

## arium® 61315

### Reverse Osmosis System



Designed to meet the demands of today's laboratories for RO treated water.

With production volumes of up to 15 l/h per hour, 99% rejection of monovalent and polyvalent ions, bacteria and viruses, automated membrane back-flushing, and with its compact design, the arium® 61315 is the ideal choice for general laboratory applications, such as dish-washing water, humidifiers and as a pretreatment step prior to ultrapure laboratory water systems.

Low-energy, high-output and thin-film composite membranes dramatically reduce the volume of water rejected to waste and increase the % of recovery.

Energy consumption is reduced by the incorporation of a uniquely designed RO module housing, which minimizes the need for solenoid valves to regulate and control permeate flow.

A single solenoid valve controls an automatic flush-fill cycle, which maintains the volume of the pressure tank while at the same time rinsing the RO modules, thereby reducing ionic concentration at the RO membrane surface. This not only conserves power, it also reduces maintenance and increases the reliability of the system.

Water storage is simplified by the use of a closed pressure tank. This significantly reduces the costs of maintenance due to airborne contamination, while reducing gas uptake by product water in storage. A pressurized storage tank provides more flexibility as to its location, without the need for an expensive distribution pump to deliver the water to a point of use.

#### Specifications

##### Unit Dimensions

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

##### System Weight

14 kg (31 lbs)

##### Operation Weight

20 kg (44 lbs)

##### Electrical Power Requirements

100-240 V AC 50-60 Hz single phase

##### Inlet Feed Water Requirements

Potable tap water feed only (feed water to meet drinking water standards of the U.S. EPA, European Union or Japan)

Minimum inlet pressure:	1 bar   15 psi
Maximum inlet pressure:	6.8 bar   100 psi
Temperature:	5°C-28°C (41°F-82.4°F)
TDS (max ppm CaCO <sub>3</sub> ):	360
Silt density index:	< 5%
Free chlorine:	0.1 ppm
Turbidity:	< 1 NTU
Iron (total as Fe):	< 0.1 ppm

##### Ambient Temperature & Humidity

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

## Laboratory Water Systems

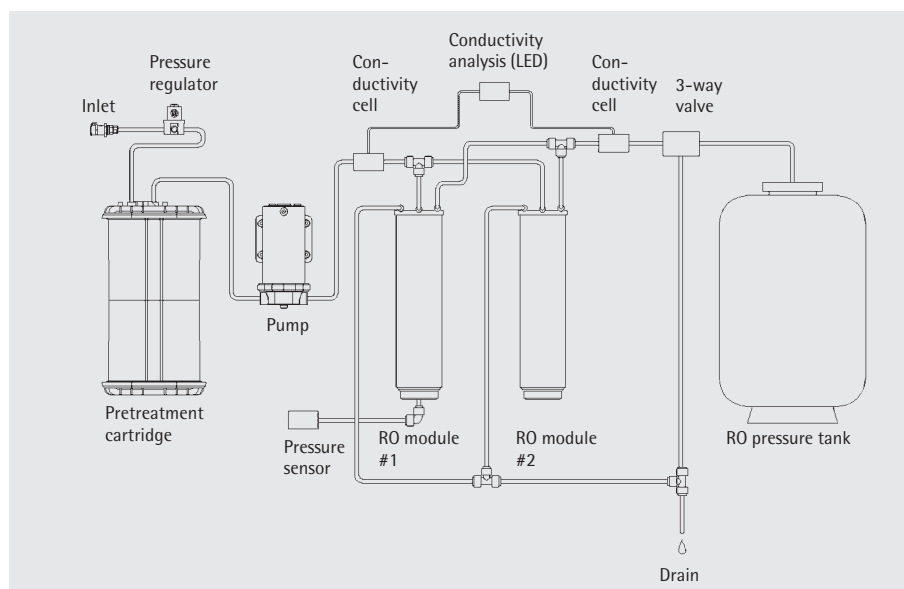
### Product Water Quality

#### Rejection rates

Monovalent ions:	Up to 98%
Polyvalent ions:	Up to 99%
Particles:	> 99%
Microorganisms:	> 99%
Dissolved organics (>300 MW):	> 99%
Reject flow [l/h]*	11-12
Product flow rate [l/h]* (± 20% at 25°C)	15
Maximum recovery rate*	Approx. 55%

\* Depends on feed water quality and temperature

### RO System Schematic



### Ordering Information

(A complete system is composed of the arium® 61315, a tank, 1 × pretreatment cartridge and 2 × RO modules)

61315	arium® RO system
613CPF05-----V	2 × pretreatment cartridges
613CPM1-----V	2 × RO modules
611CDS2	Sanitization kit for the tank (2 syringes)
613CDS2	Sanitization kit for the RO modules (2 syringes)
613AMDG1	Dispense gun for RO tank
613APV30	31-liter tank**
613APV60	76-liter tank**
613APV150	100-liter tank**
61315030F05M1A	Complete Reverse Osmosis System with a tank capacity of 31 litres** (includes arium® 61315, tank, 2 × RO modules, 2 pretreatment cartridges + sanitizing syringes for RO modules & tank)
61315060F05M1A	Complete Reverse Osmosis System with a tank capacity of 76 litres** (includes arium® 61315, tank, 2 × RO modules, 2 pretreatment cartridges + sanitizing syringes for RO modules & tank)
61315150F05M1A	Complete Reverse Osmosis System with a tank capacity of 100 litres** (includes arium® 61315, tank, 2 × RO modules, 2 pretreatment cartridges + sanitizing syringes for RO modules & tank)

\*\* at 2.5 bar inlet pressure

Sartorius AG  
Weender Landstrasse 94-108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289

[www.sartorius.com/arium](http://www.sartorius.com/arium)

USA +1.631.2544249  
UK +44.1372.737100  
France +33.1.69192100  
Italy +39.055.634041  
Spain +34.91.3588566  
Japan +81.3.33295533

Specifications subject to change without notice.  
Printed in Germany on paper that has been bleached without any use of chlorine.  
W/sart-231 · G  
Publication No.: SLG2026-e06014  
Order No.: 85030-515-38