

Preventive Maintenance and Service Contracts

Arium® Water Purification Systems



Benefits

- Consistently highest water quality
- Optimized device performance and process safety
- Trouble-free operation, reduction of downtime
- Extended device life
- Individual consumables options

Product Information

We offer regular maintenance service options to prolong the life of the Arium® water purification system and to optimize its reliability. In the long run, you are protecting your investment and guaranteeing accurate results. At the same time, you are maximizing the performance of your system as well as the efficient use of the purified water—with consistently high quality. We recommend taking out a maintenance contract tailored to your needs so you can continuously rely on your water purification system and ensure maximum productivity.

Specification

- Visual inspection
- System cleaning (for 613L and Advance L, no cleaning of the tank)
- Replacement of consumables
- Replacement of TOC module (if applicable)
- Check of feed water quality
- Calibrating the sensors
- Inspection of flow rate capacity
- Inspection of the internal system pressure
- Leakage check
- Startup test
- Issuing a maintenance report

Repairs, original spare parts, travel costs, and other expenses are not included in the scope of services and will be charged separately.

Customer Prerequisites

- Device is at the maintenance site

Optional Services

- TOC-measurement of the product water
- Consumables packages and subscription
- Annual care-free package

Remarks

Device cannot be used by staff during maintenance.

Are You 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 Goettingen
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 additional information,
visit www.sartorius.com