## S/RICH SILS

## 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 and Arium® Comfort Systems

## Simplifying Progress

