



arium® 613CPF05-----V

Reverse Osmosis Pretreatment Cartridge



Description

The Sartorius arium® 613CPF05-----V is an encapsulated cartridge, containing pretreatment media and 5 µm depth filter as a convenient, one-piece disposable. Designed for installation within the arium® 61315 RO, the pretreatment cartridge removes free chlorine and particulate matter from the system's feed water to protect the RO membranes located downstream.

Features and Benefits

Pretreatment packs utilize a patented design to encapsulate the treatment media and filter. Installation and removal are thereby simplified, reducing downtime to a minimum.

Water entering the pretreatment pack flows through the purification media bed before passing to a 5 µm depth filter. Down-flow operation over the media bed helps to remove larger particles and thus increases the service life of the particle-removing filter.

Special high-quality 20 × 50 mesh activated carbons are used to remove oxidizing agents. The small particle size carbons permit fast diffusion of contaminants and efficient adsorption or surface reaction. The small carbon particle size also allows more activated carbon mass, which results in longer service life.

A combination of proprietary catalytic media is an integral part of the pretreatment mix. The catalytic media are extremely efficient at removing free chlorine and less susceptible to the temperature and pH limits associated with activated carbon. The catalyst ensures optimum chlorine removal from feed waters of low temperature and/or high pH. Added benefits of the catalyst include limited fouling and scale prevention and bacterial control. The catalytic action reduces hardness scale (calcium and magnesium) buildup on the membrane.

Specifications

Materials of Construction

Housing:	Virgin unfilled polypropylene
Purification media:	– 20 × 50 mesh granular activated carbon – Chlorine hardness catalyst – 5 µm [nominal] PP filter cartridge
Unit Dimensions:	Width: 18 cm (7") Height: 26 cm (10.2") Depth: 11 cm (4.3")
Operation Weight:	3.5 kg

Inlet Feed Water Requirements

Potable tap water feed only	
Minimum inlet pressure:	1 bar 15 psi
Maximum inlet pressure:	6.8 bar 100 psi
TDS (max ppm CaCO ₃):	360
Silt density index:	< 5%
Turbidity:	< 1 NTU
Iron (Total as Fe):	< 0.1 ppm

Ambient Temperature and Humidity

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

Ordering Information

Order number:	613CPF05-----V
Quantity:	2 cartridges

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

USA +1.631.2544249
UK +44.1372.737100
France +33.1.69192100
Italy +39.055.505671
Spain +34.91.3586100
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-183a · G
Publication No.: SLG2035-e06022
Order No.: 85030-518-98