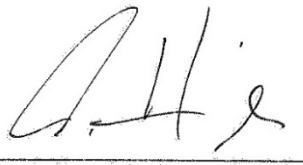
Specifications and test reports:

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

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

Frequency License (continued) Frequenzzulassung (Fortsetzung)

CETECOM ICT Services GmbH EC Identification Number 0682

authorized by the German Government



Bundesnetzagentur

BNNetzA-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 (R&TTE, Article 3.2)	EN 300 440-1 V1.4.1 (2007-08) EN 300 440-2 V1.2.1 (2007-08)	Test Report	conform

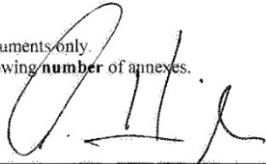
Marking: The product shall be signed with CE and our notified body number as shown right hand.

CE 0682

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

Number of annexes: **1**

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>

Technical Acceptance Certificate

Technisches Abnahmezertifikat

Technical Acceptance Certificate -



Certificate Holder: BERTHOLD TECHNOLOGIES GMBH & CO. KG
CALMBACHER STR. 22
75323 BAD WILDBAD
Germany

ISED Certification Number: 4777A-IC01X01

CTC Registration Number: 1865

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



BNetzA-CAB-03/22-51

authorized by the German
Government to act as CAB
in accordance with the MRA
EU Canada of 1st November
1998.

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.

Saarbrücken

CTC advanced GmbH

cn=Gerald Schmidt, o=CTC
advanced GmbH,
ou=SCT-141216,
email=gerald.schmidt@ctcadvan
ced.com, c=DE
2017.09.08 10:20:11 +02'00'

Certification Body



CTC advanced GmbH (formerly CETECOM ICT Services GmbH)
Untertuerkheimer Str. 66117 | German | www.ctcadvanced.com

Technical Acceptance Certificate (continued) Technisches Abnahmezertifikat (Fortsetzung)

Technical Acceptance Certificate -



Certificate Holder: BERTHOLD TECHNOLOGIES GMBH & CO. KG
CALMBACHER STR. 22
75323 BAD WILDBAD
Germany

ISED Certification Number: 4777A-IC01X12

CTC Registration Number: 1866

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



BNetzA-CAB-03/22-51

authorized by the German
Government to act as CAB
in accordance with the MRA
EU Canada of 1st November
1998.

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

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'

Certification Body

Saarbrücken

CTC advanced GmbH (formerly CETECOM ICT Services GmbH)
Untertuerkheimer Str. 66117 | German | www.ctcadvanced.com

Grant of equipment authorization

TCB

**GRANT OF EQUIPMENT
AUTHORIZATION**

TCB

Certification

Issued Under the Authority of the
Federal Communications Commission

By:

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

Date of Grant: 09/08/2017

Application Dated: 09/08/2017

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.

FCC IDENTIFIER: R9ZFCC01X01
Name of Grantee: Berthold Technologies
Equipment Class: Part 15 Field Disturbance Sensor
Notes: Concentration, Dry Mass, Water, Density
measuring system

Grant Notes

FCC Rule Parts

15.245

Frequency
Range (MHZ)

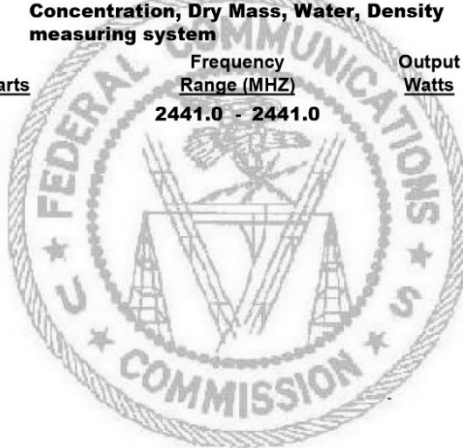
2441.0 - 2441.0

Output
Watts

Frequency
Tolerance

Emission
Designator

PC II to add additional antennas.



Grant of equipment authorization (continued)

TCB

**GRANT OF EQUIPMENT
AUTHORIZATION**

TCB

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

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

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

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.

FCC IDENTIFIER: **R9ZFCC01X12**
Name of Grantee: **Berthold Technologies**
Equipment Class: **Part 15 Field Disturbance Sensor**
Notes: **Concentration, Dry Mass, Water, Density measuring system**

Grant Notes

FCC Rule Parts
15.245

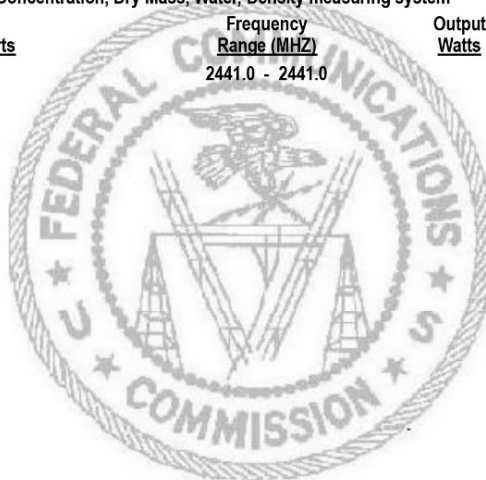
Frequency
Range (MHZ)
2441.0 - 2441.0

Output
Watts

Frequency
Tolerance

Emission
Designator

PC II to add additional antennas.



Electrical Wiring Diagram Elektrischer Anschlussplan

Line in for MicroPolar :

1. / 2. depending on configuration

1.
AC 100-240V, 45-65 Hz

2.
DC 24 V (18-36 V) or
AC 24 V -20/+5%, 40-440 Hz

Line in for MicroPolar ++:

1. / 2. depending on configuration

1.
AC 100-240V, 45-65 Hz

2.
DC 24 V (18-36 V), no reverse polarity protection

Line in für MicroPolar :

1. / 2. je nach Geräteausführung

1.
AC 100-240V, 45-65 Hz

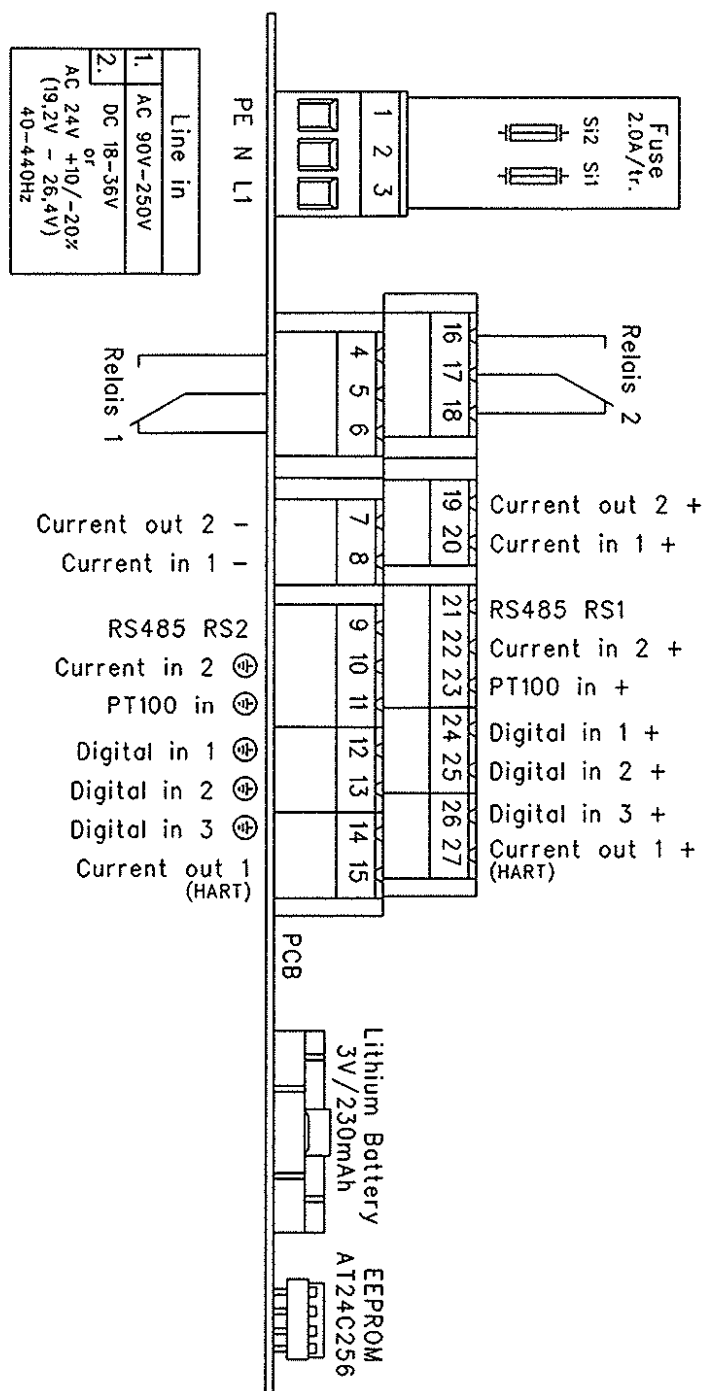
2.
DC 24 V (18-36 V) oder
AC 24 V -20/+5%, 40-440 Hz

Line in für MicroPolar ++:

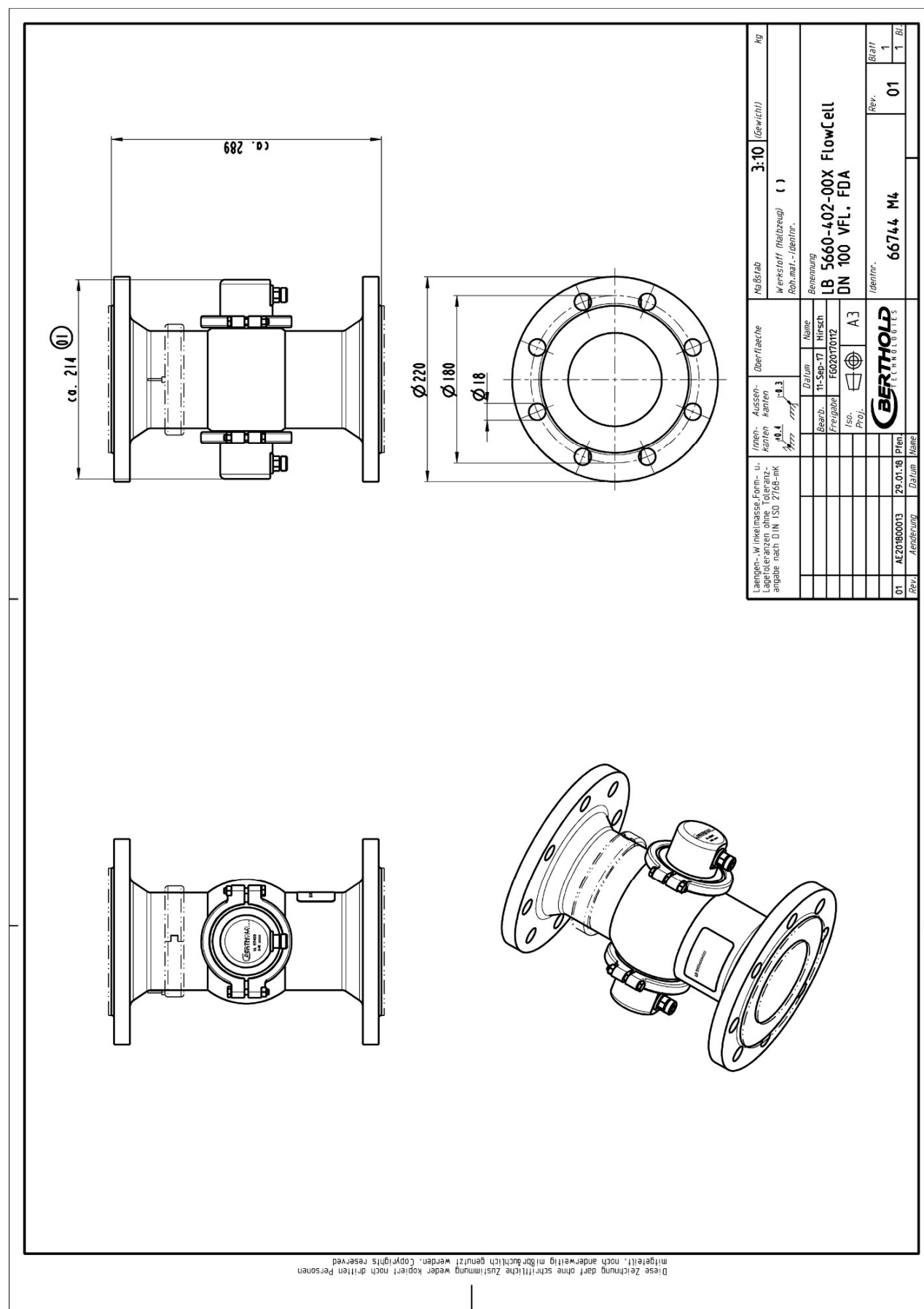
1. / 2. je nach Geräteausführung

1.
AC 100-240V, 45-65 Hz

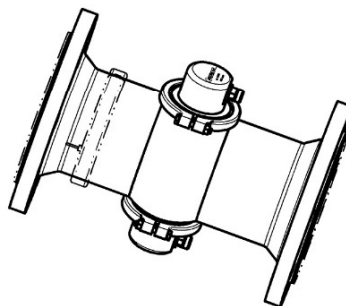
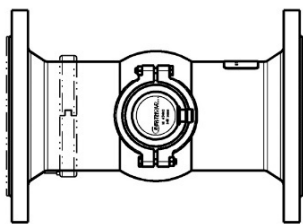
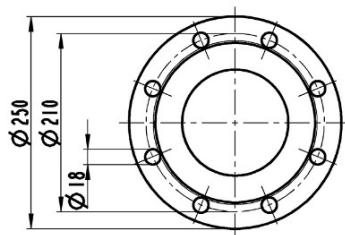
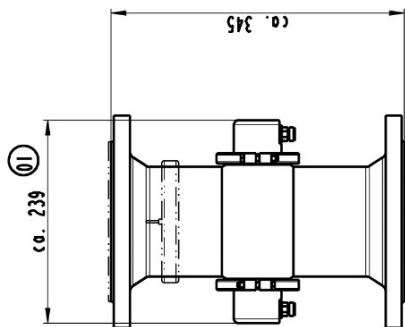
2.
DC 24 V (18-36 V), kein Verpolungsschutz



Type LB 5660-402-00X FlowCell DN 100 Flange, FDA Typ LB 5660-402-00x, Nennweite 100 mm



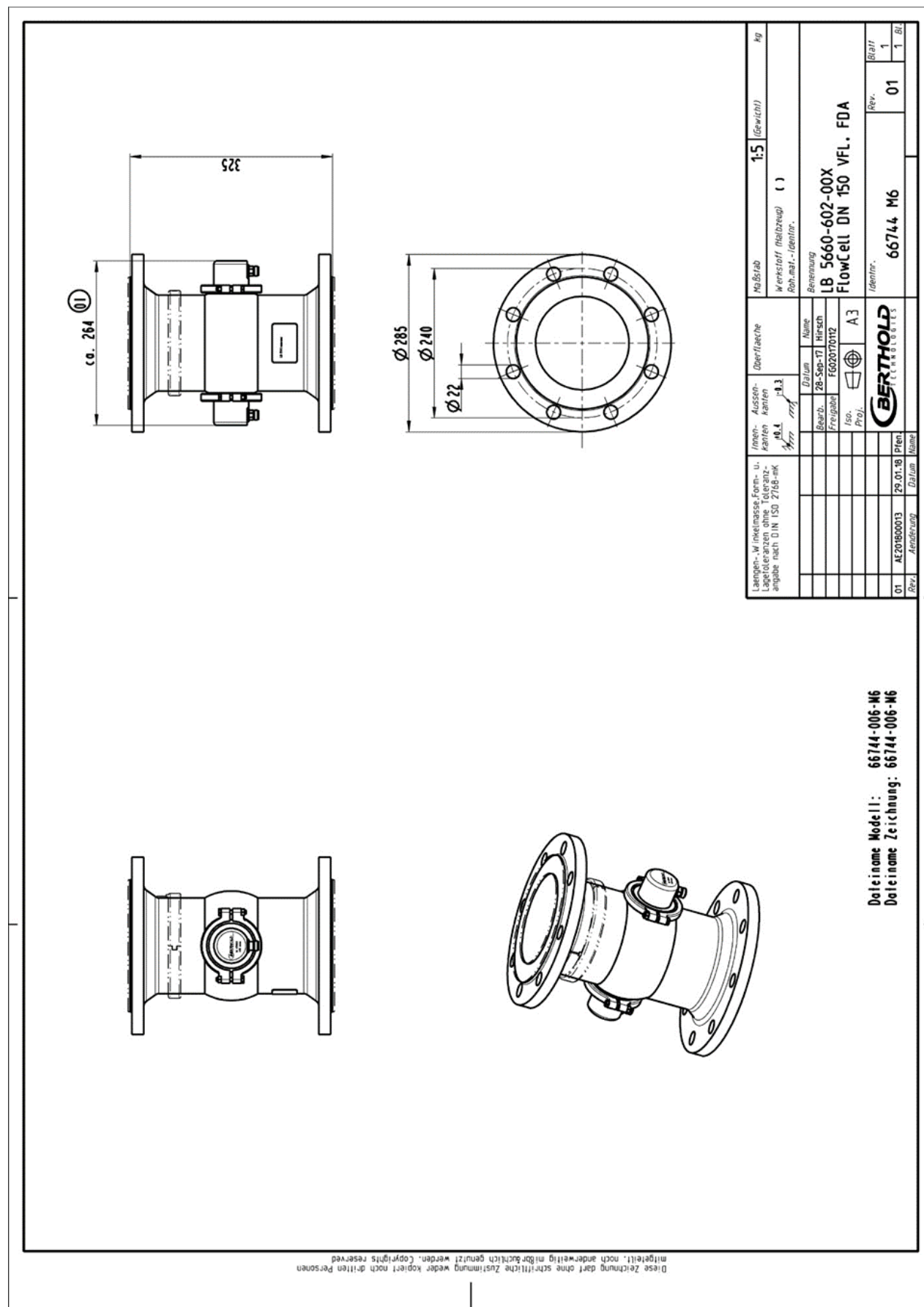
Type LB 5660-502-00X FlowCell DN 125 Flange, FDA Typ LB 5660-502-00x, Nennweite 125 mm



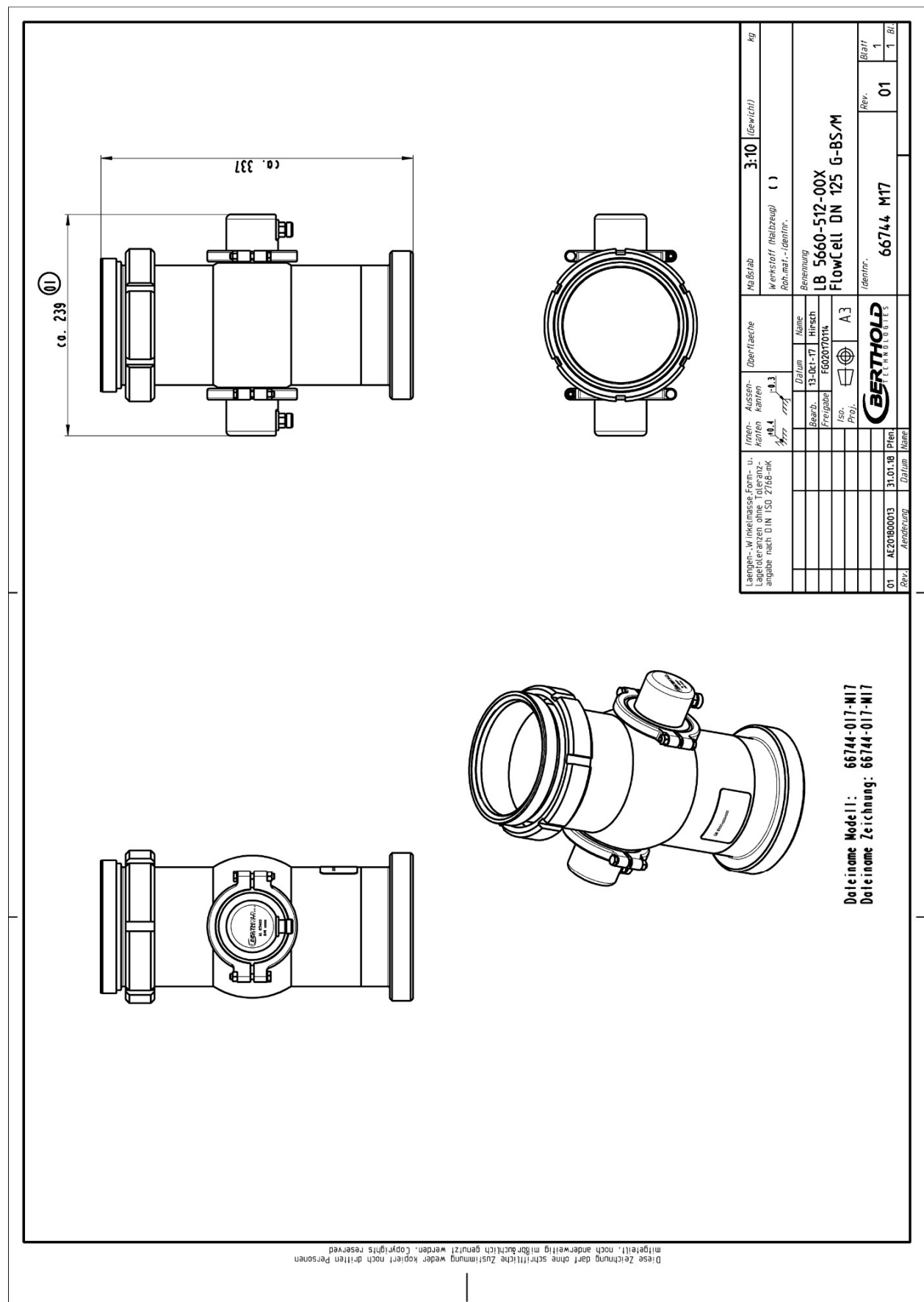
Dateiname Modell: 66744-005-M5
 Dateiname Zeichnung: 66744-005-M5

Laengen-, W inkleinasse, Form- u. Lagenanordnungen ohne Toleranzangaben nach DIN ISO 2768-mS		Innen- Ausson- kanten kanten		Ober/taetche		Maßstab	1:5 (Gewicht)	kg
AE201800013		29.01.18 Pfen.		28-Sep-17 Hirsch		LB 5660-502-00X FlowCell		
Rev.		Datum		Freigabe		DN 125 VFL. FDA		
01		29.01.18		F602017012		Identif.		
AE201800013		29.01.18		F602017012		66744 M5		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		01		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		
Rev.		Datum		Name		1		
01		29.01.18		F602017012		1		
AE201800013		29.01.18		F602017012		1		

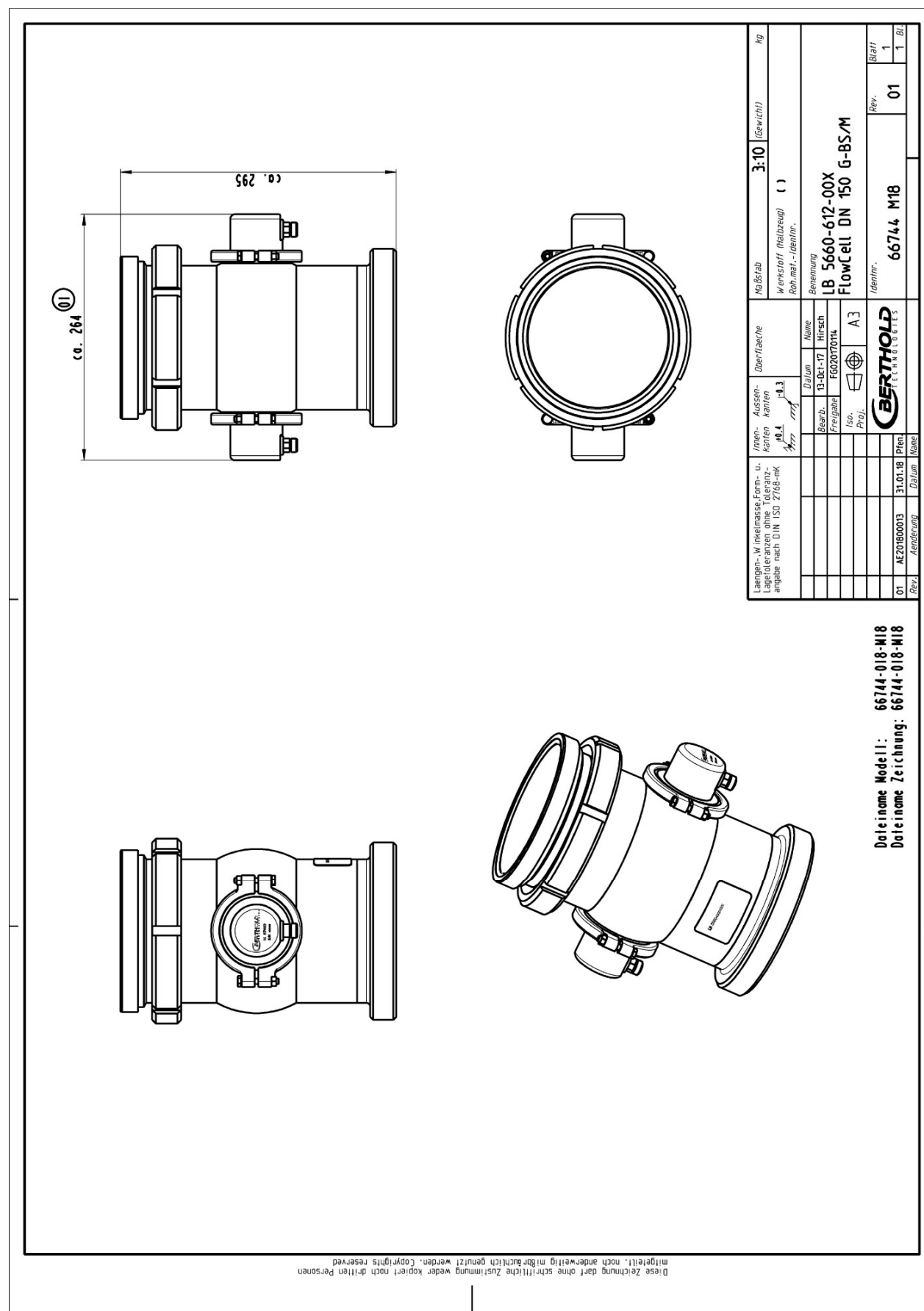
Type LB 5660-602-00X FlowCell DN 150 Flange, FDA Typ LB 5660-602-00x, Nennweite 150 mm



Type LB 5660-512-00X FlowCell DN 125 G-BS/M Typ LB 5660-512-00x, Nennweite 125 mm

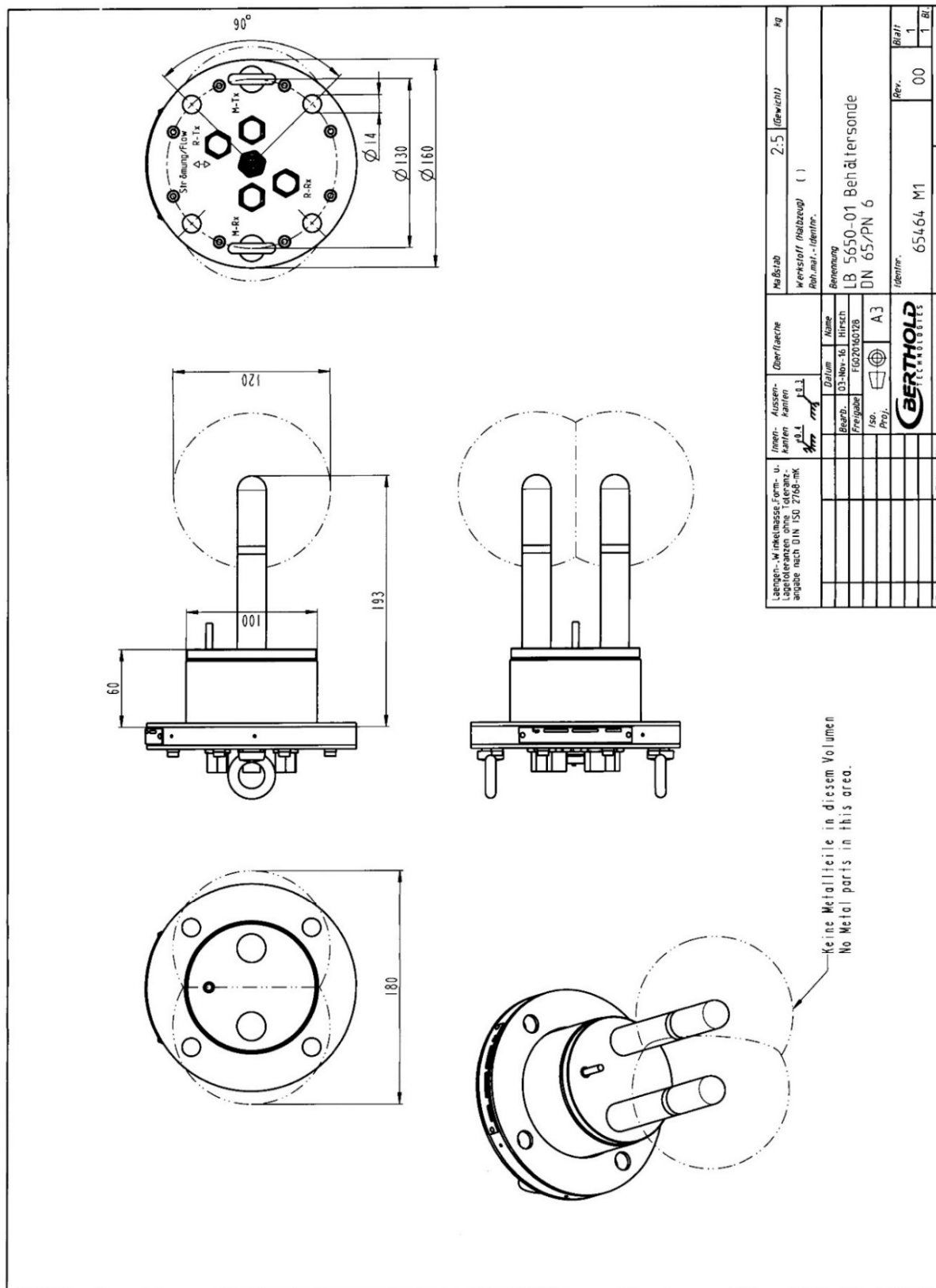


Type LB 5660-612-00X FlowCell DN 150 G-BS/M Typ LB 5660-612-00x, Nennweite 150 mm

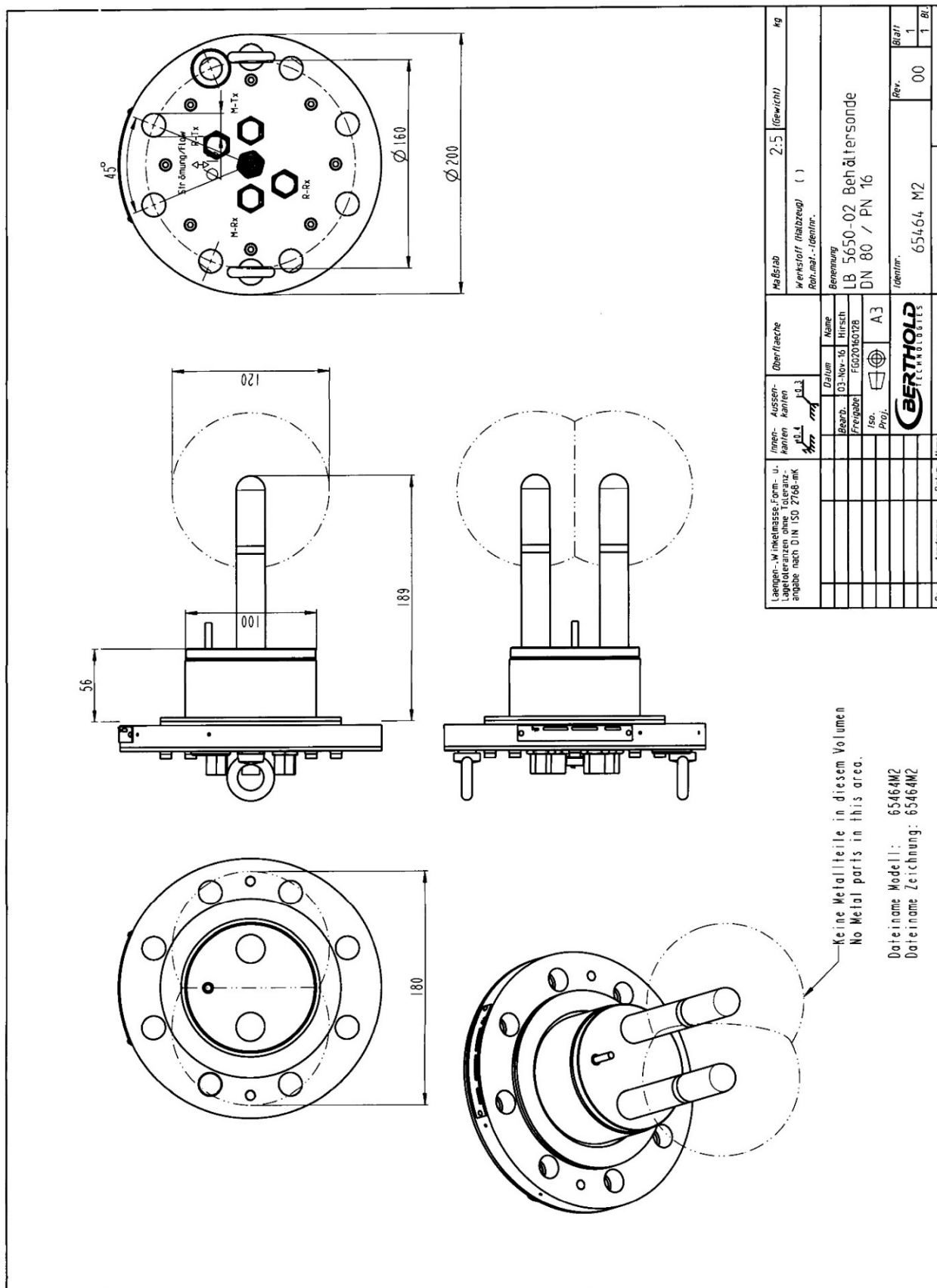


Dimensional Drawings Container Probes Maßbilder Behältersonden

Type LB 5650-01

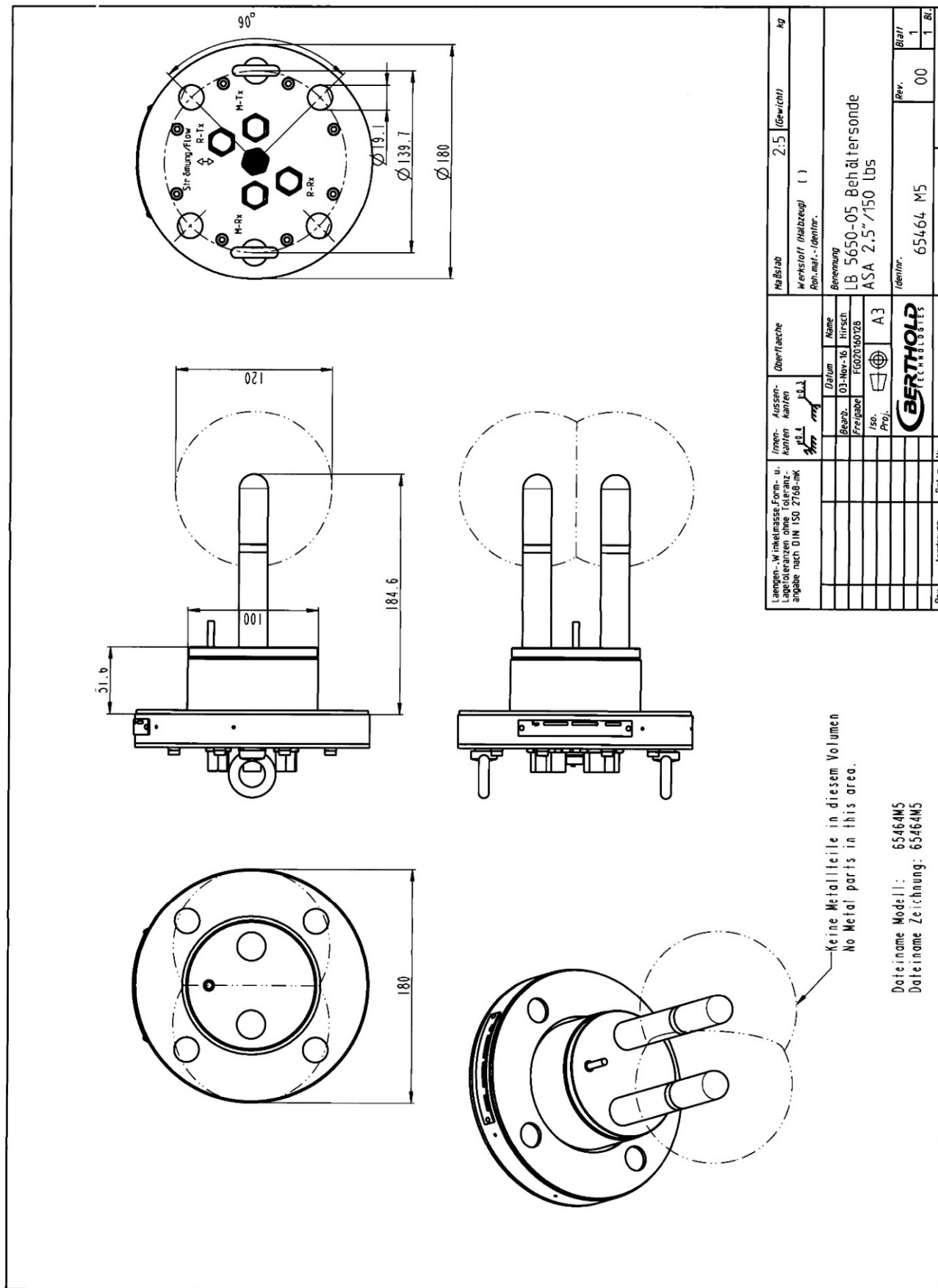


Typ LB 5650-02



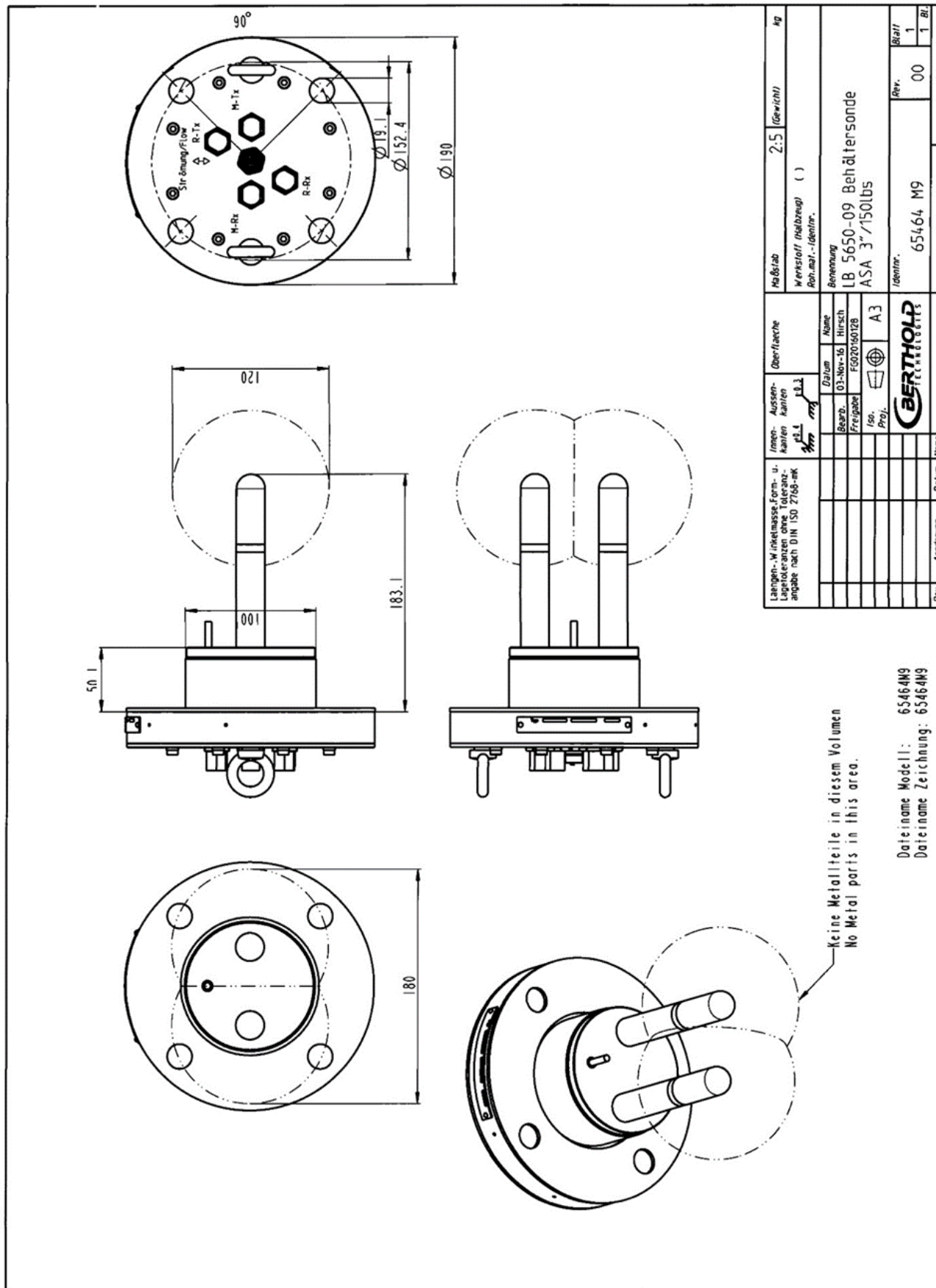
Diese Zeichnung darf ohne schriftliche Zustimmung weder kopiert noch Dritten
 mitgeteilt, noch anderweitig mündlich genutzt werden. Copyrights reserved

Type LB 5650-05



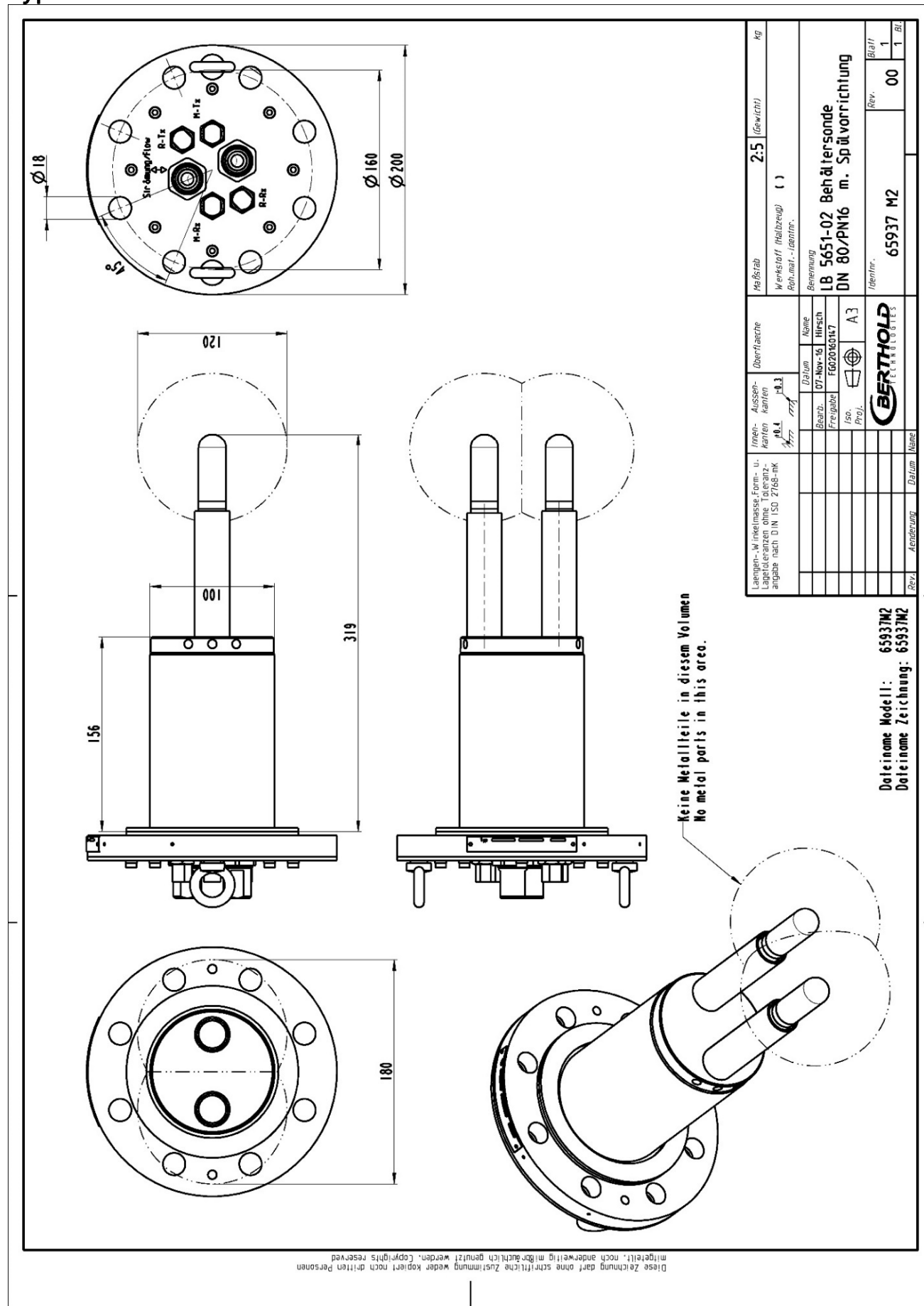
Diese Zeichnung darf ohne schriftliche Zustimmung weder kopiert noch Dritten mitgeteilt, noch anderweitig mirboräuflich genutzt werden. Copyrights reserved

Type LB 5650-09

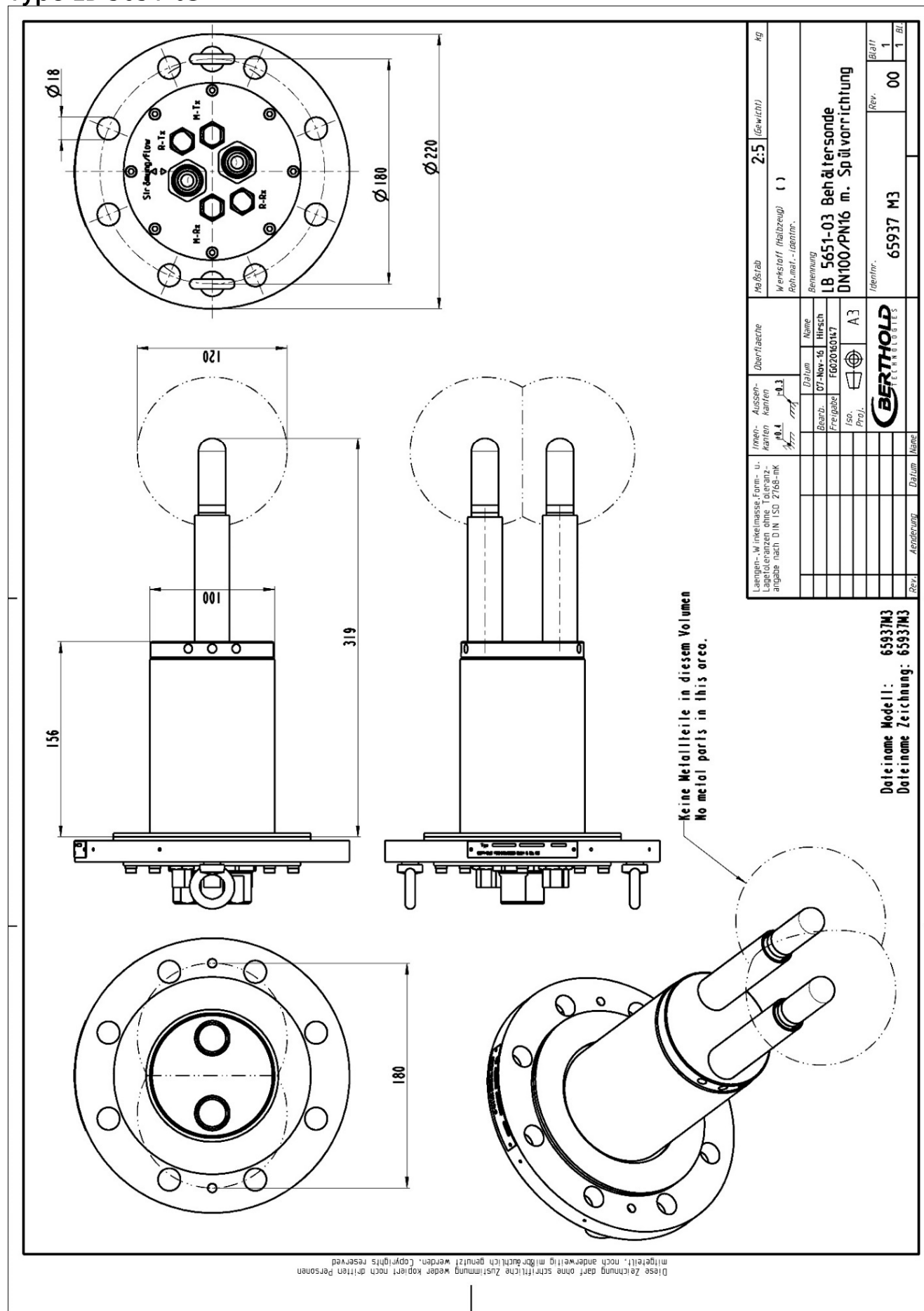


Diese Zeichnung darf ohne schriftliche Zustimmung weder kopiert noch Dritten
 mitgeteilt, noch anderweitig mirbäurlich genutzt werden. Copyrights reserved

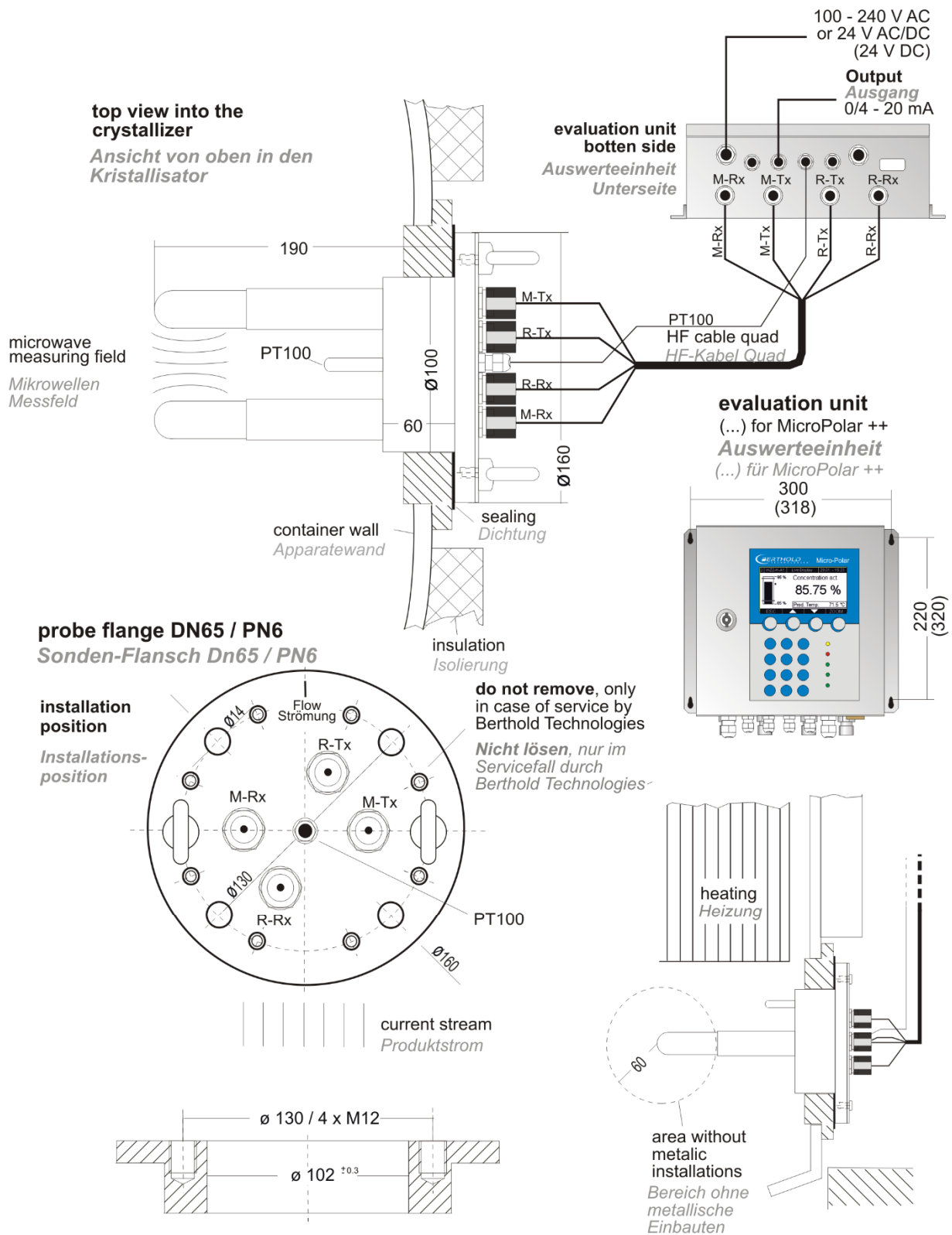
Type LB 5651-02



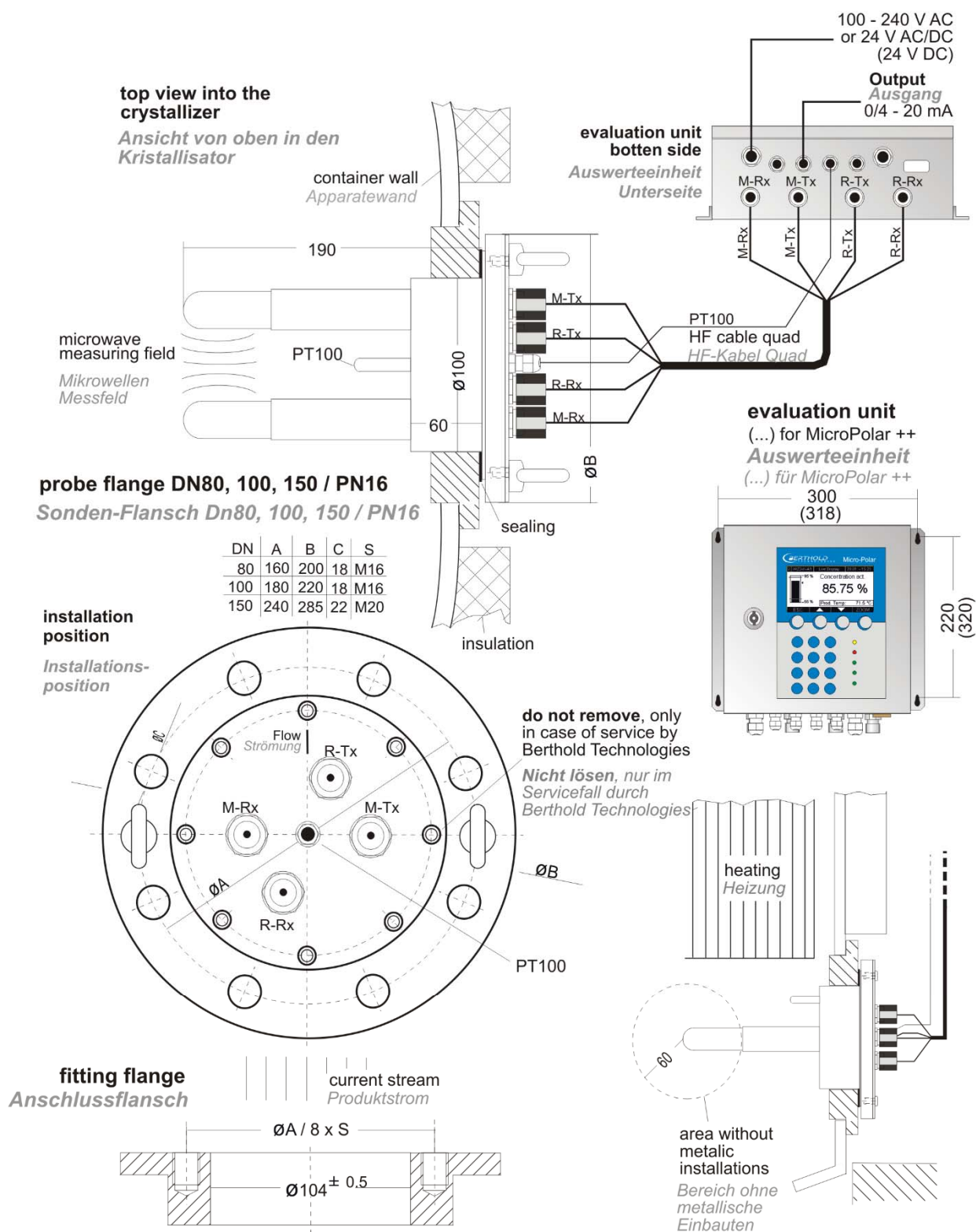
Type LB 5651-03



Installation Sheets for LB 5650 (Container Probe) Montageblätter für LB 5650 (Behältersonde)

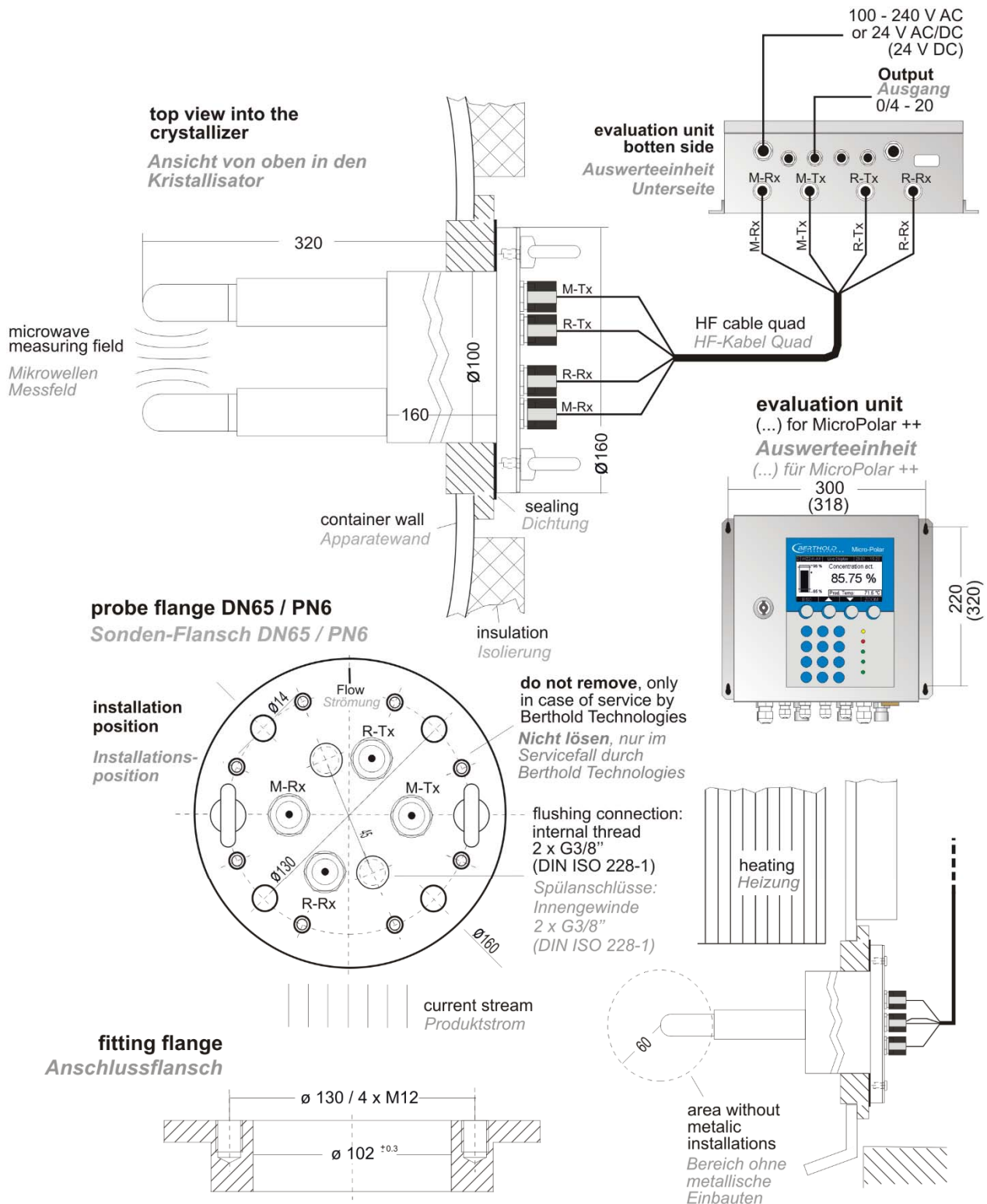


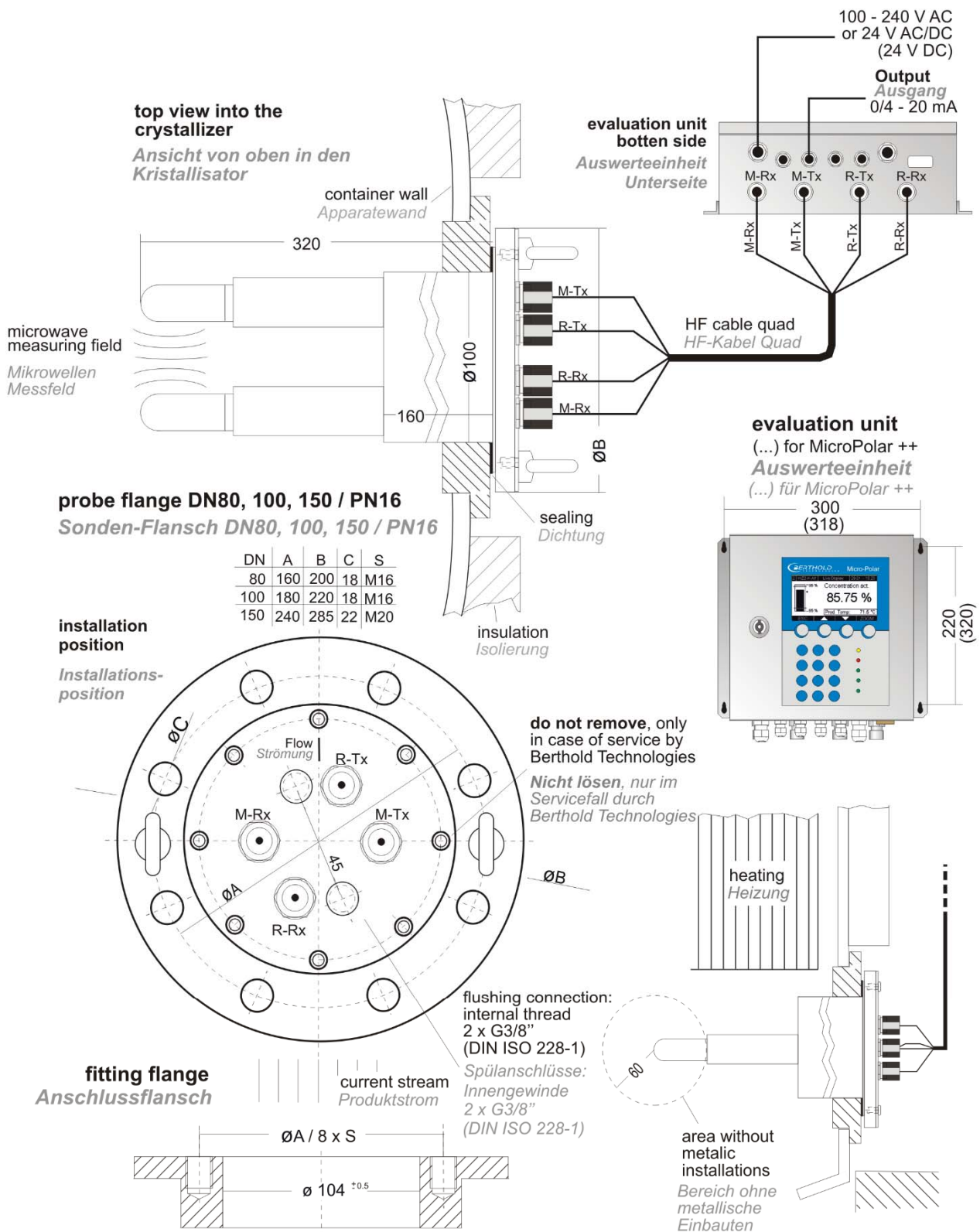
All dimensions in mm
 Alle Angaben in mm



All dimensions in mm
 Alle Angaben in mm

Installation Sheets for LB 5651 (Container Flush Probe) Montageblätter für LB 5651 (Spülsonde)





All dimensions in mm
Alle Angaben in mm