



Filtration of wine samples and measure of the wine clogging index

Product Information

When small volumes of wine need to be filtered for their presentation and tasting, MidiCaps® capsules simplify your filtration in complete safety and while respecting your wines.



From a few liters to hundreds of liters!

The filter materials of the MidiCaps® are the same as those used on the bottling lines but with surface areas 9 times smaller.

Applications

According to the filter materials chosen and their pore size:

- Clarification
- Particle retention
- Bioburden reduction
- Sterile filtration

MidiCaps® filtration capsules simplify your filtration in complete safety and while respecting your wines.

MidiCaps®: ready-to-use filtration capsules

Traceability

Each capsule is identified: type of filter, reference, pore size and batch number.

Engraving on the capsules of the conditions for use and direction of filtration.

Quality

Purge and sanitary type vent to optimize the use of the capsules. MidiCaps® have a validation guide according to FDA recommendations. Quality assurance certificates in each box.

Ease of use

Available with different connections: triclamp or flexible hose or hose barb for flexible tubing.

Sterilized by autoclaving or hot water. Contact us for advice on the regeneration and storage of capsules.

Filtration of wine samples and measurement of the clogging index of wine



5 liter vessel

Sartorius provides a stainless steel vessel for the filtration of small volumes of wine.

Application

- Filtration of wine samples with filtration capsules
- Measurement of the clogging index and Vmax

Practical and versatile

Ease of use

- Stainless steel system mounted on an height-adjustable gallow
- 0 to 6 bar pressure gauge.
- Two valves at the highest point for the intake of air and release into the atmosphere after filtration.
- A valve at the lowest point for the outlet of wine.



2 in 1!

- Equipped as standard with a support disc of 25 mm in diameter to measure the clogging index and Vmax.
- Triclamp connector for mounting a capsule filter and filtering wine samples.

Technical Specifications

Dimensions (W × D × H)	450 × 460 × 1300 mm (17 3/4 × 18 1/8 × 51 3/16 in)
Weight (approx)	17 kg (37.48 lbs)

Ordering Information

Vessel

Description	Reference
Complete 5 liter vessel	FR-INDICE5L349

Membrane for the measure of the clogging index and Vmax

Description	Reference
0.65 µm cellulose acetate membrane 25 mm in diameter (box of 100)	111 05 025N

MidiCaps® capsules for filtration of samples

Filter material	Applications	Pore size	Inlet connection	Inlet connection	References		
Positively charged fiberglass	Brightness of the wines	0.65 µm	Clamp	Olive	555 53 05 P9 SOA		
			Olive	Olive	555 53 05 P9 OOA		
		1.2 µm	Clamp	Olive	555 53 03 P9 SOA		
			Olive	Olive	555 53 03 P9 OOA		
Polypropylene	Clarification of the wines and particle retention	0.65 µm	Clamp	Olive	505 53 05 P9 SOA		
			Olive	Olive	505 53 05 P9 OOA		
		1.2 µm	Clamp	Olive	505 53 03 P9 SOA		
			Olive	Olive	505 53 03 P9 OOA		
		3 µm	Clamp	Olive	505 53 02 P9 SOA		
			Olive	Olive	505 53 02 P9 OOA		
			Clamp	Clamp	505 53 03 P9 SSA		
		5 µm	Olive	Olive	505 53 42 P9 OOA		
		Cellulose acetate	Sterilizing filtration	0.65 µm	Clamp	Clamp	5625306A9--00--A
				0.45 µm	Clamp	Olive	5625305G9--S0--A

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

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5-7
34302 Guxhagen
Phone +49.5665.407.0

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques - CS 91051
13781 Aubagne Cedex
Phone +33.442.845600

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques - CS 71058
13781 Aubagne Cedex
Phone +33.442.845600

Austria

Sartorius Stedim Austria GmbH
Modecenterstrasse 22
1030 Vienna
Phone +43.1.7965763.18

Belgium

Sartorius Stedim Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.80

Hungary

Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi
Phone +36.23.457.227

Italy

Sartorius Stedim Italy S.r.l.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)
Phone +39.055.63.40.41

Netherlands

Sartorius Stedim Netherlands B.V.
Phone +31.30.60.25.080
filtratie.nederland@sartorius-stedim.com

Poland

Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.647.38.40

Russian Federation

LLC "Sartorius Stedim RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg
Phone +7.812.327.53.27

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone +34.913.586.098

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen
Phone +41.52.354.36.36

U.K.

Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ
Phone +44.1372.737159

Ukraine

LLC "Sartorius Stedim RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918

Americas

USA

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716
Toll-Free +1.800.368.7178

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
leadsmex@sartorius.com

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800

China

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China
Phone +86.21.6878.2300

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Beijing Branch Office
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300
Phone +86.10.8042.6501

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Guangzhou Branch Office
Room 1105
Xing Guang Ying Jing Building
No. 119, Shui Yin Road
Yue Xiu District, Guangzhou 510075
Phone +86.20.3836.4193

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250

Japan

Sartorius Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan
Phone +81.3.4331.4300

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia
Phone +60.3.8996.0622

Singapore

Sartorius Stedim Singapore Pte. Ltd.
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684
Phone +65.6872.3966

South Korea

Sartorius Korea Biotech Co., Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400
Phone +82.31.622.5700



▶ www.sartorius-stedim.com