

# LoopSeries Display LB 4310

## REMOTE PROCESS MONITORING

### Remote mounting

Up to 20 m flexible display positioning for hard-to-reach measuring points.

### Rugged design

IP-certified housing for extreme environmental conditions.

### Protected

Protection at the highest level with all common Ex approvals.

### Plug & Play

Quick and easy installation/retrofitting.

### Power from the detector

No separate power supply required.

# TECHNICAL DATA & FACTS

## LoopSeries Display LB 4310

### Detector

---

|                      |                                                                                                                         |
|----------------------|-------------------------------------------------------------------------------------------------------------------------|
| Applications         | Density, Level, Level Switch                                                                                            |
| Power supply         | Via LoopSeries LB 430 detector                                                                                          |
| Housing material     | Stainless steel ISO 1.4404 / AISI 316L or Aluminium                                                                     |
| Parameters           | - Condensed status according to NE107<br>- Process value<br>- Count rates<br>- Current output<br>- Detector temperature |
| Ambient temperature  | -40 to +60 °C (-40 to +140 °F)<br>Possible temperature restrictions for explosion protection                            |
| IP protection        | IP66 for cable entry 1/2" NPT thread<br>IP68 for cable entry M20                                                        |
| Explosion protection | ATEX, IECEx, NEC/CEC, INMETRO, EAC (all pending)                                                                        |

### Special Features

---

- Supports self-diagnostics conforming to NE107
- Easy and intuitive installation, retrofit possible
- Remote wired connection to LB 430 LoopSeries detector (max. 20 m)