



# Arium® Lab Water Systems

Ensuring Reliability Through Service Contracts and Scheduled Replacement of Consumables

Simplifying Progress

**SARTORIUS**

# When It Comes to Water Quality, Play It Safe

In many laboratory processes and applications involving low detection limits, it is decisively important to consistently maintain the highest-possible water quality, as such standards are vital to achieving reliable and reproducible analysis results and high process reliability.

Every detail counts. Inadequate replacement of consumables can cause highly sensitive analyses such as high-performance liquid chromatography (HPLC) to yield non-reproducible and non-specific results.

## Arium® Lab Water Systems

Arium® Mini



Arium® Comfort



# Service Contracts and Regular Scheduled Replacement of Consumables Ensure Reliability

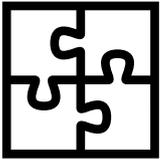
Regular replacement of system consumables and professional maintenance of your laboratory equipment by trained service specialists are essential to consistently ensuring the highest-possible water quality and preventing any risk of your Arium® lab water systems incurring unscheduled outages.

That's how you prevent any microbiological contamination of equipment and erroneous analysis results, and establish the necessary process reliability in your laboratory.



# The Right Maintenance Package for You

The choice of the optimal maintenance package depends on your individual requirements and needs in the lab. We recommend regular maintenance of your lab equipment at least once annually by specially trained service engineers so that you can always rely on your Arium® lab water systems throughout their entire service lifetime. You decide for yourself the extent to which you would like us to support you.



Several options to choose from



Maintenance tailored to your needs



Peace of mind thanks to carefree packages



Entire life cycle support

Multiple options support you in selecting the optimal maintenance package meeting your and your system's needs.

Our modular approach helps you to configure the right package. From single maintenance visits up to annual carefree package agreements, we aim to simplify your daily challenges and provide peace of mind.

The professional maintenance of your lab water system is basis and you can individually add the following:

- Consumables packages
- Consumables replacement visits by our service specialist in between the maintenance visits
- Full-service, covering breakdowns throughout the entire lifetime of your system

# Our Consumables Packages

It is essential to regularly replace the consumables of your Arium® system in order to consistently ensure untainted water quality over long-term operation.

With our consumables packages and a subscription for automatic delivery, you ensure that consumables replacement and timely reordering of supplies are not overlooked in the stressful and demanding setting of daily laboratory routines.

Besides maintenance servicing, consumables play a crucial role and are therefore part of our comprehensive package offering.

When configuring the optimal maintenance package, you decide which consumables packages you need at which point in time according to your individual requirements and needs.

You can either opt for a subscription including automatic delivery or order the consumables packages by yourself in our e-shop, in case you prefer to organize and replace the consumables by yourself.

Save precious time and money with consumables packages and automated delivery. This ensures that the right consumables are always available when you need them. In addition, you have more free storage capacity.

Easy ordering in our e-shop



Cost savings



Delivery at the right point in time



Free storage space



# Process Reliability and Consistent Water Quality



Regular  
replacement of consumables



Customized  
service contracts



Consumables packages  
and consumables subscription

## Germany

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

 For further information, visit  
[www.sartorius.com](http://www.sartorius.com)

## USA

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