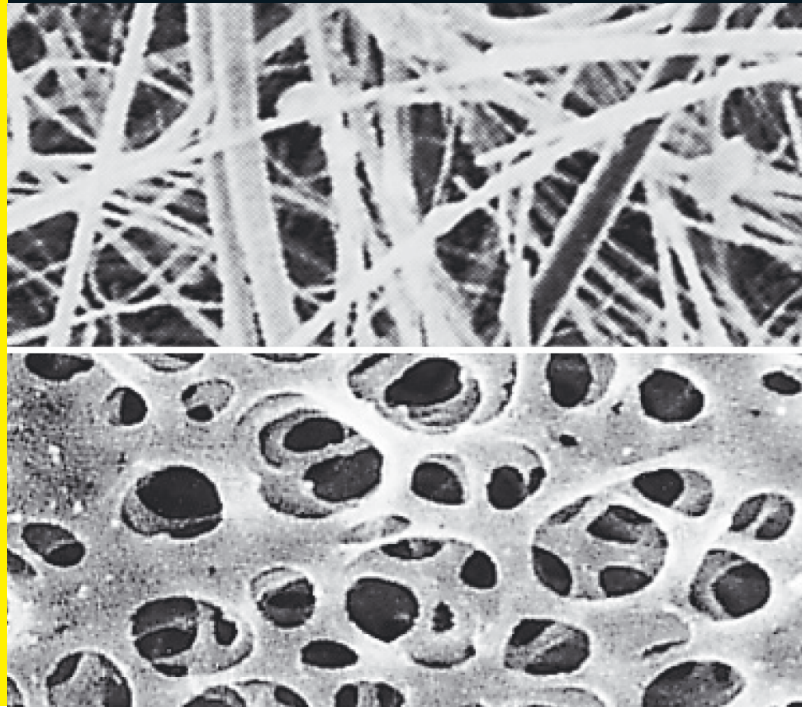


Sartoclean® GF

Adsorptive Membrane Filter
for Colloid and Bioburden
Reduction



Product Information

Sartoclean® GF filter elements combine excellent retention performance by membrane filtration with the adsorptive power of glass fiber fleeces. Therefore, Sartoclean® GF are ideally suited for removal of colloids and lipids as well as defined particle and bioburden reduction for a broad range of biopharmaceutical process fluids.

Process Contamination Control

The removal of colloidal contaminants and lipids by adsorption ensures effective downstream processing and protection of downstream equipment. In addition, bioburden reduction by membrane filtration prevents the release of pyrogens during the process and, thus, strengthens contamination control.

Performance

The combination of adsorptive glass fiber fleeces with cellulose acetate membrane filters leads to optimal total throughput performance and guarantees an economic filtration system design.

Applications

Clarification and filtration of

- Fermentation broths
- Serum
- Plasma
- Cell culture media
- Colloid and lipid containing solutions

Technical Data

Available Sizes	Filtration Area
Cartridges T-Style Maxicaps® Maxicaps®	
Size 1 (10")	0.47 m ² 5.1 ft ²
Size 2 (20")	0.94 m ² 10.2 ft ²
Size 3 (30")	1.41 m ² 15.3 ft ²
Mini Cartridges Midicaps® (Mini Cartridges only available in size 9)	
Size 7	0.05 m ² 0.54 ft ²
Size 8	0.08 m ² 0.85 ft ²
Size 9	0.16 m ² 1.7 ft ²
Size 0	0.32 m ² 3.4 ft ²

Materials

Prefilter Membrane

Cellulose Acetate

Endfilter Membrane

Cellulose Acetate

Filter Active Fleece

Glass Fiber

Support Fleece

Polypropylene

Core

Polypropylene

End Caps

Polypropylene

Capsule Housing

Polypropylene

O-Rings

Silicone

Pore Size Combination

3.0 µm + 0.8 µm (nom.)

0.8 µm + 0.65 µm (nom.)

Max. Operating Pressure

Cartridges | T-Style Maxicaps® | Maxicaps® | Mini Cartridges | Midicaps®

5 bar | 72.5 psi at 20 °C

Max. Allowable Differential Pressure

Cartridges | Mini Cartridges

5 bar | 72.5 psi at 20 °C

2 bar | 29 psi at 80 °C

T-Style Maxicaps® | Maxicaps® | Midicaps®

5 bar | 72.5 psi at 20 °C

3 bar | 43.5 psi at 50 °C

Max. Allowable Back Pressure

Cartridges | T-Style Maxicaps® | Maxicaps® | Mini Cartridges | Midicaps®

2 bar | 29 psi at 20 °C

Regulatory Compliance

- Designed, developed and manufactured in accordance with an ISO 9001 certified Quality Management System
- Meet or exceed the requirements for WFI quality standards set by the current USP
- Non-pyrogenic according to USP Bacterial Endotoxins
- Non-fiber releasing according to 21 CFR 210.3 (b) (6) and 211.72
- Pass USP Plastic Class VI Test

Sterilization

Cartridges | Mini Cartridges

- In-Line Steam Sterilization
Max. 134 °C, 0.5 bar | 7.25 psi, 20 min
Min. 25 Sterilization Cycles
- or
- Autoclaving
Max. 134 °C, 2 bar | 29 psi, 30 min
Min. 25 Sterilization Cycles

Maxicaps® | Midicaps®

- Autoclaving
Max. 134 °C, 2 bar | 29 psi, 30 min
Min. 25 Sterilization Cycles

T-Style Maxicaps®

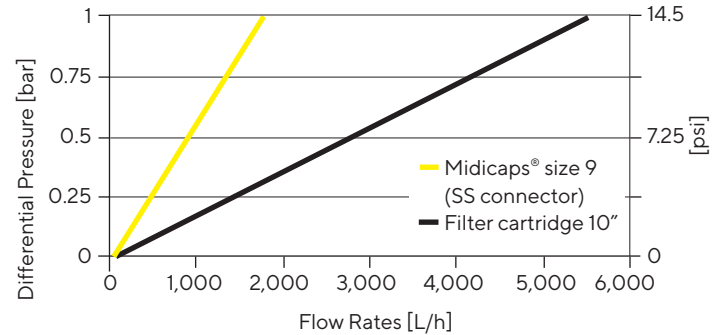
- Autoclaving
Max. 134 °C, 2 bar | 29 psi, 30 min
Min. 5 Sterilization Cycles

Technical References

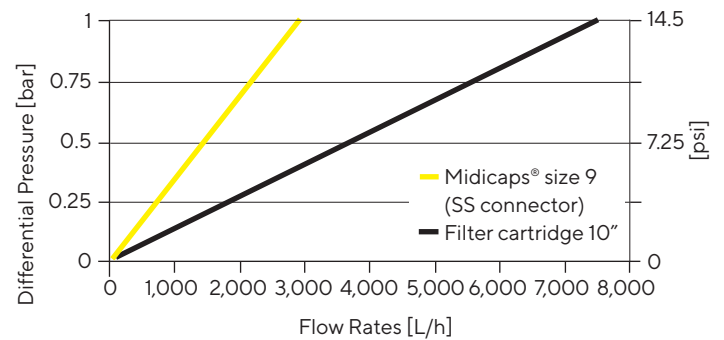
Validation Guide (SPK5788-e)
Extractables Guide (DIR 2833485)

Performance

Water Flow Rates* 0.8 µm + 0.65 µm



Water Flow Rates* 3.0 µm + 0.8 µm



* Standardized at 20 °C.



Ordering Information



Cartridge



Adapter
25: 2 flange bayonet adapter with 226 double o-ring
27: Bayonet adapter with 222 double o-ring
28: 3 flange bayonet adapter with 222 double o-ring

Pore Size
04 E: Final filter 0.8 μm , prefilter 3.0 μm
05 G: Final filter 0.65 μm , prefilter 0.8 μm

Filter Size
1: 0.47 m² | 5.1 ft² (10")
2: 0.94 m² | 10.2 ft² (20")
3: 1.41 m² | 15.3 ft² (30")



Mini Cartridge



Pore Size
04 E: Final filter 0.8 μm , prefilter 3.0 μm
05 G: Final filter 0.65 μm , prefilter 0.8 μm

Filter Size
9: 0.16 m² | 1.7 ft²

Packing Size
B: Box of 5



Maxicaps®



Pore Size
04 E: Final filter 0.8 μm , prefilter 3.0 μm
05 G: Final filter 0.65 μm , prefilter 0.8 μm

Filter Size
1: 0.47 m² | 5.1 ft²
2: 0.94 m² | 10.2 ft²
3: 1.41 m² | 15.3 ft²

Connector Inlet
S: 1½" tri-clamp 50 mm
O: ½" single stepped hose barb

Connector Outlet
S: 1½" tri-clamp 50 mm
O: ½" single stepped hose barb



T-Style Maxicaps®



Pore Size
04 E: Final filter 0.8 μm , prefilter 3.0 μm
05 G: Final filter 0.65 μm , prefilter 0.8 μm

Filter Size
1: 0.47 m² | 5.1 ft²
2: 0.94 m² | 10.2 ft²
3: 1.41 m² | 15.3 ft²

Connector Inlet
S: 1½" tri-clamp 50 mm
O: ½" single stepped hose barb

Connector Outlet
S: 1½" tri-clamp 50 mm
O: ½" single stepped hose barb



Midicaps®

560	53	■	■	■	--	■	■	--	■
Pore Size 04 E: Final filter 0.8 µm, prefilter 3.0 µm 05 G: Final filter 0.65 µm, prefilter 0.8 µm		Filter Size 7: 0.05 m ² 0.54 ft ² 8: 0.08 m ² 0.85 ft ² 9: 0.16 m ² 1.7 ft ² 0: 0.32 m ² 3.4 ft ²			Connector Inlet S: 1½" tri-clamp 50 mm O: ½" single stepped hose barb F: ¾" tri-clamp 25 mm		Connector Outlet S: 1½" tri-clamp 50 mm O: ½" single stepped hose barb F: ¾" tri-clamp 25 mm		Packing Size A: Box of 4 V: Box of 2


Check the availability of your desired configuration with your Sartorius sales representative or online at www.sartorius.com

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