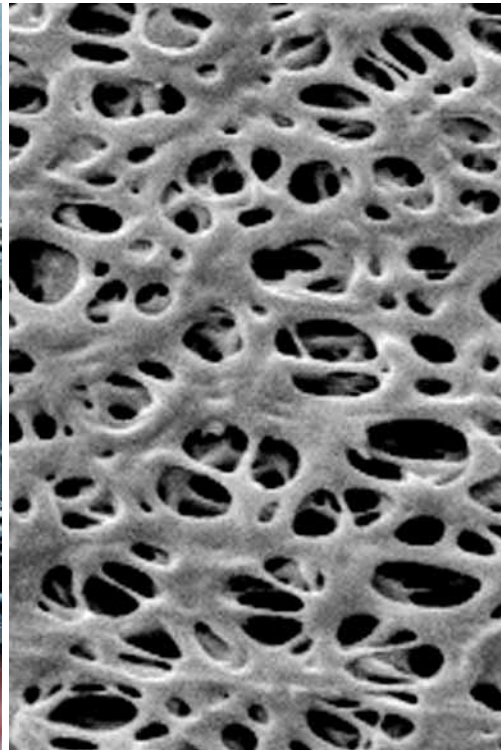




sartorius stedim
biotech

Vinosart® PS Cost efficiency at its best



turning science into solutions

Vinosart® PS at a glance

Optimal mechanical stability

Testable, maximum process
reliability for wine and sparkling
wine

Stable performance during
chemical regeneration

Long in-service lives

Validated, testable integrity

Optimized pore structure

High filtration performance



Vinosart® PS

Minimal process costs, maximum filtration performance

It's time for a new membrane technology – one that is geared towards the requirements of a highly dynamic market. The innovative Vinosart® PS filter cartridge sets new standards in technical performance and optimizes filtration costs.

Vinosart® PS has been specially designed for the filtration of wine and sparkling wines. All the practical tests conducted with this cartridge relate exclusively to wine and sparkling wine applications. This has resulted in a membrane filter that is specially optimized for the beverage filtration process.

Excellent flow rates.

These are facilitated by Vinosart® PS's high filtration performance, which is guaranteed by the new PESU (polyether sulfone) membrane filter material. This material is unique in terms of its structure and performance capability.

Broad compatibility, long in-service life.

Our new filter cartridge has another major advantage. Vinosart® PS is extremely resistant to all conventional chemical regenerating agents and processes, right across the entire pH spectrum from 1 to 14.

The cartridge's excellent chemical regeneration capability extends its service life considerably, thereby reducing the cost of filtration.

The polyether sulfone membrane provides the requisite mechanical and thermal stability. Vinosart® PS easily withstands daily 30-minute steam sterilization cycles at 121°C and thousands of pulsation cycles at a differential pressure of 2 bar (29 psi).



Summary.

Vinosart® PS is ideal for the bottling process of wine and sparkling wine, meets all customers' needs, and ensures smooth bottling operations at the customer's facilities.

Vinosart® PS's top quality provides excellent value combined with high filtration performance.

You'll find that it is well worth a trial filtration run.

Our specialists would be happy to give you a personal demonstration of the Vinosart® PS innovation along with the wide range of products available for your beverage production process.



Sartorius Stedim Biotech has the process down to a tee.

Sartorius Stedim Biotech not only develops filtration systems with optimum process reliability. It also provides equipment for the entire production process of beverages, ranging from clarification through to membrane filtration to fill level control.

Sartorius Stedim Biotech's range of products includes the following:

- Sartoflow systems for clarification or final cellar filtration
- Standard and Jumbo filter cartridges for particle-removing or polishing filtration
- Filter sheets and filter modules for depth filtration
- Membrane filter cartridges for sterile filtration of wine during bottling
- Equipment for microbiological quality testing and analytical laboratory balances
- Air and gas filter cartridges for tank venting
- Stainless steel filter cartridges for steam filtration

Technical specifications

Filter medium	Single-layer asymmetrical polyether sulfone membrane
Protective non-woven drainage layers	Polypropylene
Core outer support end caps adapters	Polypropylene
O-rings	Silicone
Filtration area	0.78 m ² / 10" cartridge
Pore sizes	0.45 µm and 0.65 µm
Heights	10" (250 mm), 20" (500 mm), 30" (750 mm), 40" (1000 mm)
Maximum differential pressure	5 bar (73 psi) at 20°C
in the direction of filtration	2 bar (29 psi) at 80°C
opposite the direction of filtration	2 bar (29 psi) at 20°C
Sterilization	See Directions for Use of the Filter Cartridges and Data Sheet
Integrity test parameters	See Directions for Use of the Filter Cartridges and Data Sheet
Recommend flow rate	Approx. 600–1000 l/h (30")

Ordering information

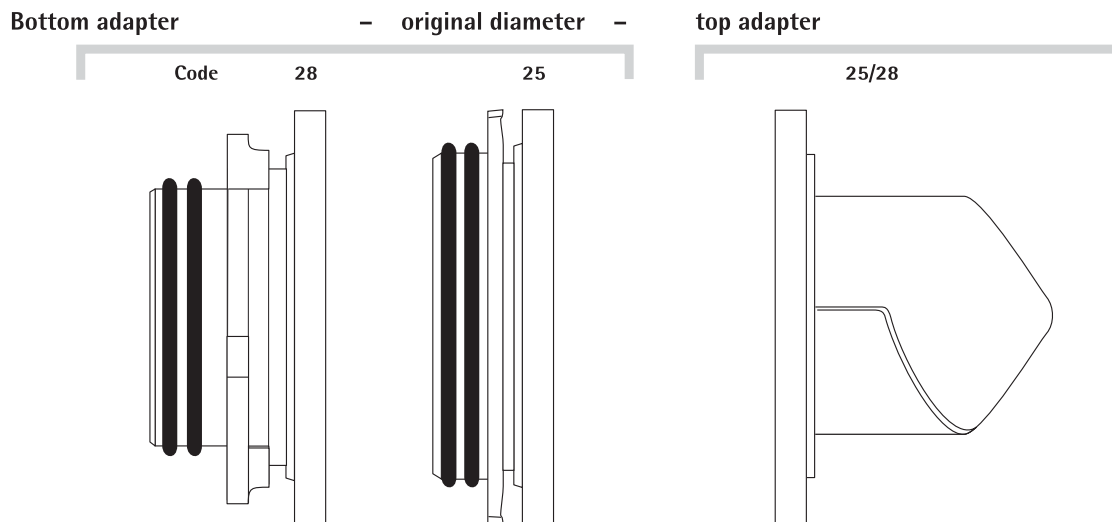
542	28	05	A	3	Vinosart® PS
					Adapter
					Pore size
					Single-layer membrane
					Height

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

Pore size

- 05 0.65 µm
- 06 0.45 µm



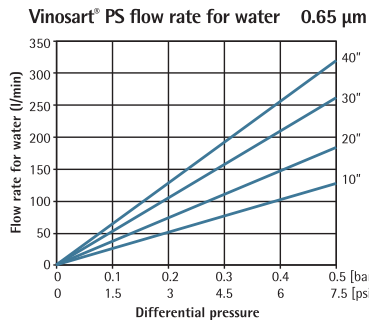
Vinosart® PS



High microbiological stability, optimal performance

Reliable retention of all wine-spoilage microorganisms at minimal filtration costs is the decisive criterion for membrane filter cartridges used in wine filtration. Vinosart® PS ideally fulfills this requirement, ensuring the highest microbiological stability of your beverages. In fact, it meets all the criteria that a membrane filter cartridge has to cover for beverage filtration:

- Integrity test capability
- Validated microbial retention
- Defined pore sizes of 0.65 µm and 0.45 µm
- Thermal sterilization
- Chemical regeneration capability



Integrity testable.

The integrity of Vinosart® PS filter cartridges can be tested daily using the Sartocheck® Mini integrity tester, for example.

The integrity test serves two important purposes. First, it determines whether the filter cartridges function properly before production begins. Second, it allows a batch to be checked immediately at the end of a day of bottling operations.

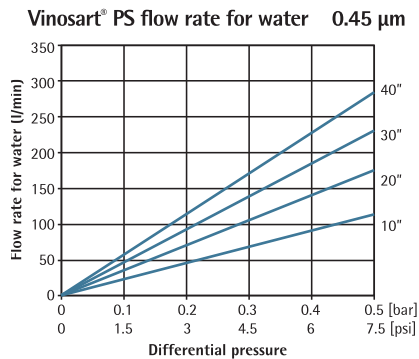
Superior membrane geometry.

The membrane inside Vinosart® PS was especially developed for the two specific F&B applications "wine- and sparkling wine filtration".

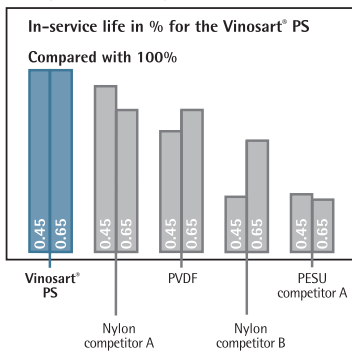
Our manufacturing process of the membrane assures an outstanding performance of Vinosart® PS since the membrane has got an asymmetrical pore geometry that ideally matches the requirements for high flow rates and total throughputs thus fractionate retention.

Summary.

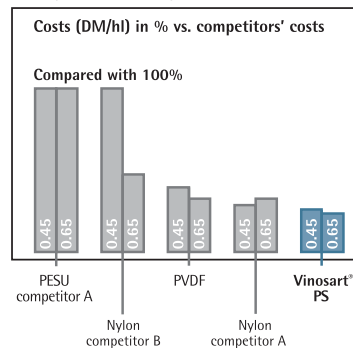
Considering the innovative benefits described above, the Vinosart® PS filter cartridge simply does more for your bottom line – for cost efficiency at its best.



In-service life comparison 0.45 µm and 0.65 µm



Comparison of filtration costs 0.45 µm und 0.65 µm



Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

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

Phone +49.551.308.0
Fax +49.551.308.3289

www.sartorius-stedim.com

Sartorius Stedim Systems GmbH
Schwarzenberger Weg 73-79
34212 Melsungen

Phone +49.5661.71.3400
Fax +49.5661.71.3702

www.sartorius-stedim.com

France

Sartorius Stedim Biotech S.A.
ZI Les Paluds
Avenue de Jouques – BP 1051
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.845619

Sartorius Stedim France SAS
ZI Les Paluds
Avenue de Jouques – CS 71058
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH
Franzosengraben 12
A-1030 Vienna

Phone +43.1.7965763.18
Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde

Phone +32.2.756.06.80
Fax +32.2.756.06.81

Denmark

Sartorius Stedim Nordic A/S
Hoerskaetten 6D, 1.
DK-2630 Taastrup

Phone +45.7023.4400
Fax +45.4630.4030

Hungary

Sartorius Stedim Hungária Kft
Kagyló u. 5
2092 Budakeszi

Phone +36.23.457.227
Fax +36.23.457.147

Italy

Sartorius Stedim Italy S.p.A.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41
Fax +39.055.63.40.526

Netherlands

Sartorius Stedim Netherlands B.V.
Edisonbaan 24
3439 MN Nieuwegein

Phone +31.30.6025080
Fax +31.30.6025099

Poland

Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.647.38.40
Fax +48.61.879.25.04

Spain

Sartorius Stedim Spain SA
C/Isabel Colbrand 10,
Oficina 70
Polígono Industrial de Fuencarral
28050 Madrid

Phone +34.90.2110935
Fax +34.91.3589623

Switzerland

Sartorius Stedim Switzerland AG
Ringstr. 24 a
8317 Tagelswangen

Phone +41.52.354.36.36
Fax +41.52.354.36.46

U.K.

Sartorius Stedim UK Limited
Longmead Business Park
Blenheim Road, Epsom
Surrey KT19 9 QQ

Phone +44.1372.737159
Fax +44.1372.726171

America

USA

Sartorius Stedim North America Inc.
5 Orville Drive
Bohemia, NY 11716

Toll-Free +1.800.368.7178
Fax +1.631.254.4253

Sartorius Stedim SUS Inc.
1910 Mark Court
Concord, CA 94520

Phone +1.925.689.6650
Toll Free +1.800.914.6644
Fax +1.925.689.6988

Sartorius Stedim Systems Inc.
201 South Ingram Mill Road
Springfield, MO 65802

Phone +1.417.873.9636
Fax +1.417.873.9275

Argentina

Sartorius Argentina S.A.
Int. A. Avalos 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Av. Dom Pedro I, 241
Bairro Vila Pires
Santo André
São Paulo
Cep 09110-001

Phone +55.11.4451.6226
Fax +55.11.4451.4369

Mexico

Sartorius de México S.A. de C.V.
Circuito Circunvalación Poniente No. 149
Ciudad Satélite
53100 Naucalpan, Estado de México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius Stedim Beijing
Representative Office
No. 33, Yu'an Road,
Airport Industrial Zone B, Shunyi District
Beijing 101300

Phone +86.10.80426516
Fax +86.10.80426580

Sartorius Stedim Shanghai
Representative Office
Room 618, Tower 1, German Centre,
Shanghai, PRC., 201203

Phone +86.21.28986393
Fax +86.21.28986392.11

Sartorius Stedim Guangzhou Office
Room 704, Broadway Plaza,
No. 233-234 Dong Feng West Road
Guangzhou 510180

Phone +86.20.8351.7921
Fax +86.20.8351.7931

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, Jakkasandra
Kunigal Road, Nelamangala Tq
Bangalore – 562 123

Phone +91.80.4350.5361
Fax +91.80.4350.5253

Japan

Sartorius Stedim Japan K.K.
KY Building, 8-11
Kita Shinagawa 1-chome
Shinagawa-ku
Tokyo 140-0001

Phone +81.3.3740.5407
Fax +81.3.3740.5406

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Stedim Singapore Pte. Ltd.
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park 2
Singapore 117528

Phone +65.6872.3966
Fax +65.6778.2494