

# LoopSeries LB 430

## ENERGY EFFICIENT, LOOP POWERED

### Remote Display

Optional display module for process values and status information.

### Different types of detectors

Optimal solution for every application and all measuring ranges

### Innovative

SiPM technology with high mechanical and electrical robustness.

### Reliable

Extensive self-diagnostics suited for preventive maintenance.

### Sustainable

Loop-powered device with low power consumption.

### Protected

Protection at the highest level with all common Ex approvals.



# TECHNICAL DATA & FACTS

## LoopSeries LB 430

### Detector

---

Applications	Level, Level Switch, Density
Measuring range point detector	50 x 50 NaI
Measuring range rod detector	500 to 2000 mm
Power supply	14 to 28 V DC current loop 4 to 20 mA
Housing material	Stainless steel ISO 1.4404/AISI 316L or Aluminium
Ambient temperature	–40 to +60 °C (–40 to +140 °F) –40 to +100 °C (–40 to +212 °F) with water cooling Possible temperature restrictions for explosion protection
Temperature stability	≤ 0.005 %/°C (–40 to +60 °C (–40 to +140 °F)) for CrystalSENS ≤ 0.02 %/°C (–40 to +60 °C (–40 to +140 °F)) for UniSENS
Max. count rate	2,000,000 cps
Water cooling for higher temperature	optional
Collimator for background suppression	optional
Analog output	4 to 20 mA potential-free, HART, max 500 Ω, passive
IP protection	IP66 (NPT 1/2"), IP68 (M20)
Explosion protection	ATEX, IECEx

### Special Features

---

- X-ray Interference Protection (XIP)
- Continuously monitored current output
- Multiple software wizards for quick and easy configuration and device operation
- Extensive self-diagnostics conforming to NE107
- Easy wiring due to loop-powered technology
- Up to 99% less energy consumption compared to externally powered devices
- Optional display module
- Software includes all applications