

Arium® Smart Station

Flexible Remote Dispenser

Advantages

- Compact: Save space integrating in your lab
- Intuitive: touch-activated color display with direct access to all important functions
- Flexible: Stepless height adjustment to fill different size containers
- Ergonomic Design: Hand-held for optimized left or right-hand operation
- Accurate: Precise volume dispense for reliable buffer and sample preparation

Product Description

The Arium® Smart Station is designed for flexible remote dispensing of pure and ultrapure water directly at the point of use.

While dispensing water into a broad range of different sized containers, the Smart Station offers constant control of all important quality parameters, at all times.

The ergonomic design supports left and right hand operation and can be easily adjusted to your need.



Compact

The Arium® Smart Station will readily fit under almost every standard laboratory shelf. This handy device will give the flexibility you need to integrate it into any lab environment and save space.

Accurate

Optimize your sample preparation and benefit from direct access to all important dispensing functions: manual, volume- controlled or predefined favorite volumes.

The various dispensing functions offers precise dispensing from 2 L/min up to manually controlled drop-by-drop* function.

Choose the Smart Station of Your Choice

The Smart Station is designed to be used for ultrapure and pure water dispense connected to your Arium® Ultrapure water system or Bagtank:

Type of System	Use with
Arium® Smart Station Ultrapure	Direct connection to Arium® Pro or Arium® Comfort supplying ultrapure Type 1 water
Arium® Smart Station Pure	Direct connection to Bagtank supplying pure water

*Depends on the connected system type

Intuitive

Easy to operate the Arium® Smart Station using the touch-activated color display, even when you are wearing laboratory gloves. The icons will guide you through the menu for intuitive, error-free operation.

Flexible

The stepless height adjustment from 8.5 – 65 cm allows you to fill different sized containers. Thanks to its angular designed stand even bulky vessels can be optimally positioned under the dispense outlet. Depending on your needs you can adjust the volume dispense from 50 mL up to 50 Liter.



Supply ultrapure water from Arium® Pro and Arium® Comfort



Supply pure water from Arium® Bagtank

Adapt the Dispense Point to Your Needs

To optimally fit the Smart Station into your lab environment there are two installation options available: bench top or wall-mounted.

Both options can be individually adjusted to left or right hand side operation, without any additional required accessories.



Flexible Dispensing

Connect up to three Smart Stations to your single ultrapure water system to serve different dispensing areas. By using the extended connection set, the distance between each Smart Stations can be set up to 4 meters. Adapted to your application you could select between different point of use filters.



Technical Specifications

Smart Station Bench-Top

Dimensions:	
Control box with stand (w × d × h)	213 × 213 × 598 mm (8.4 × 8.4 × 8.2")
Operating range fixed dispense arm (d × w × h)	428 × 476 × 835 mm (16.9 × 18.7 × 32.9")
Tubing Length:	
Distance to water system/bagtank	2 Meter
Operating range flexible hand held	0.7 Meter
Weight	Approx. 4.9 kg (10.8 lbs)

Smart Station Wall-Mounted

Dimensions:	
Control box (w × d × h)	172 × 157 × 343 mm (6.8 × 6.2 × 13.5")
Operating range dispense arm (d × w × h)	242 × 90 × 300 mm (9.5 × 3.5 × 11.8")
Tubing Length:	
Distance to water system/bagtank	2 Meter
Operating range flexible hand held	0.7 Meter
Weight	Approx. 2.4 kg (5.3 lbs)

General Specifications

Volume-controlled dispensing	50 mL - 50 Liter
Volume accuracy	± 5 % between
Power supply	100 - 240 VAC; 50 and 60 Hz, 2.5 A (max.) 2°C - 40°C
Power cord (IEC 60320-1 / C14)	Country specific

Ordering Information

Order number	Description
H2O-ARST-UP-T	Arium® Smart Station Ultrapure for benchtop installation
H2O-ARST-UP-B	Arium® Smart Station Ultrapure for wall-mounted installation
H2O-ARST-P-T	Arium® Smart Station Pure for benchtop installation
H2O-ARST-P-B	Arium® Smart Station Pure for wall-mounted installation

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

Consumables

Arium® Sterile Plus

For Sterile and Particle-Free Water Dispensing

- Excellent total throughput and flow rates
- Integrity-tested
- Validated according to HIMA and ASTM F-838-05
- Meets WFI quality standards compliant with USP, incl. USP Plastic Class VI Test
- Manufactured according to EN ISO 9001
- Easy to install
- Certified quality



Description

The Arium® Sterile Plus (Sartopore® 2 150) is a sterile, ready-to-use membrane filter capsule designed to meet the highest requirements.

The filter capsules contain a hydrophilic, heterogeneous polyethersulfone double-layer membrane, delivering excellent total throughput and flow rate performance.

The capsule is attached by a quick-connect coupling to the point of use and reliably removes particles and micro-organisms > 0.2 µm in the last water purification step.

The pleated Sterile Plus membrane filter units are validated as sterilizing-grade filters for biopharmaceutical use according to HIMA and ASTM F-838-05 guidelines (documents available). Each capsule is integrity-tested during the manufacturing process to ensure it meets the highest quality standards and safety regulations.

Technical Specifications

Materials

Membrane	Asymmetrical polyethersulfone
Filling bell assembly	Polycarbonate
Other plastics	Polypropylene

Typical Specifications for Purified Water

Flow rate (depends on the inlet pressure and the type of system)	Up to 2.0 L/min.
Bacteria	< 0.001 CFU/mL
Particle content	No particles > 0.22 µm

General Specifications

Pore sizes	0.45 µm + 0.2 µm
Filtration area	0.015 m ²
Inlet and outlet	¼" plug-in connector
Sterilization (3 cycles max.)	Autoclavable at 134°C, 1 bar (14.5 psi), 30 min.
Max. diffusion	1 mL/min. at 2.5 bar (36 psi)
Min. bubble point	3.2 bar (46.4 psi)

Ordering Information

Order Number	Description
5441307H4--CE	Arium® Sterile Plus (Sartopore® 2 150 capsule), 0.2 µm pore size; qty. per pkg.: 1

Intended Use

Device Type:

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

Arium® Cell Plus Ultrafilter

For Effective Removal of Endotoxins in Cell Culture Applications

- Effective removal of RNase | DNase
- Reliable removal of endotoxins
- High flow rate performance
- Certified quality
- Sterile-packaged



Description

The Arium® Cell Plus is a point-of-use ultrafilter for the effective removal of small molecules at 15kDa cut off.

Attached with the quick-connect coupling to the point of use, the ultrafilter provides the highest safety for your critical cell culture and life science applications by removing high efficiently RNase | DNase and Endotoxins.

A protective bell supplied with the ultrafilter additionally prevents retrograde contamination.

Moreover, the high-quality material selected for Arium® Cell Plus enables excellent total throughputs and optimal flow rates.

Technical Specifications

Materials	
Membrane	Polysulfone
Composite material	Polyurethane (PUR)
Housing	Acrylonitrile butadiene styrene (ABS)
Protective bell	Polycarbonate (PC)

Typical Specifications for Purified Water

Flow rate (depends on the inlet pressure and the type of system)	Up to 2.0 L/min.
Endotoxins	< 0.001 EU/mL
Bacteria	< 0.01 CFU/mL
RNase concentration	< 1 pg/mL
DNase concentration	< 5 pg/mL

General Specifications

Cut-off	15.000 Daltons/0.005 µm
Inlet and outlet	¼" plug-in connector
Dimensions (height × diameter)	169 mm × 50 mm (6.6" × 1.9")
Max. operating pressure	6 bar (87 psi)
Max. inlet temperature	50°C
Effective membrane area	0.5 m ² (5.4 ft ²)

Ordering Information

Order Number	Description
H2O-CUF	Arium® Cell Plus; qty. per pkg.: 1

Intended Use

Device Type:

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

Accessories

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 Smart Station Ultrapure and 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

Materials

Cable	Ethernet CAT6A
Tubing	¼" Polyethylene (PVDF PE)

Length

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

Ordering Information

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

Intended Use

Device Type:

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

Sartorius Service

We Ensure the Quality of Your Results

At Sartorius, quality products go hand in hand with professional service. With our wide service offering, we will help guarantee the safe, reliable and optimal operation of your Arium® system. Just ask us and we will even cover the entire life cycle of your laboratory water system – from commissioning to qualification to regular maintenance. Together with you, we will ensure the consistently high quality of your laboratory water purification.



Our Services at a Glance

Installation and Commissioning

Your advantage: Your system will operate reliably at peak performance from day one

Equipment Qualification (IQ | OQ)

Your advantage: You will meet all regulatory requirements (GMP | GLP)

Regular Preventative Maintenance, Including Calibration, inspection and testing of your system and exchange of consumables

Your advantages: Optimal operation of your system; reliable results; prevention of downtime or even equipment failure

Get more information now at:


www.sartorius.com/service

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
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