

Arium® Water Guard

Early detection of leaks
protects the laboratory

Benefits

- Highly sensitive optical sensor
- Audiovisual alarm signals
- Automatic water stop in the case of leakage
- High-quality material, no corrosion
- Easy to install



Product Information

Only the early detection of water leaks offers optimal protection against water damage in the laboratory. However, this is only possible to a limited extent with classic water monitors, which register a leak exclusively via the conductivity of the escaping water, as they may not detect a measurement if the conductivity is too low.

The Arium® water monitor, on the other hand, identifies water leaks using an optical sensor and is therefore able to reliably detect both pure and ultrapure water. If the optical sensor registers a leak, it automatically shuts down the water supply. To alert the user to a leak, the triggering of the sensor is accompanied by an acoustic and a visual signal.

Technical Specifications

Sensor dimensions

Diameter	5 cm
Height	2.5 cm
Cable length	2 m

Tubing connections

Input	3/4" External thread
Output	3/4" External thread
Power supply	100 – 240 VAC 50 – 60 Hz

Ordering Information

Order Number	Description
H2O-AWG-2	Arium® Water Guard, HF 1 pc

Intended Use


Device type: Arium® Advance L

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com