



Arium® Lab Water Systems

Ensuring Reliability through Service Contracts and Scheduled Replacement of Consumables

Simplifying Progress

SARTORIUS

When it comes to water quality, never make any compromises

In many laboratory processes and applications involving low detection limits, it's 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 service contract for you

Besides the necessary replacement of system consumables, 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 yourself the extent to which you would like us to support you.

Our professional maintenance encompasses the following services

- System cleaning
- Replacement of total organic carbon (TOC) analyzer module (if installed)
- Comprehensive testing of feedwater quality and product water quality
- Calibration of all sensors
- Complete check of all functions
- Leaktightness testing
- Documentation of all maintenance activities and measurement results in a test report

Decide yourself on an individual basis whether you want consumables to be replaced within the scope of our maintenance work, or whether you wish to replace them yourself independently of maintenance activities.

If you don't wish to make any compromises when it comes to regularly replacing your lab consumables, you can also have them replaced as part of a comprehensive care-free package. If you opt for such a package, our Service Engineer will visit you once each quarter and replace all consumables scheduled for replacement at that time. In such cases, all that remains for you to do yourself is replace the sterile final filter every month.

Note: In devices fitted with an integrated TOC analyzer module, this module must always be replaced with a newly calibrated module by a Service Engineer. This replacement is performed as part of maintenance work.



Replacement of consumables

It is essential to regularly replace the consumables of your Arium® system in order to consistently ensure untainted water quality over long-term operation. Our consumables packages and a consumables subscription ensure that consumables replacement and timely reordering of supplies don't get overlooked in the stressful and demanding setting of daily laboratory routines.

Consumables packages and consumables subscription

You can order all the consumables you need for your Arium® system in practical consumables packages containing all consumables that need to be replaced in the coming quarter, custom-picked and prepared. Consumables packages save you money compared to ordering items individually.

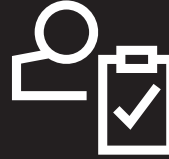
We recommend automatic delivery of your consumables packages by way of a practical subscription so that you can focus on your work and continuously rely with confidence on your Arium® system. You then automatically receive the correct package of consumables needed in the coming quarter. Our consumables subscription thereby ensures that you always have the right consumables in stock and minimum shelf-life requirements are met, while no storage space is taken up unnecessarily by overstocking inventory.



High process reliability and consistent water quality



Regular
replacement of consumables



Customized
service contracts



Consumables packages
and consumables subscription

Replacement cycles of consumables

The replacement intervals recommended below for the individual consumables are indicated in months. These intervals are based on the feedwater qualities and sampling quantities stipulated in the respective equipment specifications.

It may become necessary to replace consumables at shorter intervals, depending on the given water quality and sampling quantity.

Order numbers of consumables packages

Differing material packages result for the respective periods of use in line with the replacement cycles of the individual consumables, starting from the point in time of equipment commissioning.

Each three-month package is needed at the start of the respective quarter, requiring a corresponding order in advance. If you opt for a subscription, however, you don't have to keep reminding yourself about ordering, as we deliver all the right packages to you right on schedule.

For many equipment models, the package needed in the fourth quarter of every operating year differs to those for the year's previous quarters (see Arium® Comfort and Arium® Advance).

In every odd-numbered operating year, the "Package for 4th quarter | 1st year" is needed, while in every even-numbered operating year it's the "Package for 4th quarter | 2nd year".

Arium® Mini

Replacement cycles of consumables (in months)

Model series	Equipment model	Pretreatment cartridge	Cartridge set	Final filter	Bag	UV lamp
Mini	H2O-MM-T		6	1	6	
	H2O-MM-UV-T		6	1	6	12
Mini Plus	H2O-MA-T	6	6	1	6	
	H2O-MA-UV-T	6	6	1	6	12
Mini Essential	H2O-MU-T		6	1	6	
	H2O-MU-UV-T		6	1	6	12

Order numbers of consumables packages

Model series	Equipment model	Package for 1st quarter	Package for 2nd quarter	Package for 3rd quarter	Package for 4th quarter
Mini	H2O-MM-T	SB-12-03-0521	SB-12-03-0523	SB-12-03-0521	SB-12-03-0523
	H2O-MM-UV-T	SB-12-03-0521	SB-12-03-0523	SB-12-03-0521	SB-12-03-0526
Mini Plus	H2O-MA-T	SB-12-03-0521	SB-12-03-0524	SB-12-03-0521	SB-12-03-0524
	H2O-MA-UV-T	SB-12-03-0521	SB-12-03-0524	SB-12-03-0521	SB-12-03-0527
Mini Essential	H2O-MU-T	SB-12-03-0521	SB-12-03-0525	SB-12-03-0521	SB-12-03-0525
	H2O-MU-UV-T	SB-12-03-0521	SB-12-03-0525	SB-12-03-0521	SB-12-03-0528



Arium® Mini

Arium® Comfort

Replacement cycles of consumables (in months)

Model series	Equipment model	Cleaning set	Pretreatment cartridge	Softener cartridge	Reverse Osmosis (RO) module(s)	Cartridge set	Final filter	Bag	UV lamp
Comfort I	H2O-I-1-T/B	3	3		24	12	1		
	H2O-I-2-T/B	3	3		24	12	1		
	H2O-I-1-UV-T/B	3	3		24	12	1		12
	H2O-I-1-TOC-T/B	3	3		24	12	1		12
	H2O-I-2-UV-T/B	3	3		24	12	1		12
	H2O-I-2-TOC-T/B	3	3		24	12	1		12
Comfort II	H2O-II-1-T/B	3	3	6	24	12	1		
	H2O-II-2-T/B	3	3	6	24	12	1		
	H2O-II-1-UV-T/B	3	3	6	24	12	1		12
	H2O-II-1-TOC-T/B	3	3	6	24	12	1		12
	H2O-II-2-UV-T/B	3	3	6	24	12	1		12
	H2O-II-2-TOC-T/B	3	3	6	24	12	1		12
Bagtank	H2O-AOV-20							6	
	H2O-AOV-50							6	
	H2O-AOV-100							6	

Order numbers of consumables packages

Model series	Equipment model	Package for 1st quarter	Package for 2nd quarter	Package for 3rd quarter	Package for 4th quarter 1st year	Package for 4th quarter 2nd year
Comfort I	H2O-I-1-T/B	SB-12-03-0500	SB-12-03-0500	SB-12-03-0500	SB-12-03-0502	SB-12-03-0506
	H2O-I-2-T/B	SB-12-03-0500	SB-12-03-0500	SB-12-03-0500	SB-12-03-0502	SB-12-03-0507
	H2O-I-1-UV-T/B	SB-12-03-0500	SB-12-03-0500	SB-12-03-0500	SB-12-03-0503	SB-12-03-0508
	H2O-I-1-TOC-T/B	SB-12-03-0500	SB-12-03-0500	SB-12-03-0500	SB-12-03-0503	SB-12-03-0508
	H2O-I-2-UV-T/B	SB-12-03-0500	SB-12-03-0500	SB-12-03-0500	SB-12-03-0503	SB-12-03-0509
	H2O-I-2-TOC-T/B	SB-12-03-0500	SB-12-03-0500	SB-12-03-0500	SB-12-03-0503	SB-12-03-0509
Comfort II	H2O-II-1-T/B	SB-12-03-0500	SB-12-03-0501	SB-12-03-0500	SB-12-03-0504	SB-12-03-0510
	H2O-II-2-T/B	SB-12-03-0500	SB-12-03-0501	SB-12-03-0500	SB-12-03-0504	SB-12-03-0511
	H2O-II-1-UV-T/B	SB-12-03-0500	SB-12-03-0501	SB-12-03-0500	SB-12-03-0505	SB-12-03-0512
	H2O-II-1-TOC-T/B	SB-12-03-0500	SB-12-03-0501	SB-12-03-0500	SB-12-03-0505	SB-12-03-0512
	H2O-II-2-UV-T/B	SB-12-03-0500	SB-12-03-0501	SB-12-03-0500	SB-12-03-0505	SB-12-03-0513
	H2O-II-2-TOC-T/B	SB-12-03-0500	SB-12-03-0501	SB-12-03-0500	SB-12-03-0505	SB-12-03-0513
Bagtank	H2O-AOV-20		H2O-CBS-20		*	
	H2O-AOV-50		H2O-CBS-50		*	
	H2O-AOV-100		H2O-CBS-50		H2O-CBS-50	

* Use of 2nd bag from package for 2nd quarter



Arium® Advance



Replacement cycles of consumables (in months)

Model series	Equipment model	Cleaning set	Pretreatment cartridge	Softener cartridge	Reverse Osmosis (RO) module(s)	Bag
Advance RO	H2O-RO-1-T/B	3	3		24	
	H2O-RO-2-T/B	3	3		24	
	H2O-RO-3-T/B	3	3		24	
Advance EDI	H2O-EDI-1-T/B	3	3	6	24	
	H2O-EDI-2-T/B	3	3	6	24	
Bagtank	H2O-AOV-20					6
	H2O-AOV-50					6
	H2O-AOV-100					6

Order numbers of consumables packages

Model series	Equipment model	Package for 1st quarter	Package for 2nd quarter	Package for 3rd quarter	Package for 4th quarter 1st year	Package for 4th quarter 2nd year
Advance RO	H2O-RO-1-T/B	SB-12-03-0514	SB-12-03-0514	SB-12-03-0514	SB-12-03-0514	SB-12-03-0516
	H2O-RO-2-T/B	SB-12-03-0514	SB-12-03-0514	SB-12-03-0514	SB-12-03-0514	SB-12-03-0517
	H2O-RO-3-T/B	SB-12-03-0514	SB-12-03-0514	SB-12-03-0514	SB-12-03-0514	SB-12-03-0518
Advance EDI	H2O-EDI-1-T/B	SB-12-03-0514	SB-12-03-0515	SB-12-03-0514	SB-12-03-0515	SB-12-03-0519
	H2O-EDI-2-T/B	SB-12-03-0514	SB-12-03-0515	SB-12-03-0514	SB-12-03-0515	SB-12-03-0520
Bagtank	H2O-AOV-20		H2O-CBS-20		*	
	H2O-AOV-50		H2O-CBS-50		*	
	H2O-AOV-100		H2O-CBS-50		H2O-CBS-50	

* Use of 2nd bag from package for 2nd quarter

Arium® Pro



Replacement cycles of consumables (in months)

Model series	Equipment model	Cleaning syringe	Ultrafilter	Cartridge set	Final filter	UV lamp
Pro	H2OBASIC-T/B			12	1	
Pro DI	H2OPRO-DI-T/B	3		12	1	
Pro UF	H2OPRO-UF-T/B	3	12	12	1	
Pro UV	H2OPRO-UV-T/B	3		12	1	12
Pro UV TOC	H2OPRO-UV-T/B-TOC	3		12	1	12
Pro VF	H2OPRO-VF-T/B	3	12	12	1	12
Pro VF TOC	H2OPRO-VF-T/B-TOC	3	12	12	1	12

Order numbers of consumables packages

Model series	Equipment model	Package for 1st quarter	Package for 2nd quarter	Package for 3rd quarter	Package for 4th quarter
Pro	H2OBASIC-T/B	SB-12-03-0521	SB-12-03-0521	SB-12-03-0521	SB-12-03-0529
	If U-Pack is used	SB-12-03-0534	SB-12-03-0534	SB-12-03-0534	SB-12-03-0534
Pro DI	H2OPRO-DI-T/B	SB-12-03-0522	SB-12-03-0522	SB-12-03-0522	SB-12-03-0530
	If U-Pack is used	SB-12-03-0535	SB-12-03-0535	SB-12-03-0535	SB-12-03-0535
Pro UF	H2OPRO-UF-T/B	SB-12-03-0522	SB-12-03-0522	SB-12-03-0522	SB-12-03-0531
	If U-Pack is used	SB-12-03-0535	SB-12-03-0535	SB-12-03-0535	SB-12-03-0536
Pro UV	H2OPRO-UV-T/B	SB-12-03-0522	SB-12-03-0522	SB-12-03-0522	SB-12-03-0532
	If U-Pack is used	SB-12-03-0535	SB-12-03-0535	SB-12-03-0535	SB-12-03-0537
Pro UV TOC	H2OPRO-UV-T/B-TOC	SB-12-03-0522	SB-12-03-0522	SB-12-03-0522	SB-12-03-0532
	If U-Pack is used	SB-12-03-0535	SB-12-03-0535	SB-12-03-0535	SB-12-03-0537
Pro VF	H2OPRO-VF-T/B	SB-12-03-0522	SB-12-03-0522	SB-12-03-0522	SB-12-03-0533
	If U-Pack is used	SB-12-03-0535	SB-12-03-0535	SB-12-03-0535	SB-12-03-0538
Pro VF TOC	H2OPRO-VF-T/B-TOC	SB-12-03-0522	SB-12-03-0522	SB-12-03-0522	SB-12-03-0533
	If U-Pack is used	SB-12-03-0535	SB-12-03-0535	SB-12-03-0535	SB-12-03-0538

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Göttingen, Germany
Phone +49 (0)551 308-0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Phone (toll-free) +1 (0)800 635 2906

 Find out more at
www.sartorius.com