

arium® 611

Ultrapure Water Systems



Description

Type 1* laboratory water purification system designed to meet the requirements of reagent grade water for critical and routine analysis.

* Water exceeds ASTM, NCCLS, ISO and USP reagent water standards.

Advanced Design Features Include:

Carbon-resin Technology

A unique carbon-resin combination enables the 611 system to produce ultrapure water with a resistivity up to 18.2 MΩ × cm with Total Organic Carbon (TOC) levels as low as < 1 ppb.

Four-line Alphanumeric VDU

Continuously indicates water quality in MΩ × cm or μS/cm with compensated or non-compensated temperature.

Removable Display & Dispense Unit

Enables high quality water to be provided at point of use up to two meters from the system's installation point.

Standby Mode

Water recirculates 15 minutes every hour of inactivity to maintain purity.

Timed (Volume) Dispenser

Automated dispensing included.

Quality

Inert materials used for all wetted parts guarantee the highest water purity.

Security

The 0.2 μm PESU membrane filter capsule ensures a bacteria-free filtrate.

Specifications

Unit Dimensions

Width:	43 cm (16.85")
Height:	48 cm (19")
Depth:	33.4 cm (13.15")

Removable Display | Dispense Unit

Width:	30.2 cm (11.91")
Height:	11.3 cm (4.44")
Depth:	15.7 cm (6.21")

System Weight

611DI	15.6 kg (33 lbs)
611UF	16.3 kg (36 lbs)
611UV	16.8 kg (37 lbs)
611VF	17.4 kg (38 lbs)

Operation Weight

611DI	23.6 kg (52 lbs)
611UF	24.8 kg (55 lbs)
611UV	25.3 kg (56 lbs)
611VF	26.4 kg (58 lbs)

Clearances

Side	17 cm (7") for handling
Top	5 cm (2") for ventilation
Front	35 cm (14") for door opening

Electrical Power Requirements

100-240 V AC 50-60 Hz single phase

Interface

RS-232 serial port

Inlet Feed Water Requirements

Specific Resistivity

RO water:	20 KΩ × cm (50 μS/cm)
Distilled water:	250 KΩ × cm (4 μS/cm)
Deionized water:	50 KΩ × cm (20 μS/cm)

TOC:	< 1000 ppb
Turbidity:	< 1 N.T.U.
Silica:	< 1000 ppb

Minimum inlet pressure: Gravity

Maximum inlet pressure: 7 bar | 103 psi

Laboratory Water Systems

Ambient Temperature & Humidity

Operation:	5°C–28°C (41°F–82.4°F) 80% relative humidity non condensing
Storage:	5°C - 45°C (41°F–113°F) 80% relative humidity non-condensing

Ordering Information

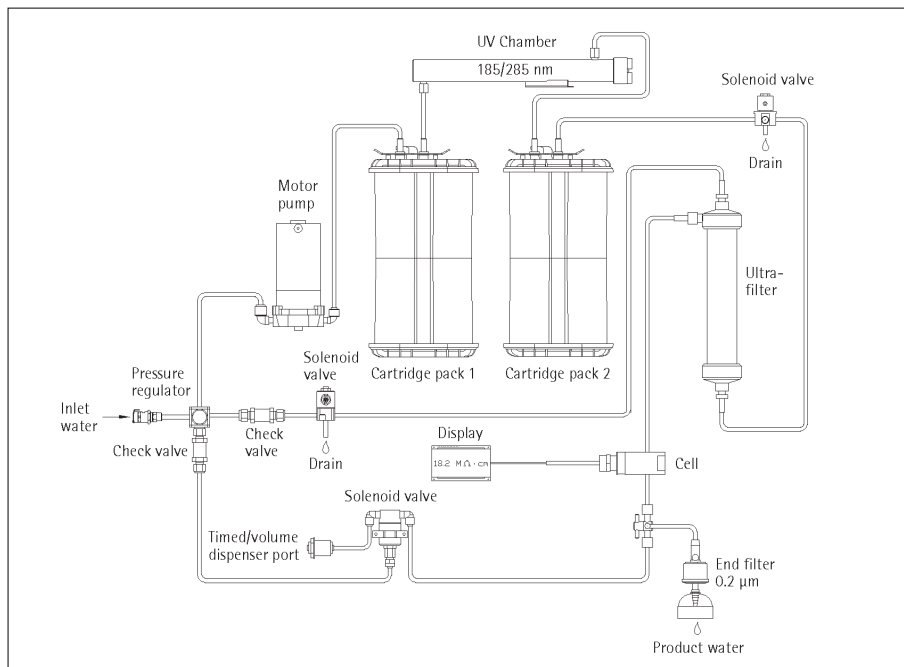
Systems

611DI	arium® basic system
611UF	arium® system with ultrafilter
611UV	arium® system with UV lamp
611VF	arium® system with ultrafilter & UV lamp

Consumables & Accessories

5441307H4-- NO--B	Sartopore 2 150 (5 final filters)
611CDS2	Sanitization kit (2 syringes)
611CDS6	Sanitization kit (6 syringes)
611CEL1	UV lamp
611CDU5	UF hollow fiber filter
611AEC1	Calibration resistor
611AMDG1	Remote Dispenser (dispense gun)
611AKD1	Removable display & dispense unit, complete with wall bracket
611APR1	Printer for data collection
Cartridge Kit	As the cartridge kit is application and feed water specific, please contact us for a complete data sheet

Flow Diagram



Product Water Quality	611DI	611UV	611UF	611VF
Resistivity (MΩ × cm)	18.2	18.2	18.2	18.2
TOC (ppb)* @ 100 ppb feed water	< 4	< 1	< 4	< 1
TOC (ppb)* @ 1000 ppb feed water	< 20	< 10	< 20	< 10
Absorbance units (AUFS @ 210 nm)	n.a.	< 0.005	n.a.	< 0.005
Endotoxin (EU/ml)	n.a.	n.a.	< 0.001	< 0.001
Particles (@ 0.2 μm)	< 1/ml	< 1/ml	< 1/ml	< 1/ml
Bacteria (CFU/1000 ml)	< 1	< 1	< 1	< 1
Flow rate (l/min)**	up to 2	up to 2	up to 1.5	up to 1.5

* depends on inlet water quality

** at a pressure of 2 bar, without final filter

Sartorius AG
Weender Landstrasse 94–108
37075 Goettingen, Germany
Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius.com

USA +1.631.2544249
UK +44.1372.737100
France +33.1.69192100
Italy +39.055.634041
Spain +34.91.3586100
Japan +81.3.33293366

Specifications subject to change
without notice.
Printed and copyrighted
by Sartorius · W/sart-236 · G
Publication No.: SLG2031-e06014
Order No.: 85030-517-58