SARTURIUS

Vinosart® PS

Cost efficiency at its best

Product Information

Vinosart® PS $0.45 \, \mu m$ | $0.65 \, \mu m$ are pleated membrane filter cartridges with a single layer asymmetrical polyether sulfone membrane inside. Vinosart® PS has been especially designed for the filtration of wine and sparkling wine. Vinosart® PS offers a very high flow rate and total throughput combined with an excellent chemical regeneration capability.



Application

Vinosart® PS is used for the final filtration prior bottling of wine and sparkling wine.

Flow rate | Total throughput

The polyether sulfone membrane has a special asymmetrical pore structure that ideally matches the requirements for the wine and champagne filtration in terms of flow rate and total throughput.

Microbiological retention

Both filters guarantee absolut retention of wine spoiling yeasts and provide highest microbiological safety for the bottling process of wine and champagne.

Chemical compatability

Vinosart® PS filter cartridges are extremely resistant to all conventional chemical regenerating agents and processes, right across the entire pH spectrum from 1 to 14.

Integrity testable

The integrity of Vinosart® PS filter cartridges can be tested daily using the Sartocheck® Mini. Integrity test parameters see technical data.

Quality

Every filter element of Vinosart® PS is integrity tested during production to guarantee highest level of security.

Technical Data

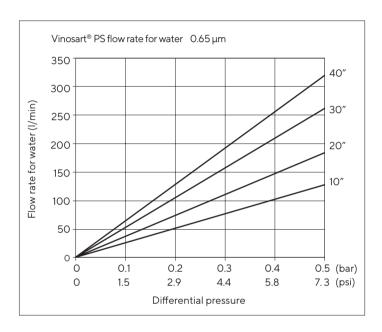
Filter Medium	Single-layer asymmetrical polyether sulfone membrane	
Protective non-woven drainage layers	Polypropylene	
Core outer support end caps adapters	Polypropylene	
O-Rings	Silicone	
Available Sizes Filtration Area	 Size 1 10" (250 mm) Size 2 20" (500 mm) Size 3 30" (750 mm) Size 4 40" (1.000 mm) 	0.78 m² 8.4 ft² 1.56 m² 16.8 ft² 2.34 m² 25.2 ft² 3.12 m² 33.6 ft²
Pore sizes	0.45 µm and 0.65 µm	
Maximum differential pressure	5 bar 72.5 psi at 20°C	
Direction of filtration	2 bar 29 psi at 80°C	
Recommend flow rate	Approx. 600–1000 l/h (30″)	

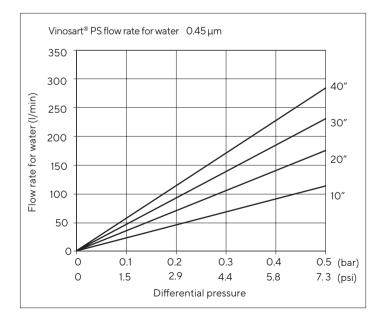
Pore size	0.45 µm	0.65 μm
Test pressure	1.5	1.0
Max. diffusion (10" element)	15 ml/mn	13 ml/mn
Test time	10 min	10 min

For more details please see the directions for use.

Available Adapters 25, 27, 28

Service life comparison of 0.45 μm | 0.65 μm membrane filters (Standard media) 100% 25% 50% 75% Vinosart® PS 0.45 0.65 Nylon 0.45 competitor A 0.65 PVDF 0.45 0.65 0.45 Nylon 0.65 competitor B PESU 0.45 competitor A 0.65





Ordering Information

Catridges	Pore Size [µm]	Pack Size [Pieces]
542**05A1	0.65	1
542**05A2	0.65	1
542**05A3	0.65	1
542**05A4	0.65	1
542**06A1	0.45	1
542**06A2	0.45	1
542**06A3	0.45	1
542**06A4	0.45	1

^{**:} Adaper Type

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