



Arium® Smart Station

The Most Flexible Solution
for Dispensing Pure and
Ultrapure Water

Simplifying Progress

SARTORIUS

Arium® Smart Station

Water. Exactly Where You Need It.



- Compact design for easy and space-saving integration in your lab environment
- Intuitive touch-activated color display with direct access to all important functions
- Flexible and stepless height adjustment to fill different size containers and beakers
- Ergonomic design can be optimized for either right- or left-hand operation
- Accurate volume dispense for fast and precise buffer and sample preparation
- Adjustable, splash-free flow rate from drop-by-drop to 2 liters per minute

In many laboratories, space is at a premium. Especially when centralized laboratory water systems are required to dispense pure and ultrapure water for your applications. To maximize your lab space capabilities, you shouldn't have to adapt your water system to the laboratory. The system should adapt to your needs.

In addition to being economical in terms of its footprint, flexibility in dispensing water is essential. Users need full control over quality parameters at the point of use and the ability to fill different sized containers at the same time across the lab.

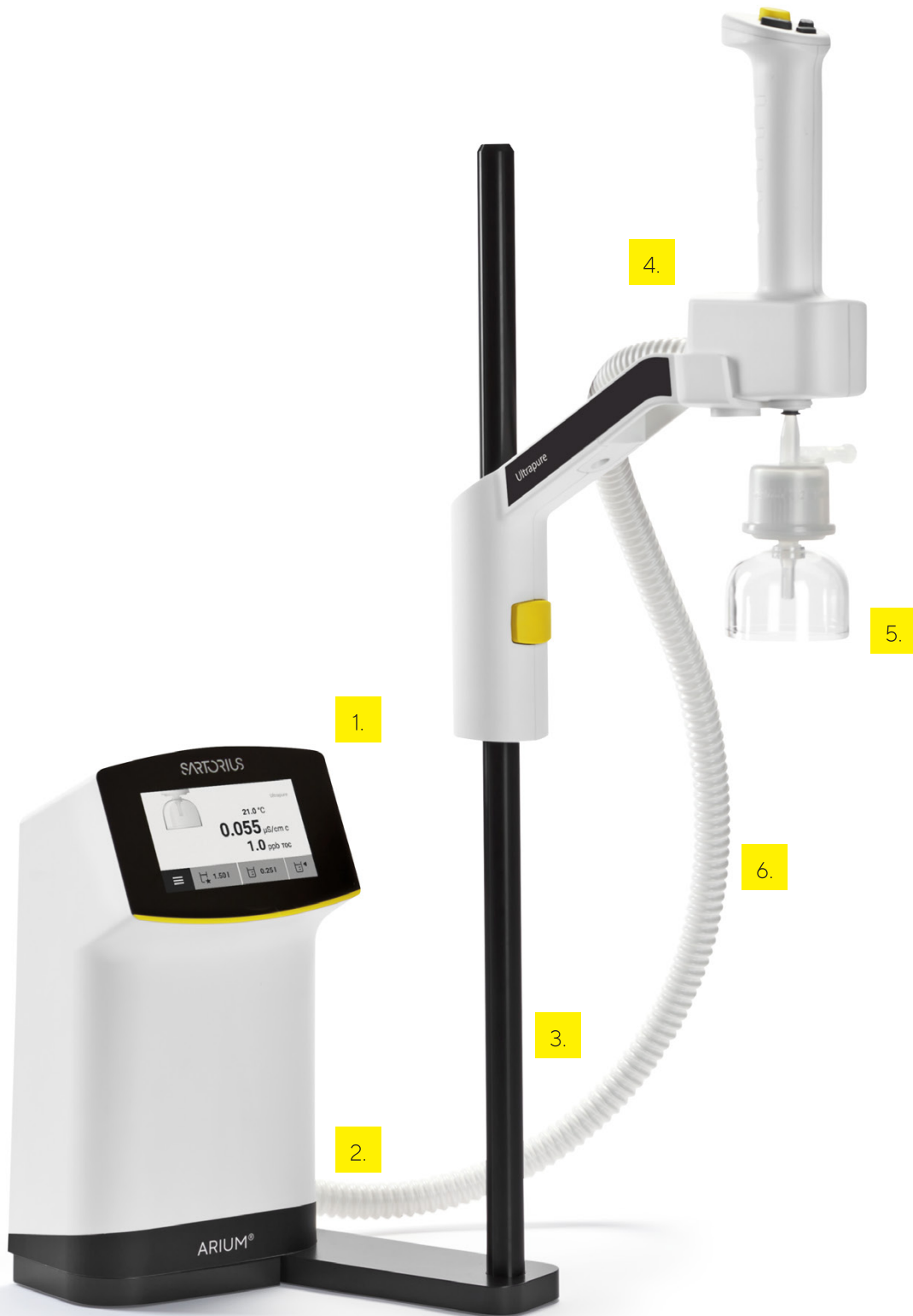
The innovative and flexible Arium® Smart Station from Sartorius addresses all of these requirements. You can dispense water, in the exact quality and quantities required for your experiments, when and where you need it.

Experience the new Arium® Smart Station and simplify your daily use of lab water.



Inspired by Your Application

The Arium® Smart Station is designed for ergonomic and flexible dispensing at the highest quality standards.





1. Full Functionality Touch Screen

- The control unit can be easily navigated using the 4.3 inch touch screen
- Displays all relevant information allowing full control of the system

2. Minimal Space Requirement

- The control unit has a small footprint with flexible positioning
- Options for bench top or wall-mounting

3. High Quality Components for Reduced Risk

- Smart recirculation ensures the highest ultrapure water quality
- Media separated valve to avoid cross contamination

4. Unbeatable Ergonomic Design for Complete Flexibility

- Designed by a team with extensive experience in ergonomics
- The dispense arm can be assembled for left and right hand operation
- Easy to fix or release the handheld from the Smart Station by magnetic connection

5. Application-Specific Filters to Meet All Your Research Needs

- Easy to exchange point of use filter due to quick connector
- Install up to 3 Smart Stations to your ultrapure water system to serve different applications.

6. Flexibility to Fill Various Vessels

- Stepless, one-hand adjustment of dispenser arm from 8.5 to 65 cm
- Dispenser arm rotates up to 360°
- Enables filling of sample tubes and bottles from 1.5 ml up to 5 L
- Easy volume dispensing including soft start | stop function and drop-by-drop

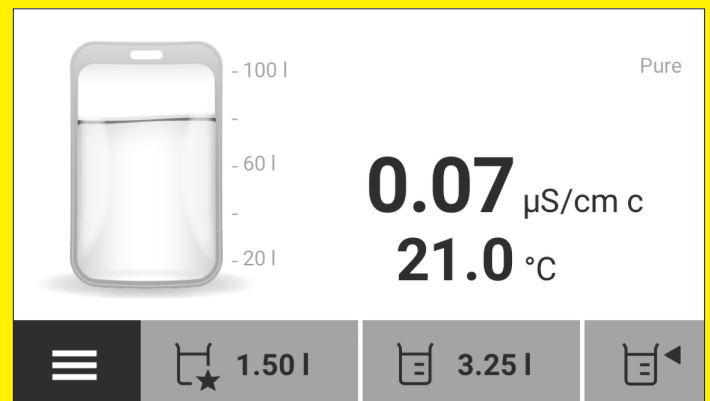
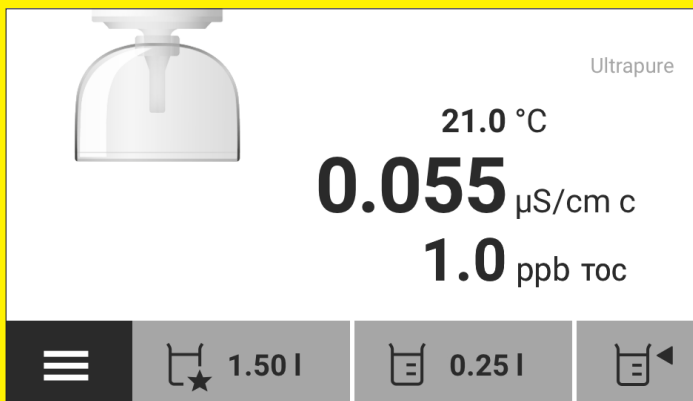
Intelligent by Design

Visual guidance of dispensing

Menu | Home function for settings and cleaning and care programs



* depends on connected system type



Click Access to All Dispense Options
Volume, favorite and manual dispense.

Designed for Pure and Ultrapure Systems
Main Water system status displayed on the screen.

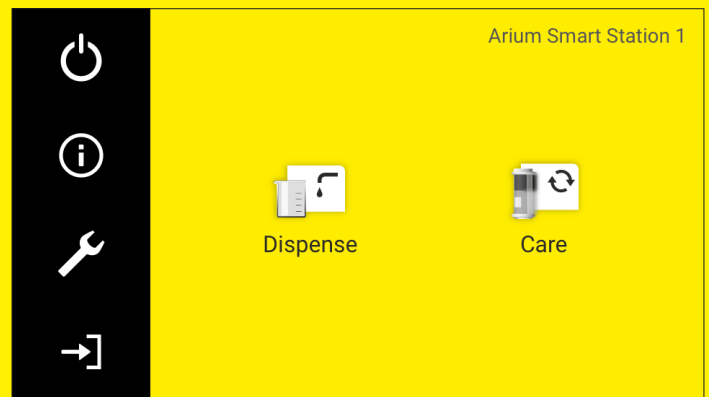


Touch-activated color display with easy-to-clean surface

Conductivity, TOC* and temperature displayed

Volume adjustment

Messages		
Service	01/31/2021	⚠
TOC Calibration	09:00 am	
Care	01/25/2021	⚠
Replace Cartridges	08:30 am	
Care	01/25/2021	⚠
Replace UV lamp	05:00 am	



Safe Operation at All Time

Important information on maintenance and calibration recommendations are clearly highlighted.

Intuitive Menu Navigation supported by Smart Icons

Works even when the user is wearing gloves.

Dispense Wherever You Need

- The Arium® Smart Station can be optimized to your lab space and fits easily into any lab environment
- Combine your Smart Station with the water system type you need to dispense pure or ultrapure water
- The Smart Station ultrapure supports multiple points of use to save costs and valuable lab space.
- Connect up to 3 Smart Stations to your ultrapure water system, the distance between each Smart Stations can be set up to 4 meters.
- Take the advantage to perform volume dispense from the Bagtank from your Smart Station pure.
- Provide water for different applications or just use the additional dispense points where its needed.



Adapted to your needs, the Smart Station has two installation options: wall-mounted or bench top.



The Smart Station can easily be adjusted to left or right handed requirements, without additional accessories.



The small footprint and adjustable arm allows you to position different vessels with varying heights and widths.

Due to its optimal height and intelligent design, the Smart Station can easily fit under a standard laboratory shelf.



SARTORIUS

ARIUM®

Ultrapur

Configure Your Smart Station

Equip your Smart Station with the best filter for your application. Select from two different final dispense filters, depending on whether you need sterile filtration or water for life science and cell culture applications.



Sterile Plus filter
with 0.2 μm pore size
removes bacteria



Cell Plus filter with 5kDa cut off
removes endotoxins,
Rnase/DNase

Choose Your Smart Station

The Smart Station is designed dispense pure and ultrapure water from your system or Bagtank.



Provide ultrapure water from Arium® Pro and
Arium® Comfort



Remote dispense unit provides pure water from
Arium® Bagtank

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 contacts, visit
www.sartorius.com