



# Sartoclean® CA

## Particle & Bioburden Reduction

### Product Information

- High product yield
- Double layer construction offers highest total throughput performance
- Mechanical and thermal resistance



Sartoclean® CA filter elements are the ideal choice for a broad range of prefiltration applications in the biopharmaceutical industry including particle removal and bioburden reduction. They offer a defined retention performance by size exclusion. The use of Sartoclean® CA prefilters avoids early blockage of downstream sterilizing-grade membrane filters and contributes significantly to an economical design of your filtration system.

#### High Product Yield

Throughout the years the cellulose acetate membranes of Sartoclean® CA filters have proven to be the membrane material with lowest nonspecific binding, ensuring high protein yield and rapid preservative recovery, enhancing your process efficiency.

#### Performance

Sartoclean® CA filters with heterogeneous double layer construction offer highest total throughput performance due to the "built-in prefiltration" to reduce filter change outs and improve economical process design.

#### Mechanical Strength

The reinforcement of the membrane results in increased mechanical and thermal resistance, which is of particular interest in applications with high differential pressure and repeated steam sterilization of filters.

#### Applications

- Plasma fractions
- Vaccines
- mAb
- Diagnostics
- Purified protein solutions
- Biological fluids
- Solutions containing preservatives

## Technical Data

Available Sizes	Filtration Area
<b>Cartridges   MaxiCaps®</b>	
Size 1 (10")	0.6 m <sup>2</sup>   6.5 ft <sup>2</sup>
Size 2 (20")	1.2 m <sup>2</sup>   12.8 ft <sup>2</sup>
Size 3 (30")	1.8 m <sup>2</sup>   19.3 ft <sup>2</sup>
<b>Mini Cartridges   MidiCaps®</b> (Mini Cartridges only available in size 9)	
Size 7	0.05 m <sup>2</sup>   0.5 ft <sup>2</sup>
Size 8	0.1 m <sup>2</sup>   1.1 ft <sup>2</sup>
Size 9	0.2 m <sup>2</sup>   2.2 ft <sup>2</sup>
Size 0	0.45 m <sup>2</sup>   4.8 ft <sup>2</sup>

### Max. Allowable Differential Pressure

**Cartridges | Mini Cartridges | MidiCaps®**  
5 bar | 72.5 psi at 20°C  
2 bar | 29 psi at 80°C

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

### Materials

**Prefilter Membrane**  
Cellulose Acetate

**Endfilter Membrane**  
Cellulose Acetate

**Support Fleece**  
Polypropylene

**Core**  
Polypropylene

**End Caps**  
Polypropylene

**Capsule Housing**  
Polypropylene

**O-Rings**  
Silicone

### Max. Allowable Back Pressure

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

### Max. Operating Pressure

**Cartridges | MaxiCaps® | Mini Cartridges | MidiCaps®**  
5 bar | 72.5 psi at 20°C

### Pore Size Combination

3.0 µm + 0.8 µm (nom.)  
0.8 µm + 0.65 µm (nom.)



## 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

#### In-Line Steam Sterilization

134°C, 0.5 bar | 7.25 psi, 20 min.

Min. 25 Sterilization Cycles

or

#### Autoclaving

134°C, 2 bar | 29 psi, 30 min.

Min. 25 Sterilization Cycles

### MaxiCaps® | MidiCaps®

#### Autoclaving

134°C, 2 bar | 29 psi, 30 min.

Min. 25 Sterilization Cycles

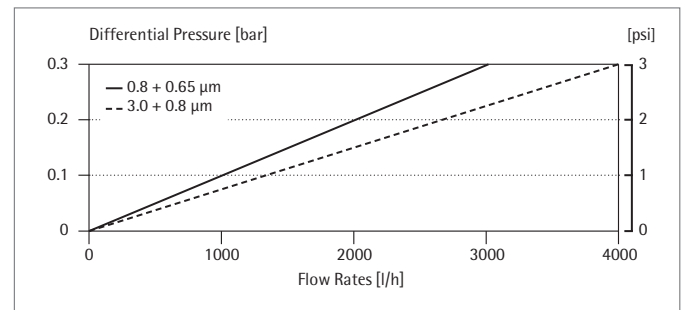
## Technical References

Validation Guide

SPK5788-e

## Performance

### Water Flow Rates for Sartoclean® CA 10" cartridges



Standardized at 20°C

## Ordering Information



### Cartridge

562				
<b>Adapter</b>	<b>Pore Size</b>	<b>Filter Size</b>		
25: 2 Flange Bayonet adapter with 226 double o-ring	04 E: Final filter 0.8 µm, prefilter 3.0 µm	1: 0.6 m <sup>2</sup>   6.5 ft <sup>2</sup> (10")		
27: Bayonet adapter with 222 double o-ring	05 G: Final filter 0.65 µm, prefilter 0.8 µm	2: 1.2 m <sup>2</sup>   12.8 ft <sup>2</sup> (20")		
28: 3 Flange Bayonet adapter with 222 double o-ring		3: 1.8 m <sup>2</sup>   19.3 ft <sup>2</sup> (30")		



### Mini Cartridge

562	15			9	--	--	--	B
<b>Pore Size</b>	<b>Filter Size</b>	<b>Packing Size</b>						
04 E: Final filter 0.8 µm, prefilter 3.0 µm	9: 0.2 m <sup>2</sup>   2.2 ft <sup>2</sup>	B: box of 5						
05 G: Final filter 0.65 µm, prefilter 0.8 µm								



### MaxiCaps®

562	73				--			
<b>Pore Size</b>	<b>Filter Size</b>	<b>Connector Inlet</b>	<b>Connector Outlet</b>					
04 E: Final filter 0.8 µm, prefilter 3.0 µm	1: 0.6 m <sup>2</sup>   6.5 ft <sup>2</sup>	S: 1½" Tri-Clamp 50 mm	S: 1½" Tri-Clamp 50 mm					
05 G: Final filter 0.65 µm, prefilter 0.8 µm	2: 1.2 m <sup>2</sup>   12.8 ft <sup>2</sup>	O: ½" single stepped hose barb	O: ½" single stepped hose barb					
	3: 1.8 m <sup>2</sup>   19.3 ft <sup>2</sup>	F: ¾" Tri-Clamp 25 mm	F: ¾" Tri-Clamp 25 mm					



### MidiCaps®

562	53				--				
<b>Pore Size</b>	<b>Filter Size</b>	<b>Connector Inlet</b>	<b>Connector Outlet</b>	<b>Packing Size</b>					
04 E: Final filter 0.8 µm, prefilter 3.0 µm	7: 0.05 m <sup>2</sup>   0.5 ft <sup>2</sup>	S: 1½" Tri-Clamp 50 mm	S: 1½" Tri-Clamp 50 mm	A: box of 4					
05 G: Final filter 0.65 µm, prefilter 0.8 µm	8: 0.1 m <sup>2</sup>   1.1 ft <sup>2</sup>	O: ½" single stepped hose barb	O: ½" single stepped hose barb	V: box of 2					
	9: 0.2 m <sup>2</sup>   2.2 ft <sup>2</sup>	F: ¾" Tri-Clamp 25 mm	F: ¾" Tri-Clamp 25 mm						
	0: 0.45 m <sup>2</sup>   4.8 ft <sup>2</sup>								

Check the availability of your desired configuration with your Sartorius sales representative or online at [www.sartorius-stedim.com](http://www.sartorius-stedim.com)

Sartorius Stedim Biotech GmbH  
 August-Spindler-Strasse 11  
 37079 Goettingen, Germany  
 Phone +49.551.308.0  
[www.sartorius-stedim.com](http://www.sartorius-stedim.com)

USA Toll-Free +1.800.368.7178  
 Argentina +54.11.4721.0505  
 Brazil +55.11.4362.8900  
 Mexico +52.55.5562.1102  
 UK +44.1372.737159  
 France +33.442.845600  
 Italy +39.055.63.40.41

Spain +34.913.586.098  
 Russian Federation  
 +7.812.327.53.27  
 Republic of South Africa  
 +27.11.315.5444  
 Japan +81.3.4331.4300  
 China +86.21.6878.2300



[www.sartorius-stedim.com](http://www.sartorius-stedim.com)