



Sartobev PS

Membrane Filter for Particle, Colloids and Bioburden Reduction in Food & Beverage Application

Membrane Filter Cartridge



Description

Sartobev PS filter cartridges feature a high-performance, single-layer pleated PES membrane. The ideal choice for the filtration in the food and beverage industry, especially developed for the prefiltration of wine and sparkling wine.

Sartobev guarantees the retention of particle and yeast cells. Enhanced total throughput performance in combination with highest product safety is achieved.

Application

Sartobev PS membrane filter cartridges are used for the final filtration prior to bottling or as an integrity testable prefilter for such beverages as the following:

- Wine
- Sparkling Wine
- Spirits

Product Performance

Sartobev PS filter cartridges incorporate a PES membrane. This new polyethersulfon membrane has been developed to meet the demanding process requirements in terms of pressure pulsation and sterilization cycles in the food and beverage industry. Sartobev PS achieves maximum total throughput to minimize the filtration costs. The membrane is chemically resistant within the pH-range of 1-14. Caustic soda and acids can be used for a chemical regeneration.

Integrity-testable and Quality

The integrity of Sartobev PS membrane filter cartridges can be easily checked by performing an integrity test before and after filtration to guarantee the highest biological safety. This membrane filter is designed and manufactured according to an ISO 9001 certified quality management system. All materials used in the construction of Sartobev PS meet FDA requirements for use in food and beverage applications.

Technical Specifications

Material

| | |
|---|---------------------------|
| Filter medium | Single-layer PES membrane |
| Protective non-woven drainage layers | Polypropylene |
| Core outer support end caps adapter | Polypropylene |
| O-rings | Silicone |

Pore Size

0.65 µm
1.0 µm

Filtration area

0.8 m² per 10" element

Available Sizes | Filtration Area

| | | |
|--------|-----|---------------------|
| Size 2 | 20" | 1.60 m ² |
| Size 3 | 30" | 2.40 m ² |
| Size 4 | 40" | 3.20 m ² |

Available Adapter

25, 28

Operating Parameters

| | |
|-------------------------------|---|
| Maximum differential pressure | 5 bar 73 psi at 20°C 2 bar 29 psi at 80°C 0.5 bar 7psi at 121°C |
| Maximum back pressure | 2 bar 29 psi at 20°C |

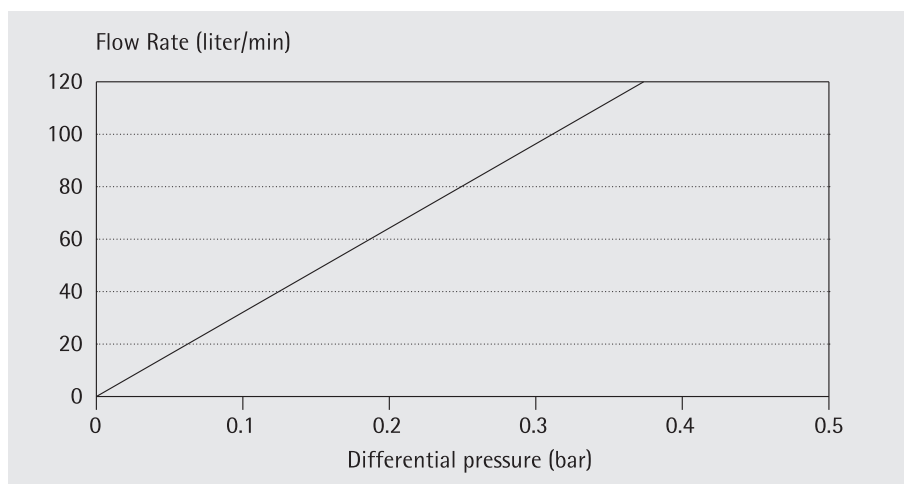
Technical Specifications

| | |
|-------------------------------|---|
| Recommended flow rate | 1.000 wine/h/30" 3.000 liter water/h/30" |
| Sterilization | Hot water (85–95°C) or Steam (110°C), at least 100 cycles |
| Microbiological recovery rate | Saccharomyces cerevisiae > 10 ⁷ /cm ² |
| ph stability | ph 1–14 |

Integrity Test Data

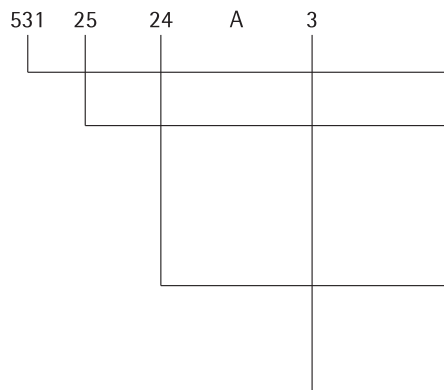
| | |
|------------------------------|------------|
| Test pressure | 0.8 bar |
| Max. diffusion (10" element) | 8 ml |
| Test time | 10 minutes |
| Bubble point | > 0.8 bar |

Flow rate for water Sartobev PS height 10"



Ordering Information

Cartridges



Sartobev PS

Adapter

- 25 S-adapter on top, 2-flange bayonet adapter with double O-ring on bottom
- 28 S-adapter on top, 3-flange bayonet adapter with double O-ring on bottom

Retention Rate

- 05 0.65 µm
- 24 1.0 µm

Height

- 2 20"
- 3 30"
- 4 40"

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.90.2110935
Russian Federation +7.812.6100821
Japan +81.3.3740.5407

Specifications subject to change without notice. Printed and copyrighted by Sartorius Stedim Biotech GmbH - W
Publication No.: SGK2011-e12083
Order No.: 85034-538-99
Ver. 08 | 2012