

SARTORIUS

Arium® Smart Station

Simplifying Water Dispensing at an Unbeatable Price

Special Launch Offer
Limited Quantity -
100 Bundles Available

**Enabling highest flexibility
for your daily application**

The Arium® Smart Station is designed for flexible remote dispensing of ultrapure water directly at the point of use. Due to its intelligent and compact design, it can be easily integrated into any lab environment.

**Don't Miss This
Special Offer!**

Get your Arium® Smart Station with an Arium® Pro UV ultrapure water system and consumables.



For just £4,685

Limited Quantity Offer

Get a fully functional Arium® Pro UV System with the new Arium® Smart Station for £4,685*

This package includes:

- Arium® Pro UV System
- Arium® Smart Station (Ultrapure)
- Arium® Pro UV Cartridge Kit
- Arium® Pro UV Sterile Filter
- Arium® Smart Station Sterile Filter

Order Information:

H2O-Action-ST-UP

* Limited quantity offer

Arium® Smart Station

Flexible remote dispenser

- Compact: Save space integrating into 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



Arium® Pro UV

Ultrapure water system

- Consistent quality: Meet or exceed ASTM Type 1 water quality to increase sensitivity of analytical results
- Optimized TOC reduction: Reduce TOC levels to a minimum and avoid background noises in chromatograms
- Reliable process stability: Variety of services (e.g. IQ/OQ), ensuring consistently reliable water quality through the entire life cycle of the system.
- Cost efficiency: High throughput cartridge for best return of investment



Please contact your Sales Representative or any of our authorized Partners for more information.

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