

Aquasart® Plus XL

Membrane Filter Cartridge for Food & Beverage Application

Product Information

- High total throughput for economic filtration
- Integrity testable by Diffusion- & Bubble Point Test
- Highly economic sizing
- Excellent filtration capabilities

Performance

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

Quality

Aquasart® Plus XL 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 XL 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.

Safety

Aquasart® Plus XL filter cartridges guarantee highest microbiological safety, which can be checked before and after every use with an integrity test (>7 LRV for B. diminuta).

Compatibility

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

Applications

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

- Non-carbonated water
- Near-water beverages

Technical Data

Available Sizes	Filtration Area	Max. Diffusion at 1.5 bar 22 psi [ml/min]	Min. Bubble Point [bar psi]
Cartridges			
Size 2	1.6 m ² 17.2 ft ²	50	2.8 40.5
Size 3	2.4 m ² 25.8 ft ²	75	2.8 40.5

Max. Allowable Differential Pressure

Cartridges

5 bar | 72.5 psi at 20°C

2 bar | 29 psi at 80°C

Max. Allowable Back Pressure

2 bar | 29 psi at 20°C

(for all elements)

Materials

Prefilter Membrane

Polyethersulfone, asymmetric

Endfilter Membrane

Polyethersulfone, asymmetric

Support Fleece

Polypropylene

(In-Line Steam sterilizable & autoclavable)

Polyester

(γ-irradiatable or γ-irradiatable | autoclavable)

Max. Operating Pressure

Cartridges

5 bar | 72.5 psi at 20°C

Pore Size Combination

0.8 μm + 0.2 μm

Regulatory Compliance

- Each individual element is tested for integrity by Diffusion test
- Fully validated as sterilizing grade filters according to ASTM current F-838-05 guidelines
- Designed, developed and manufactured in accordance with an ISO 9001 certified Quality Management System
- Meet or exceed the requirements for WFI quality standards set by the current USP
- Non pyrogenic according to USP Bacterial Endotoxins
- USP Plastic Class VI Test
- Non fiber releasing according to Code of Federal Regulations (CFR), section 210.3(b)(6) and 211.72
- Material used meet FDA requirements and fulfill EU regulation 1935/2004/EC, Regulations (EU) No. 10/2011 relating to plastic materials and article intended to come into contact with food.

Sterilization

Cartridges

In-Line Steam Sterilization

121°C, 0.3 bar, 20 min.

Min. 25 Sterilization Cycles

or

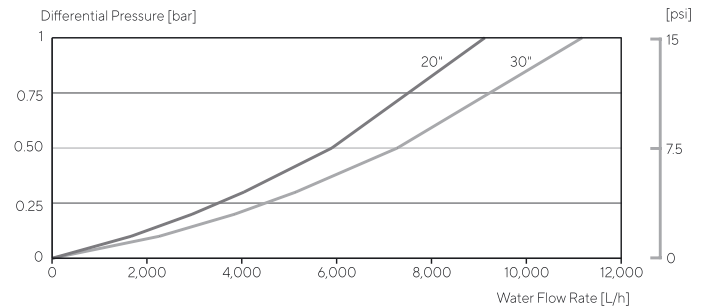
Autoclaving

121°C, 2 bar, 30 min

Min. 25 Sterilization Cycles

Performance

Water Flow Rates for Aquasart® Plus XL 20" & 30" cartridges



Technical References

Validation Guide 2552028-000

Ordering Information



54B



07

G



Adapter

25: 2 Flange Bayonet adapter
with 226 double o-ring

28: 3 Flange Bayonet adapter
with 222 double o-ring

Filter Size

2: 1.6 m² | 17.2 ft² (20")

3: 2.4 m² | 25.8 ft² (30")

Sales and Service Contacts

For further contacts, visit
sartorius.com

Germany

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

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906