

SartoScale 25

Single-Use Screening Tool for Process Development & Optimization



Minimized Product Volume for Testing

SartoScale 25 mm units feature an optimized design with 4.5 cm² effective filtration area. In combination with the extremely low hold-up volume of less than 1 ml SartoScales 25 units allow to carry out filtration trials with minimal product volume and high product recovery.

Consistency of Results

SartoScale 25 units contain the identical filter active materials as well as the support fleeces compared to process scale elements, assuring data consistency. The use of identical filter materials allows in addition to carry out adsorption studies of important formulation components at small scale.

Product Information

SartoScale 25 single-use filters are the ideal tool for screening of filter materials in bio-pharmaceutical development and process optimization. Their unique design and small surface area permits to carry out trials with lowest volumes of expensive fluid streams. Consistency of results is assured by using the identical filter materials including supportive fleeces as process scale elements.

Ease of Use

SartoScale 25 units are single use units delivered ready to use. This allows to set-up trials fast and easy to speed up bio-pharmaceutical process development. They 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.

Secure Identification & Traceability

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

Technical Data

Filter Families	Materials	Min. Bubble Point [bar psi]
Sartopore® Platinum	Polyethersulfone, surface modified	3.5 50.7
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	Nylon	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
Sartoguard NF	Polyethersulfone + PES Nanofleece	n.a.
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

Support Fleece

Polypropylene

Capsule Housing

Polypropylene

Regulatory Compliance

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

Max. Allowable Pressure

5 bar | 72.5 psi at 20°C

2 bar | 29 psi at 80°C

SartoScale 25 filter devices cannot be used in reverse direction of filtration!

Sterilization

Filter Area

4.5 cm² for all types

Autoclaving

134°C, 2 bar, 30 min
1 Sterilization Cycle



Ordering Information



SartoScale 25

	53		V	--	LX	--	C
Filter Family		Pore Size [μm]	Filter Size		Connector		Pack Size
549: Sartopore® Platinum		07H: 0.45 + 0.2	V: 4.5 cm ²		L: Female Luer Lock X: Male Luer Slip		C: box of 6 N: box of 100
544: Sartopore® 2		06G: 0.8 + 0.45					
Sartopore® 2		07H: 0.45 + 0.2					
Sartopore® 2 XLG		07G: 0.8 + 0.2					
Sartopore® 2 XLI		07I: 0.35 + 0.2					
Sartopore® 2 XLM		58M: 0.2 + 0.1					
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					
		06A: 0.45					
		07A: 0.2					
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

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