



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 work space
- 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-handed 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 should not have to adapt your water system to the laboratory, rather the system should adapt to your needs.

In addition to being space-saving 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.

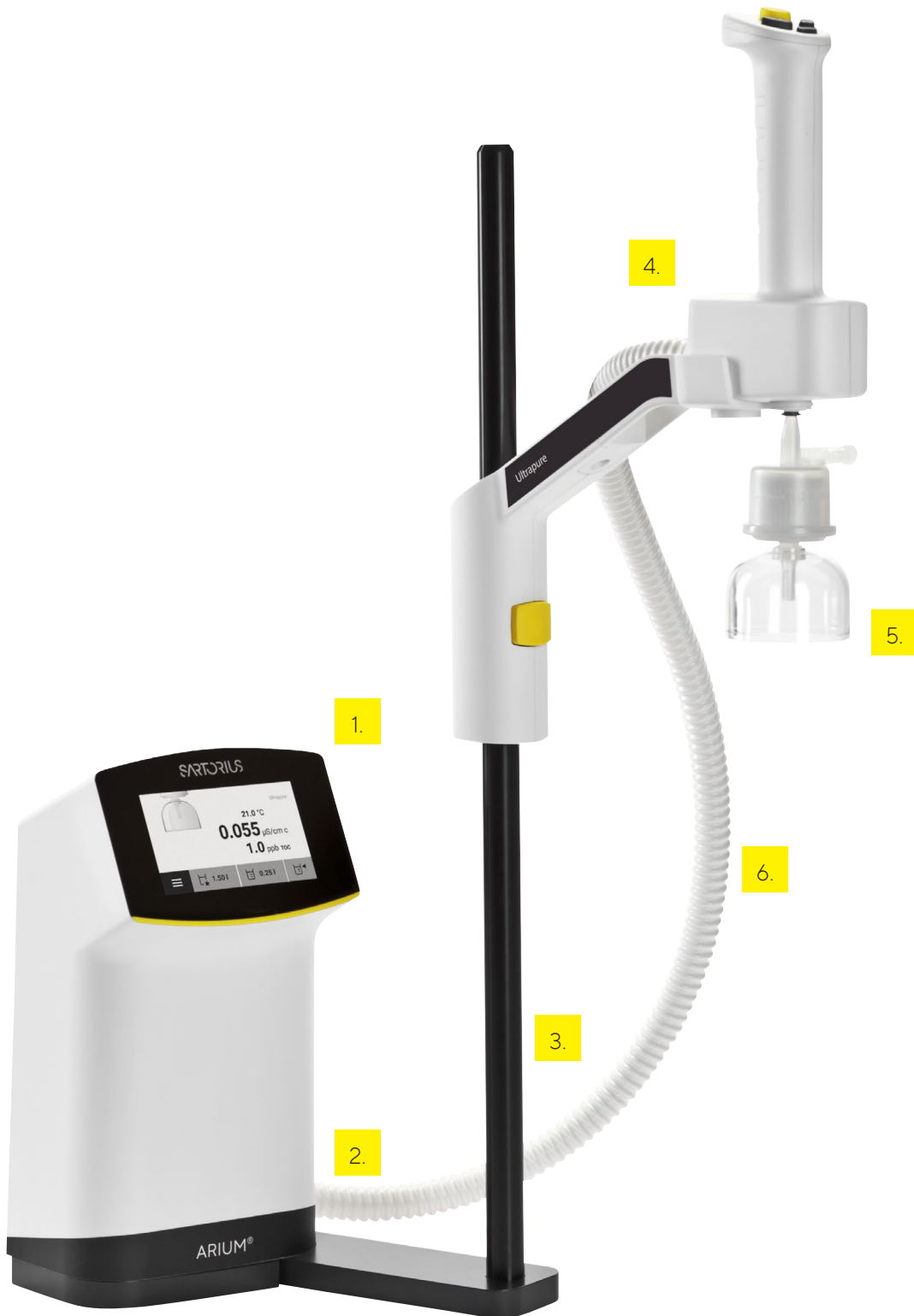
The innovative and flexible Arium® Smart Station from Sartorius addresses all of these requirements. You can dispense water, with the exact quality and quantities required for your experiments, whenever and wherever.

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 on the Arium® Smart Station Ultrapure ensures highest 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 Arium® Smart Station by magnetic connection

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

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

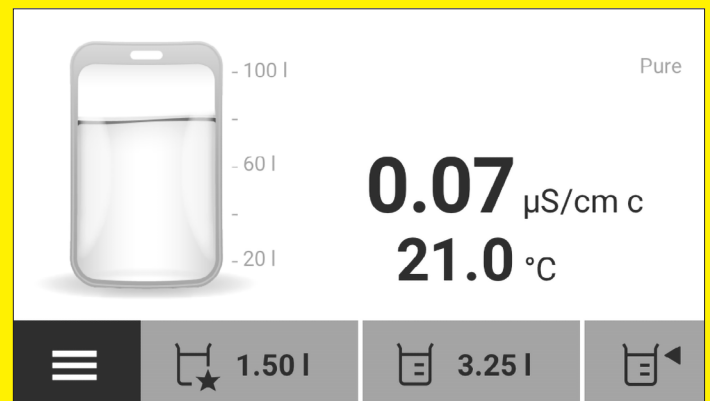
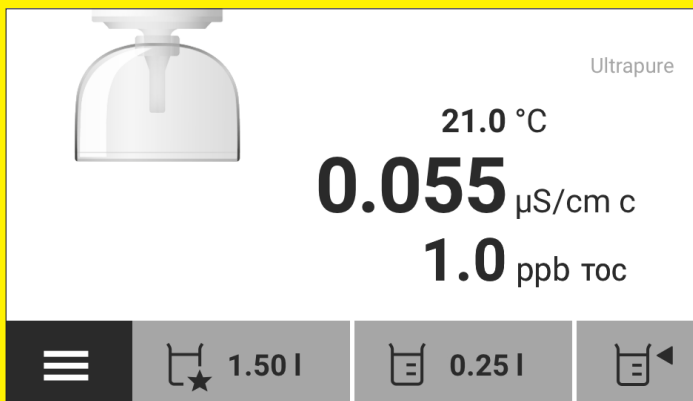
6. Flexibility to Fill Various Vessel Sizes

- Stepless, one-hand adjustment of dispenser arm from 8.5 to 65 cm
- Dispenser arm rotates up to 360°
- Enables volume dispense from 50 mL up to 50 L
- Easy volume dispensing includes soft start | stop function and drop-by-drop

Intelligent by Design

Visual guidance of dispensing

Menu | Home function for settings, cleaning and care programs



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

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



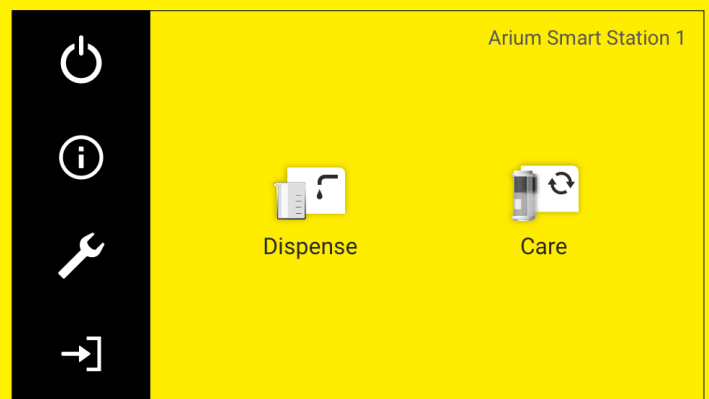
Touch-activated color display with easy-to-clean surface

Conductivity, TOC* and temperature displayed

Volume adjustment

* depends on connected system type

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 Times

Important information on maintenance and calibration recommendations are clearly highlighted.

Intuitive Menu Navigation supported by Smart Icons

Works even when the user is wearing safety gloves.

Dispense Wherever You Need

- The Arium® Smart Station can be optimized to your work space and fits easily into any lab environment
- Combine your Arium® Smart Station with the water system type you need in order to dispense pure or ultrapure water
- The Arium® Smart Station Ultrapure supports multiple points of use to save costs and valuable lab space
- Connect up to 3 Arium® Smart Stations to your ultrapure water system; the distance between each Ultrapure Smart Stations can be set up to 4 meters
- Take advantage of performing volume dispensing from the Bagtank attached to your Arium® Smart Station Pure
- Provides water for different applications and additional dispense points where needed



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



The Arium® 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 Arrium® Smart Station can easily fit under a standard laboratory shelf.



Configure Your Arium® Smart Station

Equip your Arium® Smart Station with the best filter for your application. Select from two different final dispense filters: Sterile Final Filter or CellPlus Filter - depending on your needs.



Sterile Plus filter
with 0.2 μm pore size
removes bacteria



Cell Plus filter with 15 kDa cut off
removes endotoxins,
RNase/DNase

Choose Your Arium® Smart Station

The Arium® Smart Station is designed to 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.



Data Documentation and Real-Time Monitoring

For safe and traceable operation

Data documentation in the laboratory environment are increasingly important. This functionality provides transparent tracing and documentation of product water quality parameters, as well as decentralized control of the system's operating status. With the Arium® Smart Station, you can keep track of your product water quality via one of the following three options below, to best suit your individual laboratory requirements:

Transfer Data to ELN

- Connect your device to an ELN system via Ethernet using a direct connection or an indirect connection via the local network for documentation of various parameters
- Make use of multiple available parameters, such as individual system information, product water quality data, consumable reminders, dispense reports and current messages (alarms | warnings). This enables gapless data documentation for comprehensive traceability
- Protect your data against unauthorized access by using the optionally encrypted transmission by the integrated Sartorius certificate or an individual created certificate

Request Data via Web Service

- Get remote access to your system and monitor system information, product water quality parameters, consumable reminders, alarms and dispense reports
- Connected via your local network, the system can be reached over your regular internet browser, displaying all relevant system parameters at a glance on the Web Service system
- Supported by easy order handling, including direct links to the eShop.
- The integrated print, copy and diagnosis functions offers not only easy documentation but also gives you the opportunity for data transfer to other PC applications

Transfer Data via USB-C


- Connect your Arium® Smart Station to your computer via USB-C for direct recording of the dispensed volumes, including quality data for each water dispense
- The data is automatically passed on to your computer as ASCII text and can be used by various programs for further data processing and evaluation

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