



# Aquasart® Plus

Membrane Filter Cartridge for  
Food & Beverage Application

## 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.



### 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

### 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

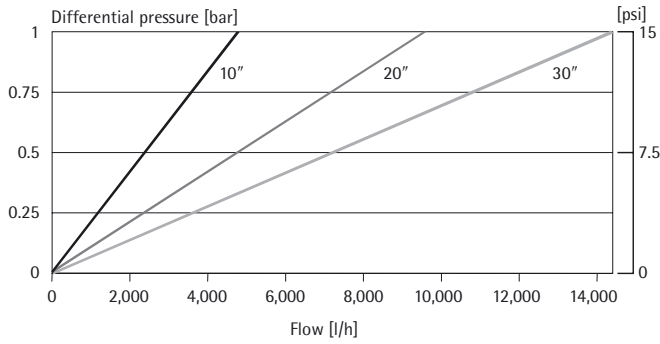
Operating Parameters	
Max. Allowable Differential Pressure	5 bar   72.5 psi at 20°C 2 bar   29 psi at 80°C
Sterilization	Hot water at 85–95°C, 30 minutes, ΔP max 0.5 bar Steam at 110°C, 30 minutes, ΔP max 0.5 bar
Max. Allowable Back Pressure	2 bar   29 psi at 20°C
Steaming cycles	min. 100

### 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

# Sales and Service Contacts

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

## Europe

### Germany

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49.551.308.0

Sartorius Stedim Systems GmbH  
Robert-Bosch-Strasse 5-7  
34302 Guxhagen  
Phone +49.5665.407.0

### France

Sartorius Stedim FMT S.A.S.  
ZI des Paluds  
Avenue de Jouques - CS 91051  
13781 Aubagne Cedex  
Phone +33.442.845600

Sartorius Stedim France SAS  
ZI des Paluds  
Avenue de Jouques - CS 71058  
13781 Aubagne Cedex  
Phone +33.442.845600

### Austria

Sartorius Stedim Austria GmbH  
Modecenterstrasse 22  
1030 Vienna  
Phone +43.1.7965763.18

### Belgium

Sartorius Stedim Belgium N.V.  
Rue Colonel Bourg 105  
1030 Bruxelles  
Phone +32.2.756.06.80

### Hungary

Sartorius Stedim Hungária Kft.  
Kagyló u. 5  
2092 Budakeszi  
Phone +36.23.457.227

### Italy

Sartorius Stedim Italy S.r.l.  
Via dell'Antella, 76/A  
50012 Antella-Bagno a Ripoli (FI)  
Phone +39.055.63.40.41

### Netherlands

Sartorius Stedim Netherlands B.V.  
Phone +31.30.60.25.080  
[filtratie.nederland@sartorius-stedim.com](mailto:filtratie.nederland@sartorius-stedim.com)

### Poland

Sartorius Stedim Poland Sp. z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn  
Phone +48.61.647.38.40

### Russian Federation

LLC "Sartorius Stedim RUS"  
Vasilyevsky Island  
5<sup>th</sup> line 70, Lit. A  
199178 St. Petersburg  
Phone +7.812.327.53.27

### Spain

Sartorius Stedim Spain, S.A.U.  
Avda. de la Industria, 32  
Edificio PAYMA  
28108 Alcobendas (Madrid)  
Phone +34.913.586.098

### Switzerland

Sartorius Stedim Switzerland AG  
Ringstrasse 24 a  
8317 Tagelswangen  
Phone +41.52.354.36.36

### U.K.

Sartorius Stedim UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9 QQ  
Phone +44.1372.737159

### Ukraine

LLC "Sartorius Stedim RUS"  
Post Box 440 "B"  
01001 Kiev, Ukraine  
Phone +380.44.411.4918

## Americas

### USA

Sartorius Stedim North America Inc.  
5 Orville Drive, Suite 200  
Bohemia, NY 11716  
Toll-Free +1.800.368.7178

### Argentina

Sartorius Argentina S.A.  
Int. A. Ávalos 4251  
B1605ECS Munro  
Buenos Aires  
Phone +54.11.4721.0505

### Brazil

Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 - SP- Brasil  
Phone +55.11.4362.8900

### Mexico

Sartorius de México, S.A. de C.V.  
Libramiento Norte de Tepotzotlan s/n,  
Colonia Barrio Tlacateco,  
Municipio de Tepotzotlan,  
Estado de México,  
C.P. 54605  
Phone +52.55.5562.1102  
[leadsmex@sartorius.com](mailto:leadsmex@sartorius.com)

## Asia | Pacific

### Australia

Sartorius Stedim Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175  
Phone +61.3.8762.1800

### China

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.  
3rd Floor, North Wing, Tower 1  
No. 4560 Jinke Road  
Zhangjiang Hi-Tech Park  
Pudong District  
Shanghai 201210, P.R. China  
Phone +86.21.6878.2300

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.  
Beijing Branch Office  
No. 33 Yu'an Road  
Airport Industrial Park Zone B  
Shunyi District, Beijing 101300  
Phone +86.10.8042.6501

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.  
Guangzhou Branch Office  
Room 1105  
Xing Guang Ying Jing Building  
No. 119, Shui Yin Road  
Yue Xiu District, Guangzhou 510075  
Phone +86.20.3836.4193

### India

Sartorius Stedim India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra  
Nelamangala Tq  
562 123 Bangalore, India  
Phone +91.80.4350.5250

### Japan

Sartorius Stedim Japan K.K.  
4th Fl., Daiwa Shinagawa North Bldg.  
8-11, Kita-Shinagawa 1-chome  
Shinagawa-ku, Tokyo, 140-0001 Japan  
Phone +81.3.4331.4300

### Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.  
Lot L3-E-3B, Enterprise 4  
Technology Park Malaysia  
Bukit Jalil  
57000 Kuala Lumpur, Malaysia  
Phone +60.3.8996.0622

### Singapore

Sartorius Stedim Singapore Pte. Ltd.  
10 Science Park Rd  
The Alpha #02-13/14  
Singapore Science Park II  
Singapore 117684  
Phone +65.6872.3966

### South Korea

Sartorius Korea Biotech Co., Ltd.  
8th Floor, Solid Space B/D,  
PanGyoYeok-Ro 220, Bundang-Gu  
SeongNam-Si, GyeongGi-Do, 463-400  
Phone +82.31.622.5700



▶ [www.sartorius-stedim.com](http://www.sartorius-stedim.com)