SVISCISVS

Product Datasheet

Arium® Advance EDI

Ultimate Reliable Electrodeionisation



Advantages

- Reliable Consistently high Type 2 water quality thanks to the latest EDI technology
- Time-saving Use of innovative Bag technology, eliminates costly tank cleaning
- Optimized water consumption Automatic with iJust
- Easy to use Display with touch function and intuitive menu

Product Description

The Arium® Advance EDI provides Type 2 water in consistently high quality. The innovative iJust automatically optimizes water consumption and ensures the conscientious use of our environmental resources. In contrast to conventional water purification systems, the unique touch display with intuitive menu navigation makes it extremely easy to use.

With a flow rate of 5 or 10 L/h, the automatic RO membrane backflush, the latest EDI technology and a constant flow rate, the Arium[®] Advance EDI is the ideal choice for general and demanding laboratory applications.

Consistently high Type 2 water quality

In addition to pretreatment and purification by reverse osmosis modules, softening and electrochemical deionization is also carried out in the third purification step. By means of this modern EDI technology, the Arium® Advance EDI safely and reliably guarantees the removal of all impurities contained in the feed water.

Innovative Bag technology

The pure water is stored in the enclosed Arium[®] Bagtank system. This guarantees optimal storage of the pure water and protects against secondary contamination. Timeconsuming tank cleaning intervals are eliminated thanks to the interchangeable bag.

iJust

iJust stands for innovative technology that optimizes water production. Amongst other things, the intelligent Arium® software controls a valve on the concentrate outlet in accordance with the data that have been entered for CaCO₃ and CO₂. iJust therefore optimizes the product water quality and water consumption.

- The highest product water quality at all times
- Optimized, economical water consumption
- Guarantees a longer life of the downstream ultrapure water systems

Display with touch function

Simply navigate intuitively in the easy-to-use and clear menu by lightly touching the display – even with gloves.

Technical Specifications

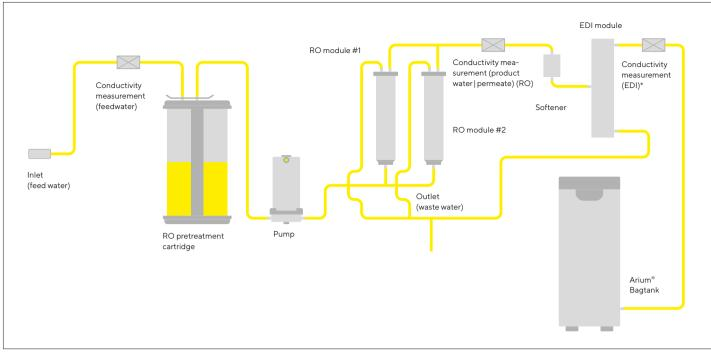
Dimensions: width × height × depth	35.0×50.1×45.1 cm
Empty weight	19 kg
Operating weight	26 kg
Power supply	100 – 240 VAC (± 10%); 50 – 60 Hz, 130 VA (max.)
Operating temperature	2°C-35°C at max. 80% relative humidity
Storage temperature	5°C-45°C at max. 80% relative humidity

Feed Water Quality

Exclusively potable tap water pursuant to the drinking water standards of the USA, the European Union, or Japan.

Input pressure'	2.0 - 6.9 bar
Temperature	2-30°C
Specific conductivity	<1,500 µS/cm compensated to 25 °C
тос	<2,000 ppb
Max. total hardness (max. CaCO₃)	360 ppm
Free chlorine	4ppm
Iron (total Fe content)	<0.1ppm
Manganese	<0.05ppm
Aluminum	<0.05ppm
CO₂ in solution	≤ 40 ppm
Fouling Index (SDI)	<5
Turbidity	<1 NTU
pH value	4-10

¹ Dynamic pressure/flow pressure 100 L/h



Flowchart Arium® Advance EDI (H2O-EDI-2-T)

Water Applications

Lab Water System Quality

	Advance EDI
Water Quality	
Type 2 Pure Water	•
Lab Water System by Daily Water Consumption	
Type 2 pure water <120 Liter/day (5 L/h)	•
Type 2 pure water <150 Liter/day (10 L/h)	•
Feed Application	
Feed ultrapure water systems	•
Feed distilled systems	•
Water for Laboratory devices (Autoclaves Washing Machine etc.)	•
General Laboratory Application	
Buffer, media and pH solutions	•
Histology	•
ELISA (Enzyme-Linked Immunosorbent Assay)	•
AAS (Atomic Absorption Spectroscopy)	•
Solutions for chemical analysis and synthesis	•

Product Water Quality

Water purification method	Particle filtration, adsorption using spherical activated carbon, catalyst, reverse osmosis, softening, electrochemical deionization, optional final particle and sterile filtration
Water type	Type 2 pure water
Production output ³	5 L/h or 10 L/h
Water dispensing flow rate⁵	Up to 3 L/min
Typical conductivity ¹	0.2-0.07 µS/cm compensated to 25 °C
Typical resistivity ¹	5–15 M Ω × cm compensated to 25 °C
Typical TOC reduction⁴	95%
Bacteria ²	<0.01 CFU/mL
Particle content ²	No particles > 0.22 µm
Particle and microorganism retention	>99%

 $^{_{1}}$ Depending on the feed water quality (CO_{^{2}} \!\leq\! 40 ppm) and temperature

² When using an Arium[®] Sterile Plus (Sartopore[®] 2 150, 0.22 μm pore size)

³ Depending on the feed water pressure, temperature, and condition of the RO module(s) ⁴ Depends on the type of organic contamination in the feed water

⁵When using an Arium[®] Bagtank design with pump, depending on hydrostatic pressure, connected accessories or end filter

Ordering Information

Arium® Advance EDI for the production of Type 2 pure water

Scope of supply: 1 Arium® Advance EDI, water guard, RO (reverse osmosis) module(s) and connection kit

Order number	Description
H2O-EDI-1-T	Arium® Advance EDI benchtop device, flow capacity 5 L/h
H2O-EDI-1-B	Arium® Advance EDI wall-mounted device, flow capacity 5 L/h
H2O-EDI-2-T	Arium® Advance EDI benchtop device, flow capacity 10 L/h
H2O-EDI-2-B	Arium® Advance EDI wall-mounted device, flow capacity 10 L/h

Accessories

Arium[®] Bagtanks

The most innovative tank system

- Integrated ventilation filter with non-return valve provides reliable protection against CO₂ pollution
- High flexibility through the 4 rollers available as an option
- Easy and fast exchange of the Arium[®] Bags
- High user safety due to the avoidance of cleaning chemicals

Description

The pure water is stored in the innovative enclosed Arium[®] Bagtank system. This system protects the prepared pure water against secondary contamination. The Sartorius Bagtank system enables consistent water quality over a prolonged period, thereby ensuring permanent, reproducible results. Unlike conventional water reservoirs, the Arium[®] Bag offers a high level of user safety and time savings, as there is no need for a complicated cleaning procedure with chemicals.

Arium[®] Bagtanks are housings which are equipped with Arium[®] Bags. The Arium[®] Bagtanks are available in 20 L, 50 L, and 100 L volumes. Their design is adaptable and saves space in any laboratory environment, and the optional rollers make this an extremely flexible system.



Integrated distributor pumps are a standard component of the 50 L and 100 L Bagtanks. A distributor pump is also available as an option for the 20 L Bagtank. In addition, a wall holder for the space-saving and user-friendly installation of this tank is also available.

Water dispensing flow rate

With pump ¹	up to 3.0 L/min
With pump, remote dispenser and Sterile Plus filter ¹	up to 2.0 L/min
Without pump ²	up to 1.5 L/min
Pump pressure	3 bar

Intended Use

Technical Specifications | Ordering Information

Materials		Order number	Description
Bagtank	Stainless steel plastic	H2O-AOV-20 ³	Arium® Bagtank 20 L, without pump, 1pc
Bag	S71 film		•
Tubing	PE silicone	H2O-AOV-50 ³	Arium® Bagtank 50 L, with pump 240 VAC, 50 Hz, 1pc
		H2O-AOV-50-US ³	Arium [®] Bagtank 50 L, with pump 115 VAC, 60 Hz, 1pc
Dimensions, excluding roller	rs and wall bracket [H×W×D]	H2O-AOV-50-W ³	Arium [®] Bagtank 50 L, without pump,
Bagtank 20	80.8×16.6×43.7 cm		1pc
Bagtank 50	85.2×25.4×58.7 cm	H2O-AOV-100 ³	Arium® Bagtank 100 L, with pump
Bagtank 100	85.2×51.4×58.7 cm		240 VAC, 50 Hz, 1pc
Bag 20 L	86.5×43.0 cm	H2O-AOV-100-US ³	Arium® Bagtank 100 L, with pump 115 VAC, 60 Hz, 1 pc
Bag 50 L	90.0×58.1cm	H2O-AOV-100-W ³	Arium [®] Bagtank 100 L, without pump, 1pc
Empty weight without Ariun Operating weight with filled		H2O-ADP-20	Pump Arium® Bagtank 20 L, 240 VAC, 50 Hz, 1pc
Bagtank 20	19 kg 40 kg	H2O-ADP-20-US	Pump Arium® Bagtank 20 L, 115 VAC, 60 Hz, 1 pc
Bagtank 50	33 kg 84 kg	H2O-ATR	Rollers for Arium [®] Bagtank 50 & bag-
Bagtank 100	47 kg 148 kg		tank 100, including fastening material, 4 pcs
Number of bags per tank		H2O-CBS-20	Arium [®] 20 L Bag for Arium [®] 20 L Bagtank, 2 pcs
Bagtank 20	1×20L	H2O-CBS-50	Arium [®] 50 L Bag for Arium [®] 50 L and
Bagtank 50	1×50 L		100 L Bagtank, 2 pcs
Bagtank 100	2×50L	H2O-ATB	Wall mount for Arium [®] Bagtank 20, 1pc
Power supply ¹	240 VAC (± 10 %), 50 Hz, 120 VA (max.)		190
Power supply US versions ¹	115 VAC (± 10 %), 60 Hz, 170 VA (max.)		
Operating temperature	2°C-35°C at max. 80% relative humidity		
	5°C-45°C at max. 80% relative		

1×¾" PLC quick-connect coupling

Water connection output		
Bagtank 20	1×¾" PLC quick-connect coupling	
Bagtank 50, Bagtank 100	2×¾" PLC quick-connect coupling	

' Bagtank 20 is supplied without a pump as standard, pump optionally

⁵ Bagtarik 20 is supplied without a particular partit particular particular particular particul

6

Arium[®] Bagtank Remote Dispenser

Ergonomic water withdrawal from the Arium $^{\otimes}$ with a working radius of up to 3.7 m

- Extended operating range of 3.7 m
- Available with height-adjustable stand or wall bracket
- Ergonomic design
- Easy-to-use
- Connection for Sterile Plus filter

Description

The Arium[®] Remote Dispenser is an ergonomically designed, easy-to-handle dosing unit which is ideally suited to the withdrawal of pure water.

Depending on the working environment, you can save space by mounting the remote dispenser on the wall or on a stand that is height-adjustable up to 70 cm. The stand enables relaxed working with optimal adaptation to the different sizes of the extraction vessels. The extended tube routing provides an operating range of 2.5 m from the Arium[®] Bagtank and a further 1.2 m from the stand.



To guarantee sterile and particle-free water removal, an Arium[®] Sterile Plus final filter with a $0.22\,\mu m$ pore size can be easily connected.

Technical Specifications | Ordering Information

Materials	
Stand	Aluminum (gray anodized)
Remote Dispenser	Plastic, white finish
Tubing	PE
Dimensions without tubing [W	/×H×D]
Remote Dispenser with stand	18.5 × 59.5 × 51.0 cm
Remote Dispenser with wall mounting bracket	9.0×10.0×28.5cm
Weight without tubing	
Remote Dispenser with stand	5.60 kg
Remote Dispenser with wall mounting bracket	0.46 kg

Order number	Description
613-AMDG'	Arium® Remote Dispenser including height-adjustable stand for connection to Arium® Bagtanks, 1pc
613-AMDG ²	Arium® Remote Dispenser including wall mounting kit for connection to Arium® Bagtanks, 1pc

Intended Use

- Arium[®] Bagtank 20*
- Arium[®] Bagtank 50
- Arium[®] Bagtank 100

* only in conjunction with an optional distributor pump

Arium[®] Water Guard

Early detection of leakages protects the laboratory

- Highly sensitive optical sensor
- Audiovisual alarm signals
- Automatic water stop in the case of leakage
- High-quality material, no corrosion
- Easy to install
- Integrated wall mounting bracket for solenoid valve

Description

Only the early detection of water leakages provides optimal protection against water damage in the laboratory. Leakages are registered by the highly sensitive optical sensor.

In contrast to conventional sensors, this sensor functions independently of conductivity measurement values as these are so low in the ultrapure water area that the activation of the guard is not guaranteed. Once a leak is detected, the water guard automatically locks the feed water inlet line. An acoustic warning is triggered immediately and the system status can be constantly controlled using the integrated LED display. With its sensitive optical sensors and high-quality materials, the Arium® Water Guard is perfect for all ultrapure and pure water systems.



Technical Specifications | Ordering Information

Sensor dimensions		
Diameter	5 cm	
Height	2.5 cm	
Cable length	2 m	

Tubing connections		
Input	%" Plug-in connector	
Output	%" Plug-in connector	
Power supply	100 - 240 VAC 50 - 60 Hz	

Order number	Description
610AWG1	Arium® Water Guard, 1pc

Intended Use

Consumables

Arium[®] Sterile Plus

Sterile and particle-free water dispensing

- Excellent service life and flow rates
- Integrity tested
- Validated according to HIMA and ASTM F-838-05
- Meets WFI quality standards pursuant to USP incl. USP plastic class VI test
- Production in accordance with DIN ISO 9001
- Easy to install
- Automatic venting
- Certified quality

Description

The Arium[®] Sterile Plus (Sartopore[®] 2 150) is a sterile, ready-to-use membrane filter capsule suitable for the most stringent requirements. Arium[®] Sterile Plus membrane filter capsules contain a hydrophilic, heterogeneous polyethersulfone double membrane. These provide an excellent service life and flow rates. The capsule is attached in the end position by means of a quick connector and reliably removes all particles and microorganisms in the last quarter purification step. A hydrophobic PTFE membrane at the

Technical Specifications | Ordering Information

Materials	
Membranes	Asym. Polyethersulfone
Bell assembly	Polycarbonate
Other plastics	Polypropylene
Pore size	0.45 µm × 0.22 µm
Filtration area	0.015 m²
Input and Output	¼" Plug-in connector
Sterilization (max. 3 cycles)	Autoclaving at 134 °C, 1 bar, 30 min.
Max. diffusion	1 mL/min @ 2.5 bar
Min. bubble point	3.2 bar

032 83 2065	

farthest point "upstream" allows for easy and clean ventilation of the capsule.

All pleated Arium[®] Sterile Plus membrane filter units are validated as sterile filters for biopharmaceutical application according to the HIMA and ASTM F-838-05 guidelines (documentation available). During the manufacturing process, every capsule is integrity-tested to meet the highest quality standards and safety regulations.

Order number	Description
5441307H4CE	Arium® Sterile Plus (Sartopore® 2 150), 0.22 µm pore size, 1 pc

Intended Use

Device type:

Arium[®] Bagtank Remote Dispenser

Arium[®] RO Pretreatment Cartridge

Safe protection of the RO module

- Fast and effective adsorption of impurities through high-grade activated carbon
- \bullet 5 μm depth filter for the retention of particles
- Highly efficient catalyst for removing free chlorine
- Patented cartridge design for easy installation

Description

The best protection for a downstream reverse osmosis (RO) membrane is the combination of spherical, catalytic effective activated carbon, a catalyst and a depth filter. It reliably removes oxidation agents, such as free chlorine, heavy-metal ions and particulate contaminants from the feed water of the system.

A special catalyst is an integral part of pre-treatment. It is particularly efficient at removing free chlorine and at a lower temperature and/or higher pH value compared to activated carbon alone.

The patented cartridge design ensures minimal time expenditure with ultra-easy installation and exchange.



Technical Specifications | Ordering Information

Materials		
Housing	High-quality polypropylene	
Cleaning media	Spherical catalytic effective activated carbon plus polypropylene filter car- tridge with nom. 5 µm separation rate	
Dimensions $[W \times H \times D]$	18×26×11cm	
Operating weight	3.5 kg	
Feed water requirements	See "Technical Specifications" page 2	

Order number	Description
H2O-CPFAD-1	Arium® RO Pretreatment Cartridge, 1 pc

Intended Use

Arium[®] RO Modules

Reverse osmosis modules with low-energy membranes

- Highly efficient reverse osmosis membranes, optimized water consumption
- Low-energy membranes for ecological and economical operation
- Backflush with product water increases the service life
- Easy replacement
- Constant flow
- Consistently high water quality

Description

The Arium[®] RO modules consist of two independent membranes whose design guarantees easy installation and reliable operation. Each of the two modules contains a low-energy reverse osmosis membrane in a polypropylene housing.

The housing has connections for feed water, permeate (product water) and concentrate (discarded water). The RO modules enable an ideal water yield, thereby optimizing the water consumption. At the same time, up to 98 % of the salts are typically retained. Thanks to the backflush



with permeate, particles and salts are removed from the surface of the membrane.

This results in a longer service life and lower system maintenance costs. In addition, this backflush function on restarting the system after a standstill allows for the immediate dispensing of high quality water.

Technical Specifications | Ordering Information

Materials	
RO membranes	Low-energy membrane made of polyamide
Housing	Polypropylene
Dimensions for each module	
Height	30.8 cm
Diameter	7.8 cm
Weight	0.468 kg
Product Water Quality	See "Technical Specifications" page 2

Order number	Description	
H2O-CRO-H-1	Arium [®] RO Module, 1pc	
H2O-CRO-H- 2	Arium [®] RO Module, 2 pcs	

Intended Use

Device type:

Arium[®] Advance EDI

Arium[®] Softener Cartridge

For maximum service life of the EDI module

- Consistently high water quality
- Long service life
- Effective CaCO₃ elimination

Description

It is sensible to soften the feed water to improve protection of the EDI module.

The cartridge reliably removes traces of alkaline earth ions from the water, thereby guaranteeing consistently high water quality and a long service life of the EDI module.



Technical Specifications | Ordering Information

Materials

Housing

High-purity polypropylene

Filling material Ultrapure ion exchange resin

Order number	Description
H2O-CSO-1	Arium® Softener Cartridge, 1pc

Intended Use

Arium[®] RO Cleaning Set

Maximum service life of the RO module

- Effective removal of scaling and metal deposits
- Elimination of organic compounds
- Dispersion of colloids
- Stable pH values
- Gentle on materials

Description

Two-stage cleaning kit for removing scaling and organic contaminants.

The alkaline substance contains non-foaming surfactants that dissolve organic compounds, disperse colloids and can be quickly removed again from the membrane surface. Cleaning efficiency depends on the pH value that is steadily maintained by buffer substances through a large temperature range.

The acidic cleaning agent to remove scaling contains chelate and reducing agents in order to dissolve metallic deposits. The ideal pH value also remains consistently low over a wide range during cleaning in this case thanks to the buffers.



Technical Specifications | Ordering Information

Ingredients	
Alkaline cleaner	HEDTA, ethanolamine, triethanol- amine
Acidic cleaner	HEDTA, phosphoric acid, citric acid

Order number	Description
H2O-CCS	Arium [®] RO Cleaning Set, 1pc

Intended Use

Device type:

Arium[®] Advance EDI

Sartorius Service

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 operation ofyour Arium[®] systems. Just ask us and we will even cover the entire life cycle of your laboratory water system – from commissioning to qualification to regular maintenance. Together with you, we will ensure the consistently high quality of your laboratory water purification.

Our 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

Get more information now at: **www.sartorius.com/service**

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 further information, visit www.sartorius.com

