

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 pure water directly at the point. Due to its intelligent design, it can easily integrate into your lab space.

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

Get your Arium® Smart Station with an Arium® Bagtank 50 and consumables.



For just 1.895€

SARTORIUS

Limited Quantity Offer

Get an Arium® Bagtank 50 with the new Arium® Smart Station for 1.895€*

This package includes:

- Arium® Bagtank 50 Liter
- Arium® Smart Station (Pure)
- Arium® Bag 50
- Arium® Smart Station Sterile Filter

Order Information:

H2O-Action-ST-P

* 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® Bagtank 50 Liter

Unique bagtank technology

- Safe: Closed system guarantees stable quality thanks to including venting filter with integrated check valve
- Time efficient: Easy bag exchange in less than 5 minutes
- Maintenance free: No need for chemical cleaning procedure
- Stable quality: Optimum water quality by prevention of a permanent biofilm thanks to special multilayer S71 bag material



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