

arium[®] mini – Compact Ultrapure Water Systems for 10 Liters per Day



arium® mini – Innovation for Efficient Research

Ultrapure water is the basis for successful analyses. Yet analytical methods are becoming more and more sensitive and thus more susceptible to interference. Consistently high water quality is decisive in ensuring the reproducibility of your results and in preventing time-consuming repeat analyses.

You can always rely on the arium® mini series of laboratory water systems to consistently deliver the highest water quality.

Regardless of the type of feed water you have, use arium® mini or arium® mini plus with our unique bagtank technology or connect arium® mini essential directly to your deionized water supply line as needed.

Type of System Feed Water*

arium® mini	Pretreated water from a supply container
arium® mini essential	Directly connects to a pretreated water supply line (RO DI EDI)
arium® mini plus	Directly connects to tap water

* For details, see the inlet water specifications.



Your arium® mini Benefits

■ Reliable

Delivers consistently high water quality for reliable and reproducible results; optionally available with UV lamp

■ Intuitive

Easy to operate using the touch-activated color display; with direct access to all important dispensing options

■ Innovative

Depending on the model type, with unique bagtank technology to save you from the hassle of time-consuming tank cleaning

■ Compact

Space-saving footprint with a width of only 28 cm to easily fit in any lab environment



Focused on Your Requirements – Convenient Operation and the Highest Quality Standards

Conveniently operate arium® mini using the touch-activated color display with self-explanatory icons – even while wearing laboratory gloves. This will not only speed up your workflows, but will also make them more accurate and reliable as a result of active error prevention – your daily lab work will become easier and more efficient.

To ensure the best ultrapure water quality and the reliability of your process, we have also implemented the highest quality standards into all functional areas. For example, you can set a limit for the purified water quality. If this limit is exceeded, further dispensing of water will be prevented or accompanied by a warning, depending on your menu configuration settings selected.

Integrated Illumination for Easy Visual Support

- Flashing light during water dispensing
 - Pulse signal emitted to accompany error messages
 - Dimmed light in standby mode
-

arium® Scientific Pack for the Highest Ultrapure Water Quality

- The highest cartridge capacity thanks to efficient purification materials
 - Optimal flow over the semiconductor-grade mixed-bed resin
 - Designed for quality standard ASTM Type 1
-

arium® SterilePlus for Effective Microbe Removal

- Validated according to HIMA and ASTM F-838-05
 - Meets quality standards in accordance with the USP
 - Reliably removes microorganisms using a 0.2 µm membrane
-

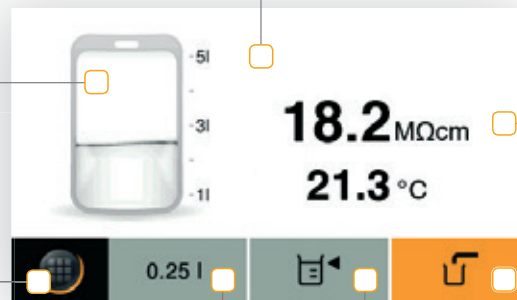


Current filling volume displayed for the bag (arium® mini & mini plus)

Touch-activated color display with easy-to-clean surface

Purified water quality and temperature displayed

Menu | Home function for settings and cleaning and care programs



Manual dispensing

Volume-controlled dispensing in 50 mL increments

Last volume dispensed is saved by the Favorites function



Intuitive Menu Navigation for Total Ease of Operation

- Icon-guided interface with touch-activated display
- Displays current measured values and messages
- Favorites function for repeated dispensing of identical volumes

arium® UV Lamp for Critical Analyses (185 | 254 nm)

- Optimal temperature gradient due to horizontal installation
- Degradation of total organic carbons (TOCs) by irradiation (≤ 5 ppb)
- Inactivates microorganisms by damaging their DNA

arium® Pretreatment Cartridge for Effective Purification

- Effective adsorption using high-quality activated carbon
- Special catalyst removes oxidants
- Impurities retained by reverse osmosis

Integrated Bagtank Technology in arium® mini and mini plus – A Revolution in Laboratory Water Purification

As a pioneer and market leader in the manufacture of single-use products for biopharmaceutical production, we have written the next chapter in our success story by implementing our innovative bagtank technology in laboratory water systems.

Originally developed for the pharmaceutical industry, our single-use bag integrated as a tank system on the side of arium® mini protects intermediately stored pure water reliably from secondary contamination.

The 5-liter bag, consisting of our specially manufactured proprietary S71 film material, is multilayer and compliant with ISO and USP standards for biocompatibility. With the quality and composition of the arium® Bag, you can optimally store your pretreated pure water for later production of Type 1 ultrapure water.

Moreover, you will enjoy the benefits of active protection against leakage provided by the integrated Aqua Stop system.

Intelligent Technology – Systematic Purification to Obtain Ultrapure Water

With intelligent technology, the arium® mini ensures continuous protection against secondary contamination and therefore the highest water quality at all times.

The closed bagtank system automatically regulates pressure equalization so that no CO₂ can penetrate. This additionally prevents contamination by ionic impurities and maintains consistently low conductivity. As an added benefit, you will profit from the especially high cartridge capacity because there is less demand placed on the ion exchangers.

Moreover, you can prevent the formation of a permanent biofilm, and therefore eliminate additional TOC loading. Depending on your requirements, you will just need to replace the bag every six months to start off fresh.



5-Liter Bag for Minimum Maintenance

- Easy bag exchange in less than 5 minutes
- No need to use chemicals for cleaning
- Practical and leak-tight connection using a quick-connect coupling

Innovative Technology to Prevent Secondary Contamination

- Optimum water quality by prevention of a permanent biofilm
- Closed system guarantees purity
- Stable quality thanks to special multilayer S71 bag material

Well-Thought-Out Functions for Reliable Performance

- Integrated pump to increase the water pressure
- Filling arium® mini with pretreated water: 2 minutes
- Filling arium® mini plus with RO permeate: 8 liters per hour

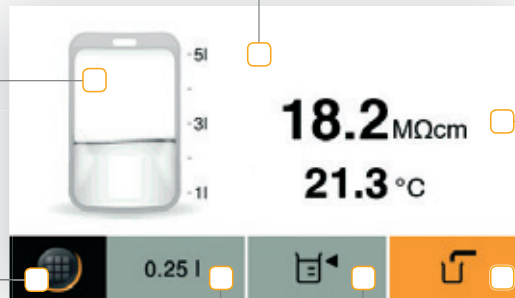


Current filling volume displayed for the bag (arium® mini & mini plus)

Touch-activated color display with easy-to-clean surface

Purified water quality and temperature displayed

Menu | Home function for settings and cleaning and care programs



Manual dispensing

Volume-controlled dispensing in 50 mL increments

Last volume dispensed is saved by the Favorites function



Intuitive Menu Navigation for Total Ease of Operation

- Icon-guided interface with touch-activated display
- Displays current measured values and messages
- Favorites function for repeated dispensing of identical volumes

arium® UV Lamp for Critical Analyses (185 | 254 nm)

- Optimal temperature gradient due to horizontal installation
- Degradation of total organic carbons (TOCs) by irradiation (≤ 5 ppb)
- Inactivates microorganisms by damaging their DNA

arium® Pretreatment Cartridge for Effective Purification

- Effective adsorption using high-quality activated carbon
- Special catalyst removes oxidants
- Impurities retained by reverse osmosis

The Right arium® mini Version for Your Daily Applications

Feed Water	arium® mini	arium® mini essential	arium® mini plus
Tap water			■
Pretreated water (RO DI line)		■	
Pretreated water (from supply container)	■		
Purified Water	arium® mini	arium® mini essential	arium® mini plus
Type 1 water	■	■	■
Type 3 water			■
Applications	arium® mini mini essential mini plus		arium® mini UV mini essential UV mini plus UV
Buffer and media preparation	■		■
Histology	■		■
Preparation of reagents, blank samples, etc.	■		■
ELISA (enzyme-linked immunosorbent assay)	■		■
AAS GF-AAS (atomic absorption spectrometry)	■		■
Photometry	■		■
SPE (solid phase extraction)			■
IC (ion chromatography)			■
ICP-MS (inductively coupled with plasma mass spectrometry)			■
GS-MS (gas chromatography coupled with mass spectrometry)			■
HPLC (high-performance liquid chromatography)			■
LC-MS (liquid chromatography coupled with mass spectrometry)			■

Sartorius Services

We Ensure the Quality of Your Results

At Sartorius, quality products go hand in hand with professional service. With our wide service offering, we will help guarantee the safe, reliable and optimal performance of your arium® mini. Just ask us and we will even cover the entire life cycle of your laboratory water system – from commissioning to equipment qualification to regular maintenance. Together with you, we will ensure the consistently high quality of your laboratory water purification.

Services at a Glance

- Installation and commissioning
Your advantage: Your system will operate reliably at peak performance from day one
- Equipment qualification (IQ | OQ)
Your advantage: You will meet all regulatory requirements (GMP | GLP)
- Regular preventative maintenance, including calibration, inspection and testing of your system and exchange of consumables
- Your advantages: Optimal operation of your system; reliable results; prevention of downtime or even equipment failure



Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49.551.308.0

France & Suisse Romande

Sartorius France
2, rue Antoine Laurent de Lavoisier
ZA de la Gaudrée
91410 Dourdan
Phone +33.1.70.62.50.00

Austria

Sartorius Austria GmbH
Modecenterstrasse 22
1030 Vienna
Phone +43.1.7965760.0

Belgium

Sartorius Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.90

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki
Phone +358.9.755.951

Hungary

Sartorius Hungária Kft.
Kagyló u. 5.
2092 Budakeszi
Phone +3623.457.227

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11
Phone +353.1.8089050

Italy

Sartorius Italy S.r.l.
Via Torino 3/5
20814 Varedo (MB)
Phone +39.0362.5557.11

Netherlands

Sartorius Netherlands B.V.
Phone +31.30.60.53.001
info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.6473830

Russian Federation

LLC "Sartorius RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg
Phone +7.812.327.53.27

Spain & Portugal

Sartorius Spain, S.A.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone Spain +34.913.586.095
Phone Portugal +351.800.855.800

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)
Phone +41.44.746.50.00

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ
Phone +44.1372.737159

Ukraine

LLS "Sartorius RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918

Americas

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1.631.254.4249
Toll-free +1.800.635.2906

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900

Canada

Sartorius Canada Inc
1173 North Service Road West, D4
Oakville, ON L6M 2V9
Phone +1.905.569.7977
Toll-Free +1.800.668.4234

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C.
Avenue Alberto del Campo 411
Floor 12 - The Office
15076 - San Isidro, Lima
Phone +51.1.441 0158

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800

China

Sartorius (Shanghai) Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China
Phone +86.21.6878.2300

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwun Tong
Kowloon, Hong Kong
Phone +852.2774.2678

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250

Japan

Sartorius Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan
Phone +81.3.3740.5408

Malaysia

Sartorius Malaysia Sdn. Bhd
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia
Phone +60.3.8996.0622

Singapore

Sartorius Singapore Pte. Ltd
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684
Phone +65.6872.3966

South Korea

Sartorius Korea Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400
Phone +82.31.622.5700

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310
Phone +66.2643.8361-6



◀ www.sartorius.com