



SartoScale

Filter Test Disposables for Use in the Biopharmaceutical Industry



Specifications

Materials

Capsule Housing:	Polypropylene
Filter Materials:	All common filter materials of Sartorius Stedim Biotech GmbH

Available Sizes | Filtration Area

Size S 17.3 cm²

Available Connectors Styles

FF, HH

F: TC Flange 25 mm (1/2")

H: 1/4" multiple stepped hose barb

Operating Parameters

Max. allowable	5 bar 72.5 psi at 20°C
Differential pressure:	2 bar 29 psi at 80°C

SartoScale filter test disposables cannot be used in reverse direction of filtration!

Description

SartoScale filter test disposables are designed to perform reliable filterability trials with 47 mm flat filter discs of original filter cartridge material. The use of disposables for filtration trials avoids time consuming preparation of filter discs in stainless steel filter holders and prevents installation mistakes of the flat filter discs.

Applications

SartoScale filter test disposables are ideally suited to perform all kind of filterability trials with the target to select the optimal membrane material for a certain application or to determine the ideal combination of prefilters and final filters with minimum product volumes.

Original Filter Material

SartoScale filter test disposables contain the original filter active material of the respective filter cartridges in order to assure reproducible test results.

Scale-up

After material selection or determination of a prefilter | final filter scheme with SartoScale filter test disposables a scale-up for flow rate and total throughput performance of the selected materials should be done using small scale pleated capsule devices (e. g. capsules of type 4).

Optimized Design

SartoScale filter test disposables feature ultra low hold up and dead volumes in order to perform filterability trials with minimized product volumes.

Reliability

SartoScale filter test disposables containing integrity testable membrane filters can be tested for integrity by a bubble-point test to assure reliable test results.

Zero-T-Test System

We recommend to use SartoScale filter test disposables together with our Zero-T Filter Test System in order to perform filtration trials effectively. The Zero-T-System consists of hardware and software modules which allow easy handling and installation of the SartoScale filter test disposables. Automatic data acquisition is achieved by the connection of a balance to a laptop. The software analyses automatically the incoming data for membrane screening and scale-up.

Availability

SartoScale filter test disposables will become available for all filter materials of Sartorius Stedim Biotech GmbH including:

- Sartopore[®] 2 544...
- Sartopore[®] 2 XLG 544...
- Sartopore[®] 2 XLI 544...
- Sartopore[®] 2 XLM 544...
- Sartopore[®] Platinum 549...
- Sartobran[®] P 523..
- Sartolon[®] 510...
- Sartoclean[®] CA 562...
- Sartoclean[®] GF 560...
- Sartopure[®] PP2 559...
- Sartopure[®] PP3 505...
- Sartopure[®] GF Plus 555...
- Sartoguard PES 547...
- Sartoguard GF 548...
- Sartoguard NF 546...

Specifications

Extractables

SartoScale filter test disposables meet or exceed the requirements for WFI quality standards set by the current USP.

Regulatory Compliance

All sterilizing grade and mycoplasma retentive SartoScale filter test disposables are randomly tested for integrity during production.

Non pyrogenic according to USP Bacterial Endotoxins

Pass USP Plastic Class VI Test

Non fiber releasing according to 21 CFR

Sterilization

Autoclaving:

134°C, 2 bar, 30 min

No In-Line Steam Sterilization

Sterilization Cycles

Autoclaving: 1 Cycle

Integrity Test Parameters

(Water wetted)

Filter type	Pore Size	Bubble-Point [bar psi]
Sartopore® 2	0.2 µm	3.2 46
Sartopore® 2	0.45 µm	2.2 32
Sartopore® 2 XLG	0.2 µm	3.2 46
Sartopore® 2 XLI	0.2 µm	3.2 46
Sartopore® Platinum	0.2 µm	3.5 50.7
Sartobran® P	0.1 µm	3.8 55
Sartobran® P	0.2 µm	3.2 46
Sartobran® P	0.45 µm	2.0 29
Sartolon®	0.2 µm	3.0 43.5
Sartoguard PES	0.1 µm nom.	2.8 40.5
Sartoguard PES	0.2 µm nom.	1.8 26
Sartoguard GF	0.1 µm nom.	2.5 36.2
Sartoguard GF	0.2 µm nom.	2.0 29

Pack Size

M = 3 Pieces per Pack

Order Information

Sartopore® 2 0.1 µm	5445358KS--**--M
Sartopore® 2 0.2 µm	5445307HS--**--M
Sartopore® 2 0.45 µm	5445306GS--**--M
Sartopore® 2 XLG 0.2 µm	5445307GS--**--M
Sartopore® 2 XLI 0.2 µm	5445307IS--**--M
Sartopore® 2 XLM 0.1 µm	5445358MS--**--M
Sartopore® Platinum 0.2 µm	5495307HS--**--M
Sartobran® P 0.1 µm	5235358HS--**--M
Sartobran® P 0.2 µm	5235307HS--**--M
Sartobran® P 0.45 µm	5235306DS--**--M
Sartolon® 0.2 µm	5105307HS--**--M
Sartoguard PES 0.1 µm nom.	5475358GS--**--M
Sartoguard PES 0.2 µm nom.	5475307FS--**--M
Sartoguard GF 0.1 µm nom.	5485358MS--**--M
Sartoguard GF 0.2 µm nom.	5485307GS--**--M
Sartoguard NF 0.1 µm nom.	5465358MS--**--M
Sartoguard NF 0.2 µm nom.	5465307GS--**--M
Sartoclean® CA 0.2 µm	5625307AS--**--M
Sartoclean® CA 0.45 µm	5625306AS--**--M
Sartoclean® CA 0.65 µm	5625305GS--**--M
Sartoclean® CA 0.8 µm	5625304ES--**--M
Sartoclean® GF 0.65 µm	5605305GS--**--M
Sartoclean® GF 0.8 µm	5605304ES--**--M
Sartopure® PP2 0.65 µm	5595305PS--**--M
Sartopure® PP2 1.2 µm	5595303PS--**--M
Sartopure® PP2 3 µm	5595302PS--**--M
Sartopure® PP2 5 µm	5595342PS--**--M
Sartopure® PP2 8 µm	5595301PS--**--M
Sartopure® PP2 20 µm	5595320PS--**--M
Sartopure® PP2 50 µm	5595350PS--**--M
Sartopure® PP3 0.45 µm	5055306PS--**--M
Sartopure® PP3 0.65 µm	5055305PS--**--M
Sartopure® PP3 1.2 µm	5055303PS--**--M
Sartopure® PP3 3 µm	5055302PS--**--M
Sartopure® PP3 5 µm	5055342PS--**--M
Sartopure® PP3 8 µm	5055301PS--**--M
Sartopure® PP3 20 µm	5055320PS--**--M
Sartopure® PP3 50 µm	5055350PS--**--M
Sartopure® PP3 100 µm	5055398PS--**--M
Sartopure® GF Plus 0.65 µm	5555305PS--**--M
Sartopure® GF Plus 1.2 µm	5555303PS--**--M

** : Connector Styles

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

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.90.2110935
Russian Federation +7.812.327.5.327
Japan +81.3.4331.4300

Specifications subject to change without notice. Printed and copyrighted by Sartorius Stedim Biotech GmbH. | W
Publication No.: SPK2069-e140203
Order No.: 85030-523-14
Ver. 02 | 2014