



Sartopore[®] 2 XLI 0.2 μm

Sterilizing Grade Filter Elements

Product Information

- Lowest filter consumption for filtration of chemically defined process fluids
- Outstanding capacity and flow performance
- Ideal solution for conventional and single-use processes
- Highly economic sizing for intermediate steps in downstream processing



30% Enlarged Effective Filtration Area

The combination of the built-in 0.35 μm pre-filter in front of a 0.2 μm final filter together with the 30% enlarged effective filtration area per filter element provide an outstanding total throughput and flow rate performance in the target applications. Thus ensuring highest process efficiency, minimized filtration costs and short filtration cycle times.

Scalability

Sartopore[®] 2 XLI 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 210 cm^2 to 2.4 m^2 per filter element guarantees greatest flexibility and most economic filter sizing.

Applications

- Ophthalmic solutions
- Chemically defined cell culture media
- High viscous large volume parenterals
- Any fully chemically defined media

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 1 | 0.8 m ² 8.6 ft ² | 21 | 3.2 46 |
| Size 2 | 1.6 m ² 17.2 ft ² | 42 | 3.2 46 |
| Size 3 | 2.4 m ² 25.8 ft ² | 63 | 3.2 46 |
| MidiCaps® Gamma MidiCaps® | | | |
| Size 7 | 0.065 m ² 0.7 ft ² | 4.5 | 3.2 46 |
| Size 8 | 0.13 m ² 1.4 ft ² | 5.5 | 3.2 46 |
| Size 9 | 0.26 m ² 2.8 ft ² | 8 | 3.2 46 |
| Size 0 | 0.52 m ² 5.6 ft ² | 16 | 3.2 46 |
| Capsules Gamma Capsules | | | |
| Size 4 | 0.021 m ² 0.22 ft ² | 1.1 | 3.2 46 |

Max. Allowable Differential Pressure

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 Size 4

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

0.35 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

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

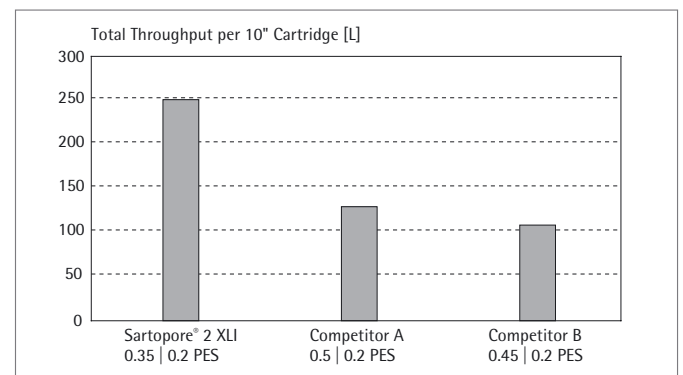
1 Sterilization Cycle

Technical References

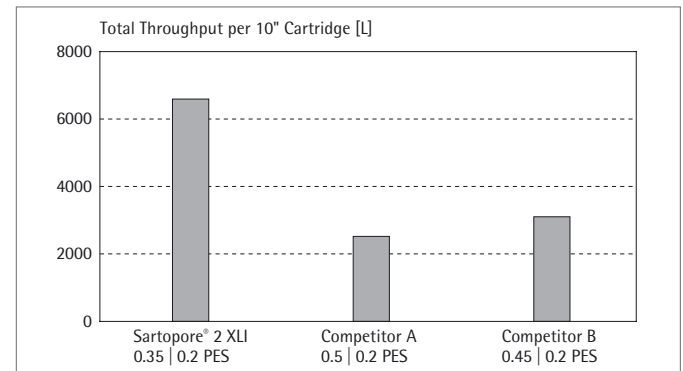
Validation Guide SPK5811-e
Extractables Guide SPK5776-e

Performance

Chemically Defined Cell Culture Media



Ophthalmic Solution



Ordering Information



Cartridge



Adapter
25: 2 Flange Bayonet adapter
with 226 double o-ring

Filter Size
1: 0.8 m² | 8.6 ft² (10")
2: 1.6 m² | 17.2 ft² (20")
3: 2.4 m² | 25.8 ft² (30")

(Standard with silicone o-ring optional with EPDM o-ring. Example: 544250711---E)



T-Style MaxiCaps®



Filter Size
1: 0.8 m² | 8.6 ft² (10")
2: 1.6 m² | 17.2 ft² (20")
3: 2.4 m² | 25.8 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®



Filter Size
1: 0.8 m² | 8.6 ft² (10")
2: 1.6 m² | 17.2 ft² (20")
3: 2.4 m² | 25.8 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: 544730711--SSIT)



MidiCaps® | Gamma MidiCaps®



Filter Size
7: 0.065 m² | 0.7 ft²
8: 0.13 m² | 1.4 ft²
9: 0.26 m² | 2.8 ft²
0: 0.52 m² | 5.6 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



Filter Size
4: 0.021 m² | 0.22 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