

# Vinosart® PS

Cost efficiency at its best

## Product Information

Vinosart® PS 0.45 µm | 0.65 µm 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 absolute retention of wine spoiling yeasts and provide highest microbiological safety for the bottling process of wine and champagne.

### Chemical compatibility

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	<ul style="list-style-type: none"><li>▪ Size 1 10" (250 mm)      0.78 m<sup>2</sup>   8.4 ft<sup>2</sup></li><li>▪ Size 2 20" (500 mm)      1.56 m<sup>2</sup>   16.8 ft<sup>2</sup></li><li>▪ Size 3 30" (750 mm)      2.34 m<sup>2</sup>   25.2 ft<sup>2</sup></li><li>▪ Size 4 40" (1.000 mm)    3.12 m<sup>2</sup>   33.6 ft<sup>2</sup></li></ul>	
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")	

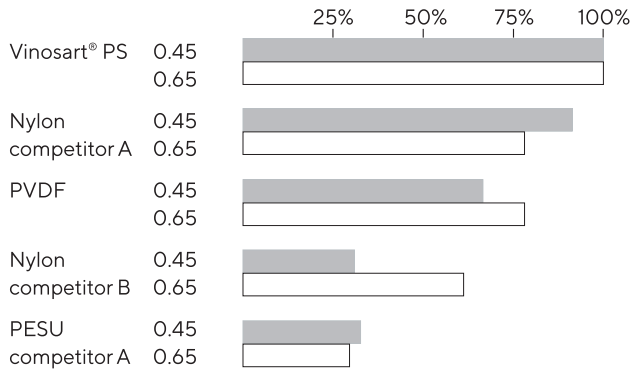
Pore size	0.45 µm	0.65 µm
Test pressure	1.5	1.0
Max. diffusion (10" element)	15 ml/mn	13 ml/mn
Test time	10 min	10 min

For more details please see the directions for use.

Available Adapters

25, 27, 28

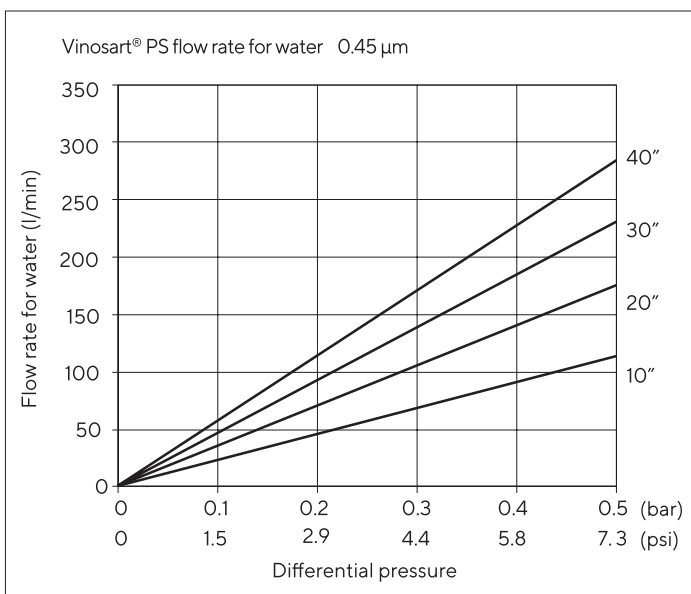
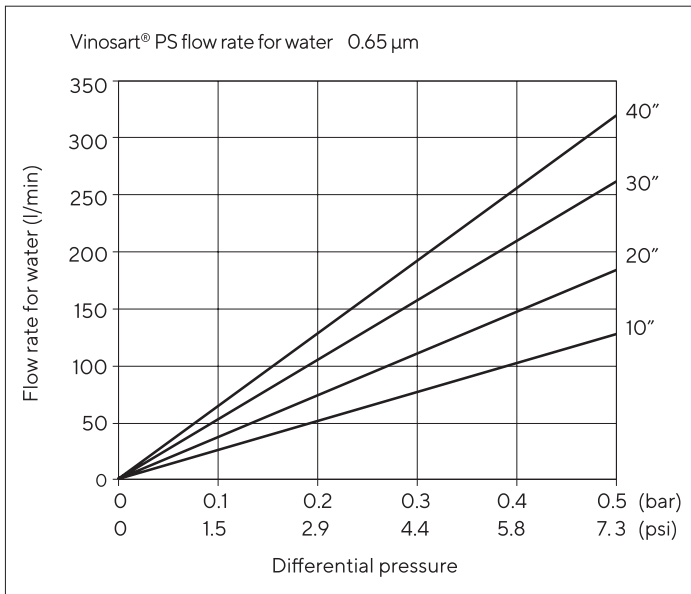
**Service life comparison of 0.45 µm | 0.65 µm membrane filters**  
(Standard media)



# Ordering Information

Cartridges	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

\*\* : Adapter 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](http://www.sartorius.com)