

Technical Data Sheet

Ionisation Chamber LB 6701M-H10

Application

Dose rate probe for photon radiation in Health Physics applications.

Measured Quantity

Ambient dose equivalent $H^*(10)$ or

Ambient dose rate equivalent $\dot{H}^*(10)$

Construction

The Ionisation Chamber is made from Aluminium with a Nitrogen gas filling at 10 bar and a radiation resistance up to 10^6 Gy. The chamber current is proportional to the dose rate, this current is converted into a +11V Norm pulse frequency in the Current/Frequency converter LB3856-23.

The Current/Frequency converter is mounted in a separate housing LB6703 which also contains the High Voltage module to operate the Ionisation Chamber.

A ^{90}Sr check source with 25 kBq activity is built in the Ionisation Chamber to continuously monitor the proper functioning of the system.

Technical Data

► **Measuring Range**

0,1 mSv/h – 100 Sv/h

► **Energy Range**

45 keV – 1,3 MeV
with regard to Cs-137 and 0°

► **Calibration Factor**

1 mSv/h per cps

► **Output pulse I-F-Converter**

Polarity:	positive
Amplitude:	+ 4,5 V in 50 Ω
Pulse Width :	2 - 5 μs
Frequency range:	0,1 Hz – 100 kHz

► **High Voltage**

1000 Volt

► **Operating Conditions**

Temperature:	0 to + 50°C
Rel. humidity:	20 to 80 %
Storage temp.:	max. 60°C

► **Protection Degree**

IP 65

► **Dimensions**

- Ionisation chamber:	80 mm Ø x 268 mm
- I-F-Converter	
Connection box:	240 x 160 x 90 mm

