



Sartopore[®] 2 0.2 μm

Sterilizing Grade Filter Elements

Product Information

- Industry standard of high-performance liquid sterile filters
- Unique hydrophilic heterogeneous double layer Polyethersulfone membrane
- Exceptional high throughput and excellent flowrates at low pressure drops
- High thermal resistance and broad chemical compatibility



High-Performance and broad compatibility

Sartopore[®] 2 filter elements feature a unique hydrophilic heterogeneous double layer design of a 0.45 μm pre-filter and 0.2 μm final filter membrane with an exceptionally high throughput and flow-rate. In addition to its outstanding performance, the Polyethersulfone membrane gives Sartopore[®] 2 0.2 μm broad chemical compatibility, including a pH-range from pH 1 to pH 14, and a high thermal resistance.

Industry standard for sterile-grade liquid filters

Sartopore[®] 2 0.2 μm filters are fully validated as sterilizing grade filter elements according to current ASTM F-838 guidelines. Each individual element is integrity tested by diffusion and bubble point test prior to release, assuring absolute reliability. Sartopore[®] 2 filter elements are designed, developed and manufactured in accordance with an ISO 9001 certified Quality Management System. The Validation and Extractables Guide are available for compliance with regulatory requirements.

Scalability

Sartopore[®] 2 0.2 μm filter elements are available in a broad range of sizes and formats to provide linear scale-up from R&D to process scale. The wide range of membrane areas from 150 cm^2 to 1.8 m^2 per filter element guarantees greatest flexibility and most economic filter sizing.

Applications

- Therapeutics
- Biological Fluids
- Ophthalmic solutions
- Injectables
- Media
- SVPs, LVPs
- Antibiotics
- WFI
- Buffers
- Chemicals
- Cleaning and sanitizing agents
- Bulk pharmaceutical products

Technical Data

Available Sizes	Filtration Area	Max. Diffusion at 2.5 bar 36 psi [ml/min]	Min. Bubble Point [bar psi]
Cartridges, T-Style MaxiCaps® MaxiCaps® Gamma MaxiCaps®			
Size 0.5 (Only Cartridge)	0.3 m ² 3.2 ft ²	10	3.2 46
Size 1	0.6 m ² 6.5 ft ²	18	3.2 46
Size 2	1.2 m ² 12.9 ft ²	36	3.2 46
Size 3	1.8 m ² 19.4 ft ²	54	3.2 46
Mini Cartridges MidiCaps® Gamma MidiCaps®			
Size 7	0.05 m ² 0.5 ft ²	4	3.2 46
Size 8	0.1 m ² 1.1 ft ²	5	3.2 46
Size 9	0.2 m ² 2.2 ft ²	7	3.2 46
Size 0 (Only MidiCaps® & Gamma MidiCaps®)	0.45 m ² 4.8 ft ²	14	3.2 46
Capsules Gamma Capsules			
Size 4	0.015 m ² 0.16 ft ²	1	3.2 46
Size 5	0.03 m ² 0.32 ft ²	2	3.2 46

Max. Allowable Differential Pressure

Mini Cartridges | Cartridges

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

T-Style MaxiCaps® | MaxiCaps® | Gamma MaxiCaps® | MidiCaps® | Gamma MidiCaps®

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

Capsules | Gamma Capsules

4 bar | 58 psi at 20°C
2 bar | 29 psi at 50°C

Max. Allowable Back Pressure

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

Materials

Prefilter Membrane
Polyethersulfone, asymmetric

Endfilter Membrane
Polyethersulfone, asymmetric

Support Fleece
Polypropylene
(In-Line Steam sterilizable & autoclavable)
Polyester
(γ-irradiatable or γ-irradiatable | autoclavable)

Core
Polypropylene

End Caps
Polypropylene

Capsule Housing
Polypropylene

O-Ring
Silicone
(other materials on request)

Filling Bell
Polycarbonate

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

Sterilization

Mini Cartridges, Cartridges

In-Line Steam Sterilization

134 °C, 0.3 bar, 20 min.

Min. 25 Sterilization Cycles

or

Autoclaving

134°C, 2 bar, 30 min

Min. 25 Sterilization Cycles

MaxiCaps®, MidiCaps® & Capsules

Autoclaving

134°C, 2 bar, 30 min

Min. 25 Sterilization Cycles (MidiCaps® & MaxiCaps®)

Min. 5 Sterilization Cycles (Capsules)

Gamma MaxiCaps®, Gamma MidiCaps® & Gamma Capsules

Gamma Irradiation

≤ 50 kGy

1 Sterilization Cycle

T-Style MaxiCaps®

Autoclaving

134°C, 2 bar, 30 min

Min. 5 Sterilization Cycles

or

Gamma Irradiation

≤ 50 kGy

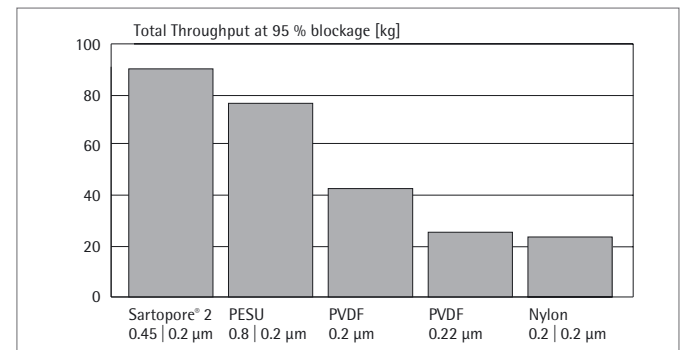
1 Sterilization Cycle

Technical References

Validation Guide	SPK5802-e
Extractables Guide	SPK5731-e

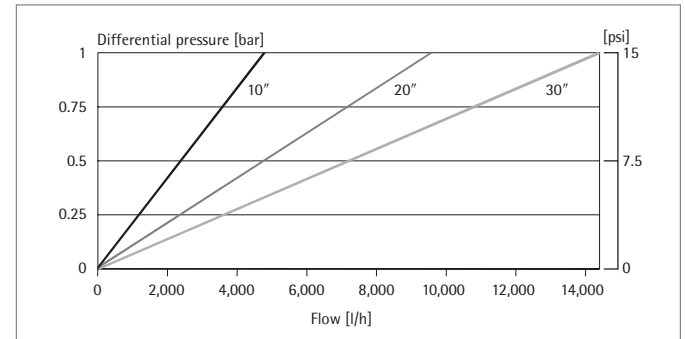
Performance

Total Throughput Comparison



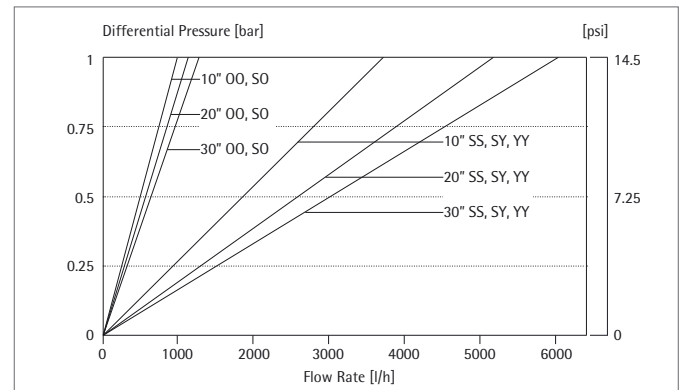
10" Cartridge format

Water Flow Rates for Standard Cartridges



Standardized at 20 °C

Water Flow Rates for T-Style MaxiCaps®



Ordering Information



Mini Cartridge

544 15 07 H B

Adapter

15: Bayonet adapter with o-ring
18: Plug Adapter with double o-ring

Filter Size

7: 0.05 m² | 0.5 ft²
8: 0.1 m² | 1.1 ft²
9: 0.2 m² | 2.2 ft²

Packing Size

B: box of 5

(Standard with silicone o-ring optional with EPDM o-ring. Example: 5441507H7---E--B)



Cartridge

544 25 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

0: 0.3 m² | 3.2 ft² (5")
1: 0.6 m² | 6.5 ft² (10")
2: 1.2 m² | 12.9 ft² (20")
3: 1.8 m² | 19.4 ft² (30")

(Standard with silicone o-ring optional with EPDM o-ring. Example: 5442507H1---E)



T-Style MaxiCaps®

544 83 07 H . G-

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

G-: γ-irradiatable and autoclavable

Connector Inlet

S: 1½" Tri-Clamp 50.5 mm
O: ½" single stepped hose barb
Y: 1" single stepped hose barb

Connector Outlet

S: 1½" Tri-Clamp 50.5 mm
O: ½" single stepped hose barb
Y: 1" single stepped hose barb

MaxiCaps® | Gamma MaxiCaps®

544 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

G-: γ-irradiatable
---: autoclavable

Connector Inlet

S: 1½" Tri-Clamp 50.5 mm
O: ½" single stepped hose barb
F: ¾" Tri-Clamp 25 mm

Connector Outlet

S: 1½" Tri-Clamp 50.5 mm
O: ½" single stepped hose barb
F: ¾" Tri-Clamp 25 mm

(Optional with vent valve design for connection of integrity tester. Example: 5447307H1--SSIT)



MidiCaps® | Gamma MidiCaps®

544 53 07 H

Filter Size

7: 0.05 m² | 0.5 ft²
8: 0.1 m² | 1.1 ft²
9: 0.2 m² | 2.2 ft²
0: 0.45 m² | 4.8 ft²

Sterilization

G-: γ-irradiatable
---: autoclavable

Connector Inlet

S: 1½" Tri-Clamp 50.5 mm
O: ½" single stepped hose barb
F: ¾" Tri-Clamp 25 mm
H: ¼" multiple stepped hose barb (only size 7 with filling bell)

Connector Outlet

S: 1½" Tri-Clamp 50.5 mm
O: ½" single stepped hose barb
F: ¾" Tri-Clamp 25 mm
H: ¼" multiple stepped hose barb (only size 7 with filling bell)

Packing Size

A: box of 4 (size 7,8,9)
V: box of 2 (size 0)

Capsules | Gamma Capsules

544 13 07 H B

Filter Size

4: 0.015 m² | 0.16 ft²
5: 0.03 m² | 0.32 ft²

Sterilization

G-: γ-irradiatable
---: autoclavable

Connector Inlet

S: ¾" Tri-Clamp 25 mm
O: ¼" multiple stepped hose barb

Connector Outlet

S: ¾" Tri-Clamp 25 mm
O: ¼" multiple stepped hose barb

Packing Size

B: box of 5

Check the availability of your desired configuration with your Sartorius sales representative or online at www.sartorius-stedim.com

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany
Phone +49.551.308.0
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