

SENSseries LB 480

COMPACT FIELD DEVICE

Various detector types

The optimal solution for every application.

Reliable & monitored

Monitored current output, continuous self-monitoring.

Easy handling

Quick Start for effective and fast commissioning. Operation via HART, PDM or AMS/DeltaV.

Safety

Safe for critical processes due to SIL2, with homogeneous redundancy SIL3.

Robust & resistant

Corrosion-resistant stainless steel housing and extremely robust design.

Explosion protection

Protection at the highest level with all common Ex approvals.



TECHNICAL DATA & FACTS

SENSseries LB 480

Detector

Applications	Density, Level, Level Switch
Power Supply	100 ... 240 VAC, 50/60 Hz, max. 12 VA 24 VDC (18 ... 32 VDC), max. 12 W
Housing Material	Stainless steel ISO 1.4301/AISI 304 or ISO 1.4404/AISI 316
Measuring range point detector	25/25 ... 150/150 mm
Measuring range rod detector	500 ... 8000 mm, cascading of up to 17 detectors possible
Ambient temperature	–40 ... +60 °C (–40 ... +140 °F) –40 ... +100 °C (–40 ... +212 °F) due to additional water cooling Observe possible temperature restrictions for explosion protection!
Temperature stability	≤ 0.002 %/°C (–40 ... +60 °C (–40 ... +140 °F)) for CrystalSENS ≤ 0.01 %/°C (–40 ... +60 °C (–40 ... +140 °F)) for SuperSENS and UniSENS ≤ 0.02 %/°C (–40 ... +60 °C (–40 ... +140 °F)) for TowerSENS
Water cooling for higher temperature	optional
Collimator for background suppression	optional
Outputs	Analog output 4 ... 20 mA potential-free, max. impedance 500 Ω, active or passive Digital output Open Collector potential-free, max. 100 mA at 5 ... 35 VDC for alarms, warning and error messages
Inputs	Analog signal input PT100 for temperature compensation (only density)
Interfaces	RS 485 for software update, cascading and compensation
IP protection	IP65, IP66, IP67, IP68, IP69K
Explosion protection	ATEX, IECEx, NEC/CEC, INMETRO, EAC
Functional safety	SIL2, SIL3

Special Features

- X-ray Interference Protection (XIP)
- Gas Property Compensation (GPC)
- Product Radiation Compensation (PRC)
- Fast measurement sample rate (Speedstar)
- Source Life Management
- Continuously monitored current output
- Quick start wizard