

Sartopure[®] PP3

Test Filter Elements

The Sartopure[®] PP3 prefilter platform is the result of decades of experience in prefiltration in pharmaceutical and biopharmaceutical processes. This development has led to a polypropylene fleece-based prefilter offering both, highest total throughput and protective abilities. Its outstanding filtration ability results in a significant reduction of the required filtration area, essentially reducing filter consumption and the overall cost of prefiltration. Sartopure[®] PP3 is the ideal choice for particle retention and protection of downstream equipment for all product contact applications.



Product Information

- High performance polypropylene fleece prefilters for particle reduction
- Advanced fleece materials with retention ratings from 0.45 – 100 µm
- Broad chemical compatibility with low extractable profile

Robust Processes

The Sartopure® PP3 series provides exceptional robustness to prefiltration applications based on the unique retention performance of its fleece material. The fleece retains particles with high efficiency even under varying process conditions, ensuring secure and reliable operation.

High Product Yield

The all-polypropylene design of Sartopure® PP3 provides low unspecific binding for highest product yield during your filtration processes.

High Flexibility

Sartopure® PP3 filter elements are available with a broad variety of retention ratings from 0.45 µm up to 100 µm making them ideally suited for numerous prefiltration applications.

Broad Compatibility and Low Extractables

Sartopure® PP3 filter elements consist entirely of polypropylene, which leads to broad chemical compatibility with a large number of solvents, acids and bases. Moreover the all-polypropylene design guarantees a small extractable profile.

Cost-Saving

The outstanding throughput performance and retention capability of Sartopure® PP3 enables downsizing of the required filtration area for pre- and final sterilizing grade filtration steps, resulting in significant cost savings.

Applications

- Protection of downstream equipment
- Clarification in vaccine production
- Ophthalmics
- Large volume parenterals
- Antibiotics & solvents
- Food & beverage

For food & beverage applications please refer to our dedicated F&B Sartopure® PP3 datasheet.

Technical Data

Available Sizes		Filtration Area [m ² ft ²]		
Cartridges T-Style Maxicaps® Maxicaps®	Retention Rate	0.45 µm	0.65 µm – 3 µm	5 µm – 100 µm
Size 1		0.4 4.3	0.45 4.8	0.65 7.0
Size 2		0.8 8.6	0.9 9.7	1.3 14
Size 3		1.2 12.9	1.35 14.5	1.95 21

Midicaps® Mini Cartridges	Retention Rate	0.45 µm	0.65 µm – 3 µm	5 µm – 100 µm
Size 7		0.04 0.43	0.05 0.54	0.065 0.7
Size 8		0.08 0.86	0.09 0.97	0.12 1.3
Size 9		0.13 1.4	0.15 1.6	0.21 2.3
Size 0*		0.26 2.8	0.30 3.2	0.42 4.5

Capsules	Retention Rate	0.45 µm	0.65 µm – 1.2 µm	3 µm – 100 µm
Size 4		0.012 0.13	0.013 0.14	0.018 0.19
Size 5		0.024 0.26	0.026 0.28	0.036 0.39

*only for Midicaps®

Max. Allowable Differential Pressure

Cartridges | Midicaps®

5 bar | 72.5 psi at 20°C
2 bar | 29 psi at 80°C

T-Style Maxicaps® | Maxicaps®

5 bar | 72.5 psi at 20°C
3 bar | 43.5 psi at 50°C

Capsules

4 bar | 58 psi at 20°C
2 bar | 29 psi at 80°C

Max. Allowable Back Pressure

2 bar | 29 psi at 20°C
(for all elements)

Max. Operating Pressure

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

5 bar | 72.5 psi at 20°C

Capsules

4 bar | 58 psi at 20°C

Materials

Filter Membrane

Polypropylene fleece

End Caps

Polypropylene

Support Fleece

Polypropylene

Housing

Polypropylene

Core

Polypropylene

O-Rings

Silicone
(optional: EPDM
or Fluoroelastomer)

Technical References

Validation Guide

85037-544-50, SPK5805-e

Extractables Analysis

EA 0.45 85037-544-51,

SPK5806-e

EA 0.65 -100 58037-544-52,

SPK5807-e

Sterilization

Cartridges

In-Line Steam Sterilization

134 °C, 20 min. at max differential pressure of 0.5 bar |
7.25 psi

Min. 25 Sterilization Cycles

or

Autoclaving

134°C, 2 bar, 30 min

Min. 25 Sterilization Cycles

Available Retention Rates

0.45 µm, 0.65 µm, 1.2 µm, 3 µm, 5 µm,
8 µm, 20 µm, 50 µm, 100 µm

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
- Pass USP Plastic Class VI Test
- Non-fiber releasing according to 21 CFR

T-Style Maxicaps® | Maxicaps® | Midicaps® |

Capsules Autoclaving

134°C, 2 bar, 30 min

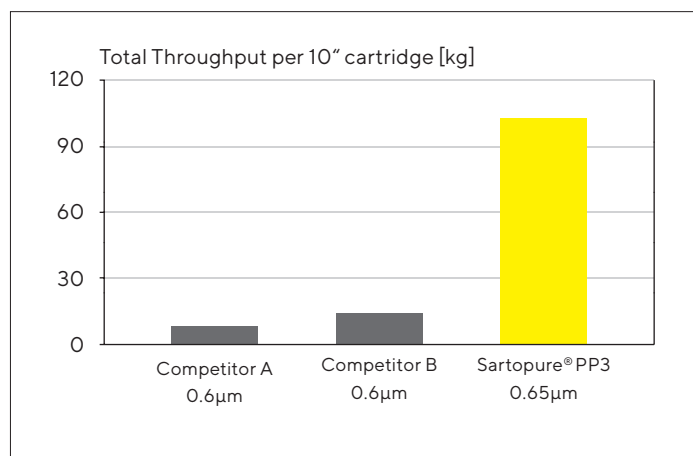
Min. 25 Sterilization Cycles

Note: T-Style Maxicaps®, Maxicaps®, Midicaps® and Capsules cannot be in-line steam sterilized

Performance

Standard Test Media 3-fold Concentrated

Total Throughput Comparison



Ordering Information



Cartridge

505	25	.	P	.
-----	----	---	---	---

Retention Rate	Filter Size
O6: 0.45 µm	1: 0.4 m ² 4.3 ft ² (10")
	2: 0.8 m ² 8.6 ft ² (20")
	3: 1.2 m ² 12.9 ft ² (30")
O5: 0.65 µm	1: 0.45 m ² 4.8 ft ² (10")
O3: 1.2 µm	2: 0.9 m ² 9.7 ft ² (20")
O2: 3 µm	3: 1.35 m ² 14.5 ft ² (30")
O4: 5 µm	1: 0.65 m ² 7 ft ² (10")
O1: 8 µm	2: 1.3 m ² 14 ft ² (20")
O20: 20 µm	3: 1.95 m ² 21 ft ² (30")
O50: 50 µm	
O98: 100 µm	



T-Style Maxicaps® | Maxicaps®

505	.	.	P	.	--	.	←
-----	---	---	---	---	----	---	---

Filter Design	Retention Rate	Filter Size	Connector Inlet	Connector Outlet
73: Maxicaps®	O6: 0.45 µm	1: 0.4 m ² 4.3 ft ² (10")	S: 1½" Tri-Clamp 50 mm	S: 1½" Tri-Clamp 50 mm
83: T-Style Maxicaps®		2: 0.8 m ² 8.6 ft ² (20")	O: ½" single stepped hose barb	O: ½" single stepped hose barb
		3: 1.2 m ² 12.9 ft ² (30")		
	O5: 0.65 µm	1: 0.45 m ² 4.8 ft ² (10")		
	O3: 1.2 µm	2: 0.9 m ² 9.7 ft ² (20")		
	O2: 3 µm	3: 1.35 m ² 14.5 ft ² (30")		
	O4: 5 µm	1: 0.65 m ² 7 ft ² (10")		
	O1: 8 µm	2: 1.3 m ² 14 ft ² (20")		
	O20: 20 µm	3: 1.95 m ² 21 ft ² (30")		
	O50: 50 µm			
	O98: 100 µm			



Midicaps®

505	53		P		--			--	
Retention Rate 06: 0.45 µm	Filter Size 7: 0.04 m ² 0.43 ft ² 8: 0.08 m ² 0.86 ft ² 9: 0.13 m ² 1.4 ft ² 0: 0.26 m ² 1.8 ft ²	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		Packing Size A: box of 4 (size 7,8,9) V: box of 2 (size 0)			
05: 0.65 µm 03: 1.2 µm 02: 3 µm	7: 0.05 m ² 0.54 ft ² 8: 0.09 m ² 0.97 ft ² 9: 0.15 m ² 1.61 ft ² 0: 0.30 m ² 3.23 ft ²								
42: 5 µm 01: 8 µm 20: 20 µm 50: 50 µm 98: 100 µm	7: 0.65 m ² 0.7 ft ² 8: 0.12 m ² 1.3 ft ² 9: 0.21 m ² 2.26 ft ² 0: 0.42 m ² 4.52 ft ²								



Mini Cartridge

505	15		P		--			--	
Retention Rate 05: 0.65 µm 03: 1.25 µm 02: 3 µm	Filter Size 7: 0.05 m ² 0.54 ft ² 8: 0.09 m ² 0.97 ft ² 9: 0.15 m ² 1.61 ft ²	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		Packing Size A: box of 4 (size 7,8,9) V: box of 2 (size 0)			
42: 5 µm 01: 8 µm 20: 20 µm 50: 50 µm 98: 100 µm	7: 0.65 m ² 0.7 ft ² 8: 0.12 m ² 1.3 ft ² 9: 0.21 m ² 2.26 ft ²								



Capsules


505	13		P		--			--	B
Retention Rate 06: 0.45 µm	Filter Size 4: 0.012 m ² 0.13 ft ² 5: 0.024 m ² 0.26 ft ²	Connector Inlet O: ¼" multiple stepped hose barb		Connector Outlet O: ¼" multiple stepped hose barb		Packing Size B: box of 5			
05: 0.65 µm 03: 1.2 µm	4: 0.013 m ² 0.14 ft ² 5: 0.026 m ² 0.28 ft ²								
02: 3 µm 42: 5 µm 01: 8 µm 20: 20 µm 50: 50 µm 98: 100 µm	4: 0.018 m ² 0.19 ft ² 5: 0.036 m ² 0.39 ft ²								

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