



## arium® SterilePlus

Sterile and particle-free water dispensing

### Advantages

- 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



### Product Description

arium® SterilePlus (Sartopore® 2 150) is a sterile, ready-to-use membrane filter capsule suitable for the most stringent requirements. Made from a hydrophilic heterogeneous polyethersulfone double membrane. The membrane filter capsules ensure excellent durability and flow rate capacities, almost free of extractables.

#### Applications

arium® SterilePlus capsules are used for the particle removal and sterile filtration of injection solutions, therapeutics, biotech products, buffers, and ultrapure water.

#### Flow Rate Capacity | Durability

A unique pleated filter design combined with the highly asymmetric pore structure of the membrane allows for excellent flow rates and great durability, especially compared to conventional flat filter systems.

#### Easy to Use

Using a ¼" quick connector, arium® SterilePlus capsules can be connected to the outlet of an arium® mini, pro, and comfort lab water system or to an arium® dispense gun.

#### Automatic Ventilation

A hydrophobic PTFE membrane at the farthest point "upstream" allows for easy ventilation of the capsule while simultaneously preventing product loss.

#### Microbiological Retention

arium® SterilePlus capsules with 0.2 µm final filter membranes are validated as sterile filters according to HIMA and ASTM F-838-05 guidelines.

#### Quality

Each and every element undergoes a diffusion test for integrity in order to meet the highest quality standards.

#### Documentation

arium® SterilePlus capsules are developed and produced in line with a DIN ISO 9001 certified quality management system. During the manufacturing process, every unit is tested for integrity in order to meet the highest quality standards and safety regulations. Validation documentation is available.

## Technical Specifications

### Design

Hydrophilic, heterogeneous double-layered polyethersulfone membrane

### Material

Final filter membrane	asymmetrical, polyethersulfone
Prefilter membrane	asymmetrical, polyethersulfone
Support fleece	polypropylene
Inner core	polypropylene
Connections	polypropylene
Housing	polypropylene
Bell assembly	polycarbonate

### Pore Size Combination

0.45  $\mu\text{m}$  + 0.2  $\mu\text{m}$

### Filtration Area

0.015  $\text{m}^2$

### Connections

Inlet	1/4 inch quick connector
Outlet	6 mm tube connector

### Max. Permitted Differential Pressure

4 bar at 20°C

2 bar at 80°C

### Max. Permitted Counter Pressure

2 bar at 20°C

### Sterilization

Autoclaving at 134°C, 1 bar, 30 min.

**Note:** no inline steam sterilization for capsules.

### Autoclaving Cycles

Max. 3

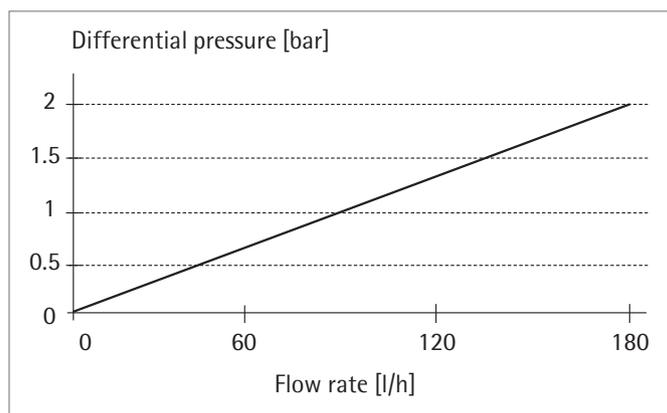
### pH Range

pH 1 to pH 14

### Integrity Test Data for arium® SterilePlus (Sartopore® 2 150 Capsules)

Pore size [ $\mu\text{m}$ ]	0.2
Test pressure [bar]	2.5
Max. diffusion [ml/min]	1
Min. B.P. [bar]	3.2

### Water Flow Rate



at 20°C

## Ordering Information

Order Number	Quantity per Package
5441307H4--CE	1

Sartorius Lab Instruments  
GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen, Germany  
Phone +49.551.308.0  
Fax +49.551.308.3289  
www.sartorius.com

USA Toll-free +1.800.635.2906  
UK +44.1372.737159  
France +33.1.70.62.50.00  
Italy +39.0362.5557.11  
Spain +34.913.586.095  
Russian Federation +7.812.327.53.27  
Japan +81.3.3740.5408

Specifications subject to change without notice.  
Copyright Sartorius Lab Instruments GmbH & Co. KG.  
Printed in the EU on paper bleached without chlorine.  
Publication No.: SLG2048-e170302  
Order No.: 85032-540-15  
Ver. 03 | 2017