

Sartopore® 2 XLM

Sterilizing-Grade & Mycoplasma Retentive Filter Elements



Product Information

- Lowest filter consumption by high filtration capacity
- Lowest risk of Mycoplasma contamination for your fermentation process
- High performance 0.1 µm Mycoplasma retentive filter
- Fully validated as sterilizing-grade filter (according to current ASTM F-838 guideline)

Background Information

It is well known that Mycoplasma can be present in commercially available cell culture media. Because Mycoplasma are able to penetrate 0.2 µm rated sterilizing-grade filters, Sartorius specifically developed the Sartopore® 2 XLM with finer filter membranes (0.1 µm) for the successful retention of Mycoplasma.

Exceptional Performance

The 0.2 µm prefilter layer is specifically designed to protect the final 0.1 µm membrane layer in cell culture media filtration applications. The optimized membrane combination together with the exceptionally high effective filtration area per filter element leads to remarkable throughput performance and Mycoplasma retention. This reduces the filtration costs per liter, the filtration time as well as the footprint of the filtration system.

Scalability

Sartopore® 2 XLM 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.

Fully Validated - Reliable Bacterial Retention

- Fully validated as sterilizing-grade filter (LRV >7/cm² filter area for *Brevundimonas diminuta* according to current ASTM F-838 guideline)
- Validated for retaining *Acholeplasma laidlawii* with an LRV7/cm² filter area using the filtration method published by the PDA Mycoplasma Task Force (PDA J Pharm Sci and Tech 2014, 68 281-296)
- This filter has also successfully been tested for *Mycoplasma orale* and *Mycoplasma fermentans*
- Sartopore® 2 XLM also retains *Leptospira licerasiae* (LRV <7/cm²)

Applications

Sterilizing-grade filtration and Mycoplasma removal from:

- Plant peptone or yeast supplemented cell culture media
- Chemically defined cell culture media
- Cell culture media components such as serum (e. g. Fetal Calf Serum)
- All other cell culture media used in biotech manufacturing

Any other application requiring below 0.2 µm filtration for enhanced sterility assurance.

Services

Sartorius Confidence® Validation Services is the perfect complement to Sartopore® 2 XLM filters.

Our services provide

- Extractables and leachables services
- Microbiological testing
- Physicochemical testing

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.

Technical Specifications

Available Sizes	Filtration Area	Max. Diffusion at 4.0 bar 58 psi [ml/min]
Cartridges, T-Style Maxicaps®, Maxicaps® Gamma Maxicaps®		
Size 0.5	0.4 m ² 4.3 ft ²	19
Size 1	0.8 m ² 8.6 ft ²	39
Size 2	1.6 m ² 17.2 ft ²	78
Size 3	2.4 m ² 25.8 ft ²	117
Midicaps® Gamma Midicaps®		
Size 7	0.065 m ² 0.7 ft ²	5
Size 8	0.13 m ² 1.4 ft ²	7
Size 9	0.26 m ² 2.8 ft ²	12
Size 0	0.52 m ² 5.6 ft ²	24
Capsules Gamma Capsules		
Size 4	0.021 m ² 0.22 ft ²	1.5



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

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

Max. Allowable Back Pressure

2 bar | 29 psi at 20°C

(for all elements)

Pore Size Combination

0.2 μm + 0.1 μm

Regulatory Compliance

- For release, 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
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® & Capsules

Autoclaving

Max. 134°C, 2 bar, 30 min
Min. 25 Sterilization Cycles (Midicaps®)
Min. 5 Sterilization Cycles (Capsules)

Gamma Maxicaps®, Gamma Midicaps® & Gamma Capsules

Gamma Irradiation

≤ 50 kGy
1 Sterilization Cycle

T-Style Maxicaps®

Autoclaving

Max. 134°C, 2 bar, 30 min
Min. 5 Sterilization Cycles

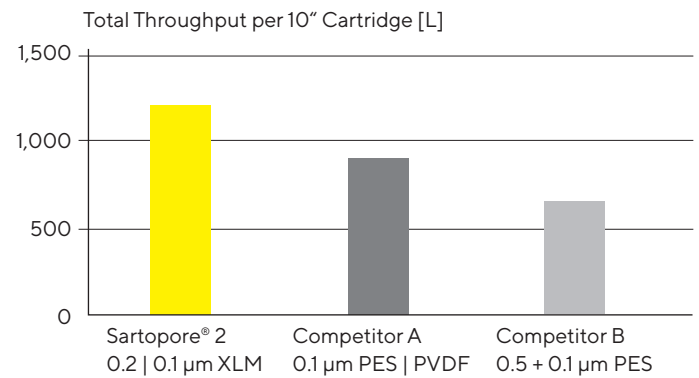
or

Gamma Irradiation

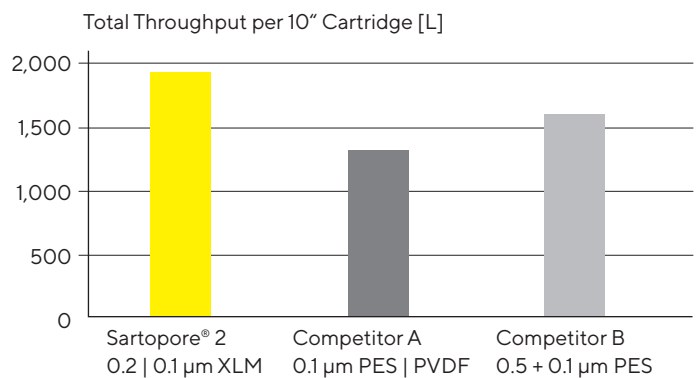
≤ 50 kGy
1 Sterilization Cycle

Performance

Plant Peptone Supplemented Cell Culture Media



Soy Supplemented Cell Culture Media



Technical References

Validation Guide 2665248
Extractables Guide SPK5823-e

Ordering Information



Cartridge

544 25 58 M ■

Adapter

25: 2 flange bayonet adapter with 226 double o-ring

Filter Size

0: 0.4 m² | 4.3 ft² (5")
 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 or Fluoroelastomer o-ring).



T-Style Maxicaps®

544 83 58 M ■ G- ■ ■

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 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 58 M ■ ■ ■ ■

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: 5447358M1G-SOIT)



Midicaps® | Gamma Midicaps®

544 53 58 M ■ ■ ■ ■ -- ■

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 58 M 4 ■ ■ ■ -- 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" single stepped hose barb

Packing Size

B: box of 5

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 **For more information, visit**
www.sartorius.com