

**New!**

# SIL Approved Radiometric Detectors

## SENSseries LB 480

The radiometric detectors SENSseries LB 480 from Berthold Technologies have been awarded SIL2 / SIL3 certification. The certificate covers the measurement applications level switch, continuous level and density. Thus Berthold Technologies is the first supplier providing SIL approved radiometric measurement systems for both level and density applications.

SIL2 certified for  
level switch,  
continuous level  
and density

intensively tested  
and SIL approved by  
TÜV Süd according  
to IEC 61508:2010

SIL3 operation in  
redundancy  
installations (1oo2)

also for cascaded  
measuring systems



[ SIL 2 ]

[ SIL 3 ]

