

# 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 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 scaffold
- 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 ¾ × 18 ⅞ × 51 ⅜ in)
Weight (approx)	117 kg (37.48 lbs)

## Ordering Information

### Vessel

Description	References
Complete 5 liter vessel	FR-INDICE5L349

### Membrane for the Measure of the Clogging Index and Vmax

Description	References
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	Hose barb	5555305P9--SO--A
			Hose barb	Hose barb	5555305P9--OO--A
		1,2 µm	Clamp	Hose barb	5555303P9--SO--A
			Hose barb	Hose barb	5555303P9--OO--A
Polypropylene	Clarification of the wines and particle retention	0,65 µm	Clamp	Hose barb	5055305P9--SO--A
			Hose barb	Hose barb	5055305P9--OO--A
		3 µm	Clamp	Hose barb	5055302P9--SO--A
			Hose barb	Hose barb	5055302P9--OO--A
		5 µm	Clamp	Clamp	5055302P9--SS--A
			Hose barb	Hose barb	5055342P9--OO--A
Cellulose acetate	Sterilizing filtration	0,65 µm	Clamp	Clamp	5625305A9--SS--A
		0,45 µm	Clamp	Hose barb	5625306G9--SO--A

**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

**France**

Sartorius Stedim France SAS  
Z.I. des Paluds  
Avenue de Jouques - CS 71058  
13781 Aubagne  
Téléphone: +33 (0)4 42 845 600

 For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)