



Sartofluor® MidiCaps® and Capsules

Membrane Filter



Description

Sartofluor® MidiCaps® and Capsules 0.2 µm rated are self contained, ready to use, sterile filter units for sterilizing grade filtration in the pharma | biotech industry. Their unique hydrophobic PTFE membrane is ideally suited for particle removal and sterilizing grade filtration of gases and for filtration of highly aggressive liquids like solvents, acids and bases.

Applications

Typical applications include sterile venting of:

- Fermenters
- Vessels
- Glass Bottles

The hydrophobic PTFE membrane is also suitable for filtration of aggressive liquids like:

- Acids | Bases
- Solvents

Easy to use

Sartofluor® MidiCaps® and Capsules are delivered as individually packed sterile units. On site, pre-use sterilization can be eliminated.

Flexibility

Sartofluor® MidiCaps® and Capsules are available with various filtration areas from 150 cm² | 0.16 ft² up to 0,45 m² | 4.8 ft² for easy adoption to any filtration process independent from the batch size.

Performance

The unique hydrophobic single layer PTFE membrane provides outstanding flow rates for gases and liquids at low differential pressure assuring most economic system design.

Cost Saving

The use of the disposable capsule design concept avoids investments into stainless steel filter housings and eliminates additional costs for cleaning of housings and cleaning validation.

Microbiological Retention

Sartofluor® MidiCaps® and Capsules 0.2 µm rated are fully validated as sterilizing grade filters according to HIMA and ASTM F-838-05 guidelines.

Quality Control

Each individual element is tested for integrity by B.-P. and Diffusion-Test prior to be released assuring absolute reliability.

Documentation

Sartofluor® MidiCaps® and Capsules are designed, developed and manufactured in accordance with a ISO 9001 certified Quality Management System. A Validation Guide is available for compliance with regulatory requirements.

Specifications

Materials

Membrane:	PTFE
Support fleece:	Polypropylene
Core:	Polypropylene
End caps:	Polypropylene
Capsule Housing:	Polypropylene
O-Rings:	EPDM
Filling Bell:	Polycarbonate

Pore Sizes

0.1 µm	(only MidiCaps®)
0.2 µm	(MidiCaps® & Capsules)
0.45 µm	(only MidiCaps®)

Available Sizes | Filtration Area

MidiCaps®

Size 7	0.05 m ² 0.54 ft ²
Size 8	0.1 m ² 1.1 ft ²
Size 9	0.2 m ² 2.2 ft ²
Size 0	0.45 m ² 4.8 ft ²

Capsules

Size 4	0.015 m ² 0.16 ft ²
Size 5	0.03 m ² 0.32 ft ²

Available Connectors MidiCaps®

SS, SO, OO, FF, FO, FH, HH (only for size 7)

Available Connectors Capsules

SS, SO, OO (Size 4)
OO (Size 5)

S:	1½" Tri-Clamp (Sanitary)
O:	½" Single stepped hose barb
F:	¾" Tri-Clamp (Sanitary)
H:	¼" Multiple stepped hose barb (with filling bell at the outlet)

Operating Parameters

Max. allowable Differential pressure:	5 bar 72.5 psi at 20°C (MidiCaps®) 4 bar 58 psi at 20°C (Capsules)
Max. allowable back Pressure:	2 bar 29 psi at 20°C

Specifications

Extractables

Sartofluor® MidiCaps® and Capsules meet, or exceed the requirements for WFI quality standards set by the current USP.

Regulatory Compliance

Individually integrity tested

Integrity test correlated to HIMA | ASTM F 838-05 Bacteria Challenge Test

Non pyrogenic according to USP Bacterial Endotoxins

Pass USP Plastic Class VI Test

Non fiber releasing according to 21 CFR

Sterilization

Autoclaving:

134°C, 2 bar, 30 min

No In-Line Steam Sterilization

Sterilization Cycles

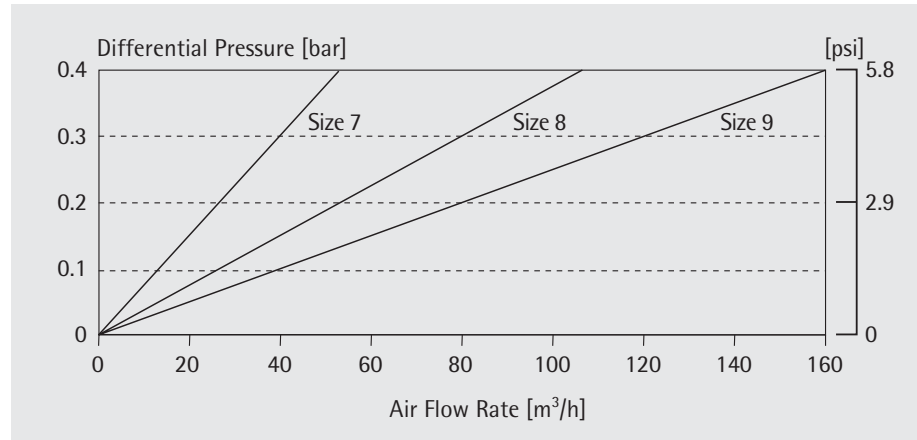
Autoclaving: Min. 25

Technical References

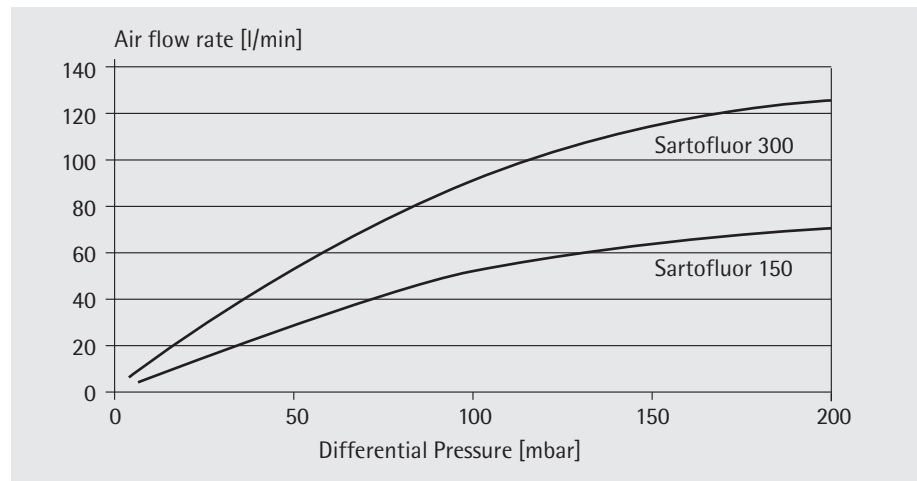
Validation Guide:

SPK5711-e (MidiCaps®)

Air Flow Rates for Sartofluor® MidiCaps®, 0.2 µm rated with SS-connector



Air Flow Rates for Sartofluor® Capsules, 0.2 µm rated with SS-connector



Order Information

Order Code	Pore size [µm]	Pack size [pieces]	Test Pressure [bar psi]	Max. Diffusion [ml min]	Min. Bubble Point [bar psi]
MidiCaps®:					
5185307T7--**--A	0.2	4	0.7 10.2	2	1.0 14.5
5185307T8--**--A	0.2	4	0.7 10.2	3	1.0 14.5
5185307T9--**--A	0.2	4	0.7 10.2	4	1.0 14.5
5185307T0--**--V	0.2	2	0.7 10.2	8	1.0 14.5
5185306T7--**--A	0.45	4	0.4 5.8	1	0.6 8.7
5185306T8--**--A	0.45	4	0.4 5.8	2	0.6 8.7
5185306T9--**--A	0.45	4	0.4 5.8	3	0.6 8.7
5185358T7--**--A	0.1	4	-	-	1.5 22
5185358T8--**--A	0.1	4	-	-	1.5 22
5185358T9--**--A	0.1	4	0.9 13.1	5	1.5 22
Capsules:					
5181307T4--**--B	0.2	4	-	-	1.0 14.5
5181307T5--00--D	0.2	10	-	-	1.0 14.5

** : Connector Styles

Sartorius Stedim Biotech GmbH

August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone +49.551.308.0

Fax +49.551.308.3289

www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178

UK +44.1372.737159

France +33.442.845600

Italy +39.055.63.40.41

Spain +34.90.2110935

Japan +81.3.3740.5407

Specifications subject to change without notice. Printed and copyrighted by Sartorius Stedim Biotech GmbH. | W
Publication No.: SPK2067-e11067

Order No.: 85030-522-62

Ver. 06 | 2011