

Mini-Switch LB 471 / DuoSeries LB 4700

2-WIRE LEVEL SWITCH MEASURING SYSTEM

Flexible operation

Switch between standard and professional mode for additional functions.

Space-saving & cost-optimised

Cost-effective and compact solution for monitoring minimum and maximum levels.



Reliable & safe

Permanent error- and self-monitoring.



Robust & resistant to interference

Optional lead collimator protects against background radiation. Corrosion-resistant stainless steel housing and extremely robust design.

TECHNICAL DATA & FACTS

Mini-Switch LB 471 / DuoSeries LB 4700

Transmitter Mini-Switch LB 471

Applications (type)	Level switch
Variants	Cassette (3U, 7HP) or plug-in card (3U, 4HP)
Power supply	18 to 30 V DC or 24 V AC +10% to -15% Option: 115/230 V AC transformer unit +10% to -15%, 47 to 65 Hz
Power consumption	max. 4 W per module
Ambient temperature	-30 to +60 °C (-22 to +140 °F)
Installation	19" rack or wall housing (IP66)
Outputs	Digital
	1 relay (SPDT) für min./max. alarm 1 relay for collective error message 1 relay for warning message, alternatively usable as min./max. alarm
	Permissible load: AC: max. 250 V, max. 1 A, max. 200 VA DC: max. 300 V, max. 1 A, max. 60 W Induction-free
Inputs	Digital
	for external empty adjust
Weight	0.3 kg per module
Detector connection	2-wire
IP protection	IP20, IP66 (depending on installation solution)

Detector Type LB 4700

Power supply	Power supply via evaluation unit
Housing material	Stainless steel 1.4301 / AISI 304 (others on request)
Measuring range	CrystalSENS: 25/25, 40/35, 50/50, 50/60 mm SuperSENS: 150/150 mm
Ambient temperature	-40 to +60 °C (-40 to +140 °F)* -40 to +100 °C (-40 to +212 °F) due to additional water cooling*
Temperature stability	≤ 0.002 %/°C (-40 to +60 °C (-40 to +140 °F)) for CrystalSENS ≤ 0.01 %/°C (-40 to +60 °C (-40 to +140 °F)) for SuperSENS
Water cooling	Optional
Collimator for background suppression	Optional
IP protection	IP66, IP67
Explosion protection	ATEX, IECEx, NEC/CEC, INMETRO, EAC, KCs

Special features

- Interference protection
- Switching between standard and professional mode
- Automatic calculation of full count rate and time constant

*Observe any temperature restrictions for explosion protection!

Berthold Technologies GmbH & Co. KG

Calmbacher Straße 22 · 75323 Bad Wildbad · Germany
+49 7081 1770 · industry@berthold.com · www.berthold.com

