

Sartolon

Sterilizing-Grade Filter Elements

- Sterile filtration of solvents and oily formulations
- Unique hydrophilic nylon membrane with broad chemical compatibility
- Heterogeneous double layer design provides outstanding throughput and exceptional flow rates

Perfectly Suited for Filtration of Solvents

Featuring a unique hydrophilic nylon (polyamide) membrane, Sartolon filters offer mechanical strength and broad chemical compatibility. They are the ideal choice for sterile filtration of solvents and oily formulations where chemical stability is needed.

Exceptional Filtration Performance

In addition to its wide range of applications, Sartolon filters provide outstanding total throughput and flow rates due to their heterogeneous double-nylon layer. The enhanced filtration performance guarantees process efficiency, reduces filtration costs and leads to short filtration cycles.



Scalability

Sartolon filter elements are available in a broad range of sizes and formats to provide linear scale-up from R&D to process scale. Membrane areas from 0.2 to 1.8 m² per filter element guarantee a high degree of flexibility and economic filter sizing.

Applications

- Solvents & Oily Formulations
- Antibiotics
- Large Volume Parenterals
- Bulk Pharmaceutical Chemicals
- ADCs

Services

Sartorius Confidence® Validation Services is the perfect complement to Sartolon.

Our services provide

- Extractables and leachables services
- Microbiological testing
- Physicochemical testing

Technical Data

Available Sizes	Filtration Area	Max. Diffusion at 2.5 bar 36 psi [ml/min]	Min. Bubble Point [bar psi]
Cartridges Maxicaps®			
Size 1	0.6 m ² 6.5 ft ²	13	3.0 43.5
Size 2	1.2 m ² 12.9 ft ²	26	3.0 43.5
Size 3	1.8 m ² 19.4 ft ²	39	3.0 43.5
Size 4 (only cartridge)	2.4 m ² 25.9 ft ²	52	3.0 43.5
Mini Cartridge Midicaps®			
Size 9	0.2 m ² 2.2 ft ²	4	3.0 43.5

Max. Allowable Differential Pressure

Cartridges | Mini Cartridges

5 bar | 72.5 psi at 20°C

2 bar | 29 psi at 80°C

Maxicaps® | Midicaps®

5 bar | 72.5 psi at 20°C

3 bar | 43.5 psi at 50°C

Max. Allowable Back Pressure

2 bar | 29 psi at 20°C

(for all elements)

in compliance with regulatory requirements. Our local teams of validation experts support you with our tailored and consultative approach to determine the most cost-effective solution and give you the confidence you need to succeed.

Materials

Prefilter Membrane

Nylon (Polyamide), reinforced with PET fleece

Endfilter Membrane

Nylon (Polyamide)

Support Fleece

Polypropylene

Core

Polypropylene

End Caps

Polypropylene

O-Rings

Silicone (optional EPDM)

Pore Size Combination

0.45 µm + 0.2 µm

Regulatory Compliance

- Each individual element is tested for integrity by Bubble Point and Diffusion test
- Fully validated as sterilizing grade filters according to ASTM current F-838 guidelines
- 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

Sterilization

Cartridges | Mini Cartridges

- In-Line Steam Sterilization
Max. 134 °C, 0.3 bar, 20 min.
Min. 25 Sterilization Cycles

or

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

Maxicaps® | Midicaps®

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

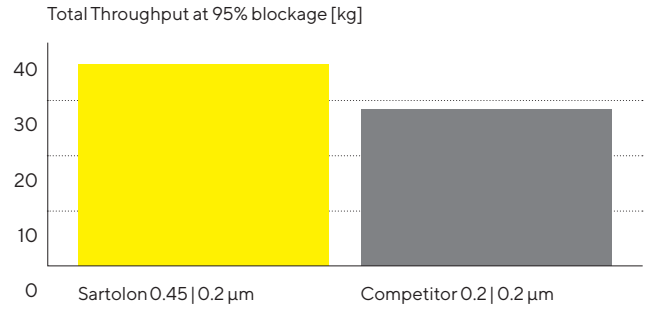
Technical References

Validation Guide SPK5716-e

Extractables Guide SPK5729-e

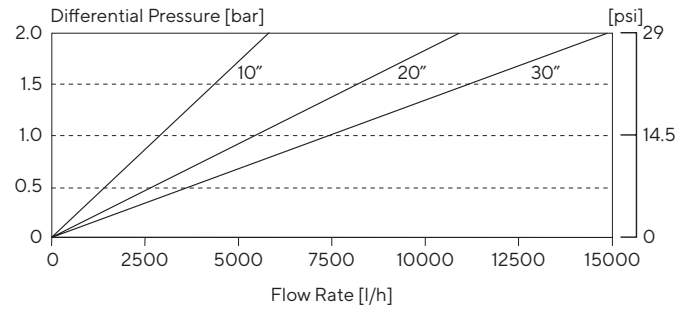
Performance

Total Throughput Comparison



10" Cartridge

Water Flow Rates Sartolon Maxicaps®



Standardized at 20°C

Ordering Information



Mini Cartridge

510 15 07 H 9 -- - -- -- B

Adapter
15: bayonet adapter with O-ring

Filter Size
9: 0.2 m² | 2.2 ft²

Packing Size
B: box of 5



Cartridge

510 ■ 07 H ■

Adapter
21: double open end with flat gaskets
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

Filter Size
1: 0.6 m² | 6.5 ft² (10")
2: 1.2 m² | 12.9 ft² (20")
3: 1.8 m² | 19.4 ft² (30")
4: 2.4 m² | 25.9 ft² (40")

(Standard with silicone O-ring, optionally with EPDM O-ring. Example 5102507H1---E)



Maxicaps®

510 73 07 H ■ -- ■ ■

Filter Size

1: 0.6 m² | 6.5 ft² (10")
2: 1.2 m² | 12.9 ft² (20")
3: 1.8 m² | 19.4 ft² (30")

Sterilization

--: autoclavable

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

(Optionally with vent valve design for connection to integrity tester. Example: 5107307H1--SSIT)



Midicaps®

510 53 07 H 9 -- ■ ■ -- A

Filter Size

9: 0.2 m² | 2.2 ft²

Sterilization

--: autoclavable

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

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