



Concentration / Moisture Measuring Systems MicroPolar Moist LB 568

*Konzentrations- / Feuchte-Messsysteme
MicroPolar Moist
LB 568*

Technical Information *Technische Information*

41990TI
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>
Transmission power <i>Sendeleistung</i>	< 10 μ W (< -20 dBm) All coaxial line power < 1 mW (< 0 dBm) <i>Alles koaxiale Leistungsleistung</i>
Applications <i>Applikationen</i>	Concentration / moisture measurement in pipelines, on conveyor belts and in chutes. <i>Konzentrations-, Feuchtemessung in Rohrleitungen, an Förderbändern und Schächten.</i>

Evaluation Unit

Auswerteinheit

Technical Data Evaluation Unit <i>Technische Daten Auswerteinheit</i>	
Housing <i>Gehäuse</i>	Wall housing made of stainless steel, see dimensional drawing MicroPolar Brix: HxWxD: 300 x 323 x 140 mm MicroPolar Brix++: HxWxD: 400 x 338 x 170 mm <i>Wandgehäuse aus Edelstahl, siehe Maßbild.</i> <i>MicroPolar Brix:HxBxT: 300 x 323 x 140 mm</i> <i>MicroPolar Brix++: HxBxT: 400 x 338 x 170 mm</i>
Protection type <i>Schutzart</i>	IP 65 <i>IP 65</i>
Weight <i>Gewicht</i>	about 8.0 kg <i>ca. 8,0 kg</i>
Ambient conditions during operation <i>Umgebungsbedingungen im Betrieb</i>	Relative humidity: max. 85 %, short-term up to 100 %, no condensation Altitude: max. 2000 m -20 ... +45 °C (253...318 K) <i>Relative Luftfeuchte: max. 85 %, kurzzeitig bis 100 %, keine Kondensation</i> <i>Höhenlage: max. 2000 m</i> <i>-20...+45 °C (253...318 K)</i>
Ambient conditions during storage <i>Umgebungsbedingung bei Lagerung</i>	-20 ... +70 °C (253 K ...343 K), Relative humidity: max. 85 %, no condensation -20...+70 °C (253...343 K) <i>Relative Luftfeuchte: max. 85 %,keine Kondensation</i>
Achievable accuracy <i>Erreichbare Genauigkeit</i>	≤ 0.1 weight % (standard deviation) depending on product and process ≤ 0,1 Gewichts % (Standardabweichung) je nach Produkt und Prozess

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>	Depending on device type: 1.) 100 ... 240 V AC, 50/60 Hz 2.) 24 V DC: 18 ... 36 V, reverse polarity protection <i>Je nach Geräteausführung: 1.) 100 ... 240 V AC, 50/60 Hz 2.) 24 V DC: 18 ... 36 V, kein Verpolungsschutz</i>
Power consumption <i>Leistungsaufnahme</i>	max. (48/60) VA, je nach Konfiguration <i>max. (48/60) VA, je nach Konfiguration</i>
Fuses <i>Sicherungen</i>	2 x 2.0 A / 250 V / T at 100 ... 240 V AC; ID no. 4403 or 2 x 6.3 A / 250 V / M at 18 ... 36 V DC; ID no. 4408 <i>2 x 2.0 A / 250 V / T bei 100 ... 240 V AC; Id.-Nr. 4403 oder 2 x 6.3 A / 250 V / M bei 18 ... 36 V DC; Id.-Nr. 4408</i>
Battery type <i>Batterietyp</i>	3 V Lithium button cell, type CR2032, ID no. 17391 <i>3 V Lithium Knopfzelle, Typ CR2032, Id.-Nr 17391</i>
Measured value <i>Messwert</i>	e.g. concentration, moisture content <i>z.B. Konzentration, Feuchtegehalt</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 Temperatúrausgabe</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> <i>DI3: Keine, Probeaufnahme, Produktauswahl</i> <i>Funktionsbeschreibung:</i> 1. <i>Messung (Start/Stopp), <u>offen</u>: Messung gestoppt, <u>geschlossen</u>: Messung gestartet bzw. Messung läuft</i> 2. <i>Messung anhalten, <u>offen</u>: Messung läuft, <u>geschlossen</u>: Messung angehalten, d.h. Mittelwerte und Stromausgang sind eingefroren</i> 3. <i>Produktauswahl über ein DI:</i> <i><u>offen</u>: Produkt 1 (P1), <u>geschlossen</u>: P2</i> <i>Produktauswahl über zwei DI's:</i> <i>DI2 & DI3 <u>offen</u>: P1</i> <i>DI2 <u>geschlossen</u> & DI3 <u>offen</u>: P2</i> <i>DI2 <u>offen</u> & DI3 <u>geschlossen</u>: P3</i> <i>DI2 & DI3 <u>geschlossen</u>: P4</i> 4. <i>Probeaufnahme: <u>offen</u>: keine Aktionen, <u>geschlossen</u>: einmalige Aufnahme startet</i>
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 <i>Konfigurationsmöglichkeiten:</i> - 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

Horn and Spiral Antennas

Horn - und Spiralantennen

Horn antenna (ID no.: 10806)

Hornantenne (Id.-Nr.: 10806)

Application <i>Anwendung</i>	Used in pairs, for example on conveyor belts and chutes for the moisture measurement in bulk goods. <i>Paarweise Anwendung z.B. an Förderbändern und Schächten zur Feuchtebestimmung in Schüttgütern.</i>
Material <i>Material</i>	Stainless steel, microwave window made of Makrolon <i>Edelstahl, Mikrowellenfenster aus Makrolon</i>
Weight <i>Gewicht</i>	1.4 kg <i>1,4 kg</i>
Temperature range <i>Temperaturbereich</i>	Ambient temperature: -20...60°C (253...333 K) Storage temperature: 10...80°C (283...353 K) <i>Umgebungstemperatur: -20...60°C (253...333 K) Lagertemperatur: 10...80°C (283...353 K)</i>
Connection <i>Anschluss</i>	1 x HF connections: N-connector, 50 Ω <i>1 x HF-Anschlüsse: N-Stecker, 50 Ω</i>

Accessories antenna fixture (ID no.: 10805)

Zubehör Antennenhalterung (Id.-Nr.: 10805)

Material <i>Material</i>	Galvanized steel <i>Verzinkter Stahl</i>
Weight <i>Gewicht</i>	3.8 kg <i>3,8 kg</i>
Dimensions <i>Maße</i>	See dimensional drawings <i>Siehe Maßzeichnungen</i>

Spiral antenna (ID no.: 15394)

Spiralantenne (Id.-Nr.: 15394)

Application <i>Anwendung</i>	Used in pairs, for example on conveyor belts and chutes for the moisture measurement in bulk goods. <i>Paarweise Anwendung z.B. an Förderbändern und Schächten zur Feuchtebestimmung in Schüttgü-tern.</i>
Material <i>Material</i>	Stainless steel, plastic <i>Edelstahl, Kunststoff</i>
Weight <i>Gewicht</i>	0.4 kg <i>0,4 kg</i>
Temperature range <i>Temperaturbereich</i>	Ambient temperature: -20...60°C (253...333 K) Storage temperature: 10...80°C (283...353 K) <i>Umgebungstemperatur: -20...60°C (253...333 K)</i>

	<i>Lagertemperatur: 10...80°C (283...353 K)</i>
Connection <i>Anschlüsse</i>	1 x HF connections: N-connector, 50 Ω 1 x HF-Anschlüsse: N-Stecker, 50 Ω
Dimensions <i>Maße</i>	See dimensional drawings <i>Siehe Maßzeichnungen</i>

Technical data Radiometric Mass per Unit Area Measurement Technische Daten radiometrische Flächengewichtsmessung

Scintillation counter Szintillationszähler	
Versions	1. With axial collimator (Id. 56942), for frontal irradiation 2. With radial collimator (Id. 56943), for radial irradiation
Varianten	1. Mit Kollimator axial (Id.-Nr. 56942), für frontale Einstrahlung 2. Mit Kollimator radial (Id.-Nr. 56943), für seitliche Einstrahlung
Crystal Kristall	CsJ 40 x 50
Material	Stainless steel Collimator: Lead, painted steel
Material	Edelstahl Kollimator: Blei, lackierter Stahl
Protection type Schutzart	IP 67
Weight	Without collimator: approx. 2 kg With axial collimator: approx. 10.6 kg With radial collimator: approx. 10 kg
Gewicht	Ohne Kollimator: ca. 2 kg Mit Kollimator axial: ca. 10.6 kg Mit Kollimator radial: ca. 10 Kg
Power supply Hilfsenergie	12...24 V DC, approx.. 1.2 W 12...24 V DC, ca. 1,2 W
Operating temperature Betriebstemperatur	-20...50 °C (253...323 K)
Storage temperature Lagertemperatur	-20 ... 60 °C (253...333 K)
Connection cable	3 m long, 7-wire, shielded (7 x 0.5 mm ²), cable connection angled 90°, temperature range: -40...70 °C (233...343 K)
Anschlusskabel	3 m lang, 7-adrig, geschirmt (7 x 0.5 mm ²), Kabelanschluss 90° gewinkelt, Temperaturbereich: -40...70 °C (233...343 K)
Maße	See dimensional drawings Siehe Maßzeichnungen
Accessory mounting options Zubehör Befestigungsmöglichkeiten	
ID-No Id.-Nr.	Description Beschreibung
56860	Scintillation counter bracket, complete Material: galvanized steel, see plastic Dimensional drawings Szintillationszähler-Halterung, komplett Material: verzinkter Stahl, siehe Kunststoff Maßzeichnungen
25668	Clamps (1 set = 2 clamps) Spannschellen (1 Satz = 2 Schellen)

Shield for pint source (LB 744x)

Abschirmung für Punktstrahler (LB 744X)

ID-No <i>Id.-Nr.</i>	Description <i>Beschreibung</i>
37624	LB 7440-D-CR, internal parts made of stainless steel LB 7440-D-CR, Innenteile aus Edelstahl
38040	LB 7440-DE-CR, stainless steel LB 7440-DE-CR, Edelstahl
38042	LB 7445-D-CR, internal parts made of stainless steel, with leakage protection LB 7445-D-CR, Innenteile aus Edelstahl, mit Auslaufschutz
38043	LB 7445-DE-CR, stainless steel, with leakage protection LB 7445-DE-CR, Edelstahl, mit Auslaufschutz

Shielding accessories

Zubehör Abschirmung

ID-No <i>Id.-Nr.</i>	Description <i>Beschreibung</i>
14716	Shielding bracket, complete for LB 7440 D <i>Abschirmungs-Halterung, komplett für LB 7440 D</i>
52752	Protective cover for shielding LB 7440/42/44 <i>Schutzhülle für Abschirmung LB 7440/42/44</i>
11213	Radiation sign (RADIOACTIVE), plastic <i>Strahlerschild (RADIOACTIVE), Kunststoff</i>
14658	Radiation sign (RADIOACTIVE), aluminum <i>Strahlerschild (RADIOACTIVE), Alu</i>

Pneumatic shutter for shielding

Pneumatischer Verschluss für Abschirmung

ID-No <i>Id.-Nr.</i>	Description <i>Beschreibung</i>
36119	Pneumatic shutter actuator with limit switch, IP 65 <i>Pneumatischer Verschlussantrieb mit Endschalter, IP 65</i>

Daten für pneumatischen Verschlussantrieb

Compressed air:	min. 4 x 4 bar max. 4 x 7 bar Connection: G 1/8
<i>Druckluft:</i>	<i>min. 4 x 4 bar max. 4 x 7 bar Anschluss: G 1/8</i>
Air quality:	Clean as usual for pneumatic tools, oil-free
<i>Luftqualität:</i>	<i>Sauber wie für Druckluft-Werkzeuge üblich, ölfrei</i>
Temperature range:	-20...80 °C
<i>Temperaturbereich:</i>	

Limit switch unit, signaling options for OPEN / CLOSE

Endschaltereinheit, Optionen für Signalisierung AUF / ZU

Option	IP 65 2 contacts (OPEN/CLOSE) 48 V DC, 1A <i>IP 65 2 Kontakte (AUF/ZU) 48 V DC, 1A</i>
--------	---

Point source Cs-137
Punktstrahler Cs-137

ID no.: <i>Id.-Nr.:</i>	Made of stainless steel <i>Aus Edelstahl</i>
54712-08	Pkt.-Str. Cs-137 370 MBq (10 mCi) – SSC-200
54712-12	Pkt.-Str. Cs-137 1110 MBq (30 mCi) – SSC-200
54712-13	Pkt.-Str. Cs-137 1850 MBq (50 mCi) – SSC-200
ID no.: <i>Id.-Nr.:</i>	Description <i>Beschreibung</i>
54281-08	Pkt.-Str. Cs-137 370 MBq (10 mCi) – SSC-100
54281-12	Pkt.-Str. Cs-137 1110 MBq (30 mCi) – SSC-100
54281-13	Pkt.-Str. Cs-137 1850 MBq (50 mCi) – SSC-100

Technical Data Measuring Chute

Technische Daten Messschacht

Measuring chute, complete Messschacht, komplett	
Application <i>Anwendung</i>	For moisture and concentration determination in bulk material. <i>Zur Feuchte- und Konzentrationsbestimmung in Schüttgütern.</i>
Variants / chute material <i>Varianten / Schachtmaterial</i>	1. Polypropylen homo polymer (PP-H) ID no. 57957 2. Polyvinylidene fluoride (PVDF), ID no. on request 1. Polypropylen-Homopolymer (PP-H), Id.-Nr. 57957 2. Polyvinylidenfluorid (PVDF), Id.-Nr. auf Anfrage
Components <i>Komponenten</i>	- Chute - Assembly plate - 4 brackets - two RF angle connectors - Fastening material - Schacht - Montageplatte - 4 Halterungen - zwei RF-Winkelverbinder - Befestigungsmaterial
Weights <i>Gewichte</i>	Only the chute: Version 1: approx. 10 Kg Version 2: on request Measuring chute, complete: Version 1: approx. 33 Kg Version 2: on request <i>Nur der Schacht:</i> <i>Variante 1: ca. 10 Kg</i> <i>Variante 2: auf Anfrage</i> <i>Messschacht komplett:</i> <i>Variante 1: ca. 33 Kg</i> <i>Variante 2: auf Anfrage</i>
Temperature range <i>Temperaturbereich</i>	Environment: 0...50 °C (273...323 K) Storage: 10...80 °C (283...353 K) Product temperature: Version 1: > 0...90 °C (273...363 K) Version 2: > 0...140 °C (273 ... 413 K) <i>Umgebung: 0...50 °C (273...323 K)</i> <i>Lagerung: 10...80 °C (283...353 K)</i> <i>Produkttemperatur:</i> <i>Variante 1: > 0...90 °C (273...363 K)</i> <i>Variante 2: > 0...140 °C (273 ... 413 K)</i>

Assembly plate, brackets <i>Montageplatte, Halterungen</i>	Material: Stainless steel, galvanized steel <i>Material: Edelstahl, verzinkter Stahl</i>
Dimensions <i>Maße</i>	See dimensional drawings <i>Siehe Maßzeichnungen</i>

Technical Data HF Cable

Technische Daten HF Kabel

HF cable Quad <i>HF-Kabel Quad</i>	
Material/Material	Cable sheath: Polyethylene (PE) <i>Kabelmantel: Polyethylen (PE)</i>
Protection type <i>Schutzart</i>	IP 68 in the screwed-on state <i>IP 68 im aufgeschraubten Zustand</i>
Temperature <i>Temperatur</i>	In operation: -40 ... +85 °C When installing: -40 ... +85 °C <i>Bei Betrieb: -40 ... +85 °C</i> <i>Bei Installation: -40 ... +85 °C</i>
Dämpfungsbelaag <i>Dämpfungsbelaag</i>	approx. 0.3 dB/m <i>ca. 0,3 dB/m</i>
Cable length [m] <i>Kabellänge [m]</i>	ID no. <i>Ident.-Nr.</i>
0.5	11473
1.0	11474
1.5	11475
2.0	11476
2.5	11477
3.0	11478
3.5	11479
4.0	11480

Format of Serial Data Output RS232

Headline

Date·Time→State→Status→DetectorStatus→Synchronizer→Product→Att→Phi→Phi(f=0)→

R2→Corr→Tint→IN1→IN2→PT100→C→Cm→C2→C2m→CpsM→DetTemp→Mqua→Density¶

Following lines

01.01.2005·00:00:00→0000→0→0→5→1→0.43→5.30→0.07→0.98→

1 2 3 4 5 6 7 8 9 10

0.0→0.0→0.0→0.0→75.36→75.00→0.00→0.00→3653→35.8→10.25→0.8612¶

11 12 13 14 15 16 17 18 19 20 21 22

Column no.	Description	Format
1	Date and time	DD.MM.YY·HH:MM:SS
2	State	4 digits, HEX
3	Status: Information on the quality of the last measurement	0 : Measurement OK < 0 : Error
4	Detector status: Information on the quality of the last measurement	0 : Measurement OK < 0 : Error
5	Product synchronization	5 : not active 1: still asynchronous 0: all values synchronous -1: Error -2: Time too short for syn. -3: Speed outside range
6	Product number	X (1 to 4)
7	Attenuation [dB]	X.XX
8	Phase [°/GHz]	X.XX
9	Dispersion of phase regression	X.XX
10	Correlation of phase regression	X.XX
11	Device temperature [temperature unit]	X.X
12	Current input 1 [unit of current input]	X.X
13	Current input 2 [unit of current input]	X.X
14	PT100 temperature [temperature unit]	X.X
15	Concentration 1 live	X.XX
16	Concentration 1 averaged	X.XX
17	Concentration 2 live	X.XX
18	Concentration 2 averaged	X.XX
19	Averaged count rate [counts/s]	X
20	Detector temperature [°C]	X.X
21	Mass per unit area [g/cm²]	X.XX
22	Density [g/cm³]	X.XXXX

Special characters

“→” Tabulation

“¶” Carriage return + Line feed

“.” Spaces

Format der seriellen Datenausgabe RS232

Titelzeile

Date·Time→State→Status→DetectorStatus→Synchronizer→Product→Att→Phi→
 R2→Corr→Tint→IN1→IN2→PT100→C→Cm→C2→C2m→CpsM→DetTemp→Mqua→Density¶

Folgezeilen

01.01.2005:00:00:00→0000→0→0→5→1→0.43→5.30→0.07→0.98→
 1 2 3 4 5 6 7 8 9 10

0.0→0.0→0.0→0.0→75.36→75.00→0.00→0.00→3653→35.8→10.25→0.8612→¶
 11 12 13 14 15 16 17 18 19 20 21 22

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

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@Berthold.com
www.Berthold.com

EC-Declaration

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

Description: **Concentration- and Moisture-Measuring Systems
Micro-Polar 2, Micro-Polar 2 ++,
Micro-Polar Moist, and Micro-Polar Moist ++**

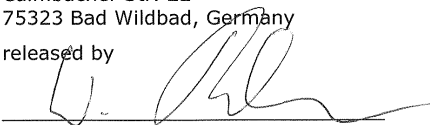
Type: **LB 567-XX and LB 568-XX**

	Richtlinie und Änderungen	angewendete Normen
EMC	2004/108/EC	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
LVD	2006/95/EC	EN 61010 part 1 2002-08

This declaration is issued by the manufacturer

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

released by


Dr. Wilfried Reuter – Technical Director
Bad Wildbad, 28th of April 2010

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

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

Sparkasse PF-CW	75323 Bad Wildbad	Konto/Account No. 8 045 003 (BLZ 666 500 85)	SWIFT-BIC PZHSDE66	IBAN: DE37 6665 0085 0008 0450 03
Volksbank	75119 Pforzheim	Konto/Account No. 957 004 (BLZ 666 900 00)	SWIFT-BIC VBPFDE66	IBAN: DE85 6669 0000 0000 9570 04
Commerzbank	75105 Pforzheim	Konto/Account No. 6 511 120 (BLZ 666 800 13)	SWIFT-BIC DRESDEFF66	IBAN: DE05 6668 0013 0651 1120 00

EG-Konformitätserklärung



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 der nachfolgend bezeichneten Geräte / Systeme / Anlagen in der von uns in den Verkehr gebrachten Ausführung den unten genannten einschlägigen EG-Richtlinien entspricht.

Produktbezeichnung: **Konzentrations- und Feuchte-Messsysteme
Micro-Polar 2, Micro-Polar 2 ++,
Micro-Polar Moist und Micro-Polar Moist ++**

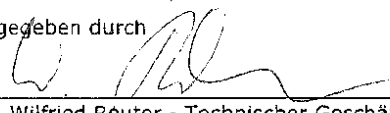
Typenbezeichnung: **LB 567-XX und LB 568-XX**

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 EN 61000-4-11 1994-08 +A1:2001-02 Namur NE21 2004
Niederspg.	2006/95/EG	EN 61010 Teil 1 2002-08

Diese Erklärung wird verantwortlich für den Hersteller

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

abgegeben durch


Dr. Wilfried Reuter - Technischer Geschäftsführer
Bad Wildbad, den 28. April 2010

Registered office / Firm of the manufacturer
Personlich haftende Gesellschaft / Fully liable Partnership
"Berthold Technologies" / "Berthold Technologies"
Berthold Technologies / Fully liable Partnership
Oftersheim / Germany
Deutsche Firmennummer / German Tax No.
VAT-Reg. No.

Stichtag / Date of issue
BERTHOLD TECHNOLOGIES
Calmbacher Str. 22
75323 Bad Wildbad, Germany
Dr. Wilfried Reuter
DE 33056111
49718708100
01 976666-0

Spezialvers. PC / Software
Vollversion
Calmbacher

75323 Bad Wildbad
75323 Bad Wildbad
75323 Bad Wildbad

Kontaktdaten No. 41990 TI (Rev. 00/06/2018)
Kontaktdaten No. 957 000 (Rev. 00/06/2018)
Kontaktdaten No. 957 000 (Rev. 00/06/2018)

SWP-PC-PC-PC-PC-PC-PC
SWP-PC-PC-PC-PC-PC-PC
SWP-PC-PC-PC-PC-PC-PC

Barcode: 6666 0000 0000 0000 00
Barcode: 6666 0000 0000 0000 00
Barcode: 6666 0000 0000 0000 00

Frequency License *Frequenzzulassung*

TCB

GRANT OF EQUIPMENT
AUTHORIZATION

TCB

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

CETECOM ICT Services GmbH
Untertuerkheimer Strasse 6-10
66117 Saarbruecken,
Germany

Date of Grant: 02/10/2016

Application Dated: 10/07/2015

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: R9ZFCC02X03
Name of Grantee: Berthold Technologies
Equipment Class: Ultra Wideband Transmitter
Notes: Concentration / Moisture / Dry Mass /
Density Measuring System

Grant Notes

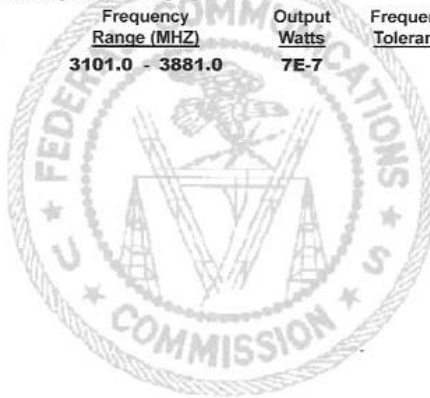
FCC Rule Parts
15F

Frequency
Range (MHZ)
3101.0 - 3881.0

Output
Watts
7E-7

Frequency
Tolerance

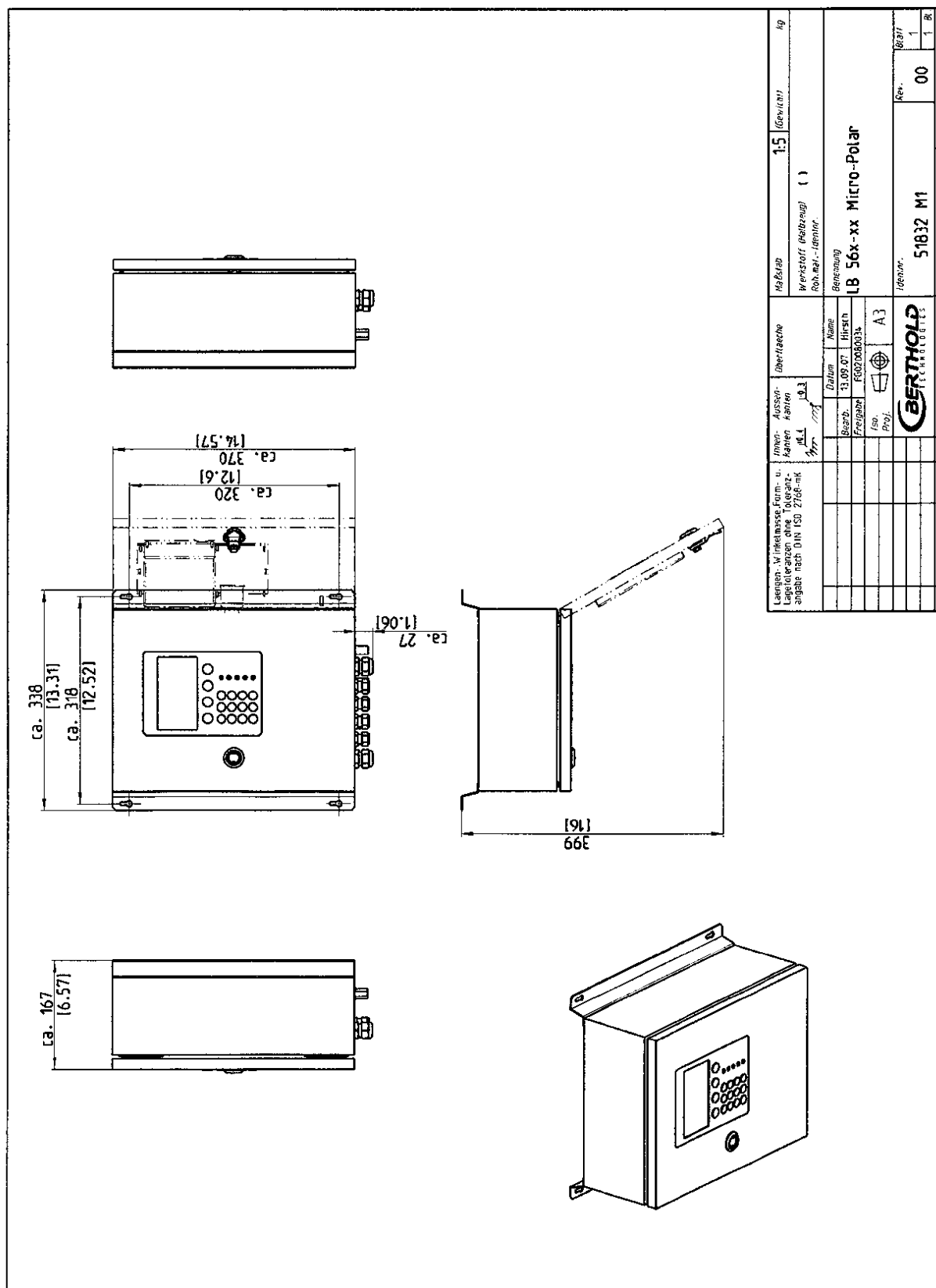
Emission
Designator



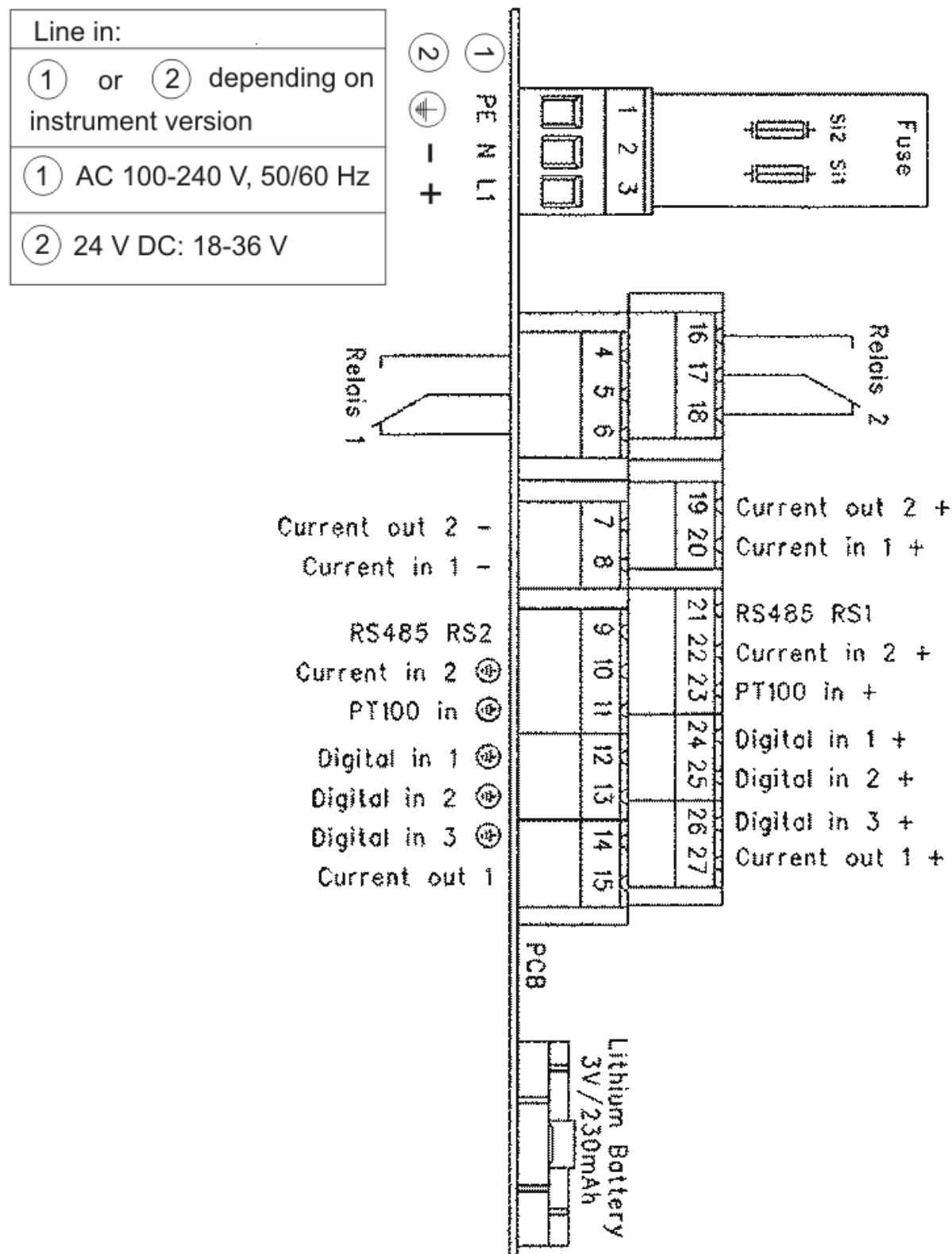
Dimensional Drawings

Maßzeichnungen

Evaluation Unit
Auswertereinheit

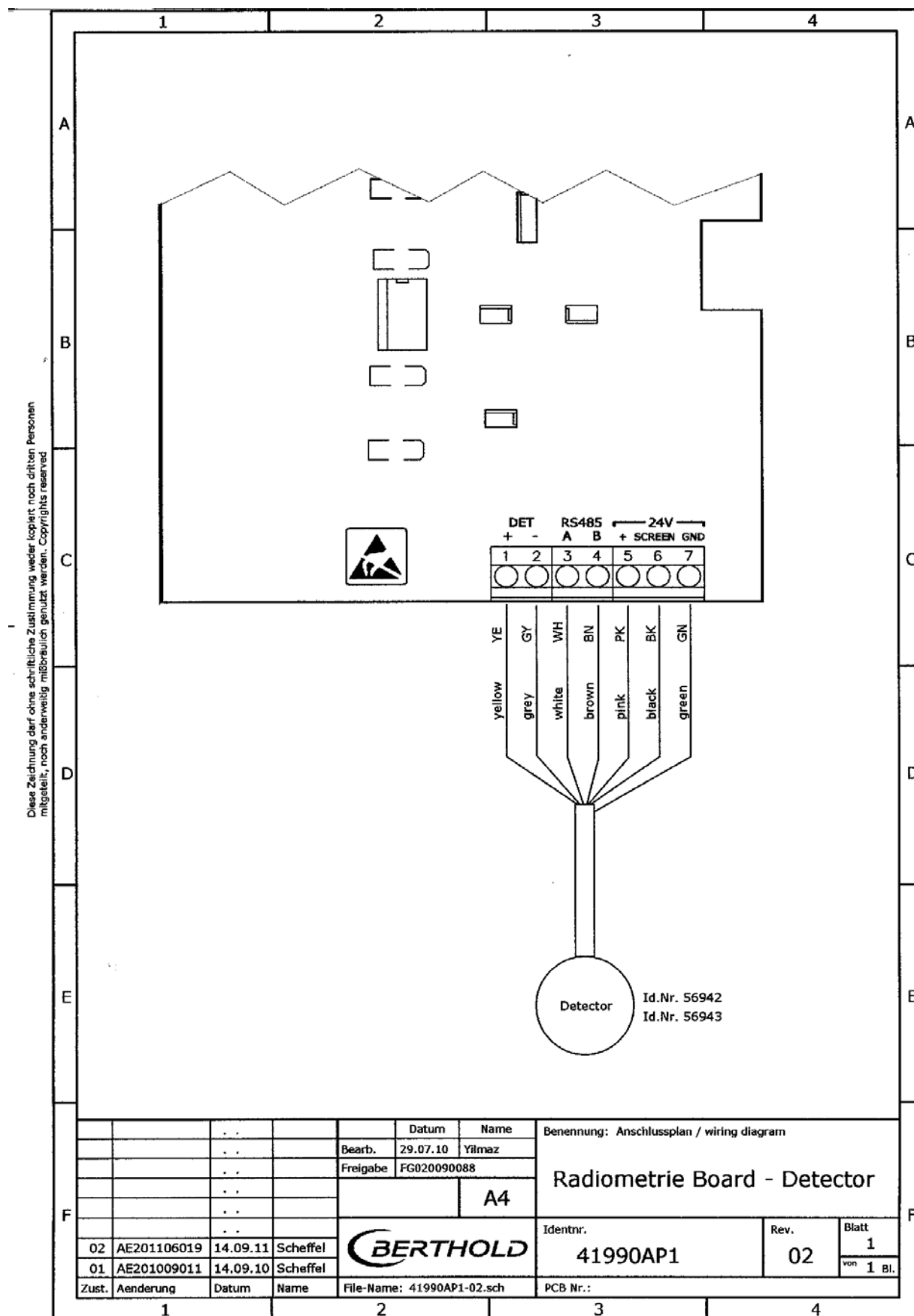


Electrical Wiring Diagram *Elektrischer Anschlussplan*



Electrical Wiring Diagram Scintillation Counter

Elektrischer Anschlussplan Szintillationszähler

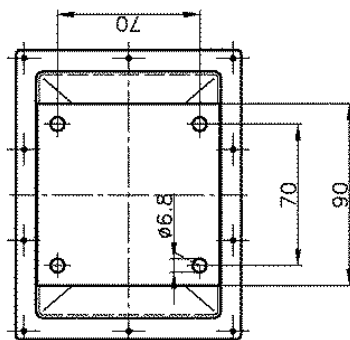
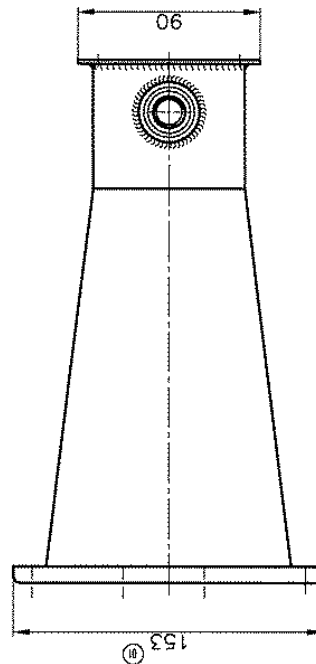
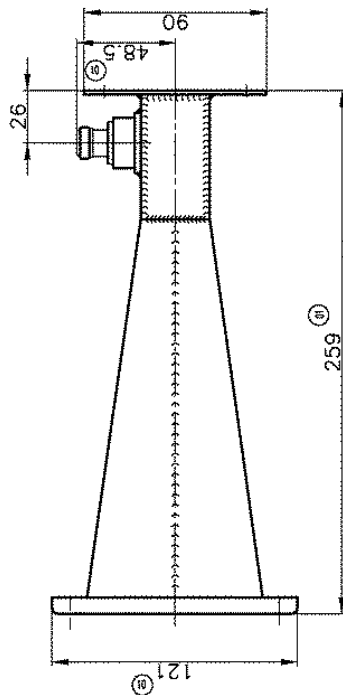


Dimensional Drawings Horn and Spiral Antennae

Maßbilder Horn- und Spiralantennen

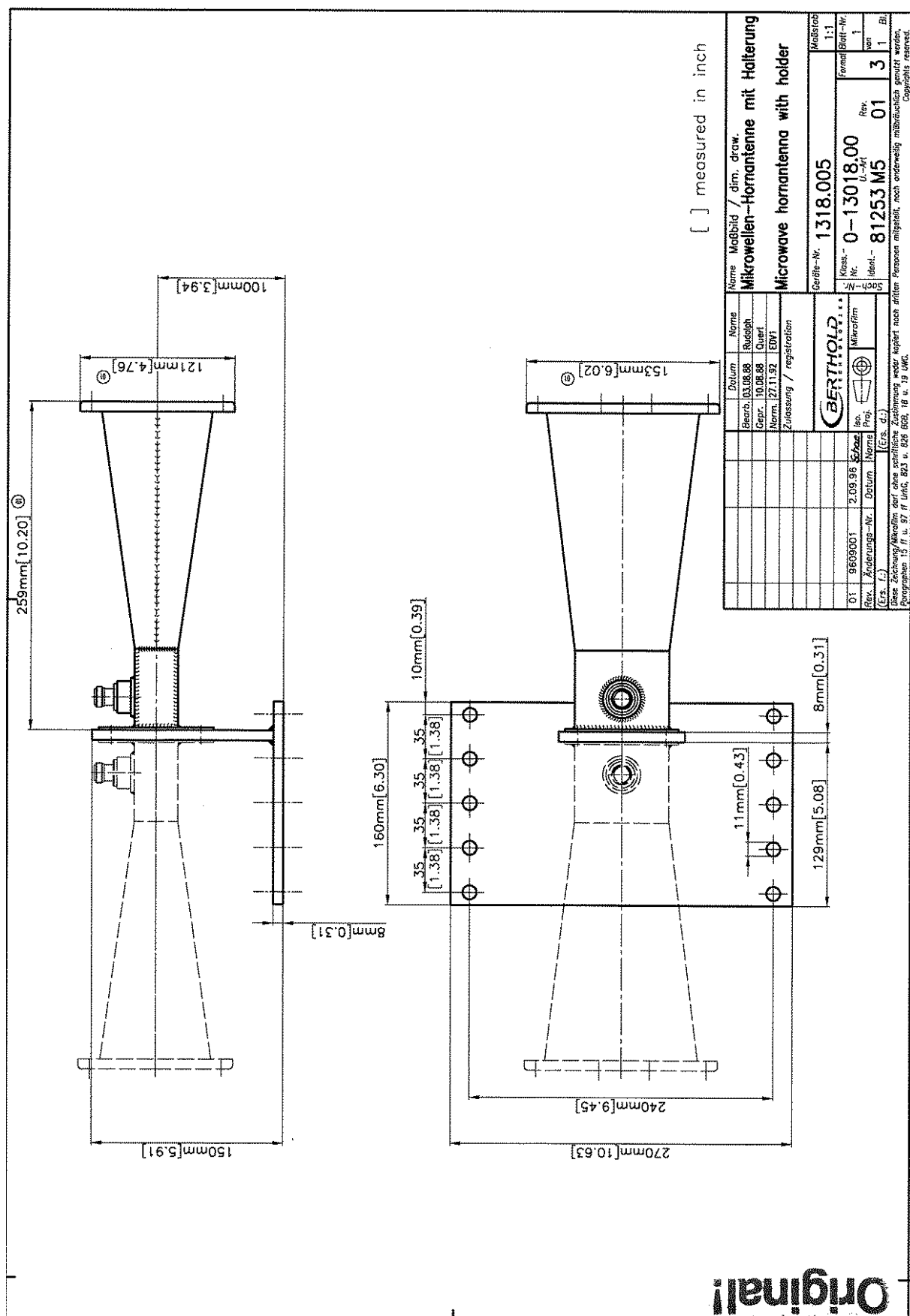
Horn Antenna and Horn Antenna Holder

Hornantenne und Hornantennenhalter



Name		Maßbild / dim. draw.	Name		Maßstab
Mikrowellen Hornantenne (Micro Moist)			Mikrowellen hornantenna (Micro Moist)		1:1
Geräte-Nr. 1318.004			Klass.-Nr. 0-13018.00		1
Revis.-Nr. 81253 M4			Revis.-Nr. 01		3
Zeich.-Nr. 01			Zeich.-Nr. 01		1
Datum 24.11.87			Datum 24.11.87		
Gepr. 28.06.88			Gepr. 28.06.88		
Norm. 27.11.92			Norm. 27.11.92		
Zulassung / registration			Zulassung / registration		
Berthold			Berthold		
01 9603001 2.09.96			01 9603001 2.09.96		
Revis.-Nr. 01			Revis.-Nr. 01		
Datum 24.11.87			Datum 24.11.87		
Gepr. 28.06.88			Gepr. 28.06.88		
Norm. 27.11.92			Norm. 27.11.92		
Zulassung / registration			Zulassung / registration		

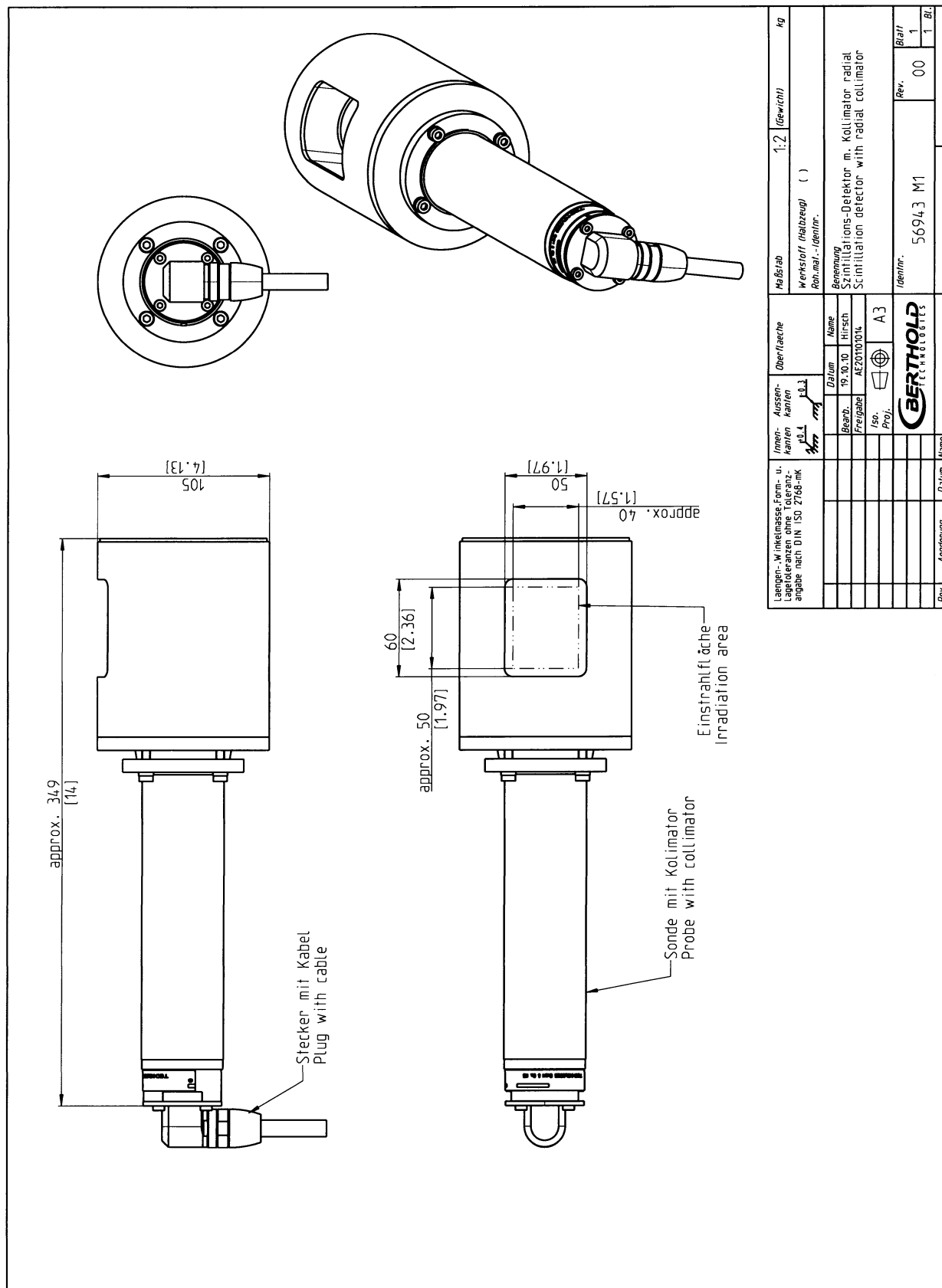
Original!



Original!

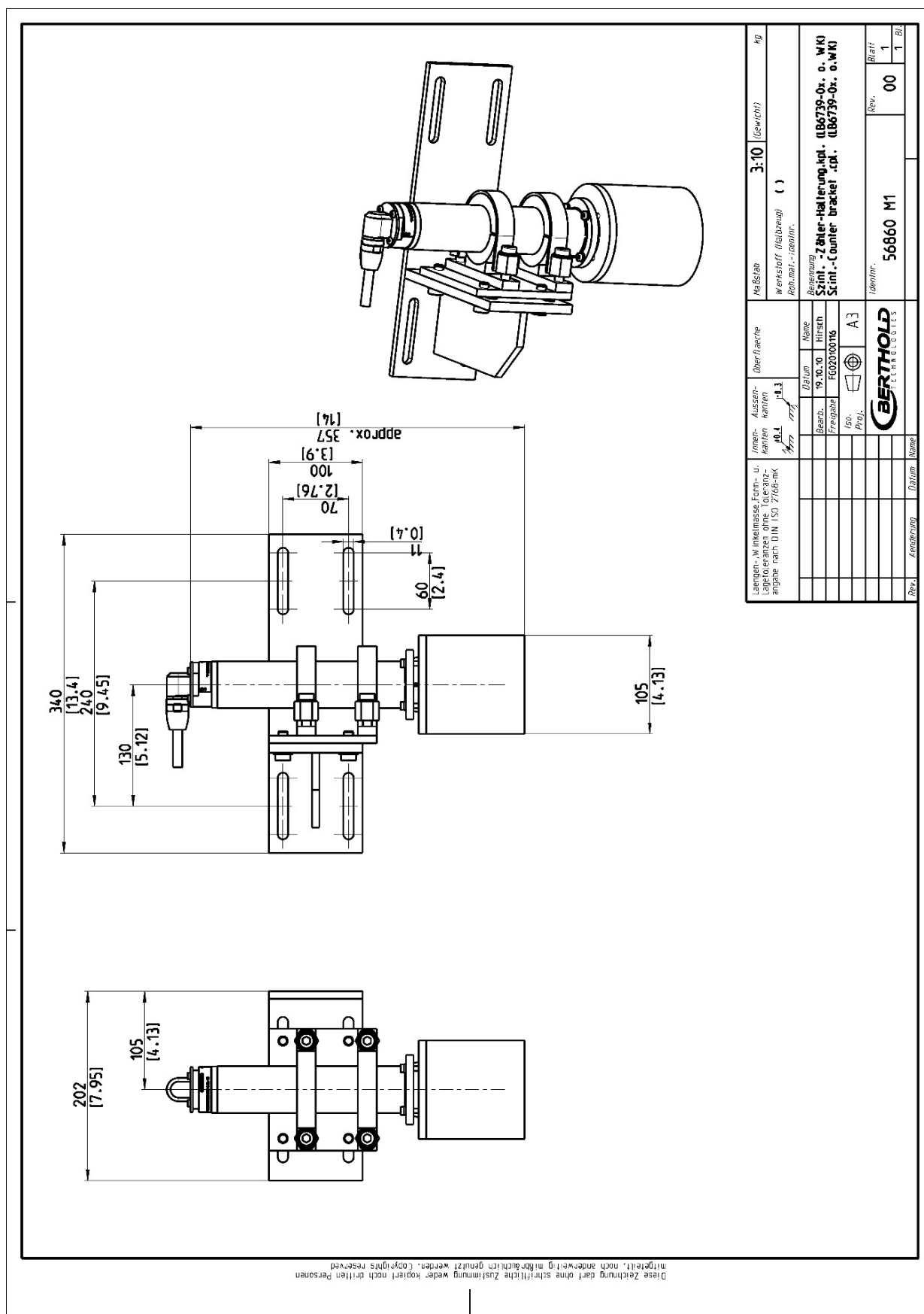


Scintillation Counter with Radial Collimator Szintillationszähler, Kollimator radial



Diese Zeichnung darf ohne schriftliche Zustimmung weder geändert noch kopiert werden. Copyrights reserved

Scintillation Counter with Bracket *Szintillationszähler mit Halterung*



Installation Proposal at the Conveyor Belt Einbauvorschlag am Förderband

