

Arium® Smart Station

Water. Exactly where you need it.

The Arium® Smart Station is designed for flexible remote dispensing of pure and ultrapure water directly at the point of use.

While dispensing water into a broad range of different sized containers, the Smart Station offers constant control of all important quality parameters, at all times. The ergonomic design supports left and right hand operation and can be easily adjusted to your need.

Features and Benefits

- 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
- Adjustable, splash-free flow rate from drop-by-drop to 2 liters per minute



Discover more about
Arium® Smart Station



Please scan!