



Sartopore® 2 XLG 0.2 µm

Sterilizing Grade Filter Elements

Product Information

- Lowest filter consumption for filtration of cell culture media
- Outstanding capacity and flow performance
- Ideal solution for conventional and single-use processes
- Highly economic sizing for intermediate steps in downstream processing



Reduce 50–90% Filtration Costs

Sartopore® 2 XLG uses a coarser pre-filter layer (0.8 µm) in front of the sterilizing 0.2 µm rated final membrane. The coarser pre-filter layer avoids early blocking in many biological liquids. The unique filtration capacity can reduce filtration costs by 50–90% for many processes.

Scalability

Sartopore® 2 XLG 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 cm² to 2.4 m² per filter element guarantees greatest flexibility and most economic filter sizing.

Applications

- Filtration of cell culture media such as:
 - Plant peptone or yeast supplemented cell culture media
 - Serum containing cell culture media
 - Other cell culture media used in biotech manufacturing
- Post cell harvest filtration
- Downstream Intermediates (before and after UF | DF and chromatography steps)

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	23	3.2 46
Size 2	1.6 m ² 17.2 ft ²	46	3.2 46
Size 3	2.4 m ² 25.8 ft ²	69	3.2 46
MidiCaps® Gamma MidiCaps®			
Size 7	0.065 m ² 0.7 ft ²	5	3.2 46
Size 8	0.13 m ² 1.4 ft ²	6	3.2 46
Size 9	0.26 m ² 2.8 ft ²	9	3.2 46
Size 0	0.52 m ² 5.6 ft ²	18	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

0.8 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

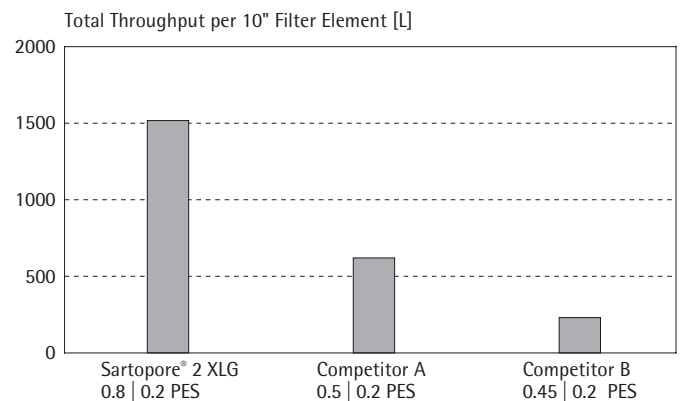
SPK5772-e

Extractables Guide

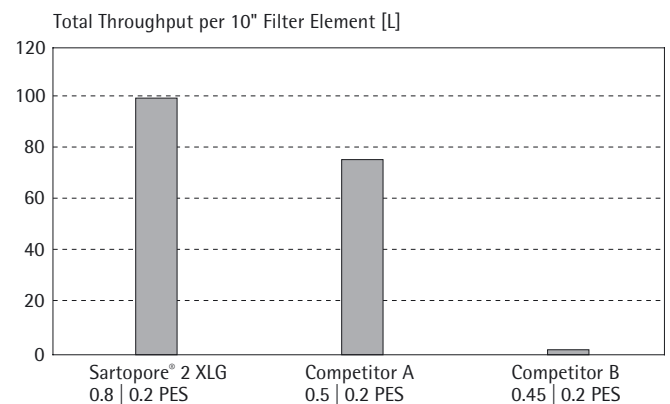
SPK5775-e

Performance

Soy Peptone Supplemented Cell Culture Media



Monoclonal Antibody Pool



Antibody Concentration: 47.5 mg/ml

Ordering Information



Cartridge

544 25 07 G 1

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: 5442507G1---E)



T-Style MaxiCaps®

544 83 07 G 1 G- 1 1

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®

544 73 07 G 1 1 1 1

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: 5447307G1--SOIT)



MidiCaps® | Gamma MidiCaps®

544 53 07 G 1 1 1 1 -- 1

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

544 13 07 G 4 1 1 1 -- B

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