

LoopSeries LB 430

ENERGY EFFICIENT, LOOP POWERED

Innovative

SiPM technology with high mechanical and electrical robustness.

Sustainable

Current loop powered device with low power consumption.

User friendly

Quick and easy configuration/operation with software wizards.

Remote Display

Optional display module for process values and status information.

Reliable

Extensive self-diagnosis suited for preventive maintenance.

Protected

Protection at the highest level with all common Ex approvals.



TECHNICAL DATA & FACTS

LoopSeries LB 430

Detector

Applications	Density, Level, Level Switch
Power supply	14 ... 28 V DC current loop 4 ... 20 mA
Housing material	Stainless steel ISO 1.4404/AISI 316L or Aluminium
Ambient temperature	–40 ... +60 °C (–40 ... +140 °F) –40 ... +100 °C (–40 ... +212 °F) with additional water cooling Observe possible temperature restrictions for explosion protection!
Temperature stability	≤ 0.5 %/°C (–40 ... +60 °C (–40 ... +140 °F))
Max. count rate	2 000 000 cps
Water cooling for higher temperature	optional
Collimator for background suppression	optional
Analog output	4 ... 20 mA potential-free, HART, max 500 Ω, passive
IP protection	IP65, IP66, IP67, IP68
Explosion protection	ATEX, IECEx, NEC/CEC, INMETRO

Special Features

- X-ray Interference Protection (XIP)
- Continuously monitored current output
- Multiple software wizards for easy configuration and device operation
- Extensive self-diagnosis in conformity to NE107
- Easy wiring due to loop-powered technology
- Power savings up to 99% compared to external powered devices
- Optional display module
- All applications included in one software