

# Aquasart® Plus

Membrane Filter Cartridge for  
Food & Beverage Application



## Applications

Aquasart® Plus membrane filter cartridges are used for final filtration prior to bottling for products, such as the following:

- Non-carbonated water
- Near-water beverages

## Product Information

Aquasart® Plus filter cartridges feature a high-performance, double-layer pleated hydrophilic polyethersulfone membrane. Specially developed for the filtration of water and non-carbonated drinks. Aquasart® Plus filter cartridges guarantee the highest biological product safety in combination with enhanced total throughput performance.

## Performance

Aquasart® Plus filter cartridges incorporate a special double layer membrane. This new combination, coarse prefilter membrane followed by a final membrane, has been developed to meet the demanding process requirements in terms of high total throughput performance. In combination with an optimally designed prefilter, Aquasart® Plus achieves maximum service life performance and minimizes filtration costs.

## Safety

Aquasart® Plus filter cartridges guarantee highest microbiological safety, which can be checked before and after every use with an integrity test.

## Quality

Aquasart® Plus membrane filters are designed and manufactured according to an ISO 9001 certified quality management system. All raw materials which are used to produce Aquasart® Plus are in accordance with EU regulation 1935/2004/EC or are listed in the Regulations (EU) No. 10/2011 relating to plastic materials and article intended to come into contact with food. All used materials meet FDA requirements.

## Technical Data

### Specifications

Materials	
Prefilter Membrane	Polyethersulfone, asymmetric
Endfilter Membrane	Polyethersulfone, asymmetric
Support Fleece	Polypropylene
Core	Polypropylene
End Caps	Polypropylene
O-Rings	Silicone (optional EPDM or Fluoroelastomer)

Pore Size
0.45 µm + 0.2 µm
0.20 µm + 0.1 µm

### Available Sizes | Filtration Area

Cartridges		
Size 2	20"	1.2 m <sup>2</sup>  12.9 ft <sup>2</sup>
Size 3	30"	1.8 m <sup>2</sup>  19.4 ft <sup>2</sup>
Size 4	40"	2.4 m <sup>2</sup>  25.9 ft <sup>2</sup>

### Available Adapters

### Cartridges

25, 28

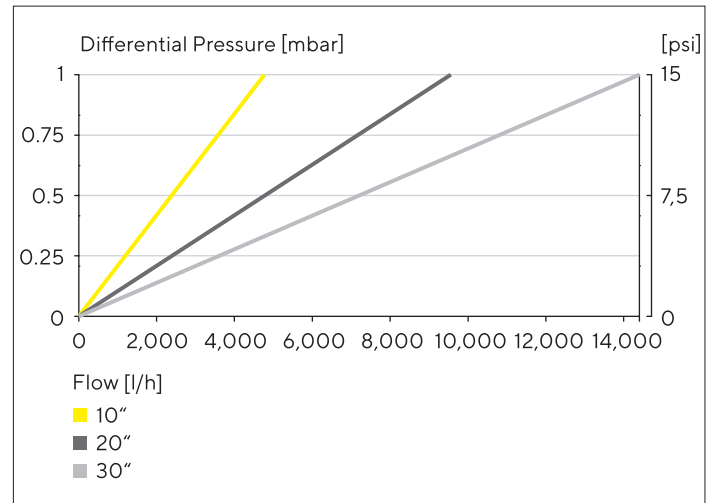
Operating Parameters	
Max. Allowable Differential Pressure	5 bar   72.5 psi at 20°C 2 bar   29 psi at 80°C
Sterilization	Hot water at 85–95°C, 30 minutes, ΔP max 0.5 bar Steam at 110°C, 30 minutes, ΔP max 0.5 bar
Max. Allowable Back Pressure	2 bar   29 psi at 20°C
Steaming cycles	min. 100

## Compatibility

Aquasart® Plus filter cartridges fulfill all requirements of the current USP for “Sterile Water for Injection”. This includes testing for Toxicity according to USP Plastic Class VI Test.

Aquasart® Plus filter cartridges are fully suitable for use in the food & beverage industry pH-compatibility: pH 1 to pH 14.

## Water Flow Rates for 10", 20" and 30" Cartridges 0.1 µm and 0.2 µm



## Ordering Information

Order Code	Pore Size [µm]	Pack Size [Pieces]	Test Pressure [bar   psi]	Max. Diffusion [ml/min]	Min. Bubble Point [bar   psi]
<b>Standard Cartridges</b>					
538**07H2	0.2	1	2.5   36	36	3.2   46
538**07H3	0.2	1	2.5   36	54	3.2   46
538**07H4	0.2	1	2.5   36	72	3.2   46
538**58K3	0.1	1	4.0   58	72	n.a
538**58K4	0.1	1	4.0   58	96	n.a

\*\* Adapter 25, 28

**Germany**

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

**USA**

Sartorius Stedim North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 368 7178

 For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)