



# 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.



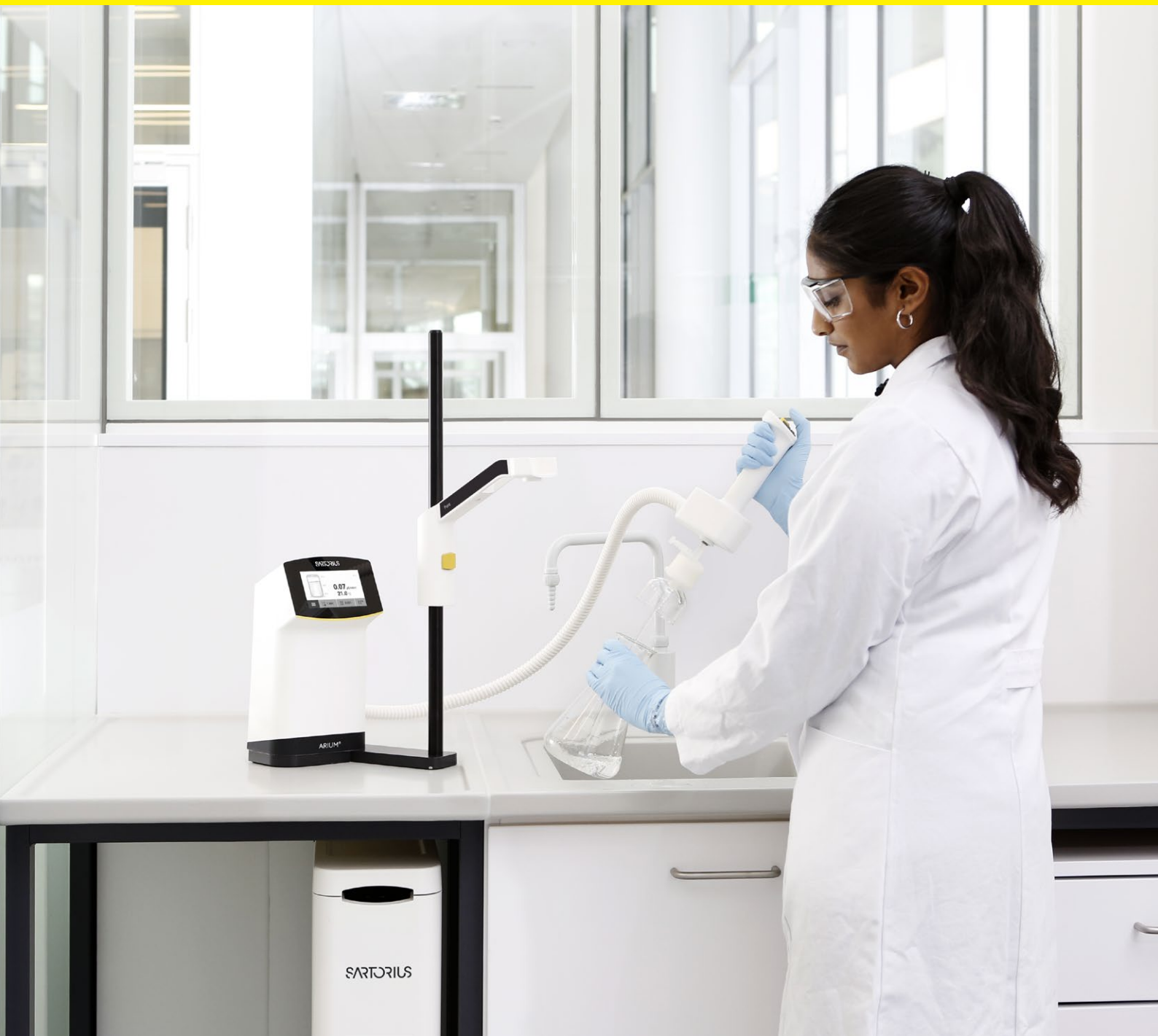
- Compact design for easy and space-saving integration in your lab environment
- Intuitive touch-activated color display with direct access to all important functions
- Flexible and stepless height adjustment to fill different size containers and beakers
- Ergonomic design can be optimized for either 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.

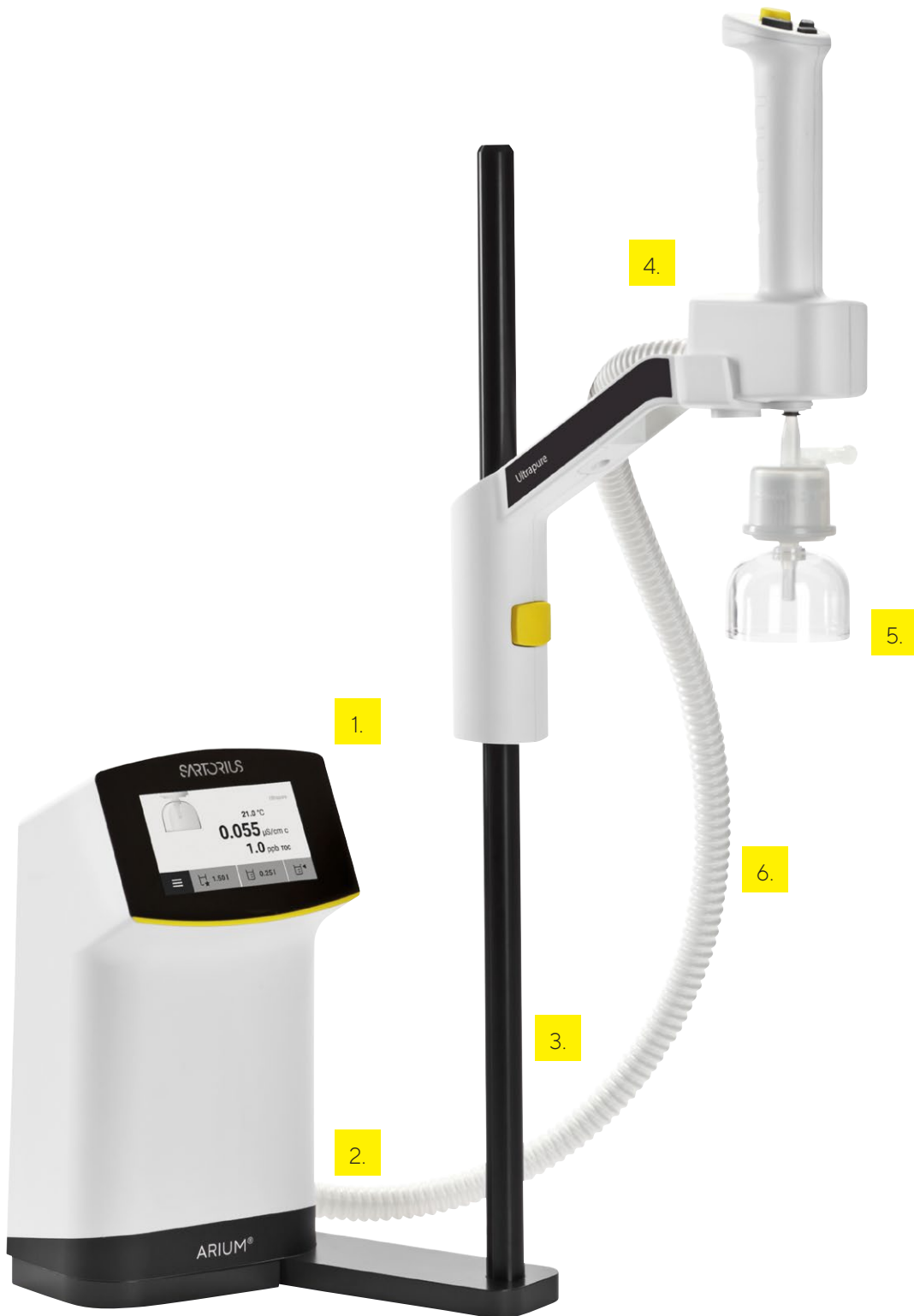
The innovative and flexible Arium® Smart Station from Sartorius addresses all of these requirements. You can dispense water, in the exact quality and quantities required for your experiments, when and where you need it.

**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.

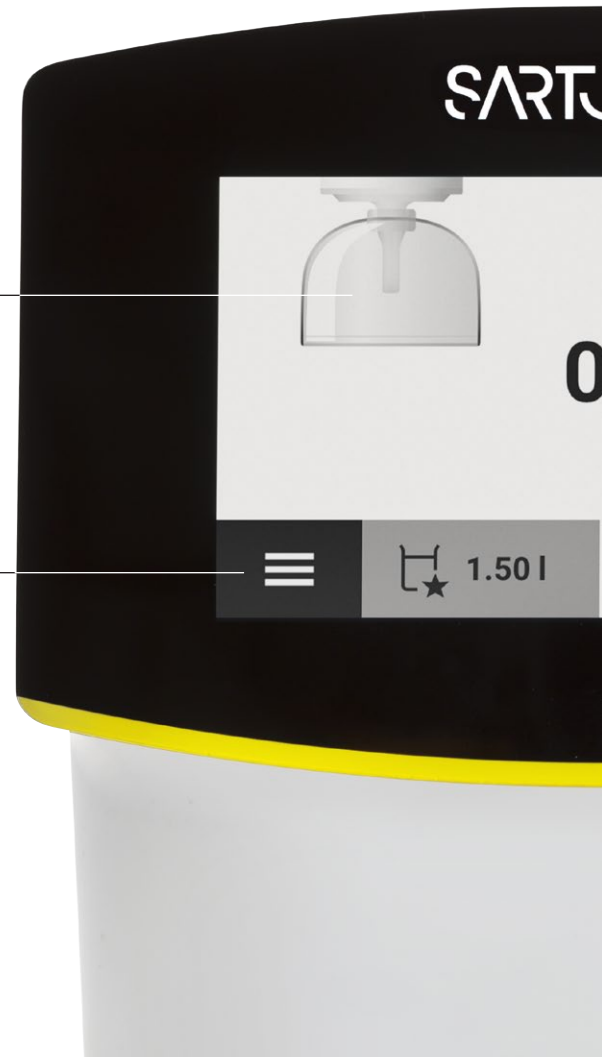
### 6. Flexibility to Fill Various Vessels

- 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

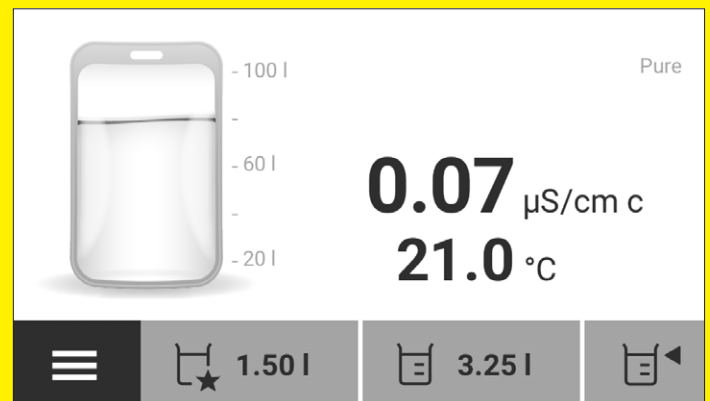
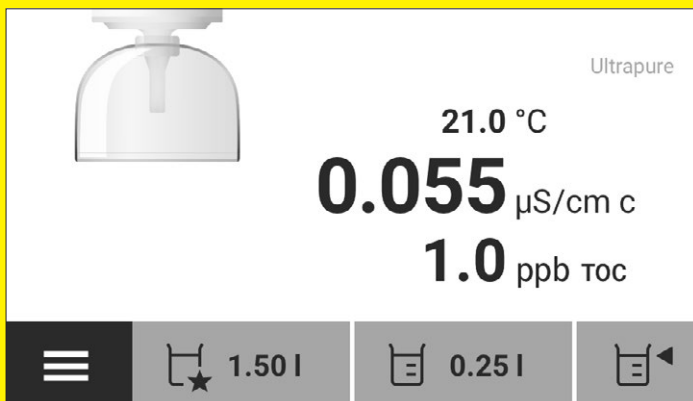
# Intelligent by Design

Visual guidance of dispensing

Menu | Home function for settings and cleaning and care programs



\* depends on connected system type



**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.

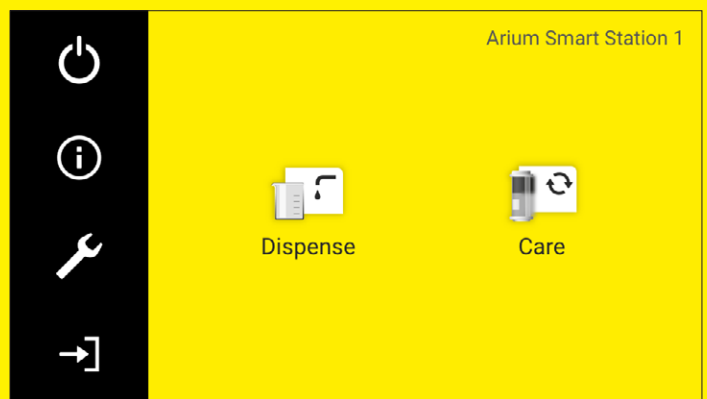


Touch-activated color display with easy-to-clean surface

Conductivity, TOC\* and temperature displayed

Volume adjustment

Messages		
<b>Service</b>	01/31/2021	▲
TOC Calibration	09:00 am	
<b>Care</b>	01/25/2021	▲
Replace Cartridges	08:30 am	
<b>Care</b>	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 it's 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.



# 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  $\mu\text{m}$  pore size  
removes bacteria



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

# Choose Your Smart Station

The 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

### For safe and traceable operation

Data documentation in the laboratory environment are becoming increasingly important. This functionality provides transparent tracing and documentation of product water quality parameters, as well as decentralized control of the systems operating status. With the Arium® Smart Station, you can always keep track of your product water quality via one of the following three options, 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.
- The large number of available parameters, such as individual system information, product water quality data, consumable reminder, dispense reports and current messages (alarms | warnings), enable gapless data documentation for comprehensive traceability.
- For secure transfer, all data can be protected against unauthorized access by encryption using a Sartorius or an individual 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 system Web Service.
- Supported by easy order handling, including direct links to the eShop. With print, copy and diagnosis functions this 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, which 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](http://www.sartorius.com)