

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 m² 12.9 ft²
Size 3	30"	1.8 m² 19.4 ft²
Size 4	40"	2.4 m² 25.9 ft²

Available Adapters

Cartridges

25, 28

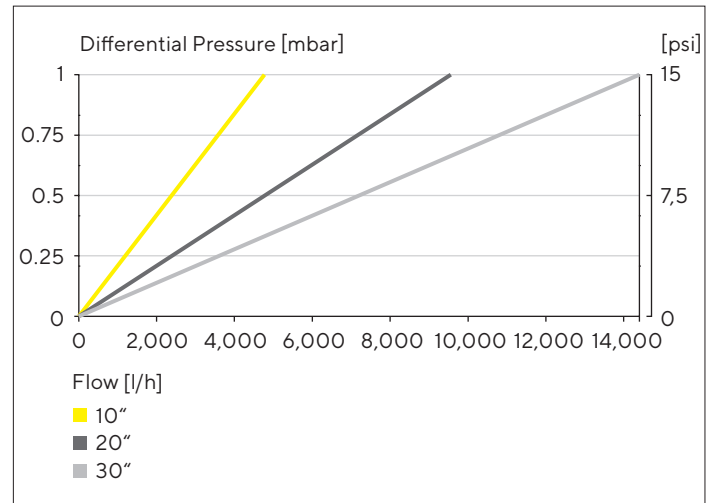
Operating Parameters	
Max. Allowable Differential Pressure	5 bar 72.5 psi at 20°C 2 bar 29 psi at 80°C
Sterilization	Hot water at 85 – 95°C, 30 minutes, ΔP max 0.5 bar Steam at 110°C, 30 minutes, ΔP max 0.5 bar
Max. Allowable Back Pressure	2 bar 29 psi at 20°C
Steaming cycles	min. 100

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]
Standard Cartridges					
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