

Installation

Arium® Water Purification Systems



Benefits

- Perfectly functioning devices
- Easy and rapid deployment
- Long-term highest water quality
- Professionally trained operating personnel
- Extended device life

Product Information

Professional installation of your Arium® system by our Sartorius Service is the easiest and safest way to ensure flawless startup of your new water purification system. We set up your system for you and guarantee its correct operation. So you can rely on the highest water quality and optimal results from your Arium® system right from the start.

Specification

- Checking the equipment delivery for completeness and any transport damage
- Installation of the device at the place of use (no wall mounting)
- Connection to the supply line
- Initial startup
- Leakage check
- Inspection of flow rate capacity
- Inspection of conductivity values
- Issuing an installation report
- Instruction of operating personnel (no application training)
- Issuing an installation report
- Consultation on device maintenance and maintenance cycles

Customer Prerequisites

- Device is at the installation site and has been sufficiently acclimatized
- The installation place is freely accessible
- To wall-mount the device, the included wall bracket must be pre-installed by the customer
- Supply lines available at the installation site, max. 1 m away from the location of the system
- Feed water pressure according to device specifications
- Atmospherically vented drain with odor trap (max. 1 m away from the location of the system)
- Splash-proof wall outlet (240 V, 50 Hz), protected with 16 A fuse, max. 1 m away from the location of the system

Optional Services

- GLP | GMP compliant equipment qualification

You are interested?


Get in touch with us right away. We will be happy to provide you with a non-binding service offer. You can also find further information on our website: www.sartorius.com/en/services

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further contacts, visit
www.sartorius.com