



Sartofluor[®] LG MaxiCaps[®] & Cartridges

Membrane Filtration of aggressive media

Product Information

The hydrophobic ePTFE membrane material is widely used for air|gas filters. PTFE is chemically extremely resistant and is a powerful membrane material also for solvent filtration.



Description

It is known that the requirements on membrane construction are somewhat different for air and liquid filtration. Therefore, the design of Sartofluor[®] LG (LG = liquid & gases) cartridges and MaxiCaps[®] was specially adapted to the needs of liquid filtration. This consequently resulted in a modified construction with a different pleating structure by which the effective filtration area is reduced compared to Sartofluor[®] GA (GA = gases & air). However, the modified construction is beneficial with regard to hydrodynamic flow and thus Sartofluor[®] LG filters are especially recommended for solvent filtration but can also be used as an air|gas filter.

Applications

Sartofluor[®] LG filters are available in cartridge and MaxiCaps[®] format. Sartofluor LG MaxiCaps[®] improve the process security of sterile filtration of aggressive media (acids and bases) and solvents. There is no need to open the filter housing after filtration. The capsule design allows filtration of such media without any handling of the contaminated filter cartridge post use.

Chemical Resistance

Sartofluor[®] LG filters show a high mechanical and chemical robustness. A compatibility list for standard solvents is part of the Validation Guide. The compatibility for other substances can be proven in specific validation studies.

Design

The high quality of the used PTFE membrane allows the construction as a single-layer design with compromising safety. The mechanical robustness of the membrane does not require a double-layer construction and therefore maximizes the flow rate and avoids potential problems during sterilization.

Documentation

Sartofluor[®] LG Cartridges and MaxiCaps[®] are designed, developed and manufactured in accordance with an ISO 90001 certified Quality Management System. A Validation Guide is available for compliance with regulatory requirements.

Cost

Sartofluor[®] LG MaxiCaps[®] remove the need for investment in stainless steel or PVDF filter housings and an inventory of spare parts such as valves and o-rings.

Technical Data

Specifications

Materials	
Filter Membrane	PTFE (single layer)
Support Fleece	Polypropylene
Core	Polypropylene
End Caps	Polypropylene
O-Rings	MaxiCaps®: EPDM (Fluoroelastomer as accessory in the package) Cartridges: Silicone (optional EPDM or Fluoroelastomer)

Pore Size

0.45 µm
0.2 µm
0.1 µm (only Cartridges)

Available Sizes | Filtration Area

MaxiCaps®		
Size 1	10"	0.5 m ² 5.4 ft ²
Size 2	20"	1.0 m ² 10.8 ft ²
Size 3	30"	1.5 m ² 16.1 ft ²

Available Connectors MaxiCaps®

S: Tri-Clamp 50 mm (1½")
F: Tri-Clamp 25 mm (¾")

Available Adapters Cartridges

25: 2 Flange Bayonet adapter with 226 double o-ring
(others on request)

Operating Parameters

Max. Allowable Differential Pressure	5 bar 72.5 psi at 20°C (Cartridges, MaxiCaps®) 2 bar 29 psi at 80°C (Cartridges) 3 bar 43.5 psi at 50°C (MaxiCaps®)
Max. Allowable Back Pressure	2 bar 29 psi at 20°C

Extractables

Sartofluor® LG MaxiCaps® and Cartridges meet, or exceed the requirements for WFI quality standards set by the current USP.

Regulatory Compliance

Individually integrity tested

Integrity test correlated to HIMA/ASTM
F-838-05 Bacteria Challenge Test

Non-pyrogenic according to USP Bacterial Endotoxins

Meets USP Plastics Class VI biological
reactivity test, in vivo

Non-fiber releasing according to 21 CFR

Sterilization

Autoclaving

134°C, 2 bar | 29 psi, 30 min

No In-Line Steam Sterilization

In-Line Steam Sterilization

134°C, 20 min. at max differential pressure
of 0.5 bar | 7.25 psi (only Cartridges)

Note: No In-Line Steam Sterilization for MaxiCaps®

Sterilization Cycles

Autoclaving: Min. 25

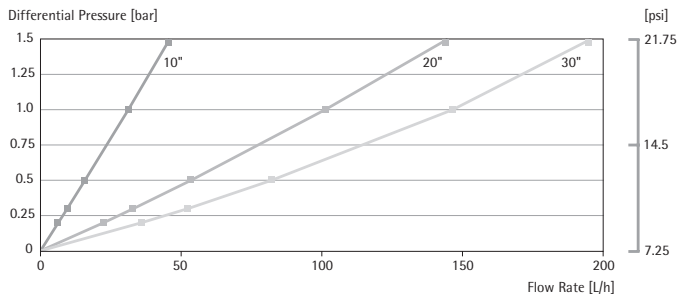
In-Line Steam Sterilization

min. 25 (only Cartridges)

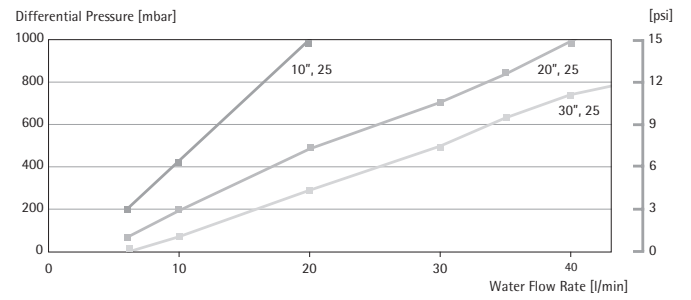
Technical References

Validation Guide SPK5708-e

Liquid Flow Rates for Sartofluor® LG MaxiCaps® 0.2 µm



Liquid Flow Rates for Sartofluor® LG Filter Cartridges 0.2 µm



Ordering Information

Order Code	Pore Size [µm]	Pack Size [Pieces]
MaxiCaps®		
5187306T1--**	0.45	1
5187307T1--**	0.2	1
5187307T2--**	0.2	1
5187307T3--**	0.2	1
Cartridges		
5182506T1----LG	0.45	1
5182506T2----LG	0.45	1
5182506T3----LG	0.45	1
5182507T1----LG	0.2	1
5182507T2----LG	0.2	1
5182507T3----LG	0.2	1
5182558T1----LG	0.1	1
5182558T2----LG	0.1	1
5182558T3----LG	0.1	1

** Connector Styles
(first code letter represents the inlet, the second the outlet)

Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5 - 7
34302 Guxhagen

Phone +49.5665.407.0
Fax +49.5665.407.2200

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques - CS 91051
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.845619

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques - CS 71058
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH
Franzosengraben 12
1030 Vienna

Phone +43.1.7965763.18
Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde

Phone +32.2.756.06.80
Fax +32.2.756.06.81

Hungary

Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi

Phone +36.23.457.227
Fax +36.23.457.147

Italy

Sartorius Stedim Italy S.p.A.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41
Fax +39.055.63.40.526

Netherlands

Sartorius Stedim Netherlands B.V.

Phone +31.30.60.25.080
Fax +31.30.60.25.099

filtratie.nederland@sartorius-stedim.com

Poland

Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.647.38.40
Fax +48.61.879.25.04

Russian Federation

LLC "Sartorius ICR"
Uralskaya str. 4, Lit. B
199155, Saint-Petersburg

Phone +7.812.327.5.327
Fax +7.812.327.5.323

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone +34.902.110.935
Fax +34.91.358.96.23

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen

Phone +41.52.354.36.36
Fax +41.52.354.36.46

U.K.

Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ

Phone +44.1372.737159
Fax +44.1372.726171

America

USA

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716

Toll-Free +1.800.368.7178
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900
Fax + 55.11.4362.8901

Mexico

Sartorius de México S.A. de C.V.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius Stedim Biotech (Beijing) Co. Ltd.
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300

Phone +86.10.80426516
Fax +86.10.80426580

Sartorius Stedim Biotech (Beijing) Co. Ltd.
Shanghai Branch Office
3rd Floor, North Wing, Tower 1
No. 4560 Jin Ke Road
Pudong District, Shanghai 201210

Phone +86.21.68782300
Fax +86.21.68782332 | 68782882

Sartorius Stedim Biotech (Beijing) Co. Ltd.
Guangzhou Representative Office
Unit K, Building 23
Huihua Commerce & Trade Building
No. 80 Xianlie Middle Road
Guangzhou 510070

Phone +86.20.37618687 | 37618651
Fax +86.20.37619051

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India

Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.4331.4300
Fax +81.3.4331.4301

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Stedim Singapore Pte. Ltd.
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528

Phone +65.6872.3966
Fax +65.6778.2494

South Korea

Sartorius Korea Biotech Co., Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700
Fax +82.31.622.5799



◀▶ www.sartorius-stedim.com