

Sartoclean[®] CA

Particle & Bioburden Reduction



Product Information

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

High Product Yield

Throughout the years the cellulose acetate membranes of Sartoclean[®] CA filters have proven to be a membrane material with very low non-specific binding, ensuring high protein yield and preservative recovery.

Performance

The Sartoclean® CA filter heterogeneous, double layer construction offers high total throughput performance due to the “built-in prefiltration”, which reduces filter change outs and improves 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
Cartridges T-Style Maxicaps® Maxicaps®	
Size 1 (10")	0.6 m ² 6.5 ft ²
Size 2 (20")	1.2 m ² 12.8 ft ²
Size 3 (30")	1.8 m ² 19.3 ft ²

Mini Cartridges Midicaps® (Mini Cartridges only available in size 9)	
Size 7	0.05 m ² 0.5 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

Materials

Prefilter Membrane

Cellulose Acetate

Endfilter Membrane

Cellulose Acetate

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 2822632)

Performance

Water Flow Rates

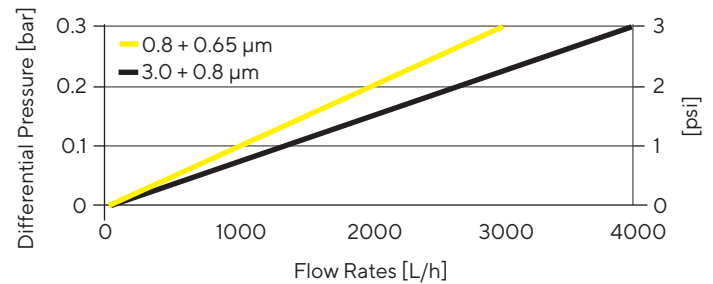


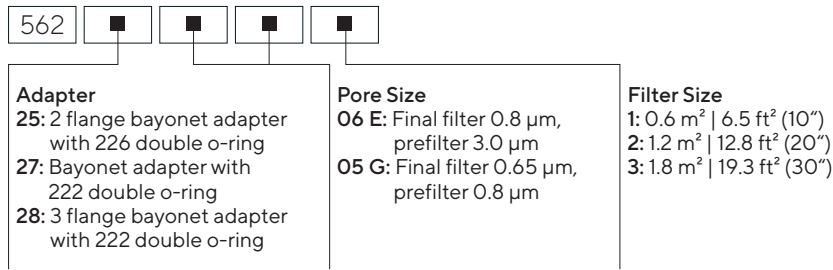
Figure 1: Water flow rates for 10" Sartoclean® CA cartridges. Standardized at 20 °C.



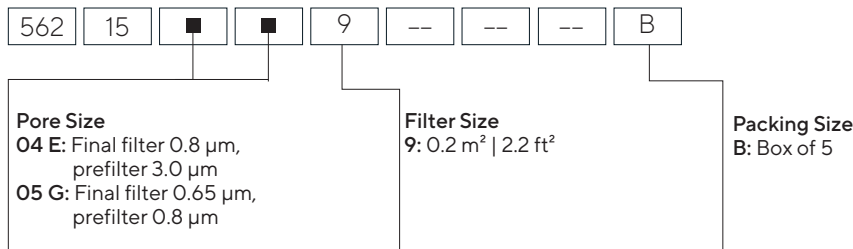
Ordering Information



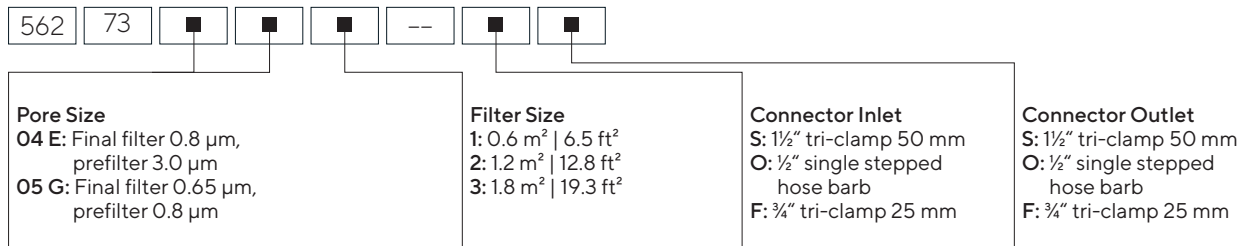
Cartridge



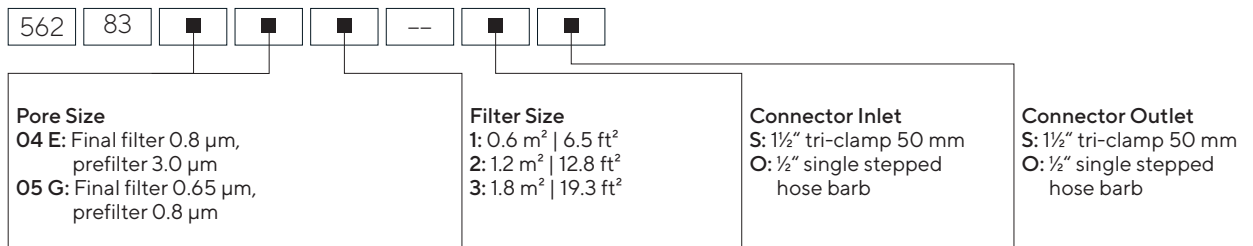
Mini Cartridge



Maxicaps®

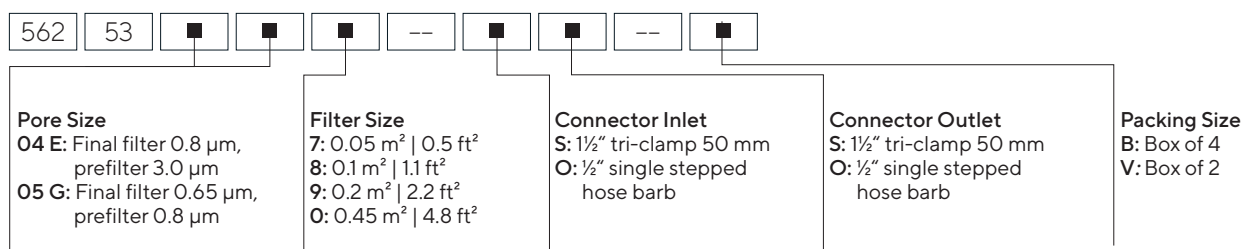


T-Style Maxicaps®





Midicaps®




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 Göttingen
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