SVISCISAS

Product Datasheet

Sartoclean® GF

Colloid & Bioburden Reduction Maxicaps[®] & Midicaps[®]



Product Information

Sartoclean[®] GF Maxicaps[®] and Midicaps[®] are selfcontained, ready to use filter units for a broad range of prefiltration applications in the biopharmaceutical industry. Sartoclean[®] GF Maxicaps[®] and Midicaps[®] combine absolute retention performance by membrane filtration with high adsorptive power by glass fiber fleeces. Therefore the filters are ideally suited for removal of colloids, lipids, defined particle retention and bioburden reduction.

Applications

Sartoclean[®] GF Maxicaps[®] and Midicaps[®] are widely used for prefiltration in biotech manufacturing processes to protect subsequent downstream processing equipments. Sartoclean[®] GF Maxicaps[®] and Midicaps[®] are ideally suited for the biodurden reduction as well as effective colloid and lipid removal from:

- Fermentation Broth
- Serum
- Plasma
- Cell Culture Media
- Colloid and Lipid containing solutions

Performance

Due to the combination of high adsorptive glass fiber fleeces with membrane filters Sartoclean® GF Maxicaps® and Midicaps® assure optimal total throughput performance. Therefore the Midicaps® & Maxicaps® allows a more economical filtration system design.

Process Safety

The removal of collodial contaminations and lipids by adsorption allows an effective downstream processing and bioburden reduction by membrane filtration prevents the formation of pyrogenes during the process – resulting in an increased process safety, especially for biotech derived fluids.

Flexibility

Sartoclean® GF Maxicaps® and Midicaps® are available with various filtration areas from 500 cm² | 0.54 ft² up to 1.8 m² | 19.4 ft² for easy adoption to any filtration process independent from the batch size

Technical Data

Specifications

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
Filling Bell	Polycarbonate

Pore Size Combination

0.8 + 0.65 μm, 3.0 + 0.8 μm

Available Sizes | Filtration Area

Maxicaps®			
Size 1	10″	0.6 m ² 6.5 ft ²	
Size 2	20″	1.2 m ² 12.9 ft ²	
Size 3	30″	1.8 m² 19.4 ft²	

Scalability

Consistent and predictable scale-up and down trials can reliably be performed as all Sartoclean[®] GF Maxicaps[®] and Midicaps[®] are produced with the same type of membrane and identical materials of construction.

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.

Documentation

Sartoclean® GF Maxicaps® and Midicaps® 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.

Midicaps®		
0.05 m² 0.54 ft²		
0.1 m ² 1.1 ft ²		
0.2 m ² 2.2 ft ²		
0.45 m ² 4.8 ft ²		

Available Connectors Maxicaps®, Midicaps®

- S: Tri-Clamp 50 mm (1½")
- O: $\frac{1}{2}$ single stepped hose barb
- F: Tri-Clamp 25 mm (¾")
- H: ¼" multiple stepped hose barb

(only Midicaps[®] size 7 with filling bell)

Operating Parameters	
Max. Allowable Differential Presure	5 bar 72.5 psi at 20°C (Maxicaps®, Midicaps®) 2 bar 29 psi at 80°C (Midicaps®) 3 bar 43.5 psi at 50°C (Maxicaps®)
Max. Allowable Back Pressure	2 bar 29 psi at 20°C

Extractables

Sartoclean[®] GF Maxicaps[®] and Midicaps[®] meet or exceed the requirements for WFI quality standards set by the current USP.

Regulatory Compliance

Non pyrogenic according to USP Bacterial Endotoxins

Pass USP Plastic Class VI Test

 $(0.65 \,\mu\text{m})$ with SS Connector

Non fiber releasing according to 21 CFR

Water Flow Rates for Sartoclean® GF Midicaps®

Sterilization

Autoclaving 134°C, 2 bar|29 psi, 30 min

No In-Line Steam Sterilization

Sterilization Cycles

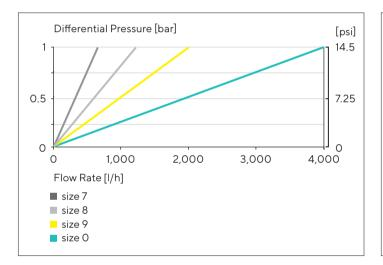
Autoclaving: Min. 25

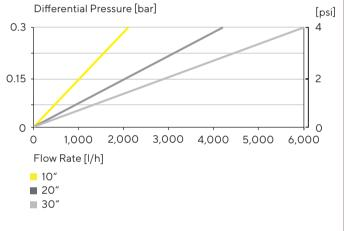
Technical References

Validation Guide

SPK5788-e

Water Flow Rates for Sartoclean $^{\circ}$ GF Maxicaps $^{\circ}$ (0.65 $\mu m)$ with SS Connector





Ordering Information

Order Code	Pore Size [µm]	Pack Size [Pieces]
Maxicaps®		
5607304E1**	0.8	1
5607304E2**	0.8	1
5607304E3**	0.8	1
5607305G1**	0.65	1
5607305G2**	0.65	1
5607305G3**	0.65	1

Midicaps®		
5605304E7**A	0.8	4
5605304E8**A	0.8	4
5605304E9**A	0.8	4
5605304E0**V	0.8	2
5605305G7**A	0.65	4
5605305G8**A	0.65	4
5605305G9**A	0.65	4
5605305G0**V	0.65	2

** Connector Styles

(first code letter represents the inlet, the second the outlet)

Germany

USA

Sartorius Stedim Biotech GmbH August-Spindler-Strasse 11 37079 Goettingen Phone +49 551 308 0 Sartorius Stedim North America Inc. 565 Johnson Avenue Bohemia, NY 11716 Toll-Free +1 800 368 7178

For further contacts, visit www.sartorius.com

Specifications subject to change without notice. Copyright Sartorius Stedim Biotech GmbH. DIR No.: 2648627-000-00 Order No.: 85030-531-66 Status: 10 | 2020