

Do you want to perform more tests with SartoScale 25 or scale-up your process?

SartoScale 25

Sartopore® 2 XLM: 5445358MV--LX--C
Sartoguard PES 0.1 µm: 5475358GV--LX--C
Sartoguard GF 0.1 µm: 5485358MV--LX--C
Sartoguard NF 0.1 µm: 5465358MV--LX--C

Capsules size 4 for scale-up

Sartopore® 2 XLM: 5441358M4--**--B
Sartoguard PES 0.1 µm: 5471358G4--**--B
Sartoguard GF 0.1 µm: 5481358M4--**--B
Sartoguard NF 0.1 µm: 5461358M4--**--B

** : available with SS, SO, OO connection

Cartridges for process scale

Sartopore® 2 XLM: 5442558M#
Sartoguard PES 0.1 µm: 5472558G#
Sartoguard GF 0.1 µm: 5482558M#
Sartoguard NF 0.1 µm: 5462558M#

: available in 10" (1), 20" (2), 30" (3) size

For other formats, please contact your local sales representative.

Contact & Information

Sartorius Stedim Biotech GmbH
August-Spindler-Straße 11
37079 Göttingen, Germany
Phone +49.551.308.2281
Email AS_Filtration@sartorius.com

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1.800.368.8200
Email techsupport@sartorius-stedim.com

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

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

Mycoplasma Removal Kit

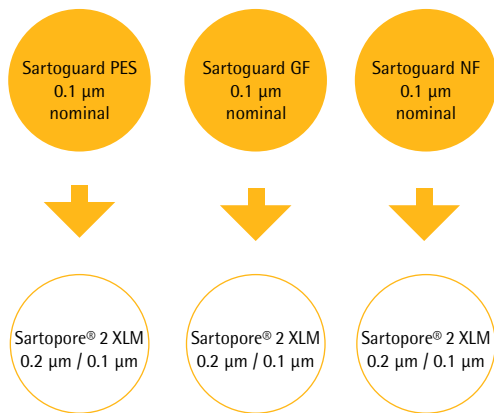
Find the ultimate choice for secure and cost effective Mycoplasma removal



Specifications subject to change without notice. Printed in Germany on paper that has been bleached without any use of chlorine.
Publication No.: SPKI1004-r166601 - Order No.: 85037-555-08 - Ver. 06 | 2016 - KT

How to use this kit?

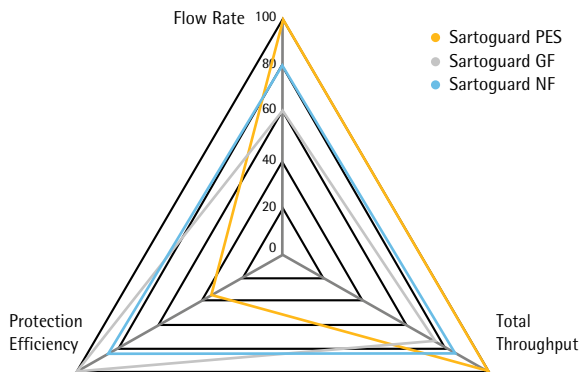
Check out the different combinations of Sartoguard prefilters and Sartopore® 2 XLM 0.1 µm sterile filter for your media:



For more information please visit www.sartorius-stedim.com

Sartoguard Family

- Optimal protection of Sartopore® 2 XLM
- 3–4 times higher capacity of Sartopore® 2 XLM
- Flexible adaptation with different filter materials for your individual media:
 - Sartoguard PES
Optimized for highest flow rate and total throughput
 - Sartoguard GF
Optimized for maximum turbidity reduction and protection efficiency
 - Sartoguard NF
Best combination of flow rate, total throughput and clarification



Sartopore® 2 XLM

- High security due to LRV of 7/cm² filtration area (TR >11/cartridge) for Mycoplasma
- Perfect cell growth and productivity due to minimized adsorption of growth factors
- Highest total throughput performance reduces filtration costs

