

Arium® Water Purification Systems

Arium® Laboratory-Grade Water Purification Systems With Added Value

Simplifying Progress



Arium® Pro

Application-orientated and flexible to meet the highest demands

- Modular System selection specifically for your application
- Flexible Perfect integration into any laboratory
- Easy to use Display with touch function and intuitive menu
- Fast Favorites function with direct access for recurring volumes
- Compatible with Arium® Smart Station



Arium® Mini

Compact Ultrapure Water Systems for 10L per Day

- Constant low conductivity for reliable results
- Integrated Bagtank technology for best economic water storage
- Easy-to-use touch display for intuitively menu navigation
- Space-saving unit with a width of only 28 cm (11 inches)
- Highest quality standards made in Germany



Arium® Smart Station

The Most Flexible Solution for Dispensing Pure and Ultrapure Water

- Compact: Save space integrating in 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
- Compatible with Arium® Pro