



Sartopore® Platinum

Sterilizing Grade Filter Elements

Product Information

- Smaller filters, less filters by high filtration capacity
- Smaller dimensioning of assemblies – smaller bags, less tubing, hardware
- Faultless integrity testing - no risk of wetting issues
- Highest product yield by lowest adsorption



Unique Surface Modification

A patented membrane hydrophilization process is used to permanently modify the membrane surface.

This technology provides those membrane surface properties that are responsible for the outstanding wettability and low protein binding character of the Sartopore® Platinum membrane, even after extreme thermal and chemical stress, without affecting wettability and integrity testing.

Perfect Solution for Single-Use Processes

The perfect wettability of the PES membrane leads to drastic reduction of the required flushing volumes (up to 95% less water needed) which is highly beneficial especially for single-use processes. Due to the low flushing volume the required waste bags can be significantly downsized by which costs are reduced and handling is optimized.

Maximum Yield for High-Value Products

Besides high filtration capacity the surface modification leads to lowest protein adsorption of any PES membrane in the market by which product yield can be maximized.

Reliable Integrity Testing

Imperfect wetting is the most frequent reason for failed integrity tests. Especially for single-use processes re-wetting possibility is limited.

The extraordinary wetting behavior of Sartopore® Platinum thus helps to eliminate this important risk factor. Using Sartopore® Platinum the process of integrity testing meets a yet unmet degree of reliability.

Flexible Cartridge Formats

Sartopore® Platinum filter cartridges are available in standard (0.6 m²/10") and large membrane version (1 m²/10"). This allows highly economic filter sizing and also capacity adaptation without replacing stainless steel housings.

Applications

- Production of high-value biologicals like monoclonal antibodies
- Final conjugated bulk
- Point of fill filtration
- Especially beneficial for single-use filtration processes
- Blood & plasma processes
- Ophthalmics

Technical Data

Available Sizes	Filtration Area	Max. Diffusion at 2.5 bar 36 psi [ml/min]	Min. Bubble Point [bar psi]
Cartridges			
Size 1	0.6 m ² 6.5 ft ²	15	3.5 51
Size 2	1.2 m ² 12.9 ft ²	30	3.5 51
Size 3	1.8 m ² 19.4 ft ²	45	3.5 51
Cartridges, T-Style MaxiCaps[®], MaxiCaps[®]			
Size 1	1.0 m ² 10.8 ft ²	25	3.5 51
Size 2	2.0 m ² 21.5 ft ²	50	3.5 51
Size 3	3.0 m ² 32.3 ft ²	75	3.5 51
MidiCaps[®] Gamma MidiCaps[®]			
Size 7	0.065 m ² 0.7 ft ²	4	3.5 51
Size 8	0.13 m ² 1.4 ft ²	5	3.5 51
Size 9	0.26 m ² 2.8 ft ²	7	3.5 51
Size 0	0.52 m ² 5.6 ft ²	14	3.5 51
Capsules Gamma Capsules			
Size 4	0.021 m ² 0.22 ft ²	1.1	3.5 51

Max. Allowable Differential Pressure

Cartridges

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

T-Style MaxiCaps[®], 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
(optional EPDM or
Fluoroelastomer)

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

Cartridges

In-Line Steam Sterilization (dry or wet steaming)

134°C, 0.3 bar, 20 min.

Min. 25 Sterilization Cycles

or

Autoclaving

134°C, 2 bar, 30 min

Min. 25 Sterilization Cycles

MidiCaps® & Capsules

Autoclaving

134°C, 2 bar, 30 min

Min. 25 Sterilization Cycles (MidiCaps®)

Min. 5 Sterilization Cycles (Capsules)

Gamma MidiCaps® & Gamma Capsules

Gamma Irradiation

≤ 50 kGy

1 Sterilization Cycle

T-Style MaxiCaps® & MaxiCaps®

Autoclaving

134°C, 2 bar, 30 min

Min. 5 Sterilization Cycles

or

Gamma Irradiation

≤ 50 kGy

1 Sterilization Cycle

Technical References

Validation Guide

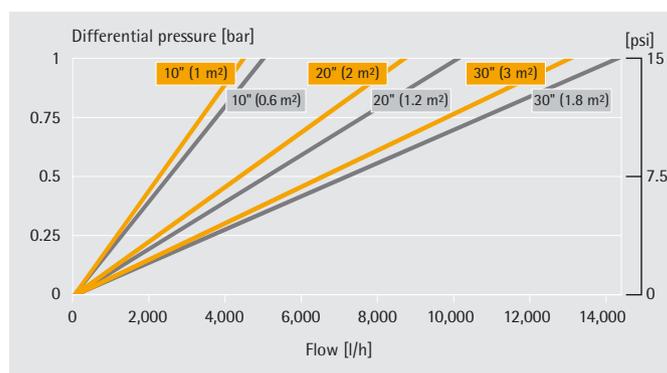
SPK5795-e

Extractables Guide

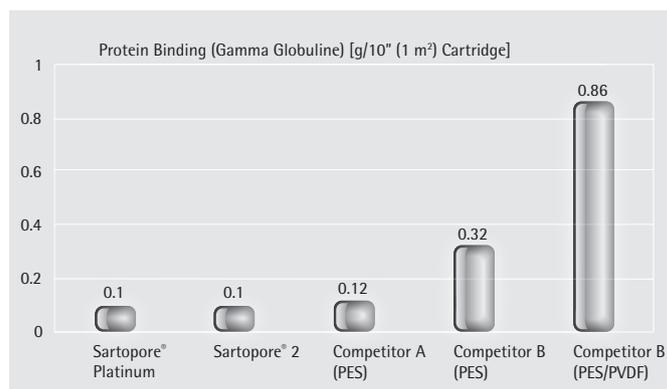
SPK5799-e

Performance

Water Flow Rates 10", 20", 30"



Unspecific Protein Binding



Ordering Information



Cartridge

549 25 07 H

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

Filter Size
1: 1.0 m² | 10.8 ft² (10")
2: 2.0 m² | 21.5 ft² (20")
3: 3.0 m² | 32.3 ft² (30")
1-S: 0.6 m² | 6.5 ft² (10")
2-S: 1.2 m² | 12.9 ft² (20")
3-S: 1.8 m² | 19.4 ft² (30")

(Standard with silicone o-ring optional with EPDM o-ring. Example: 5492507H1---E)



T-Style MaxiCaps®

549 83 07 H G-

Filter Size
1: 1.0 m² | 10.8 ft² (10")
2: 2.0 m² | 21.5 ft² (20")
3: 3.0 m² | 32.3 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®

549 73 07 H G-

Filter Size
1: 1.0 m² | 10.8 ft² (10")
2: 2.0 m² | 21.5 ft² (20")
3: 3.0 m² | 32.3 ft² (30")

Sterilization
G-: γ-irradiatable and 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: 5497307H1--SOIT)



MidiCaps® | Gamma MidiCaps®

549 53 07 H

Filter Size
7: 0.065 m² | 0.7 ft²
8: 0.13 m² | 1.4 ft²
9: 0.26 m² | 2.8 ft²
O: 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

549 13 07 H 4

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