

# LB 147 / LB 148

Portable Contamination Monitors



 **BERTHOLD**

# PERSONAL CONTAMINATION MONITORS

## Equipment Concept

The Personal Contamination Monitors LB 147 and LB 148 are contamination monitors designed for use in the radiation protection field to detect contamination caused by alpha and beta/gamma radiation. Low-maintenance thin-layer ZnS (Ag) scintillators are used as detectors. The user can interact with the system via a touch panel on a graphical display.

The LB 148 version has four automatically tracked hand detectors that allow both the palms and the back of the hands to be measured simultaneously. With the LB 147 version, an automatic prompt for back of hand measurement can be activated. The LB 147 is also available in a narrow version LB 147 Slim with a different positioning of the hand detectors. There are two variants: the detectors are facing the user (ID 45356-15) or the detectors are facing away from the user (ID 45356-18). The LB 147 and LB 148 are also available as an Alpha Sense version with special modification of the protective grid for lower  $\alpha$ -detection limits. All monitor versions have a detachable probe for clothing measurements.

Each detector has its own calibration factor for each nuclide and its own spill-over factor for each alpha nuclide. More than 50 nuclides with calibration factors according to DIN ISO 7503-1 and A-100 are already stored in the software.

The devices can be operated as monitors for simultaneous  $\alpha$ - and  $\beta/\gamma$ -measurement or only as pure  $\alpha$ - or pure  $\beta/\gamma$ -monitors, whereby the measurement results can be displayed in the unit Bq/cm<sup>2</sup> or cps. The detection limits and characteristic limits are calculated for each measured value in accordance with DIN ISO 11929 and saved together with the measured values.

The ambient radiation is continuously checked separately for each detector and compensated for as a long-term average value during the contamination measurement. The monitors also check for fluctuations in the background immediately before starting the measurement.

The menu navigation is clearly laid out and easy to operate. The built-in power supply unit automatically adapts to the various AC voltage supplies.



LB 147 with signal tower



LB 147 SLIM with averted detectors

## Functions and Accessories LB 147 and LB 148

Numerous service functions permit easy measurements for necessary periodical tests. This includes automatic calibration routines for all detectors and a fast system test. All in- and outputs can be checked very quickly by means of further service functions.

Users can be registered using an optional transponder or magnetic card reader. The HFK programme (optional) can be used to evaluate the personal measurement data and record it in a database.

An optional signal tower can be connected via one of the USB (host) interfaces or via the three relays built into the devices. A door control unit can also be connected here with configuration options in the software.

A remote parameter programme is available for configuring and documenting the monitors. In addition to direct communication, the application can also be used offline to create or edit setting files. The monitors have the following interfaces: 5 USB

(host), 1 USB (device), 1 Ethernet 10/100/1000 Mb, 1 RS232 and 1 RS485. In addition to signal towers, transponders, magnetic card readers or printers can also be connected to the USB (host) interfaces (optional).

Up to 10.000 measurement data can be stored and transferred to external computers via the USB or Ethernet interface. Alternatively, the serial interface can also be used. A comprehensive communication protocol allows integration into a measurement network.

Reading out the measurement data and carrying out software updates on the monitors is quick and easy via one of the five USB (host) interfaces.



LB 148 with detectors for back of hands

# TECHNICAL SPECIFICATIONS

## Technical data LB 147 / LB 148

Detectors		A: LB 147/LB 148, B: Alpha Sense	
Radiation detector	ZnS(Ag) scintillator		
Material entrance window	2 µm x 3 µm metallized, plastic (0.4 mg/cm²)		
Hand Detectors			
Dimensions entrance window	150 mm x 230 mm		
Sensitive area	345 cm²		
Zero effect	α-channel about 0.1cps β-γ-channel about 15cps		
Typ. Efficiencies (according to ISO 7503-1)		A	B
	<sup>241</sup> Am	33%	43%
	<sup>14</sup> C	20%	25%
	<sup>36</sup> Cl	49%	70%
	<sup>90</sup> Sr/ <sup>90</sup> Y	52%	66%
Typ. MDA (according to ISO 11929)	<sup>241</sup> Am	0.026	0.019
	<sup>14</sup> C	0.649	0.460
with 10s measuring	<sup>36</sup> Cl	0.135	0.078
time in Bq/cm²	<sup>90</sup> Sr/ <sup>90</sup> Y	0.126	0.088
Foot Detectors			
Dimensions entrance window	150 mm x 370 mm		
Sensitive area	555 cm²		
Background	α-channel ca. 0.2 cps β-γ-channel ca. 40 cps		
Typ. Efficiencies (according to ISO 7503-1)		A	B
	<sup>241</sup> Am	19%	40%
	<sup>14</sup> C	21%	21%
	<sup>36</sup> Cl	54%	71%
	<sup>90</sup> Sr/ <sup>90</sup> Y	43%	64%
Typ. MDA (according to ISO 11929)	<sup>241</sup> Am	0.055	0.026
	<sup>14</sup> C	0.954	0.961
with 10s measuring	<sup>36</sup> Cl	0.191	0.137
time in Bq/cm²	<sup>90</sup> Sr/ <sup>90</sup> Y	0.254	0.153
Ambient Conditions			
Temperature range	- 5 °C bis + 40 °C		
Rel. humidity	0% bis 90%, non condensing		

Mechanical Data	
LB 147/Alpha Sense	about 65 x 125 x 80 cm³ (W x H x D)
LB 147 Slim	about 45 x 125 x 90 cm³ (W x H x D)
LB 148	about 85 x 125 x 82 cm³ (W x H x D)
LB 147/Slim/Alpha Sense	approx. 35 kg
LB 148/Alpha Sense	approx. 45 kg
Electronics	
Display	Colour 7" Touch Panel, 800 x 480 Pixel
Ports	5 USB(Host), 1 USB(Device), 1 Ethernet 10/100/1000 Mb, 1 RS 232, 1 RS 485
External signal transmitters	3 relays max. 50 V, 5A
Power supply	100 - 240 VAC, wide-range input
Power consumption	approx. 14.5 W, Fuse 2A
Systems	
LB 147	45356-10
LB 147 Slim	45356-15 (with facing detectors)
LB 147 Slim	45356-18 (with averted detectors)
LB 147 Alpha Sense	45356-16
Accessories (optional)	
	Ident. No.
Mobile set (only for LB 147)	52874
Transponder system (RFID)	66164
Chip/Card (RFID)	59495/59503
Magnetic card reader	66442
Magnetic card	34481
Signal tower for LB 147 (red/green)	65252
Signal tower for LB 148 (red/green)	65252
Calibration source <sup>36</sup> Cl/ <sup>241</sup> Am	29336/25509
HFM program	42849-042
Stomach and thyroid probe*	49950
Stomach probe*	53516
Thyroid probe*	53517
*only for LB 147, only for export	
Subject changes without prior notice.	

## TRANSFORMING SCIENCE INTO SOLUTIONS



Experience and expertise are of great importance to be able to ensure safety-relevant measurements properly and reliably. With more than 70 years of experience in planning and design, installation and commissioning, calibration, documentation and service of radiation protection measurement systems, we continue to support our customers in their task to continuously optimize their work processes and to ensure the safety of the environment and personnel.

### Berthold Technologies GmbH & Co. KG

Calmbacher Straße 22 · 75323 Bad Wildbad · Germany  
+49 7081 1770 · nuclear@berthold.com · [www.berthold.com/rp](http://www.berthold.com/rp)

© Berthold Technologies. All rights reserved. All trademarks are the property of Berthold Technologies or their respective owners. Berthold Technologies reserves the right to implement technical improvements and/or design changes without prior notice.