

Arium® Mini

Designed to deliver pure performance

Compact Arium® Mini laboratory water systems have been designed for Type 1 ultrapure water requirements of 10 liters per day and are ideal for use in preparation of buffers, media and samples, both in life sciences applications and in analytical laboratory procedures. A large, touch-activated screen and intuitive menu navigation ensure exceptionally easy operation. Regardless of your type of feed water available, use Arium® Mini or Arium® Mini plus featuring our unique bagtank technology or connect Arium® Mini essential directly to your deionized water supply line in the lab.

Features

- Reliable: Delivers consistently high water quality for reliable and reproducible results
- Intuitive: Touch-activated color display with direct access to all important dispensing options
- Innovative: With unique bagtank technology, depending on the type of system; saves time-intensive tank cleaning
- Compact: Space-saving, with a width of only 28 cm



Discover more about
Arium® Mini Lab Water System



Please scan!