

Arium® Pro

Tailored to Your Specific Requirements

As a reliable source of ultrapure water, the Arium® Pro series offers a flexible and modular system which, compared to conventional devices, demonstrates excellent added value. All systems meet and exceed the ASTM Type 1 water quality standards and ensure the best reproducible results in their class. Up to 2 liters of consistently high-quality ultrapure water with a conductivity of $0.055 \mu\text{S}/\text{cm}$ can be dispensed per minute. The ultrapure water is virtually microorganism-free when an Arium® SterilePlus final filter is used. The patented Sartorius technology, the SD card slot, the long service life and low maintenance requirement distinguish the Arium® Pro systems as an easy-to-use, economical and reliable Type 1 ultrapure water system.

Features

- 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



Discover more about
Arium® Pro Lab Water System



Please scan!