

# Preventive Maintenance

## Arium® Water Purification Systems

### Consumables Included



### Benefits

- Consistently highest water quality
- Optimized device performance
- Trouble-free operation, reduction of downtime
- Extended device life
- Consumables included

### Product Information

We offer a regular maintenance service 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 that you take out a maintenance contract so that you can fully 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)
- Consumables and their replacement
- 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

## Remarks

Device cannot be used by staff during maintenance.

## 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](http://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](http://www.sartorius.com)