

Arium® Level Sensor

Practical filling of separate tanks

- Flexible water transport to any location
- Any tank system can be filled

Description

The level sensor makes it easy to connect an external water storage tank and subsequently fill a tank with ultra-pure water.



Technical Specifications | Ordering Information

Level sensor length	88 mm
Connection diameter	2.03 cm (max.)
Drill hole	1.65 cm
Cable length	3 m

Order number	Description
H2O-ALS1	Arium® Level Sensor, 1 pc

Intended Use

Device type:

- Arium® Pro DI, Pro UF, Pro UV, and Pro VF