

arium® advance RO

Reverse Osmosis System with iJust



Reverse osmosis system for water purification

Optimized water consumption – automatic with iJust

Easy-to-use – unique glass display with touch function

Easy handling – intuitive menu navigation

Description

The arium® advance RO provides type 3 reverse osmosis water of the highest quality. The innovative iJust makes the best use of environmental resources by automatically optimizing water consumption. Unlike conventional water purification systems, the unique display with touch functions in combination with the intuitive menu navigation creates an easy to use system.

With a flow rate of up to 8, 16 or 24 l/h, the automatic RO membrane backflush and a constant flow rate, the arium® advance RO is the optimal choice for daily laboratory applications.

Applications

- Feed water for ultrapure water systems
- Rinsing of glass and laboratory containers
- Feed water for various laboratory devices, such as humidifiers, autoclaves and lab glassware washers

iJust

iJust is based on an innovative technology which optimizes the water production. The intelligent arium® software controls a valve at the concentrate outlet according to measurement data for CaCO_3 and CO_2 . iJust thereby optimizes product water quality and water consumption

- Premium product water quality at all times
- Optimized, economic water consumption
- Guaranteed longer life of down-stream 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.

High system flexibility

The space-saving installation of the device as bench-top, wall mounted or built-in unit integrates it perfectly into any laboratory environment.

Technical Specifications

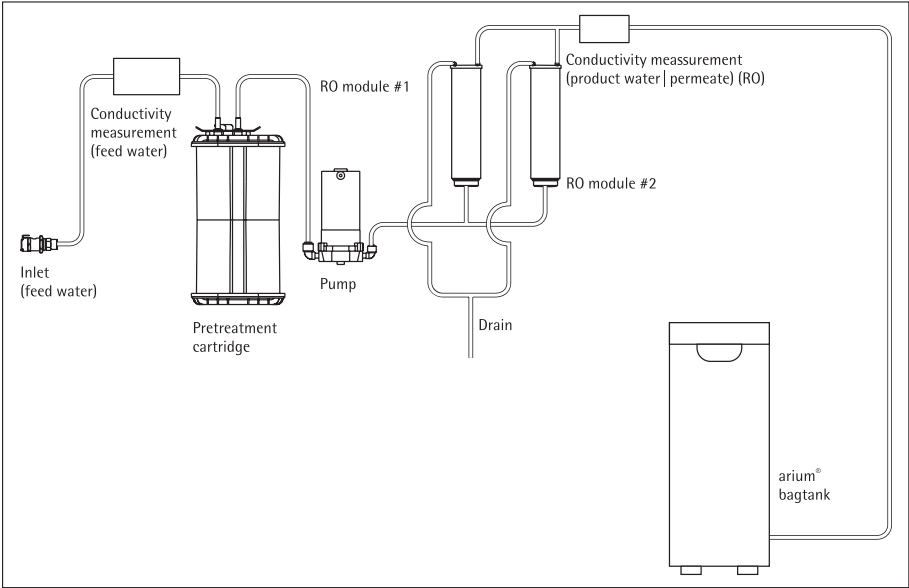
Technical Specifications

| | |
|--------------------------------|--|
| Water purification methods | Particle filtration, spherical activated carbon, catalyst, reverse osmosis, optional final particle and sterile filtration |
| Dimensions: Width×height×depth | 35.0×50.1×45.1 cm |
| Empty weight | approx. 15 kg |
| Operating weight | approx. 22 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 |

arium® advance RO reverse osmosis system for producing type 3 pure water

Equipment supplied: 1 arium® advance H2O-RO, RO (reverse osmosis) Module(s) and Connection Set

| Order no. | Description |
|------------|---|
| H2O-RO-1-T | arium® advance RO Bench-Top system in a compact design for every laboratory workstation, product water flow rate 8 l/h |
| H2O-RO-1-B | arium® advance RO space-saving Wall-Mounted system with integrated wall bracket, product water flow rate 8 l/h |
| H2O-RO-2-T | arium® advance RO Bench-Top system in a compact design for every laboratory workstation, product water flow rate 16 l/h |
| H2O-RO-2-B | arium® advance RO space-saving Wall-Mounted system with integrated wall bracket, product water flow rate 16 l/h |
| H2O-RO-3-T | arium® advance RO Bench-Top system in a compact design for every laboratory workstation, product water flow rate 24 l/h |
| H2O-RO-3-B | arium® advance RO space-saving Wall-Mounted system with integrated wall bracket, product water flow rate 24 l/h |



Flowchart arium® advance RO (H2O-RO-2-T)

Product water quality

| | |
|---|-------------------|
| Water type | Pure water Type 3 |
| Product water flow rate ³ | 8, 16 or 24 l/h |
| Water dispensing flow rate ⁴ | 3 l/min |
| Typical conductivity ¹ | < 20 µS/cm |
| Typical resistivity ¹ | < 0.05 MΩ × cm |
| Microorganism content ² | < 1 CFU/1000 ml |
| Particle content ² | < 1/ml |
| Typical ion retention | up to 98% |
| Retention of dissolved organic substances (MW > 300 Dalton) | > 99 % |
| Particle and microorganism retention | > 99 % |

Feed Water Quality

Exclusively potable water according to the drinking water regulations of the USA, the European Union, or Japan.

| | |
|---|------------------------------------|
| Input pressure | 0.5 – 6.9 bar, recommended > 2 bar |
| Temperature | 2 – 30 °C |
| Specific conductivity | < 1500 µS/cm compensated to 25 °C |
| TOC | < 2000 ppb |
| Max. permanent hardness (max. CaCO ₃) | 360 ppm |
| Free chlorine | < 4 ppm |
| Iron (total Fe content) | < 0.1 ppm |
| Fouling Index (SDI) | < 5 |
| Turbidity | < 1 NTU |
| pH value | 4 – 10 |

¹ Depending on the feed water and temperature

² When using an arium® Sterile Filter (Sartopore® 2 150)

³ Depending on the feed water pressure, temperature, and condition of the RO Modules

⁴ Depending on the arium® bagtank design, hydrostatic pressure, connected accessories and end filter

arium® bagtanks

The most innovative tank system



- Integrated venting filter with check valve securely protects against CO₂ contamination
- Optional rollers provide highest flexibility
- Easy replacement of arium® bags, with no time-consuming sanitization
- High operation safety, no use of cleaning agents

Description

Pure water is stored in the innovative, closed arium® bagtank system. Here, the prepared pure water is securely protected from secondary contaminations. The arium® bagtank system enables high water quality over an extended period, thereby providing for long-term, reproducible results. In contrast to typical water tanks, the arium® bag provides a high degree of user safety and saves time by the elimination of complex sanitization processes with chemicals.

arium® bagtanks are housings that are equipped with arium® bags. The arium® bagtanks are available in 20 l, 50 l, and 100 l volumes. Their space-saving design is adaptable to any laboratory environment, and the optional rollers make the system extremely flexible.

Integrated distribution pumps are a standard components of the 50 l and 100 l bagtanks. An optional additional pump is also available for a 20 l bagtank. Also available is a wall bracket for a space-saving and user-friendly installation of the arium® bagtank 20.

Water dispensing flow rate

| | |
|---|-----------------|
| with pump ¹ | up to 3.0 l/min |
| with pump, dispense gun and sterile-grade filter ² | up to 2.0 l/min |
| without pump ² | up to 1.5 l/min |

Intended Use

System type:

arium® comfort I and comfort II,
arium® advance RO and advance EDI

¹ bagtank 20 is supplied without pump, pump optionally available

² Value only applies for bagtank 20 dispensing site at the same height or lower than the tank connection

³ Note: The arium® bag is not included in the equipment supplied with the arium® bagtank

Technical Specifications

Materials

| | |
|---------|----------------------------|
| bagtank | Stainless steel plastics |
| bag | S71 Film |
| Tubing | PE silicon |

Dimensions, without rollers and wall brackets [H × W × D]

| | |
|-------------|-----------------------|
| bagtank 20 | 80.8 × 16.6 × 43.7 cm |
| bagtank 50 | 85.2 × 25.4 × 58.7 cm |
| bagtank 100 | 85.2 × 51.4 × 58.7 cm |
| bag 20 l | 86.5 × 43.0 cm |
| bag 50 l | 90.0 × 58.1 cm |

Empty weight without arium® bag | Operating weight with filled arium® bag

| | |
|-------------|----------------|
| bagtank 20 | 19 kg 40 kg |
| bagtank 50 | 33 kg 84 kg |
| bagtank 100 | 47 kg 148 kg |

Number of bags per tank

| | |
|-------------|--------------|
| bagtank 20 | 1 × 20 liter |
| bagtank 50 | 1 × 50 liter |
| bagtank 100 | 2 × 50 liter |

| | |
|---------------------------|---------------------------------------|
| Power supply ¹ | 240 VAC (± 10%), 50 Hz, 120 VA (max.) |
|---------------------------|---------------------------------------|

| | |
|--|---------------------------------------|
| Power supply, US versions ¹ | 115 VAC (± 10%), 60 Hz, 170 VA (max.) |
|--|---------------------------------------|

| | |
|-----------------------|--|
| Operating temperature | 2°C–35°C at max. 80% relative humidity |
|-----------------------|--|

| | |
|---------------------|--|
| Storage temperature | 5°C–45°C at max. 80% relative humidity |
|---------------------|--|

| | |
|------------------------|-----------------------------|
| Water connection input | 1 × 3/8" PLC Quick Coupling |
|------------------------|-----------------------------|

Water connection output

| | |
|-------------------------|-----------------------------|
| bagtank 20 ¹ | 1 × 3/8" PLC Quick Coupling |
| bagtank 50, bagtank 100 | 2 × 3/8" PLC Quick Coupling |

Order Number Description

| | |
|-----------------------------|---|
| H20-AOV-20 ³ | arium® bagtank 20 l, without pump, qty. 1 unit |
| H20-AOV-50 ³ | arium® bagtank 50 l, with pump 240 VAC 50 Hz, qty. 1 unit |
| H20-AOV-50-US ³ | arium® bagtank 50 l, with pump 115 VAC 60 Hz, qty. 1 unit |
| H20-AOV-50-W ³ | arium® bagtank 50 l, without pump, qty. 1 unit |
| H20-AOV-100 ³ | arium® bagtank 100 l, with pump 240 VAC 50 Hz, qty. 1 unit |
| H20-AOV-100-US ³ | arium® bagtank 100 l, with pump 115 VAC 60 Hz, qty. 1 unit |
| H20-AOV-100-W ³ | arium® bagtank 50 l, without pump, qty. 1 unit |
| H20-ADP-20 | arium pump® bagtank 20 l, 240 VAC 50 Hz, qty. 1 unit |
| H20-ADP-20-US | arium pump® bagtank 20 l, 115 VAC 60 Hz, qty. 1 unit |
| H20-ATR | Rollers for arium® bagtank 50 & bagtank 100, incl. fittings, qty. 4 units |
| H20-CBS-20 | arium® 20 l bag for arium® 20 l bagtank, qty. 2 units |
| H20-CBS-50 | arium® 50 l bag for arium® 50 l and 100 l bagtank, qty. 2 units |
| H20-ATB | Wall bracket for arium® bagtank 20, qty. 1 unit |

arium® bagtank Dispense Gun

Ergonomic water dispensing from the arium® bagtank with a working radius of up to 3.7 m



- Extended work area of 3.7 m
- Available with height-adjustable stand or with wall brackets
- Ergonomic design
- Easy-to-use
- Sterile-grade filter can be connected

Description

The arium® dispense gun is an ergonomically designed, easy-to-use dispensing unit and is perfectly suited for dispensing pure water.

Depending on the work environment, either position the dispense gun on the wall to save space or on the 70 cm height-adjustable stand. Moreover, the stand lets you work in a relaxed position and allows optimal adjustment to different sized sampling vessels. The extended tube guide ensures a work area 2.5 m away from the arium® bagtank or pressure tank and a further 1.2 m from the stand.

A sterile-grade Sartopore® 2 150 capsule with 0.2 µm pore size can be easily connected for guaranteed sterile and particle-free water dispensing.

| Order Number | Description |
|--------------|--|
| 613-AMDG1 | arium® Dispense Gun inclusive height-adjustable Stand, qty. 1 unit |
| 613-AMDG2 | arium® Dispense Gun inclusive wall mounting kit, qty. 1 unit |

Technical Specifications

Materials

| | |
|--------------|--------------------------|
| Stand | Aluminum (gray anodized) |
| Dispense gun | Plastic, white finish |
| Tubing | PE |

Dimensions without tubing [W × H × D]

| | |
|---|-----------------------|
| Dispense gun with stand | 18.5 × 59.5 × 51.0 cm |
| Dispense gun with wall mounting bracket | 9.0 × 10.0 × 28.5 cm |

Weight without tubing

| | |
|---|---------|
| Dispense gun with stand | 5.60 kg |
| Dispense gun with wall mounting bracket | 0.46 kg |

Intended Use for arium bagtanks:

arium® bagtank 20*
arium® bagtank 50
arium® bagtank 100
arium® pressure tank 30
arium® pressure tank 50
arium® pressure tank 70
arium® pressure tank 100

* suitable only with optionally available distribution pump

arium® Water Guard

Early Detection of Leakages Protects the Laboratory



- Highly sensitive optical sensor
- Audiovisual alarm signals
- Automatic water-stop in case of leakage
- High-grade material, non-corrosive
- Easy to install
- Integrated wall mounting bracket for magnetic 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 pure and ultrapure water area that the activation of the guard is not guaranteed. Once a leakage 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. The sensitive, optical sensors and high-grade materials mean that the arium® water guard is perfect for all ultrapure and pure water systems.

| Order Number | Description |
|--------------|---------------------------------|
| 610AWG1 | arium® Water Guard, qty. 1 unit |

Technical Specifications

Sensor Dimensions

| | |
|--------------|--------|
| Diameter | 5 cm |
| Height | 2.5 cm |
| Cable length | 2 m |

Tubing Connections

| | |
|--------------|----------------------|
| Input | 3/8" plug connection |
| Output | 3/8" plug connection |
| Power supply | 100–240 VAC 50–60 Hz |

Intended Use

System type:
arium® comfort I and comfort II
arium® pro, pro DI, pro UF, pro UV and pro VF
arium® advance RO and EDI
arium® 611, 612 and 613

arium® Sterile Filter

Sterile and Particle-free Water Dispensing



- Excellent service lifetime and flow rate
- Integrity tested
- Validated acc. to HIMA and ASTM F-838-05
- Meets WFI quality standards acc. to USP incl. USP plastic class VI test
- Manufacture acc. to DIN ISO 9001
- Easy to install
- Automatic venting
- Certified quality

Description

The arium® Sterile Filter (Sartopore® 2 150) is a sterile, ready-to-use membrane filter capsule for the most stringent requirements. Sartopore® 2 150 membrane filter capsules contain a hydrophilic, heterogeneous polyethersulfone double membrane. It enables an excellent service life and output. The capsule is attached, by a quick connector, at the final position and reliably removes all particles and micro-organisms in the last water purification step. A hydrophobic PTFE membrane at the farthest point „upstream“ allows for easy and clean venting of the capsule.

All pleated Sartopore® 2 membrane filter units are validated as sterile filters for bio-pharmaceutical applications according to the HIMA and ASTM F-838-05 guidelines (documentation available). During the manufacturing process, every unit is tested for integrity in order to meet the highest quality standards and safety regulations.

Technical Specifications

Materials

| | |
|----------------------------------|---|
| Membranes | Asym. Polyethersulfone |
| Filling bell | Polycarbonate |
| Other plastics | Polypropylene |
| Pore size | 0.45 µm + 0.2 µm |
| Filtration area | 0.015 m ² |
| Input and output | 1/4" plug connection |
| 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 |

Intended Use

On Dispense Guns and Display-Dispense Unit for system type:
arium® comfort I and comfort II
arium® pro, pro DI, pro UF, pro UV and pro VF
arium® 611
arium® bagtank Dispense Gun
arium® Dispense Gun

Order Number

Description

| | |
|------------------|--|
| 5441307H4--CE--B | arium® Sterile Filter (Sterile-grade Sartopore® 2 150 capsules), 0.2 µm pore size, qty. 5 units |
|------------------|--|

arium® RO Pretreatment Cartridge

Safe Protection of the RO Modules



- Quick and effective adsorption of impurities by high-grade activated carbon
- 5 µm depth filter for the retention of particles
- Highly efficient catalyst for removing free chlorine
- Easy to install by patented cartridge design

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 oxidizing agents, such as free chlorine, heavy metal ions and particles, from the feed water.

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

Apart from preventing deposits, the catalyst reduces fouling processes and inhibits micro-biological growth. The patented cartridge design saves time with its easy installation and replacement.

Technical Specifications

Materials

| | |
|----------------------------|---|
| Housing | High-grade polypropylene |
| Cleaning media | Spherical catalytic effective activated carbon plus polypropylene filter cartridge with nom. 5 µm separation rate |
| Dimensions [W × H × D] | 18 × 26 × 11 cm |
| Operating weight | 3.5 kg |
| Feed water requirements | see "Technical specifications" page 2 |

Intended Use

System type:
arium® advance RO and advance EDI
61316, 61215

Order Number

Description

| | |
|----------------|--|
| 613CPF05-----V | arium® RO Pretreatment Cartridge Set, qty. 2 units |
|----------------|--|

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 contain a low-energy reverse osmosis membrane in a polypropylene housing. The housing has connections for feed water, permeate (product water) and concentrate (waste water). The RO Modules typically enable high recovery rates. This optimizes the water consumption while still retaining at least up to 98% of the ions. 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 allows for the immediate dispensing of high quality water.

Technical Specifications

Materialien

| | |
|--------------|---------------------------------------|
| RO membranes | Low-energy membrane made of polyamide |
| Housing | Polypropylene |

Dimension of each Module

| | |
|-----------------------|---------------------------------------|
| Height | 30.8 cm |
| Diameter | 7.8 cm |
| Weight | 0.468 kg |
| Product water quality | see "Technical specifications" page 2 |

Intended Use

System type:
arium® comfort I
arium® advance RO
arium® 61316 and 61215

Order Number

Description

| | |
|---------------|---------------------------------|
| 613CPM4-----V | arium® RO Modules, qty. 2 units |
| 613CPM4 | arium® RO Module, qty. 1 unit |

arium® RO Module Cleaning Set

Maximum Service Life of the RO Module



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

Description

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

The alkaline substance contains non-foaming tensides that dissolve organic compounds, disperse colloids and can be quickly removed from the membrane surface. The cleaning efficiency depends on the pH value that can be safely maintained over a wide temperature range by the contained buffer substances.

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

Technical Specifications

Ingredients

| | |
|------------------|--------------------------------------|
| Alkaline cleaner | HEDTA, ethanolamine, triethanolamine |
| Acidic cleaner | HEDTA, phosphoric acid, citric acid |

Intended Use

System type:
arium® comfort I and comfort II
arium® advance RO and advance EDI
arium® 612 and 613

Order Number

Description

| | |
|---------|--|
| H20-CCS | arium® RO Module Cleaning Set, qty. 1 unit |
|---------|--|

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Weighing Technology GmbH
Weender Landstrasse 94-108
37075 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

France & Suisse Romande

Sartorius France S.A.S.
ZI Les Paluds
Avenue de Jouques – CS 31090
13781 Aubagne Cedex
Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Austria GmbH
Franzosengraben 12
1030 Vienna
Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde
Phone +32.2.756.06.71
Fax +32.2.253.45.95

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki
Phone +358.9.755.951
Fax +358.9.755.95.292

Hungary

Sartorius Hungária Kft.
Kagyló u. 5.
2092 Budakeszi
Phone +3623.457.227
Fax +3623.457.147

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11
Phone +353.1.8089050
Fax +353.1.8089388

Italy

Sartorius Italy S.r.l.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)
Phone +39.055.63.40.41
Fax +39.055.63.40.526

Netherlands

Sartorius Netherlands B.V.
Edisonbaan 24
3439 MN Nieuwegein
Phone +31.30.6053001
Fax +31.30.6052917

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.6473830
Fax +48.61.6473839

Russian Federation

LLC "Sartorius ICR"
and LLC "Biohit"
Uralskaya str. 4, Lit. B
199155, Saint-Petersburg
Phone +7.812.327.5.327
Fax +7.812.327.5.323

Scandinavia

Sartorius Nordic A/S
Hoerskaetten 6D
2630 Taastrup, Denmark
Phone +45.7023.4400
Fax +45.4630.4030

Spain & Portugal

Sartorius Spain S.A.
Offices in Madrid:
C/ Isabel Colbrand, 10-12, of. 70
28050 Madrid
Phone Spain +34.902.123.367
Phone Portugal +351.800.855.800
Fax Spain +34.91.358.84.85
Fax Portugal +351.800.855.799

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)
Phone +41.44.746.50.00
Fax +41.44.746.50.50

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ
Phone +44.1372.737159
Fax +44.1372.729927

America

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1.631.254.4249
Toll-free +1.800.635.2906
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Av. Dom Pedro I, 241
Bairro Vila Pires
Santo André
São Paulo
Cep 09110-001
Phone +55.11.4451.6226
Fax +55.11.4451.4369

Canada

Sartorius Canada Inc.
2179 Dunwin Drive #4
Mississauga, ON L5L 1X2
Phone +1.905.569.7977
Toll-Free +1.800.668.4234
Fax +1.905.569.7021

Mexico

Sartorius de México S.A. de C.V.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México
Phone +52.5555.62.1102
Fax +52.5555.62.2942

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius Scientific
Instruments (Beijing) Co., Ltd.
33 Yu An Road, Airport Industrial Park
Zone B,
Shunyi District, Beijing 101300,
P.R.China
Phone +86.10.8042.6300
Fax +86.10.8042.6486

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwung Tong
Kowloon, Hong Kong
Phone +852.2774.2678
Fax +852.2766.3526

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Stedim Japan K.K.
4th Fl., Diawa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan
Phone +81.3.4331.4300
Fax +81.3.4331.4301

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
Fax +60.3.8996.0755

Singapore

Sartorius Singapore Pte. Ltd
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528
Phone +65.6872.3966
Fax +65.6778.2494

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
Fax +82.31.622.5799

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310
Phone +66.2643.8361-6
Fax +66.2643.8367

