

Product Information

LB 126

portable

Dose and Doserate Monitor



Handheld monitor for fast and reliable measurements of ambient dose equivalent $H^*(10)$ in photon radiation fields

Technical Data:

- Detector: energy compensated proportional counter tube LB 6006-H10
- Measuring range: 50 nSv/h to 50 mSv/h
- Energy range, polar response: 30 keV to 1,3 MeV, $\pm 45^\circ$
- Selectable Units: $\mu\text{Sv/h}$, μSv or mrem/h mrem
- free selectable alarm thresholds for dose and dose rate
- Microprocessor electronics with charging circuit for 4 rechargeable batteries (AA cells)
- Data storage for 1000 measurements
- Real time clock, battery-buffered
- Signal and alarm indication
- RS485 networking for up to 32 devices
- Temperature range: -10 bis +40 °C

