

# Arium® Extended Connection Set

## Flexible enhance your working space

- Extend the operating range up to 4 meter
- Fast and reliable connection

## Description

The extended tubing set is used for the Arium® Smart Station Pure for uncomplicated extension of the standard supplied tubing set.

With the extension set the operating range can be easily increased from 2 up to 4 meter.



## Technical Specifications | Ordering Information

### Materials

Cable	Ethernet CAT6A
Tubing	¼" Polyethylene (PVDF   PE)

### Length

Operating range	4 meter
-----------------	---------

### Order number

H2O-ATES-UP	Arium® Tubing Extension Set Ultrapure
H2O-ATES-P	Arium® Tubing Extension Set Pure

### Description

Benchtop and wall-mounted edition can be assembled for left or right hand side, without additional equipment required.

### Intended Use

Device type:

- Arium® Smart Station Ultrapure
- Arium® Smart Station Pure