

Sartobev PS

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



Product Information

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.

Description

Sartobev PS guarantees the retention of particles 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
- Water

Product Performance

Sartobev PS filter cartridges incorporate a PES membrane. This new polyethersulfone 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 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 the raw materials used to produce the Sartobev PS cartridge comply with the European Union Regulation (EC) No. 1935/2004 as well as the Regulation (EU) No. 10/2011 concerning plastic materials and articles intended to come into contact with foodstuffs.

These guidelines for plastics allow the use in food and beverage applications.

All materials used meet the requirements of the FDA.

Technical Data

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 Adapters

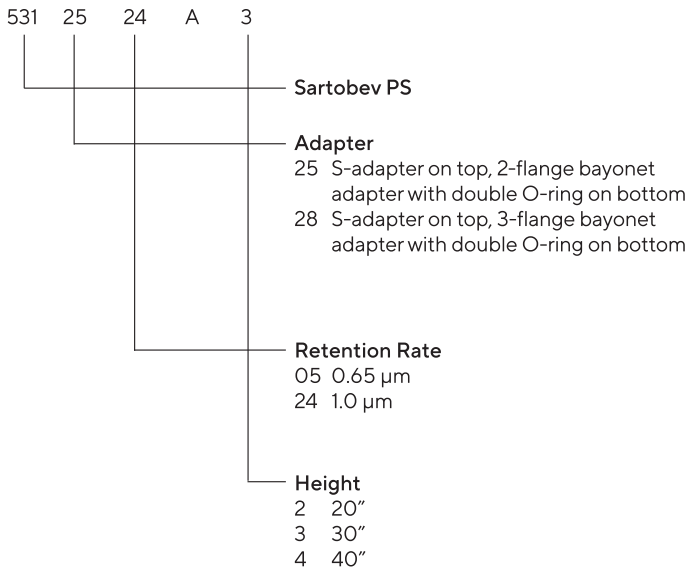
25, 28

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 |

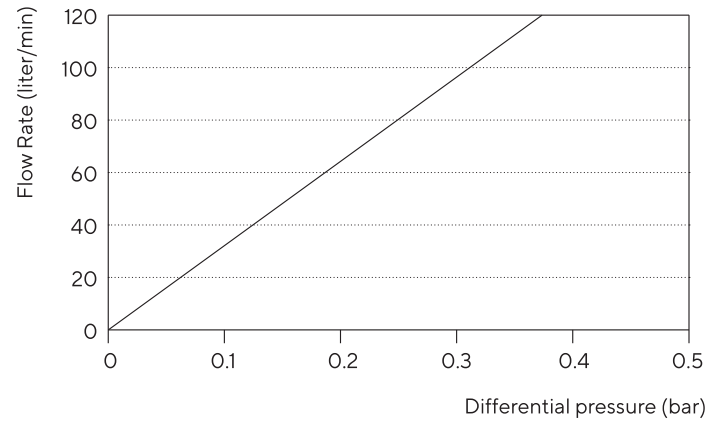
Ordering Information

Cartridges



Flow rate for water

Sartobev PS height 10"



Technical Specifications

| | |
|---|---|
| Recommended flow rate for 30" cartridge | Wine: 15 to 20 hl/h Water: 3m ³ /h |
| Sterilization | Hot water (85-95°C) or Steam (110°C), at least 100 cycles |
| Microbiological recovery rate | Saccharomcyes 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 |

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