

Sartoscale

Single-Use Screening Tool for Process Development & Optimization



Minimal Product Use for Testing

Sartoscales are available in two sizes: Sartoscale 25 with 4.5 cm² effective filtration area (EFA) and Sartoscale 47 that offers an EFA of 17.3 cm². These compact device constructions offer extremely low hold-up volumes (Sartoscale 25: ≤ 1 mL) and enable filtration trials with minimal product volume and high product recovery.

Consistency of Results

Sartoscale single-use filters are the ideal tool for screening of filter materials in biopharmaceutical development and process optimization. They contain the identical filter active materials as well as the support fleeces as their process filter counterparts to assure data consistency for scale-up. In addition, the use of identical filter materials allows to carry out adsorption studies with formulation components or high-value product streams at a small scale.

Ease of Use

Sartoscale units are single-use units delivered ready-to-use. This facilitates trial set-up to speed up biopharmaceutical process development. Sartoscale 25 feature a female luer lock inlet connection and venting port as well as a male luer slip outlet connection for easy integration into every testing setup. Inlet, outlet and venting port are closed with caps to protect the units before use. Sartoscale 47 feature tri-clamp or hose barb connectors on inlet and outlet as well as an upstream venting port.

Secure Identification & Traceability

Each Sartoscale unit is laser-labeled with the product name, pore size or retention rate, lot number and piece number assuring reliable filter identification and complete traceability. In addition the Sartoscale 25 carries a 2-d barcode for automatic data transfer.

Technical Data

Filter Families	Filter Material	Min. Bubble Point [bar psi]
Sartopore® Platinum 0.2 µm	Polyethersulfone, surface modified	3.5 50.7
Sartopore® Platinum HB 0.2 µm	Polyethersulfone, surface modified	4.1 59.5
Sartopore® 2 0.45 µm	Polyethersulfone	2.2 32
Sartopore® 2 0.2 µm		3.2 46
Sartopore® 2 XLG		3.2 46
Sartopore® 2 XLI		3.2 46
Sartopore® 2 XLM		n.a.
Sartobran® P 0.45 µm	Cellulose Acetate	2.0 29
Sartobran® P 0.2 µm		3.2 46
Sartolon 0.2 µm	Polyamide	3.0 43.5
Sartoguard PES 0.1 µm nom.	Polyethersulfone	2.8 40.5
Sartoguard PES 0.2 µm nom.		1.8 26
Sartoguard GF 0.1 µm nom.	Polyethersulfone + Glass Fiber	n.a.
Sartoguard GF 0.2 µm nom.		2.3 33
Sartoclean® CA	Cellulose Acetate	n.a.
Sartoclean® GF	Cellulose Acetate + Glass Fiber	n.a.
Sartopure® PP3	Polypropylene	n.a.
Sartopure® GF Plus	Glass Fiber	n.a.

Materials

Filter Material

Depending on product family (see table)

Support Fleece

Polypropylene

Capsule Housing

Polypropylene

Max. Allowable Pressure

5 bar | 72.5 psi at 20°C

2 bar | 29 psi at 80°C

Sartoscale filter devices cannot be used in reverse direction of filtration!

Regulatory Compliance

- Non-pyrogenic according to USP Bacterial Endotoxins
- Pass USP Plastic Class VI Test
- Non-fiber releasing according to 21 CFR



Available Sizes | Filtration Area

Sartoscale 25 | 4.5 cm²
Sartoscale 47 | 17.3 cm²

Sterilization

Autoclaving

Max. 134 °C, 2 bar, 30 min
Min. 1 Sterilization Cycle

Ordering Information

Sartoscale



	53		V	--	LX	--	C
Filter Family	Pore Size [µm]	Filter Size	Connectors Sartoscale 25	Special Type -- : Standard version	Pack Size Sartoscale 25		
549: Sartopore® Platinum	07H: 0.45 + 0.2	V: 4.5 cm² S: 17.3 cm²	L: Female Luer lock (inlet) X: Male Luer slip (outlet)	HB: High Bubble Point	C: box of 6 N: box of 100		
544: Sartopore® 2 Sartopore® 2 Sartopore® 2 XLG Sartopore® 2 XLI Sartopore® 2 XLM	06G: 0.8 + 0.45 07H: 0.45 + 0.2 07G: 0.8 + 0.2 07I: 0.35 + 0.2 58M: 0.2 + 0.1		Connectors Sartoscale 47 FF: tri-clamp 25 mm (inlet & outlet) HH: ¼" multiple stepped hose barb (inlet & outlet)		Pack Size Sartoscale 47 M: box of 3		
523: Sartobran® P	06D: 0.65 + 0.45 07H: 0.45 + 0.2						
510: Sartolon	07H: 0.45 + 0.2						
547: Sartoguard PES	58G: 0.1 nom. 07F: 0.2 nom.						
548: Sartoguard GF	58M: 0.1 nom. 07G: 0.2 nom.						
546: Sartoguard NF	58M: 0.1 nom. 07G: 0.2 nom.						
562: Sartoclean® CA	04E: 3.0 + 0.8 05G: 0.8 + 0.65						
560: Sartoclean® GF	04E: 3.0 + 0.8 05G: 0.8 + 0.65						
	Retention Rate [µm]						
505: Sartopure® PP3	06P: 0.45 05P: 0.65 03P: 1.2 02P: 3.0 42P: 5.0 01P: 8.0 20P: 20.0 50P: 50.0 98P: 100.0						
555: Sartopure® GF Plus	05P: 0.65 03P: 1.2						

Example

Sartopore® Platinum: 5495307HV-LX-C

Check the availability of your desired configuration with your Sartorius sales representative or online at www.sartorius.com

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 further contacts, visit
www.sartorius.com