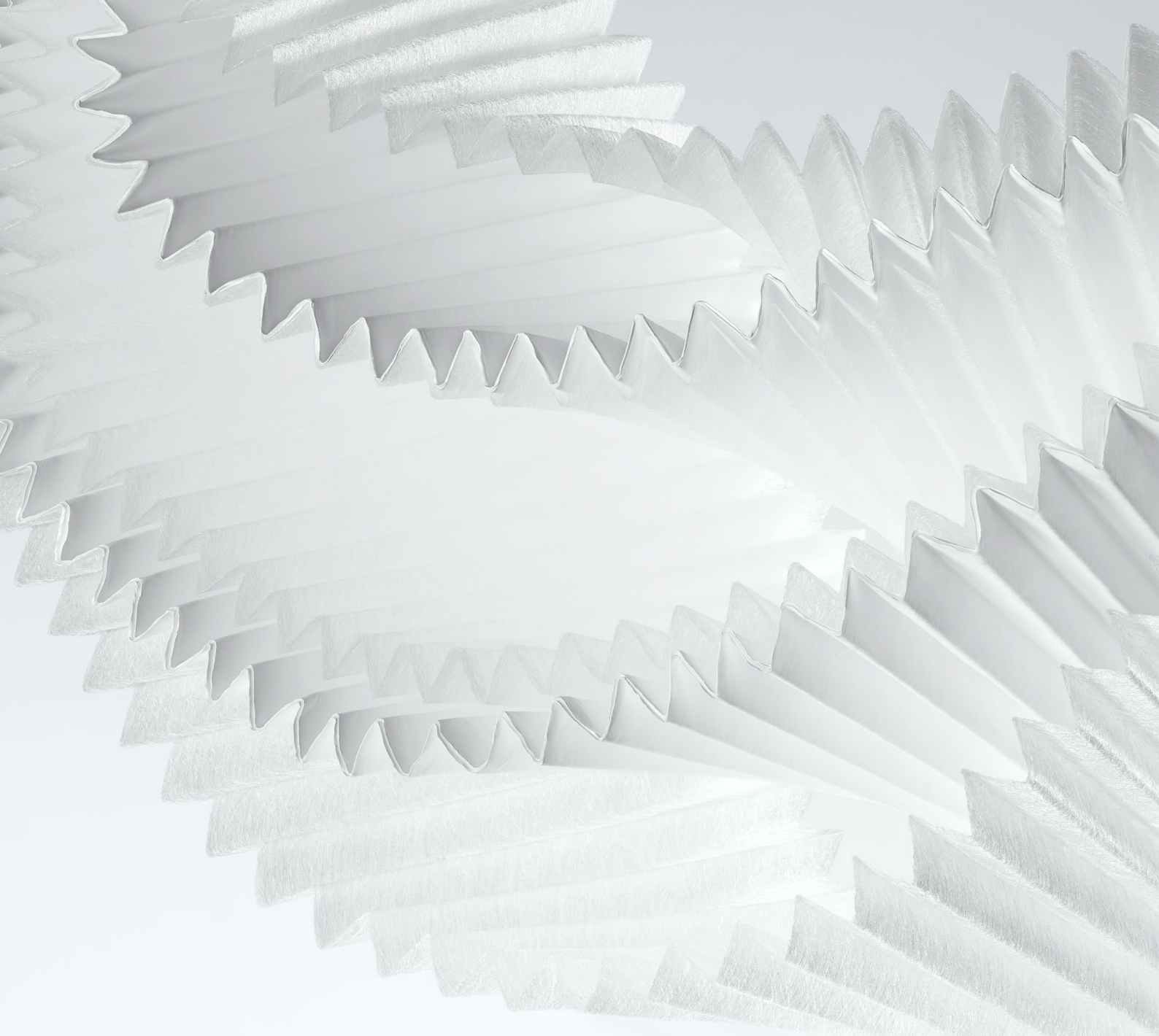




Filter Portfolio Overview

Simplifying Progress

SARTORIUS

Table of Contents

Application Filter Matrix	4
Our Portfolio	10
Flexibility in Design	18
How to Order	21
Sartopore® Family	22
Sartoguard Family	30
Sartobran® P Family	32
Sartoclean® Family	33
Sartolon Family	35
Sartopure® Family	36
Sartofluor® Family	40
Midisart®	41
Virosart®	42
Filter Transfer Sets	46
Maxicaps® MR	48





Application | Filter Matrix

This matrix provides a guideline for selection of the right filter in a given application backed up by decades of experience in the biopharmaceutical industry. However, it is recommended carrying out small scale filtration trials to identify the optimal filter combination based on the actual product and process conditions.

Biotech

		Particle Reduction			Prefilters Bioburden Reduction			
		Sartopure®		Sartoclear®	Sartoguard		Sartoclean®	
		PP3	GF Plus		PES	GF	GF	CA
		PP	GF		PES	PES + GF	CA + GF	CA
Unit Operation	Application							
Media Filtration	Serum-containing		■		■	■		
	Serum-free				■			
	Protein-free							
	Chemically defined							
Buffer & Auxiliaries Filtration	pH adjustment for Cell Culture Fermentation	■						
	Buffer for Cell Removal & Downstream	■			■			
	ADC Solvent Filtration							
	Formulation Buffer							
	Adjuvants				■			
Harvest Cell Removal	mAbs, rec. Proteins			■		■		
	Viral Vaccines	■	■			■	■	
	Viral Vectors	■	■			■	■	■
	Plasmid DNA	■	■	■		■		
Downstream Intermediates	mAbs, rec. Proteins				■			
	Viral Vaccines				■		■	■
	Viral Vectors				■			■
	Plasmid DNA				■			■
	mRNA				■			
Filtration after Virus Inactivation	mAbs, rec. Proteins	■	■			■		
Form & Fill	mAbs, rec. Proteins							
	Viral Vaccines							■
	Viral Vectors							■
	Plasmid DNA				■			
	Lipid Nano Particle				■			

■ recommended
 ■ alternatively recommended

Mycoplasma Retentive Filters		Sterile Filters							
Sartopore 2®		Sartopore 2			Sartopore Evo®	Sartopore® Platinum		Sartobran® P	Sartolon
XLM		0.2 µm	XLG	XLI	0.2 µm	0.2 µm	HB	0.2 µm	
PES		PES	PES	PES	mod. PES	mod. PES	mod. PES	CA	PA
	■		■						
	■		■						
	■		■						
	■		■	■					
		■	■						
		■	■						
									■
		■	■						
			■			■	■		
								■	
								■	
			■		■				
			■					■	
			■					■	
			■						
			■					■	
					■	■	■		
				■	■	■		■	
					■	■	■	■	
					■	■	■		
					■	■	■		

Application | Filter Matrix

Pharma

		Particle Reduction			Prefilters Bioburden Reduction			
		Sartopure®		Sartoclear®	Sartoguard		Sartoclean®	
		PP3	GF Plus		PES	GF	GF	CA
		PP	GF		PES	PES + GF	CA + GF	CA
Material ▶								
Classical Pharma								
SVP		■	■		■			
LVP		■			■			
API	Antibiotics (solvents)	■						
	Water-based	■			■			
Opthalmics		■	■		■			
Blood & Plasma								
Albumin & Globulines	Intermediate Process Filtration		■	■		■		
	Filtration after Virus Inactivation	■	■			■		
	Form & Fill				■			
Clotting Factors	Intermediate Process Filtration						■	■
	Form & Fill							■
Others								
	Water	■						
	Oily formulations	■						
	Solvents	■						

Professional Support When You Need It

Make the choice that helps you bring your products to market rapidly.

Our global expert teams are available to support you with solutions for process design, application development as well as product and process specific Validation Services including Extractables| Leachables studies, physicochemical and microbiological testing.

Visit us at: www.sartorius.com/validation-service

■ recommended
 ■ alternatively recommended

Mycoplasma Retentive Filters		Sterile Filters								
Sartopore® 2		Sartopore® 2			Sartopore Evo®	Sartopore® Platinum		Sartobran® P	Sartolon	Sartofluor®
XLM	0.2 µm	XLG	XLI	0.2 µm	0.2 µm	HB	0.2 µm			LG
PES	PES	PES	PES	mod. PES	mod. PES	mod. PES	CA	PA		PTFE
	■	■			■					
	■	■								
								■		■
	■							■		
	■		■		■		■			
		■		■						
		■		■						
		■		■		■	■			
		■								
		■					■			
	■	■								
		■						■		
								■		■

Application | Filter Matrix

Air & Gas Filtration

		Midisart®		Sartopore® Air
				2000
		PES	PTFE	PES
Material ▶				
Compressed Gas				
Venting of tanks	Pressure resistant			
	Non-pressure resistant			
Glas & plastic bottles			■	
Venting, inflating & leak testing of SU bags assemblies		■		
Venting of WFI loops (high temperature)				
Bioreactors Fermentors	Single-use	■		■
	Stainless steel		■	
Protection of integrity test devices				■
Autoclave				
Lyophilization				
PUPSIT operation	Single-use			■
	Stainless steel			

Virus Retention

		Prefilters		Virus Retentive Filters
		Virosart® Max	Sartopore® 2 XLM	Virosart® HF
		Polyamide	PES	mod. PES
Material ▶				
Biotech				
mAb, rec. Proteins	Virus Retention DSP	■		■
	Virus Retention for Media Filtration		■	
Viral Vaccines, Viral Vectors and Cell Therapy	Virus Retention for Media Filtration		■	
Blood & Plasma				
Plasma Derivates	Virus Retention DSP	■		

■ recommended
 ■ alternatively recommended

Sartopure®	Sartofluor®	
GA ¹	GA	HR
Hydrophob. GF	PTFE	PTFE
	■	
	■	
■		
	■	
		■
	■	
	■	
	■	■
	■	
	■	

¹ Non-sterilizing grade

Virosart® CPV	Virosart® Media	Virosart® HC
PES	PES	mod. PES
■		
	■	
	■	
		■

Our Portfolio

Sartopore Evo®

Sterilizing Grade 0.8 | 0.2 µm

Filter Material		Filter Designs	Sterilization Methods
Prefilter Membrane:	modified Polyethersulfone	Mini Cartridges	In-line steamable or autoclavable
Final Membrane:	modified Polyethersulfone	Cartridges	In-line steamable or autoclavable [S]
		T-Style Maxicaps®	Autoclavable or γ-irradiatable (hybrid version) [S]
		Maxicaps®	Autoclavable or γ-irradiatable (hybrid version) [S]
		Midicaps®	Autoclavable or γ-irradiatable (hybrid version) [S]
		Capsules	Autoclavable or γ-irradiatable (hybrid version) [S]
		Sartoscales	Autoclavable

Sartopore® Platinum | Platinum HB

Sterilizing Grade 0.45 | 0.2 µm

Filter Material		Filter Designs	Sterilization Methods
Prefilter Membrane:	modified Polyethersulfone	Cartridge	In-line steamable or autoclavable
Final Membrane:	modified Polyethersulfone	T-Style Maxicaps®	Autoclavable or γ-irradiatable (hybrid version) [S]
		Maxicaps®	Autoclavable or γ-irradiatable (hybrid version) [S]
		Midicaps®	Autoclavable [S]
			γ-irradiatable
		Capsules	Autoclavable [S]
			γ-irradiatable
		Sartoscales	Autoclavable

Sartopore® 2

Bioburden Reduction 0.8 | 0.45 µm
Sterilizing Grade 0.45 | 0.2 µm

Filter Material		Filter Designs	Sterilization Methods
Prefilter Membrane:	Polyethersulfone	Mini Cartridges	In-line steamable or autoclavable
Final Membrane:	Polyethersulfone	Cartridges	In-line steamable or autoclavable
		T-Style Maxicaps®	Autoclavable or γ-irradiatable (hybrid version)
		Maxicaps®	Autoclavable
			γ-irradiatable
		Midicaps®	Autoclavable [S]
			γ-irradiatable
		Capsules	Autoclavable [S]
			γ-irradiatable
		Sartoscales	Autoclavable

Sartopore® 2 XLG

Sterilizing Grade 0.8 | 0.2 µm

Filter Material		Filter Designs	Sterilization Methods
Prefilter Membrane:	Polyethersulfone	Cartridges	In-line steamable or autoclavable
Final Membrane:	Polyethersulfone	T-Style Maxicaps®	Autoclavable or γ-irradiatable (hybrid version)
		Maxicaps®	Autoclavable γ-irradiatable
		Midicaps®	Autoclavable [S] γ-irradiatable
		Capsules	Autoclavable [S] γ-irradiatable
		Sartoscales	Autoclavable

Sartopore® 2 XLI

Sterilizing Grade 0.35 | 0.2 µm

Filter Material		Filter Designs	Sterilization Methods
Prefilter Membrane:	Polyethersulfone	Cartridges	In-line steamable or autoclavable
Final Membrane:	Polyethersulfone	T-Style Maxicaps®	Autoclavable or γ-irradiatable (hybrid version)
		Maxicaps®	Autoclavable γ-irradiatable
		Midicaps®	Autoclavable [S] γ-irradiatable
		Capsules	Autoclavable [S] γ-irradiatable
		Sartoscales	Autoclavable

Sartopore® 2 XLM

Sterilizing Grade |
Mycoplasma Retention 0.2 | 0.1 µm

Filter Material		Filter Designs	Sterilization Methods
Prefilter Membrane:	Polyethersulfone	Cartridges	In-line steamable or autoclavable
Final Membrane:	Polyethersulfone	T-Style Maxicaps®	Autoclavable or γ-irradiatable (hybrid version)
		Maxicaps®	Autoclavable γ-irradiatable
		Midicaps®	Autoclavable [S] γ-irradiatable
		Capsules	Autoclavable [S] γ-irradiatable
		Sartoscales	Autoclavable

[S] delivered in sterile condition (autoclaved)

Sartoguard PES

Bioburden Reduction | 0.8 | 0.1 µm (nom.)
Particle Retention 1.2 | 0.2 µm (nom.)

Filter Material	Filter Designs	Sterilization Methods
Prefilter Membrane: Polyethersulfone	Cartridges	In-line steamable or autoclavable
Final Membrane: Polyethersulfone	T-Style Maxicaps®	Autoclavable or γ-irradiatable (hybrid version)
	Maxicaps®	Autoclavable
		γ-irradiatable
	Midicaps®	Autoclavable [S]
		γ-irradiatable
	Capsules	Autoclavable [S]
		γ-irradiatable
	Sartoscales	Autoclavable

Sartoguard GF

Bioburden Reduction | 0.2 | 0.1 µm (nom.)
Particle Retention 0.8 | 0.2 µm (nom.)

Filter Material	Filter Designs	Sterilization Methods
Filter Active Fleece: Glass Fiber	Cartridges	In-line steamable or autoclavable
Prefilter Membrane: Polyethersulfone	T-Style Maxicaps®	Autoclavable
Final Membrane: Polyethersulfone	Maxicaps®	Autoclavable
	Midicaps®	Autoclavable
	Capsules	Autoclavable
	Sartoscales	Autoclavable

Sartobran® P

Bioburden Reduction 0.65 | 0.45 µm
Sterilizing Grade 0.45 | 0.2 µm

Filter Material	Filter Designs	Sterilization Methods
Prefilter Membrane: Cellulose Acetate	Mini Cartridges	In-line steamable or autoclavable
Final Membrane: Cellulose Acetate	Cartridges	In-line steamable or autoclavable
	T-Style Maxicaps®	Autoclavable [S]
	Maxicaps®	Autoclavable [S]
	Midicaps®	Autoclavable [S]
	Capsules	Autoclavable [S]
	Sartoscales	Autoclavable

[S] delivered in sterile condition (autoclaved)

Sartoclean® CA

Particle Retention 3 | 0.8 µm (nom.)
0.8 | 0.65 µm (nom.)

Filter Material	
Prefilter Membrane:	Cellulose Acetate
Final Membrane:	Cellulose Acetate

Filter Designs	Sterilization Methods
Mini Cartridges	In-line steamable or autoclavable
Cartridges	In-line steamable or autoclavable
T-Style Maxicaps®	Autoclavable S
Maxicaps®	Autoclavable S
Midicaps®	Autoclavable S
Sartoscales	Autoclavable

Sartoclean® GF

Particle Retention 3 | 0.8 µm (nom.)
0.8 | 0.65 µm (nom.)

Filter Material	
Filter Active Fleece:	Glass Fiber
Prefilter Membrane:	Cellulose Acetate
Final Membrane:	Cellulose Acetate

Filter Designs	Sterilization Methods
Mini Cartridges	In-line steamable or autoclavable
Cartridges	In-line steamable or autoclavable
T-Style Maxicaps®	Autoclavable S
Maxicaps®	Autoclavable S
Midicaps®	Autoclavable S
Sartoscales	Autoclavable

Sartolon

Sterilizing Grade 0.45 | 0.2 µm

Filter Material	
Prefilter Membrane:	Polyamide (Nylon)
Final Membrane:	Polyamide (Nylon)

Filter Designs	Sterilization Methods
Mini Cartridges	In-line steamable or autoclavable
Cartridges	In-line steamable or autoclavable
Maxicaps®	Autoclavable S
Midicaps®	Autoclavable S
Sartoscales	Autoclavable

Sartopure® PP3

Particle Retention ¹ (for all)	50 µm	3 µm
	20 µm	1.2 µm
	8 µm	0.65 µm
	5 µm	0.45 µm

Filter Material		Filter Designs	Sterilization Methods
Filter Active Fleece:	Polypropylene	Mini Cartridges	In-line steamable or autoclavable
		Cartridges	In-line steamable or autoclavable
		T-Style Maxicaps®	Autoclavable
		Maxicaps®	Autoclavable
		Midicaps®	Autoclavable
		Capsules	Autoclavable
		Sartoscales	Autoclavable

Sartopure® GF Plus

Particle Retention ¹ (for all)	1.2 µm
	0.65 µm

Filter Material		Filter Designs	Sterilization Methods
Filter Active Fleece:	Glass Fiber	Mini Cartridges	In-line steamable or autoclavable
		Cartridges	In-line steamable or autoclavable
		Maxicaps®	Autoclavable
		Midicaps®	Autoclavable
		Sartoscales	Autoclavable

¹ nom.



Sartopore® Air

Sterilizing Grade 0.2 µm

Filter Material		Filter Designs	Sterilization Methods
Membrane:	hydrophobic Polyethersulfone	Maxicaps®	γ-irradiatable
		Midicaps®	γ-irradiatable
		Capsules	γ-irradiatable

Sartopure® GA

Bioburden Reduction | Particle Retention 0.2 µm (nom.)

Filter Material		Filter Designs	Sterilization Methods
Filter Active Fleece:	hydrophobic Glass Fiber	Mini Cartridges	In-line steamable or autoclavable
		Cartridges	In-line steamable or autoclavable

Sartofluor® GA | LG

Sterilizing Grade 0.2 µm (GA)
Sterilizing Grade 0.2 µm (LG)

Filter Material		Filter Designs	Sterilization Methods
Membrane:	PTFE	Mini Cartridges	In-line steamable or autoclavable
		Junior	In-line steamable or autoclavable
		Cartridges	In-line steamable or autoclavable
		Maxicaps®	Autoclavable S
		Midicaps®	Autoclavable S
		Capsules	Autoclavable S

Sartofluor® HR

Sterilizing Grade 0.2 µm

Filter Material		Filter Designs	Sterilization Methods
Membrane:	PTFE	Cartridges	In-line steamable or autoclavable

S delivered in sterile condition (autoclaved)

Virosart® HF

Virus Retention 20 nm nom.

Filter Material		Filter Designs	Sterilization Methods
Membrane:	Modified Polyethersulfone	Process modules	γ-irradiation
Format:	Hollow fiber	Mid-scale modules	γ-irradiation
		Lab modules	γ-irradiation [S]

Virosart® Media

Virus Retention 20 nm nom.

Filter Material		Filter Designs	Sterilization Methods
Membrane:	Polyethersulfone	Process modules	γ-irradiation
Format:	Hollow fiber	Mid-scale modules	γ-irradiation
		Lab modules	γ-irradiation [S]

Virosart® CPV

Virus Retention 20 nm nom.

Filter Material		Filter Designs	Sterilization Methods
Membrane:	Polyethersulfone	Cartridges	In-line steamable or autoclavable
		Maxicaps®	Autoclavable
		Midicaps®	Autoclavable
		Capsules Size 4	Autoclavable

Virosart® HC

Virus Retention 20 nm nom.

Filter Material		Filter Designs	Sterilization Methods
Membrane:	Modified Polyethersulfone	Cartridges	In-line steamable or autoclavable
		T-Style Maxicaps®	Autoclavable
		Midicaps®	Autoclavable
		Capsules Size 4	Autoclavable

Virosart® Max

Retention Rate 0.1 μm

Filter Material		Filter Designs	Sterilization Methods
Membrane:	Polyamide	Cartridges	In-line steamable or autoclavable
		T-Style Maxicaps®	Autoclavable [S] *
		Midicaps®	Autoclavable [S] *
		Capsules Size 4	Autoclavable [S] *

[S] delivered in sterile condition (γ-irradiated)

* [S] delivered in sterile condition (autoclaved)



SARTORIUS
VIROSART® MAX
V-STYLE MAXICAPS®
Platform for VIROSART® CPV/HC

Model: ViroSart-MAX
Max. Working Volume: 100 L
Max. Operating Pressure: 1.5 bar
Max. Operating Temperature: 37°C
Max. Stirring Speed: 120 rpm
Max. Stirring Power: 1.5 kW
Max. Stirring Torque: 15 Nm
Max. Stirring Time: 120 min
Max. Stirring Frequency: 120 Hz
Max. Stirring Amplitude: 120 mm
Max. Stirring Diameter: 120 mm
Max. Stirring Length: 120 mm
Max. Stirring Width: 120 mm
Max. Stirring Height: 120 mm
Max. Stirring Depth: 120 mm
Max. Stirring Area: 120 m²
Max. Stirring Volume: 120 L
Max. Stirring Weight: 120 kg
Max. Stirring Length: 120 m
Max. Stirring Width: 120 m
Max. Stirring Height: 120 m
Max. Stirring Depth: 120 m
Max. Stirring Area: 120 m²
Max. Stirring Volume: 120 L
Max. Stirring Weight: 120 kg

Max. Working Volume: 100 L
Max. Operating Pressure: 1.5 bar
Max. Operating Temperature: 37°C
Max. Stirring Speed: 120 rpm
Max. Stirring Power: 1.5 kW
Max. Stirring Torque: 15 Nm
Max. Stirring Time: 120 min
Max. Stirring Frequency: 120 Hz
Max. Stirring Amplitude: 120 mm
Max. Stirring Diameter: 120 mm
Max. Stirring Length: 120 mm
Max. Stirring Width: 120 mm
Max. Stirring Height: 120 mm
Max. Stirring Depth: 120 mm
Max. Stirring Area: 120 m²
Max. Stirring Volume: 120 L
Max. Stirring Weight: 120 kg

Flexibility in Design

Capsules



Connector Type O
¼" multiple stepped hose barb



Connector Type S
¾" Tri-Clamp
25 mm



Hollow Fiber Lab Modules



Connector Inlet Type M
Male Luer Lock

Connector Outlet Type L
Female Luer Lock

Midicaps®



Connector Type S
1½" Tri-Clamp
50.5 mm



Connector Type O
½" single stepped hose barb



Connector Type F*
¾" Tri-Clamp
25 mm



Hollow Fiber Mid Scale Modules



Connector Inlet Type F
Sanitary ¾" Tri-Clamp
25 mm

Connector Outlet Type O
½" single stepped hose barb



T-Style Maxicaps®



Connector Type S
1½" Tri-Clamp
50.5 mm



Connector Type O
½" single stepped hose barb



Hollow Fiber Process Modules



Connector Inlet Type S
Sanitary 1 ½" Tri-Clamp
50.5 mm

Connector Outlet Type F
Sanitary ¾" Tri-Clamp
25 mm

Maxicaps®



Connector Type S
1½" Tri-Clamp
50.5 mm



Connector Type O
½" single stepped hose barb



Connector Type F*
¾" Tri-Clamp
25 mm



■ Multi-use

■ Single-use

* On request

Sartoscale 25



Connector Inlet Type L
Female Luer Lock for
venting | bleeding

Connector Outlet Type X
Male Luer Slip

Sartoscale 47



Connector Type F
¾" Tri-Clamp 25 mm

Mini Cartridges



Adapter 15
Bayonet adapter
with o-ring



Adapter 18 *
Plug adapter with
double o-ring



Cartridges



Adapter 25
2 Flange Bayonet adapter
with 226 double o-ring



Adapter 27 *
Adapter with
222 double o-ring



Adapter 28 *
3 Flange Bayonet adapter
with 222 double o-ring



Small Filters for Special Cases

Midisart®



**Standard Stepped
Hose Barb**
6 – 12 mm



**Small Stepped
Hose Barb**
4 – 12 mm



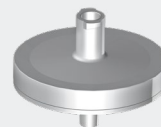
Thread
⅛" NPT



Tri-Clamp
¾" Tri-Clamp
25 mm



Minisart®



Connector Inlet
Female Luer Lock

Connector Outlet
Male Luer Slip

In this brochure small scale virus filters in the Minisart® format are described. Minisart® pre-filters, sterilizing grade filters and air filter can be found in the brochure "Minisart® Syringe Filter Family".



How to Order

Filter Design		Pore Size Prefilter		Sterilization		Special Additions	
83	T-Style Maxicaps®	A	No prefilter	(T-Style) Maxicaps®		P	Pharmaceutical grade
53	Midicaps®	P	Polypropylene layer	Midicaps® Capsules		GA	Gas Filtration
13	Capsule Size 4 & 5	T	Single layer PTFE membrane	G - γ-irradiatable hybrid*		LG	Liquid Filtration
73	Maxicaps®	E	3 μm	-- autoclavable		IT	Vent valve for connecting Sartocheck® (only Maxicaps®)
--	Hollow Fiber Modules	F	1.2 μm	Hollow Fiber Modules		HB	High Bubble Point
Cartridge Adapter Type		G	0.8 μm	C - prepared for γ-irradiation			
25 27 28		D	0.65 μm	CG γ-irradiated			
Mini Cartridge Adapter Type		H	0.45 μm				
15 18		I	0.35 μm				
14 (only Sartofluor® Junior)		M	0.2 μm				
		V	Virosart® CPV/HC				
		N	Virosart® Max				
549	73	07	H	1	G -	SS	IT
Filter Type		Pore Size Final Filter		Filter Size		Connectors	
505	Sartopure® PP3	50	= 50 μm	Cartridges (T-Style) Maxicaps®		Mini Standard Cartridges	
510	Sartolon	20	= 20 μm	0	5" (only cartridges)	--	Silicone O-Rings
517	Sartofluor® HR	01	= 8 μm	1	10"	- E	EPDM O-Rings
518	Sartofluor®	42	= 5 μm	2	20"	- V	Fluoroelastomer O-Rings
519	Sartopore® Air	02	= 3 μm	3	30"	(T-Style) Maxicaps® Midicaps® Hollow Fiber Process + mid scale Modules	
523	Sartobran® P	03	= 1.2 μm	4	40" (only cartridges)	S	Tri-Clamp 50.5 mm (1½")
539	Virosart® HC	04	= 0.8 μm	Capsules		O	½" single stepped hose barb
544	Sartopore® 2	05	= 0.65 μm	4		F	Tri-Clamp 25 mm (¾")
545	Virosart® CPV	06	= 0.45 μm	5		Capsules	
547	Sartoguard PES	07	= 0.2 μm	Mini Cartridges Midicaps®		O	¼" multiple stepped hose barb
548	Sartoguard GF	58	= 0.1 μm	7		S	Tri-Clamp 25 mm (¾")
549	Sartopore® Platinum	28	= 20 nm	8		Sartoscale 47	
555	Sartopure® GF Plus			9		F	Tri-Clamp 25mm (¾")
559	Sartopure® GA			Sartoscale		Sartoscale 25	
560	Sartoclean® GF			S	Sartoscale 47	L	Female Luer Lock
562	Sartoclean® CA			V	Sartoscale 25	X	Male Luer Slip
599	Sartopore Evo®			Hollow Fiber Modules		Hollow Fiber Lab Modules	
54A	Virosart® Max			T	1.7 cm²	L	Female Luer Lock
3VI	Virosart® HF			B	5 cm²	M	Male Luer Slip
3V2	Virosart® Media			D	200 cm²		
				F	0.2 m²		
				G	0.3 m²		
				H	0.8 m²		
				I	1 m²		
				M	2.4 m²		
						Note: the first code letter represents the inlet, the second the outlet.	

* hybrid = autoclavable or γ-irradiatable

Sartopore® Family

Sartopore® Evo

		Size 0 (5")	Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.4 m ² 4.3 ft ²	0.8 m ² 8.6 ft ²	1.6 m ² 17.2 ft ²	2.4 m ² 25.8 ft ²
Cartridges	0.8 + 0.2	5992507G0	5992507G1	5992507G2	5992507G3
T-Style Maxicaps® γ-irradiatable or autoclavable	0.8 + 0.2		5998307G1G-**-	5998307G2G-**-	5998307G3G-**-
Maxicaps® γ-irradiatable or autoclavable	0.8 + 0.2		5997307G1G-**-	5997307G2G-**-	5997307G3G-**-

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.065 m ² 0.7 ft ²	0.13 m ² 1.4 ft ²	0.26 m ² 2.8 ft ²	0.52 m ² 5.6 ft ²
Mini Cartridges	0.8 + 0.2	5991507G7-----B	5991507G8-----B	5991507G9-----B	
Midicaps® γ-irradiatable or autoclavable	0.8 + 0.2	5995307G7G-**-A	5995307G8G-**-A	5995307G9G-**-A	5995307G0G-**-V

		Size 4
	Pore Size	0.021 m ² 0.22 ft ²
Capsules® γ-irradiatable or autoclavable	0.8 + 0.2	5991307G4G-**-B

		25	47
	Pore Size	4.5 cm ²	17.3 cm ²
Sartoscale	0.8 + 0.2	5995307GV--LX--C	5995307GS--**-M

** = description of available filter design or connector | adapter types see page 18 | 19
Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

Sartopore® Platinum

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.6 m ² 6.5 ft ²	1.2 m ² 12.9 ft ²	1.8 m ² 19.4 ft ²
Cartridges standard pleated	0.45 + 0.2	5492507H1-S	5492507H2-S	5492507H3-S

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	1.0 m ² 10.8 ft ²	2.0 m ² 21.5 ft ²	3.0 m ² 32.3 ft ²
Cartridges	0.45 + 0.2	5492507H1	5492507H2	5492507H3
T-Style Maxicaps® γ-irradiatable or autoclavable	0.45 + 0.2	5498307H1G-**-	5498307H2G-**-	5498307H3G-**-
Maxicaps® γ-irradiatable or autoclavable	0.45 + 0.2	5497307H1G-**-xx	5497307H2G-**-xx	5497307H3G-**-xx

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.065 m ² 0.7 ft ²	0.13 m ² 1.4 ft ²	0.26 m ² 2.8 ft ²	0.52 m ² 5.6 ft ²
Midicaps®	0.45 + 0.2	5495307H7-**- -A	5495307H8-**- -A	5495307H9-**- -A	5495307H0-**- -V
Gamma Midicaps®	0.45 + 0.2	5495307H7G-**- -A	5495307H8G-**- -A	5495307H9G-**- -A	5495307H0G-**- -V

		Size 4
	Pore Size	0.021 m ² 0.22 ft ²
Capsules	0.45 + 0.2	5491307H4-**- -B
Gamma Capsules	0.45 + 0.2	5491307H4G-**- -B

		25	47
	Pore Size	4.5 cm ²	17.3 cm ²
Sartoscale	0.45 + 0.2	5495307HV- -LX- -C	5495307HS- -FF- -M

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

xx = IT vent valve design suitable for connection of Sartocheck® integrity tester

Sartopore® Platinum HB

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	1.0 m ² 10.8 ft ²	2.0 m ² 21.5 ft ²	3.0 m ² 32.3 ft ²
Cartridges	0.45 + 0.2	5492507H1- - - -HB	5492507H2- - - -HB	5492507H3- - - -HB
T-Style Maxicaps® γ-irradiatable or autoclavable	0.45 + 0.2	5498307H1G- **HB	5498307H2G- **HB	5498307H3G- **HB
Maxicaps® γ-irradiatable or autoclavable	0.45 + 0.2	5497307H1G- **HB	5497307H2G- **HB	5497307H3G- **HB

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.065 m ² 0.7 ft ²	0.13 m ² 1.4 ft ²	0.26 m ² 2.8 ft ²	0.52 m ² 5.6 ft ²
Midicaps®	0.45 + 0.2	5495307H7- - **HBA	5495307H8- - **HBA	5495307H9- - **HBA	5495307H0- - **HBV
Gamma Midicaps®	0.45 + 0.2	5495307H7G- **HBA	5495307H8G- **HBA	5495307H9G- **HBA	5495307H8G- **HBV

		Size 4
	Pore Size	0.021 m ² 0.22 ft ²
Capsules	0.45 + 0.2	5491307H4- - **HBB
Gamma Capsules	0.45 + 0.2	5491307H4G- **HBB

		25	47
	Pore Size	4.5 cm ²	17.3 cm ²
Sartoscale	0.45 + 0.2	5495307HV- -LXHBC	5495307HS- -FFHBM

** = description of available filter design or connector | adapter types see page 18 | 19
 Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

Sartopore® 2

		Size 0 (5")	Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.3 m ² 3.2 ft ²	0.6 m ² 6.5 ft ²	1.2 m ² 12.9 ft ²	1.8 m ² 19.4 ft ²
Cartridges	0.8 + 0.45	544**06G0	544**06G1	544**06G2	544**06G3
	0.45 + 0.2	544**07H0	544**07H1	544**07H2	544**07H3
Maxicaps®	0.8 + 0.45		5447306G1- -**	5447306G2- -**	5447306G3- -**
	0.45 + 0.2		5447307H1- -** ^{xx}	5447307H2- -** ^{xx}	5447307H3- -** ^{xx}
Gamma Maxicaps®	0.8 + 0.45		5447306G1G- -** ^{xx}	5447306G2G- -** ^{xx}	5447306G3G- -** ^{xx}
	0.45 + 0.2		5447307H1G- -** ^{xx}	5447307H2G- -** ^{xx}	5447307H3G- -** ^{xx}
T-Style Maxicaps® γ-irradiatable or autoclavable	0.8 + 0.45		5448306G1G- -**	5448306G2G- -**	5448306G3G- -**
	0.45 + 0.2		5448307H1G- -**	5448307H2G- -**	5448307H3G- -**

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.05 m ² 0.5 ft ²	0.1 m ² 1.1 ft ²	0.2 m ² 2.2 ft ²	0.45 m ² 4.8 ft ²
Mini Cartridges	0.8 + 0.45	544**06G7- ---- -B	544**06G8- ---- -B	544**06G9- ---- -B	
	0.45 + 0.2	544**07H7- ---- -B	544**07H8- ---- -B	544**07H9- ---- -B	
Midicaps®	0.8 + 0.45	5445306G7- -** -A	5445306G8- -** -A	5445306G9- -** -A	5445306G0- -** -V
	0.45 + 0.2	5445307H7- -** -A	5445307H8- -** -A	5445307H9- -** -A	5445307H0- -** -V
Gamma Midicaps®	0.8 + 0.45	5445306G7G- -** -A	5445306G8G- -** -A	5445306G9G- -** -A	5445306G0G- -** -V
	0.45 + 0.2	5445307H7G- -** -A	5445307H8G- -** -A	5445307H9G- -** -A	5445307H0G- -** -V

		Size 4	Size 5
	Pore Size	0.015 m ² 0.16 ft ²	0.03 m ² 0.32 ft ²
Capsules	0.8 + 0.45		5441306G5- -OO- -B
	0.45 + 0.2	5441307H4- -** -B	5441307H5- -OO- -B
Gamma Capsules	0.45 + 0.2	5441307H4G- -** -B	5441307H5G- -OO- -B

		25	47
	Pore Size	4.5 cm ²	17.3 cm ²
Sartoscale	0.8 + 0.45	5445306GV- -LX- -C	5445306GS- -FF- -M
	0.45 + 0.2	5445307HV- -LX- -C	5445307HS- -FF- -M

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

^{xx} = IT vent valve design suitable for connection of Sartocheck® integrity tester

Sartopore® 2 XLG

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.8 m ² 8.6 ft ²	1.6 m ² 17.2 ft ²	2.4 m ² 25.8 ft ²
Cartridges	0.8 + 0.2	5442507G1	5442507G2	5442507G3
Maxicaps®	0.8 + 0.2	5447307G1- **xx	5447307G2- **xx	5447307G3- **xx
Gamma Maxicaps®	0.8 + 0.2	5447307G1G- **xx	5447307G2G- **xx	5447307G3G- **xx
T-Style Maxicaps® γ-irradiatable or autoclavable	0.8 + 0.2	5448307G1G- **	5448307G2G- **	5448307G3G- **

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.065 m ² 0.7 ft ²	0.13 m ² 1.4 ft ²	0.26 m ² 2.8 ft ²	0.52 m ² 5.6 ft ²
Midicaps®	0.8 + 0.2	5445307G7- **- -A	5445307G8- **- -A	5445307G9- **- -A	5445307G0- **- -V
Gamma Midicaps®	0.8 + 0.2	5445307G7G- **- -A	5445307G8G- **- -A	5445307G9G- **- -A	5445307G0G- **- -V

		Size 4
	Pore Size	0.021 m ² 0.22 ft ²
Capsules	0.8 + 0.2	5441307G4- **- -B
Gamma Capsules	0.8 + 0.2	5441307G4G- **- -B

		25	47
	Pore Size	4.5 cm ²	17.3 cm ²
Sartoscale	0.8 + 0.2	5445307GV- -LX- -C	5445307GS- -FF- -M

** = description of available filter design or connector | adapter types see page 18 | 19
 Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

Sartopore® 2 XLI

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.8 m ² 8.6 ft ²	1.6 m ² 17.2 ft ²	2.4 m ² 25.8 ft ²
Cartridges	0.35 + 0.2	5442507I1	5442507I2	5442507I3
Maxicaps®	0.35 + 0.2	5447307I1- -** ^{xx}	5447307I2- -** ^{xx}	5447307I3- -** ^{xx}
Gamma Maxicaps®	0.35 + 0.2	5447307I1G-**- ^{xx}	5447307I2G-**- ^{xx}	5447307I3G-**- ^{xx}
T-Style Maxicaps® γ-irradiatable or autoclavable	0.35 + 0.2	5448307I1G-**-	5448307I2G-**-	5448307I3G-**-

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.065 m ² 0.7 ft ²	0.13 m ² 1.4 ft ²	0.26 m ² 2.8 ft ²	0.52 m ² 5.6 ft ²
Midicaps®	0.35 + 0.2	5445307I7- -**- -A	5445307I8- -**- -A	5445307I9- -**- -A	5445307I0- -**- -V
Gamma Midicaps®	0.35 + 0.2	5445307I7G-**- -A	5445307I8G-**- -A	5445307I9G-**- -A	5445307I0G-**- -V

		Size 4
	Pore Size	0.021 m ² 0.22 ft ²
Capsules	0.35 + 0.2	5441307I4- -**- -B
Gamma Capsules	0.35 + 0.2	5441307I4G-**- -B

		25	47
	Pore Size	4.5 cm ²	17.3 cm ²
Sartoscale	0.35 + 0.2	5445307IV- -LX- -C	5445307IS- -FF- -M

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

^{xx} = IT vent valve design suitable for connection of Sartocheck® integrity tester

Sartopore® 2 XLM

		Size 0 (5")	Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.4 m ² 4.3 ft ²	0.8 m ² 8.6 ft ²	1.6 m ² 17.2 ft ²	2.4 m ² 25.8 ft ²
Cartridges	0.2 + 0.1	5442558M0	5442558M1	5442558M2	5442558M3
Maxicaps®	0.2 + 0.1		5447358M1- -** xx	5447358M2- -** xx	5447358M3- -** xx
Gamma Maxicaps®	0.2 + 0.1		5447358M1G- ** xx	5447358M2G- ** xx	5447358M3G- ** xx
T-Style Maxicaps® γ-irradiatable or autoclavable	0.2 + 0.1		5448358M1G- **	5448358M2G- **	5448358M3G- **

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.065 m ² 0.7 ft ²	0.13 m ² 1.4 ft ²	0.26 m ² 2.8 ft ²	0.52 m ² 5.6 ft ²
Midicaps®	0.2 + 0.1	5445358M7- -**- -A	5445358M8- -**- -A	5445358M9- -**- -A	5445358M0- -**- -V
Gamma Midicaps®	0.2 + 0.1	5445358M7G- **- -A	5445358M8G- **- -A	5445358M9G- **- -A	5445358M0G- **- -V

		Size 4
	Pore Size	0.021 m ² 0.22 ft ²
Capsules	0.2 + 0.1	5441358M4- -**- -B
Gamma Capsules	0.2 + 0.1	5441358M4G- **- -B

		25	47
	Pore Size	4.5 cm ²	17.3 cm ²
Sartoscale	0.2 + 0.1	5445358MV- -LX- -C	5445358MS- -FF- -M

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

xx = IT vent valve design suitable for connection of Sartocheck® integrity tester

Sartopore® Air

		Size 1 (10")	Size 2 (20")
	Pore Size	0.53 m ² 5.7 ft ²	1.06 m ² 11.4 ft ²
Maxicaps® γ-irradiatable	0.2	5197307A1G-**-IT	5197307A2G-**-IT

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.05 m ² 0.54 ft ²	0.1 m ² 1.1 ft ²	0.17 m ² 1.83 ft ²	0.34 m ² 3.66 ft ²
Midicaps® γ-irradiatable	0.2	5195307A7G-**-A	5195307A8G-**-A	5195307A9G-**-A	5195307A0G-**-V

		Size 4
	Pore Size	0.016 m ² 0.17 ft ²
Capsules γ-irradiatable	0.2	5191307A4G-**-B

** = description of available filter design or connector | adapter types see page 18 | 19
Units per package: A = 4 pieces | B = 5 pieces | V = 2 pieces

Sartoguard Family

Sartoguard PES

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size ¹	0.8 m ² 8.6 ft ²	1.6 m ² 17.2 ft ²	2.4 m ² 25.8 ft ²
Cartridges	0.8 + 0.1	5472558G1	5472558G2	5472558G3
	1.2 + 0.2	5472507F1	5472507F2	5472507F3
T-Style Maxicaps® γ-irradiatable or autoclavable	0.8 + 0.1	5478358G1G-**-	5478358G2G-**-	5478358G3G-**-
	1.2 + 0.2	5478307F1G-**-	5478307F2G-**-	5478307F3G-**-
Maxicaps®	0.8 + 0.1	5477358G1-**-	5477358G2-**-	5477358G3-**-
	1.2 + 0.2	5477307F1-**-	5477307F2-**-	5477307F3-**-
Gamma Maxicaps®	0.8 + 0.1	5477358G1G-**-	5477358G2G-**-	5477358G3G-**-
	1.2 + 0.2	5477307F1G-**-	5477307F2G-**-	5477307F3G-**-

		Size 7	Size 8	Size 9	Size 0
	Pore Size ¹	0.065 m ² 0.7 ft ²	0.13 m ² 1.4 ft ²	0.26 m ² 2.8 ft ²	0.52 m ² 5.6 ft ²
Midicaps®	0.8 + 0.1	5475358G7-**- -A	5475358G8-**- -A	5475358G9-**- -A	5475358G0-**- -V
	1.2 + 0.2	5475307F7-**- -A	5475307F8-**- -A	5475307F9-**- -A	5475307F0-**- -V
Gamma Midicaps®	0.8 + 0.1	5475358G7G-**- -A	5475358G8G-**- -A	5475358G9G-**- -A	5475358G0G-**- -V
	1.2 + 0.2	5475307F7G-**- -A	5475307F8G-**- -A	5475307F9G-**- -A	5475307F0G-**- -V

		Size 4
	Pore Size ¹	0.021 m ² 0.22 ft ²
Capsules	0.8 + 0.1	5471358G4-**- -B
	1.2 + 0.2	5471307F4-**- -B
Gamma Capsules	0.8 + 0.1	5471358G4G-**- -B
	1.2 + 0.2	5471307F4G-**- -B

		25	47
	Pore Size ¹	4.5 cm ²	17.3 cm ²
Sartoscale	0.8 + 0.1	5475358GV- -LX- -C	5475358GS- -FF- -M
	1.2 + 0.2	5475307FV- -LX- -C	5475307FS- -FF- -M

¹ nom.

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

Sartoguard GF

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size ¹	0.45 m ² 4.8 ft ²	0.9 m ² 9.6 ft ²	1.35 m ² 14.5 ft ²
Cartridges	0.2 + 0.1	5482558M1	5482558M2	5482558M3
	0.8 + 0.2	5482507G1	5482507G2	5482507G3
T-Style Maxicaps®	0.2 + 0.1	5488358M1- -**	5488358M2- -**	5488358M3- -**
	0.8 + 0.2	5488307G1- -**	5488307G2- -**	5488307G3- -**
Maxicaps®	0.2 + 0.1	5487358M1- -**	5487358M2- -**	5487358M3- -**
	0.8 + 0.2	5487307G1- -**	5487307G2- -**	5487307G3- -**

		Size 8	Size 9	Size 0
	Pore Size ¹	0.08 m ² 0.9 ft ²	0.15 m ² 1.6 ft ²	0.3 m ² 3.2 ft ²
Midicaps®	0.2 + 0.1	5485358M8- -** -A	5485358M9- -** -A	5485358M0- -** -V
	0.8 + 0.2	5485307G8- -** -A	5485307G9- -** -A	5485307G0- -** -V

		Size 4
	Pore Size ¹	0.012 m ² 0.13 ft ²
Capsules	0.2 + 0.1	5481358M4- -** -B
	0.8 + 0.2	5481307G4- -** -B

		25	47
	Pore Size ¹	4.5 cm ²	17.3 cm ²
Sartoscale	0.2 + 0.1	5485358MV- -LX- -C	5485358MS- -FF- -M
	0.8 + 0.2	5485307GV- -LX- -C	5485307GS- -FF- -M

¹ nom.

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

Sartobran® P Family

Sartobran® P

		Size 0 (5")	Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.3 m ² 3.2 ft ²	0.6 m ² 6.5 ft ²	1.2 m ² 12.9 ft ²	1.8 m ² 19.4 ft ²
Cartridges	0.65 + 0.45	523**06D0- - - -P	523**06D1- - - -P	523**06D2- - - -P	523**06D3- - - -P
	0.45 + 0.2	523**07H0- - - -P	523**07H1- - - -P	523**07H2- - - -P	523**07H3- - - -P
Maxicaps®	0.65 + 0.45		5237306D1- -** xx	5237306D2- -** xx	5237306D3- -** xx
	0.45 + 0.2		5237307H1- -** xx	5237307H2- -** xx	5237307H3- -** xx
T-Style Maxicaps®	0.65 + 0.45		5238306D1- -**	5238306D2- -**	5238306D3- -**
	0.45 + 0.2		5238307H1- -**	5238307H2- -**	5238307H3- -**

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.05 m ² 0.54 ft ²	0.1 m ² 1.1 ft ²	0.2 m ² 2.2 ft ²	0.45 m ² 4.8 ft ²
Mini Cartridges	0.65 + 0.45	5231506D7- - - - -B	5231506D8- - - - -B	5231506D9- - - - -B	
	0.45 + 0.2	523**07H7- - - - -B	523**07H8- - - - -B	523**07H9- - - - -B	
Midicaps®	0.65 + 0.45	5235306D7- -** -A	5235306D8- -** -A	5235306D9- -** -A	5235306D0- -** -V
	0.45 + 0.2	5235307H7- -** -A	5235307H8- -** -A	5235307H9- -** -A	5235307H0- -** -V

		Size 4	Size 5
	Pore Size	0.015 m ² 0.16 ft ²	0.03 m ² 0.32 ft ²
Capsules	0.65 + 0.45		5231306D5- -OO- -B
	0.45 + 0.2	5231307H4- -** -B	5231307H5- -OO- -B

		25	47
	Pore Size	4.5 cm ²	17.3 cm ²
Sartoscale	0.65 + 0.45	5235306DV- -LX- -C	5235306DS- -FF- -M
	0.45 + 0.2	5235307HV- -LX- -C	5235307HS- -FF- -M

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

xx = IT vent valve design suitable for connection of Sartocheck® integrity tester

Sartoclean® Family

Sartoclean® CA

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size ¹	0.6 m ² 6.5 ft ²	1.2 m ² 12.8 ft ²	1.8 m ² 19.3 ft ²
Cartridges	3.0 + 0.8	562**04E1	562**04E2	562**04E3
	0.8 + 0.65	562**05G1	562**05G2	562**05G3
T-Style Maxicaps®	3.0 + 0.8	5628304E1- -**	5628304E2- -**	5628304E3- -**
	0.8 + 0.65	5628305G1- -**	5628305G2- -**	5628305G3- -**
Maxicaps®	3.0 + 0.8	5627304E1- -**	5627304E2- -**	5627304E3- -**
	0.8 + 0.65	5627305G1- -**	5627305G2- -**	5627305G3- -**

		Size 7	Size 8	Size 9	Size 0
	Pore Size ¹	0.05 m ² 0.5 ft ²	0.1 m ² 1.1 ft ²	0.2 m ² 2.2 ft ²	0.45 m ² 4.8 ft ²
Mini Cartridges	3.0 + 0.8			5621504E9- ---- -B	
	0.8 + 0.65			5621505G9- ---- -B	
Midicaps®	3.0 + 0.8	5625304E7- ***- -A	5625304E8- ***- -A	5625304E9- ***- -A	5625304E0- ***- -V
	0.8 + 0.65	5625305G7- ***- -A	5625305G8- ***- -A	5625305G9- ***- -A	5625305G0- ***- -V

		25	47
	Pore Size ¹	4.5 cm ²	17.3 cm ²
Sartoscale	3.0 + 0.8	5625304EV- -LX- -C	5625304ES- -FF- -M
	0.8 + 0.65	5625305GV- -LX- -C	5625305GS- -FF- -M

¹ nom.

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

Sartoclean® GF

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size ¹	0.47 m ² 5.1 ft ²	0.94 m ² 10.2 ft ²	1.41 m ² 15.3 ft ²
Cartridges	3.0 + 0.8	560**04E1	560**04E2	560**04E3
	0.8 + 0.65	560**05G1	560**05G2	560**05G3
T-Style Maxicaps®	3.0 + 0.8	5608304E1- -**	5608304E2- -**	5608304E3- -**
	0.8 + 0.65	5608305G1- -**	5608305G2- -**	5608305G3- -**
Maxicaps®	3.0 + 0.8	5607304E1- -**	5607304E2- -**	5607304E3- -**
	0.8 + 0.65	5607305G1- -**	5607305G2- -**	5607305G3- -**

		Size 7	Size 8	Size 9	Size 0
	Pore Size ¹	0.05 m ² 0.54 ft ²	0.08 m ² 0.85 ft ²	0.16 m ² 1.7 ft ²	0.32 m ² 3.4 ft ²
Mini Cartridges	3.0 + 0.8			5601504E9- ---- -B	
	0.8 + 0.65			5601505G9- ---- -B	
Midicaps®	3.0 + 0.8	5605304E7- -** -A	5605304E8- -** -A	5605304E9- -** -A	5605304E0- -** -V
	0.8 + 0.65	5605305G7- -** -A	5605305G8- -** -A	5605305G9- -** -A	5605305G0- -** -V

		25	47
	Pore Size ¹	4.5 cm ²	17.3 cm ²
Sartoscale	3.0 + 0.8	5605304EV- -LX- -C	5605304ES- -FF- -M
	0.8 + 0.65	5605305GV- -LX- -C	5605305GS- -FF- -M

¹ nom.

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

Sartolon Family

Sartolon

		Size 1 (10")	Size 2 (20")	Size 3 (30")	Size 4 (40")
	Pore Size	0.6 m² 6.5 ft²	1.2 m² 12.9 ft²	1.8 m² 19.4 ft²	2.4 m² 25.9 ft²
Cartridges	0.45 + 0.2	510**07H1	510**07H2	510**07H3	510**07H4
Maxicaps®	0.45 + 0.2	5107307H1- -** xx	5107307H2- -** xx	5107307H3- -** xx	

		Size 9
	Pore Size	0.2 m² 2.2 ft²
Mini Cartridges	0.45 + 0.2	5101507H9- - - - -B
Midicaps®	0.45 + 0.2	5105307H9- -** -A

		25	47
	Pore Size	4.5 cm²	17.3 cm²
Sartoscale	0.45 + 0.2	5105307HV- -LX- -C	5105307HS- -FF- -M

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | M = 3 pieces

xx = IT vent valve design suitable for connection of Sartocheck® integrity tester

Sartopure® Family

Sartopure® PP3 0.45 µm

		Size 1 (10")	Size 2 (20")	Size 3 (30")	Size 4 (40")
	Retention Rate ¹	0.4 m² 4.3 ft²	0.8 m² 8.6 ft²	1.2 m² 12.9 ft²	1.6 m² 17.2 ft²
Cartridges	0.45	505**06P1	505**06P2	505**06P3	505**06P4
T-Style Maxicaps®	0.45	5058306P1- -**	5058306P2- -**	5058306P3- -**	
Maxicaps®	0.45	5057306P1- -**	5057306P2- -**	5057306P3- -**	

		Size 7	Size 8	Size 9	Size 0
	Retention Rate ¹	0.04 m² 0.43 ft²	0.08 m² 0.86 ft²	0.13 m² 1.4 ft²	0.26 m² 2.8 ft²
Midicaps®	0.45	5055306P7- -** -A	5055306P8- -** -A	5055306P9- -** -A	5055306P0- -** -V

Sartopure® PP3 0.65 – 3 µm

		Size 1 (10")	Size 2 (20")	Size 3 (30")	Size 4 (40")
	Retention Rate ¹	0.45 m² 4.8 ft²	0.9 m² 9.7 ft²	1.35 m² 14.5 ft²	1.8 m² 19.4 ft²
Cartridges	3	505**02P1	505**02P2	505**02P3	
	1.2	505**03P1	505**03P2	505**03P3	505**03P4
	0.65	505**05P1	505**05P2	505**05P3	505**05P4
T-Style Maxicaps®	3	5058302P1- -**	5058302P2- -**	5058302P3- -**	
	1.2	5058303P1- -**	5058303P2- -**	5058303P3- -**	
	0.65	5058305P1- -**	5058305P2- -**	5058305P3- -**	
Maxicaps®	3	5057302P1- -**	5057302P2- -**	5057302P3- -**	
	1.2	5057303P1- -**	5057303P2- -**	5057303P3- -**	
	0.65	5057305P1- -**	5057305P2- -**	5057305P3- -**	

		Size 7	Size 8	Size 9	Size 0
	Retention Rate ¹	0.05 m² 0.54 ft²	0.09 m² 0.97 ft²	0.15 m² 1.61 ft²	0.30 m² 3.23 ft²
Mini Cartridges	3	5051502P7- ---- -B	5051502P8- ---- -B	5051502P9- ---- -B	
	1.2	5051503P7- ---- -B	5051503P8- ---- -B	5051503P9- ---- -B	
	0.65	5051505P7- ---- -B	5051505P8- ---- -B	5051505P9- ---- -B	
Midicaps®	3	5055302P7- -** -A	5055302P8- -** -A	5055302P9- -** -A	5055302P0- -** -V
	1.2	5055303P7- -** -A	5055303P8- -** -A	5055303P9- -** -A	5055303P0- -** -V
	0.65	5055305P7- -** -A	5055305P8- -** -A	5055305P9- -** -A	5055305P0- -** -V

¹ nom.

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | V = 2 pieces

Sartopure® PP3 5 – 50 µm

		Size 1 (10")	Size 2 (20")	Size 3 (30")	Size 4 (40")
Retention Rate ¹		0.65 m ² 7 ft ²	1.3 m ² 14 ft ²	1.95 m ² 21 ft ²	2.6 m ² 28 ft ²
Cartridges	50	505**50P1	505**50P2	505**50P3	
	20	505**20P1	505**20P2	505**20P3	
	8	505**01P1	505**01P2	505**01P3	505**01P4
	5	505**42P1	505**42P2	505**42P3	505**42P4
T-Style Maxicaps®	50	5058350P1- -**	5058350P2- -**	5058350P3- -**	
	20	5058320P1- -**	5058320P2- -**	5058320P3- -**	
	8	5058301P1- -**	5058301P2- -**	5058301P3- -**	
	5	5058342P1- -**	5058342P2- -**	5058342P3- -**	
Maxicaps®	50	5057350P1- -**	5057350P2- -**	5057350P3- -**	
	20	5057320P1- -**	5057320P2- -**	5057320P3- -**	
	8	5057301P1- -**	5057301P2- -**	5057301P3- -**	
	5	5057342P1- -**	5057342P2- -**	5057342P3- -**	

		Size 7	Size 8	Size 9	Size 0
Retention Rate ¹		0.065 m ² 0.7 ft ²	0.12 m ² 1.3 ft ²	0.21 m ² 2.26 ft ²	0.42 m ² 4.52 ft ²
Mini Cartridges	20			5051520P9- ---- -B	
	8			5051501P9- ---- -B	
	5	5051542P7- ---- -B	5051542P8- ---- -B	5051542P9- ---- -B	
Midicaps®	50	5055350P7- **- -A	5055350P8- **- -A	5055350P9- **- -A	5055350P0- **- -V
	20	5055320P7- **- -A	5055320P8- **- -A	5055320P9- **- -A	5055320P0- **- -V
	8	5055301P7- **- -A	5055301P8- **- -A	5055301P9- **- -A	5055301P0- **- -V
	5	5055342P7- **- -A	5055342P8- **- -A	5055342P9- **- -A	5055342P0- **- -V

¹ nom.

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | V = 2 pieces

Sartopure® PP3 Capsules & Sartoscale

		Size 4	Size 5
	Retention Rate ¹	0.012 m ² 0.13 ft ²	0.024 m ² 0.26 ft ²
Capsules	0.45	5051306P4- **- -B	5051306P5- **- -B

		Size 4	Size 5
	Retention Rate ¹	0.013 m ² 0.14 ft ²	0.026 m ² 0.28 ft ²
Capsules	1.2	5051303P4- **- -B	5051303P5- **- -B
	0.65	5051305P4- **- -B	5051305P5- **- -B

		Size 4	Size 5
	Retention Rate ¹	0.018 m ² 0.19 ft ²	0.036 m ² 0.39 ft ²
Capsules	50	5051350P4- -SS- -B	
	20	5051320P4- **- -B	5051320P5- -OO- -B
	8	5051301P4- **- -B	5051301P5- -OO- -B
	5	5051342P4- **- -B	5051342P5- -OO- -B
	3	5051302P4- **- -B	5051302P5- -OO- -B

		25	47
	Retention Rate ¹	4.5 cm ²	17.3 cm ²
Sartoscale	50	5055350PV- -LX- -C	5055350PS- -FF- -M
	20	5055320PV- -LX- -C	5055320PS- -FF- -M
	8	5055301PV- -LX- -C	5055301PS- -FF- -M
	5	5055342PV- -LX- -C	5055342PS- -FF- -M
	3	5055302PV- -LX- -C	5055302PS- -FF- -M
	1.2	5055303PV- -LX- -C	5055303PS- -FF- -M
	0.65	5055305PV- -LX- -C	5055305PS- -FF- -M
	0.45	5055306PV- -LX- -C	5055306PS- -FF- -M

¹ nom.

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: B = 5 pieces | C = 6 pieces | M = 3 pieces

Sartopure® GF Plus

		Size 1 (10")	Size 2 (20")	Size 3 (30")	Size 4 (40")
	Retention Rate ¹	0.4 m ² 4.3 ft ²	0.8 m ² 8.6 ft ²	1.2 m ² 12.9 ft ²	1.6 m ² 17.2 ft ²
Cartridges	1.2	555**03P1	555**03P2	555**03P3	555**03P4
	0.65	555**05P1	555**05P2	555**05P3	555**05P4
T-Style Maxicaps®	1.2	5558303P1- -**	5558303P2- -**	5558303P3- -**	
	0.65	5558305P1- -**	5558305P2- -**	5558305P3- -**	
Maxicaps®	1.2	5557303P1- -**	5557303P2- -**	5557303P3- -**	
	0.65	5557305P1- -**	5557305P2- -**	5557305P3- -**	

		Size 7	Size 8	Size 9	Size 0
	Retention Rate ¹	0.04 m ² 0.4 ft ²	0.08 m ² 0.9 ft ²	0.12 m ² 1.3 ft ²	0.25 m ² 2.7 ft ²
Mini Cartridges	1.2	5551503P7- ----B	5551503P8- ----B	5551503P9- ----B	
	0.65	5551505P7- ----B	5551505P8- ----B	5551505P9- ----B	
Midicaps®	1.2	5555303P7- ***-A	5555303P8- ***-A	5555303P9- ***-A	5555303P0- ***-V
	0.65	5555305P7- ***-A	5555305P8- ***-A	5555305P9- ***-A	5555305P0- ***-V

		25	47
	Retention Rate ¹	4.5 cm ²	17.3 cm ²
Sartoscale	1.2	5555303PV- -LX- -C	5555303PS- -FF- -M
	0.65	5555305PV- -LX- -C	5555305PS- -FF- -M

Sartopure® GA

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Retention Rate ¹	0.7 m ² 7.5 ft ²	1.4 m ² 15.1 ft ²	2.1 m ² 22.6 ft ²
Cartridges	0.2	559**07P1- ---GA	559**07P2- ---GA	559**07P3- ---GA

		Size 9
	Retention Rate ¹	0.2 m ² 2.2 ft ²
Mini Cartridges	0.2	5591507P9- ---GAB

¹ nom.

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces

Sartofluor® Family

Sartofluor®

		Size 0 (5")	Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.375 m² 4.05 ft²	0.75 m² 8.1 ft²	1.5 m² 16.1 ft²	2.25 m² 24.2 ft²
Cartridges GA	0.2	518**07T0- ---GA	518**07T1- ---GA	518**07T2- ---GA	518**07T3- ---GA
	Pore Size		0.5 m² 5.4 ft²	1.0 m² 10.8 ft²	1.5 m² 16.1 ft²
Cartridges LG	0.2		518**07T1- ---LG	518**07T2- ---LG	518**07T3- ---LG
Maxicaps®	0.2		5187307T1- -**	5187307T2- -**	

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.05 m² 0.5 ft²	0.1 m² 1.1 ft²	0.2 m² 2.2 ft²	0.45 m² 4.8 ft²
Mini Cartridges	0.2	518**07T7- ----B	518**07T8- ----B	518**07T9- ----B	
Junior	0.2	5181407T7- ----B			
Midicaps®	0.2	5185307T7- **- -A	5185307T8- **- -A	5185307T9- **- -A	5185307T0- **- -V

		Size 4	Size 5
	Pore Size	0.015 m² 0.16 ft²	0.03 m² 0.32 ft²
Capsules	0.2	5181307T4- **- -B	5181307T5- -OO- -D

Sartofluor® HR

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.56 m² 6.7 ft²	1.12 m² 12.1 ft²	1.68 m² 18.1 ft²
Cartridges	0.2	5172507T1- ---HR	5172507T2- ---HR	5172507T3- ---HR

** = description of available filter design or connector | adapter types see page 18 | 19

Units per package: A = 4 pieces | B = 5 pieces | C = 6 pieces | V = 2 pieces | M = 3 pieces | D = 10 pieces

Midisart®

Midisart® 2000

Pore Size	Connectors	Box of 12 (sterile)	Box of 25 (sterile)	Box of 100 (unsterile)	Plastic bucket* of 100 (unsterile)	Box of 500 (unsterile)
0.2 µm	Stepped hose barb	17805-----E	17805-----G	17805-----UPN	17805-----UQN	17805-----UPQ
0.2 µm	½ NPT	17805-----NPE	17805-----NPG			
0.2 µm	Small hose barb (4–12 mm)			17877-----UPN		
0.2 µm	TC 25			17805-----TCN		
0.45 µm	Stepped hose barb	17804-----E	17804-----G			

Sartopore® Air Midisart®

Pore Size	Connectors	Box of 12 (unsterile)	Box of 100 (unsterile)	Plastic bucket* of 100 (unsterile)	Box of 500 (unsterile)
0.2 µm	Stepped hose barb	17805-----AIE	17805-----AIN	17805-----AQN	17805-----AIQ

* = Plastic bag in plastic bucket (double packaging) for transfer into clean room. Memory discs enclosed.

Virosart®

Virosart® HF

	Pore Size	0.8 m ² 8.6 ft ²	2.4 m ² 25.8 ft ²
Process module Sterile (γ-irradiated)	20 nm	3VI- -28-HCGSF	3VI- -28-MCGSF
Process module γ-irradiatable	20 nm	3VI- -28-HC-SF	3VI- -28-MC-SF

	Pore Size	200 cm ² 0.22 ft ²	0.2 m ² 2.15 ft ²
Mid-scale module Sterile (γ-irradiated)	20 nm	3VI- -28-DCGFO	3VI- -28-FCGFO
Mid-scale module γ-irradiatable	20 nm	3VI- -28-DC-FO	3VI- -28-FC-FO

	Pore Size	1.7 cm ²	5 cm ²
Lab module Sterile (γ-irradiated)	20 nm	3VI- -28-TCGML- -V	3VI- -28-BCGML- -V
Lab module Integrity tested; Sterile (γ-irradiated)	20 nm	3VI- -28-TCGMLITV	3VI- -28-BCGMLITV

Virosart® Media

	Pore Size	1 m ² 10.8 ft ²
Process module sterile (γ-irradiated)	20 nm	3V2- -28-IVGSF
Process module γ-irradiatable	20 nm	3V2- -28-IV-SF

	Pore Size	0.3 m ² 3.2 ft ²
Mid-scale module Sterile (γ-irradiated)	20 nm	3V2- -28-GVGFO
Mid-scale module γ-irradiatable	20 nm	3V2- -28-GV-FO

	Pore Size	5 cm ²
Lab module Sterile (γ-irradiated)	20 nm	3V2- -28-BVGML- -V
Lab module Integrity tested; Sterile (γ-irradiated)	20 nm	3V2- -28-BVGMLITV

Virosart® CPV

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.7 m ² 7.5 ft ²	1.4 m ² 15 ft ²	2.1 m ² 22.6 ft ²
Cartridge (Autoclavable Steamable)	20 nm	5452528V1	5452528V2	5452528V3
Maxicaps® (Autoclavable)	20 nm	5457328V1- -SS	5457328V2- -SS	5457328V3- -SS

		Size 9
	Pore Size	0.2 m ² 2.15 ft ²
Midicaps® (Autoclavable)	20 nm	5455328V9- -FF- -V

		Size 4
	Pore Size	180 cm ² 0.19 ft ²
Capsule (Autoclavable)	20 nm	5451328V4- -OO- -B 5451328V4- -FF- -B

		5 cm ²
Minisart® (Autoclavable)	20 nm	545VM- - - - -B
Minisart® Integrity tested Sterile (by ETO)	20 nm	545VM- - - - -ITA

Virosart® HC

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.9 m ² 9.7 ft ²	1.8 m ² 19.4 ft ²	2.7 m ² 29 ft ²
Cartridge (Autoclavable Steamable)	20 nm	5392528V1	5392528V2	5392528V3
T-Style Maxicaps® (Autoclavable)	20 nm	5398328V1- -SS	5398328V2- -SS	5398328V3- -SS

		Size 9
	Pore Size	0.27 m ² 2.9 ft ²
Midicaps® (Autoclavable)	20 nm	5395328V9- -FF- -V

		Size 4
	Pore Size	240 cm ² 0.27 ft ²
Capsule (Autoclavable)	20 nm	5391328V4- -FF- -B

		5 cm ²
Minisart® (Autoclavable)	20 nm	539VM-----B
Minisart® Integrity tested Sterile (by ETO)	20 nm	539VM-----ITA

Virosart® Max

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.7 m ² 7.5 ft ²	1.4 m ² 15 ft ²	2.1 m ² 22.6 ft ²
Cartridge (Autoclavable Steamable)	0.1 µm	54A2558N1	54A2558N2	54A2558N3
T-Style Maxicaps® (Autoclavable)	0.1 µm	54A8358N1- -SS	54A8358N2- -SS	54A8358N3- -SS

		Size 9
	Pore Size	0.21 m ² 2.26 ft ²
Midicaps® Sterile (Autoclaved)	0.1 µm	54A5358N9- -FF- -V

		Size 4
	Pore Size	220 cm ² 0.24 ft ²
Capsule Sterile (Autoclaved)	0.1 µm	54A1358N4- -FF- -B

	Pore Size	5 cm ²	21 cm ²
Minisart® Sartolab (Autoclavable)	0.1 µm	54AMI-----B	54ASA-----V

Filter Transfer Sets

The Fast Solution

Standardized basic Transfer Sets that are quickly available:
Pre-Designed Solutions (PDS)

Sartopore®
Platinum

Sartopore® 2,
XLM and XLG

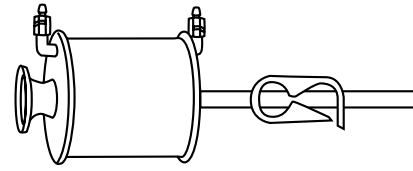


500 mm
TPE tubing



Opta®

MPX Connector



Sealing & Welding
C-Flex® tube for sterile
weld with Biowelder® TC

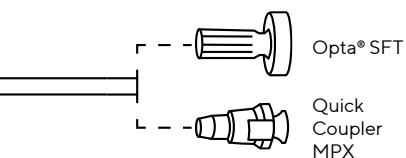
Sartopore® Platinum 0.45 | 0.2 µm

Part Number	Tubing Outlet Connector	Filter	Lot size*
FFU500235	½" ID Clear C-Flex® 500 mm Opta® male ½"	Midicaps® size 7	4
FFU500236		Midicaps® size 8	4
FFU500237		Midicaps® size 9	4
FFU500238		Midicaps® size 0	4
FFU500239		Maxicaps® T-Style size 1	4
FFU500240		Maxicaps® T-Style size 2	4
FFU500241		Maxicaps® T-Style size 3	4
FFU500242	½" ID Clear C-Flex® 500 mm MPX male + sealing cap	Midicaps® size 7	4
FFU500243		Midicaps® size 8	4
FFU500244		Midicaps® size 9	4
FFU500245		Midicaps® size 0	4
FFU500246		Maxicaps® T-Style size 1	4
FFU500247		Maxicaps® T-Style size 2	4
FFU500248		Maxicaps® T-Style size 3	4

Sartopore® 2 0.45 | 0.2 µm

Part Number	Tubing Outlet Connector	Filter	Lot size*
FFU500214	½" ID Clear C-Flex® 500 mm Opta® male ½"	Midicaps® size 7	4
FFU500215		Midicaps® size 8	4
FFU500216		Midicaps® size 9	4
FFU500217		Midicaps® size 0	4
FFU500249		Maxicaps® T-Style size 1	4
FFU500250		Maxicaps® T-Style size 2	4
FFU500251		Maxicaps® T-Style size 3	4
FFU500252	½" ID Clear C-Flex® 500 mm MPX male + sealing cap	Midicaps® size 7	4
FFU500253		Midicaps® size 8	4
FFU500254		Midicaps® size 9	4
FFU500255		Midicaps® size 0	4
FFU500256		Maxicaps® T-Style size 1	4
FFU500257		Maxicaps® T-Style size 2	4
FFU500258		Maxicaps® T-Style size 3	4

C-Flex® is a registered trademark of Saint-Gobain Performance Plastics Corporation.



The Flexible Solution

Configurable filter and sensor Transfer Sets with a broad range of tubing and connector options: Configure-To-Order (CTO)

Sartorius Filters

BioPAT® Flow

BioPAT® Pressure

The Special Solution

Customized solutions for complex assemblies that need to be designed from scratch: Engineer-To-Order (ETO)

[Click here to download our Single-Use Filtration and Sensor Solutions brochure and learn more about our CTO & ETO Solutions](#)

Sartopore® 2 XLG 0.8 | 0.2 µm

Part Number	Tubing Outlet Connector	Filter	Lot size*
FFU500220	½" ID Clear C-Flex® 500 mm	Midicaps® size 7	4
FFU500221		Midicaps® size 8	4
FFU500222		Midicaps® size 9	4
FFU500223		Midicaps® size 0	4
FFU500259		Maxicaps® T-Style size 1	4
FFU500260		Maxicaps® T-Style size 2	4
FFU500261		Maxicaps® T-Style size 3	4
FFU500262	½" ID Clear C-Flex® 500 mm	Midicaps® size 7	4
FFU500263		Midicaps® size 8	4
FFU500264		Midicaps® size 9	4
FFU500172		Midicaps® size 0	4
FFU500265		Maxicaps® T-Style size 1	4
FFU500266		Maxicaps® T-Style size 2	4
FFU500267		Maxicaps® T-Style size 3	4

Sartopore® 2 XLM 0.2 | 0.1 µm

Part Number	Tubing Outlet Connector	Filter	Lot size*
FFU500268	½" ID Clear C-Flex® 500 mm	Midicaps® size 7	4
FFU500269		Midicaps® size 8	4
FFU500270		Midicaps® size 9	4
FFU500271		Midicaps® size 0	4
FFU500272		Maxicaps® T-Style size 1	4
FFU500273		Maxicaps® T-Style size 2	4
FFU500274		Maxicaps® T-Style size 3	4
FFU500275	½" ID Clear C-Flex® 500 mm	Midicaps® size 7	4
FFU500276		Midicaps® size 8	4
FFU500277		Midicaps® size 9	4
FFU500278		Midicaps® size 0	4
FFU500226		Maxicaps® T-Style size 1	4
FFU500279		Maxicaps® T-Style size 2	4
FFU500225		Maxicaps® T-Style size 3	4

* = It can be only ordered the minimum order quantity or a multiple of it.

Maxicaps® MR

Large-Scale Single-Use
Filtration Device:
Maxicaps® MR

Ready-to-use, pre-assembled,
pre-sterilized

Configuration

Filter



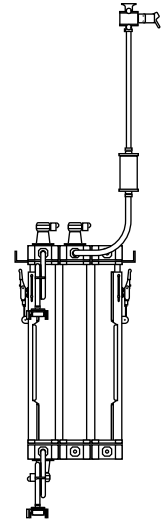
Venting



Inlet | Outlet
Tubing



Inlet | Outlet
Connectors



Configuration Options

Filter	Retention Rate Pore Size	Material	Delivery Condition	Number of 30" Cartridges	Filtration Area (Min.)	Filtration Area (Max.)
Sartoguard PES	0.1 µm 0.2 µm	Polyetersulfone	Sterile	2 – 9	4.8 m² 51.6 ft²	21.6 m² 232.2 ft²
Sartoguard GF	0.1 µm 0.2 µm	Polyetersulfone	Non-Sterile	2 – 9	2.7 m² 29 ft²	12.15 m² 130.5 ft²
Sartopure® GF Plus	0.65 µm 1.2 µm	Glass Fiber	Non-Sterile Sanitary	2 – 9	2.4 m² 25.8 ft²	10.8 m² 116.1 ft²
Sartopure® PP3	0.45 µm 0.65 µm 1.2 µm 3 µm 5 µm 8 µm 20 µm 50 µm 100 µm	Polypropylene	Non-Sterile Sanitary	2 – 9	2.4 m² 25.8 ft²	17.55 m² 189 ft²
Sartoclean® CA	0.65 µm 0.8 µm	Cellulose Acetate	Sanitary	2 – 9	3.6 m² 38.8 ft²	16.2 m² 174.6 ft²
Sartoclean® GF	0.65 µm 0.8 µm	Cellulose Acetate	Sanitary	2 – 9	2.82 m² 30.6 ft²	12.69 m² 137.7 ft²
Sartopore® Platinum	0.2 µm	Modified Polyethersulfone	Sterile	2 – 9	6 m² 64.6 ft²	27 m² 290.7 ft²
Sartopore® 2	0.2 µm 0.45 µm	Polyetersulfone	Sterile	2 – 9	3.6 m² 38.8 ft²	16.2 m² 174.6 ft²
Sartopore® 2 XLG	0.2 µm	Polyetersulfone	Sterile	2 – 9	4.8 m² 51.6 ft²	21.6 m² 232.2 ft²
Sartopore® 2 XLI	0.2 µm	Polyetersulfone	Sterile	2 – 9	4.8 m² 51.6 ft²	21.6 m² 232.2 ft²
Sartopore® 2 XLM	0.1 µm	Polyetersulfone	Sterile	2 – 9	4.8 m² 51.6 ft²	21.6 m² 232.2 ft²
Sartobran® P	0.2 µm 0.45 µm	Cellulose Acetate	Sanitary	2 – 9	3.6 m² 38.8 ft²	16.2 m² 174.6 ft²
Virosart® Max	0.1 µm	Polyamide	Sanitary	2 – 9	4.2 m² 45.2 ft²	18.9 m² 203.4 ft²

Venting

Basic

Advanced

Inlet | Outlet Tubing

Tuflux® SIL | 750 mm

Sani-Tech® STHT-R® | 750 mm

Pharma-50 | 750 mm

Tuflux® TPE | 1,500 mm

C-Flex® 374 | 1,500 mm

Inlet | Outlet Connectors

AseptiQuik® G

AseptiQuik® L

Opta® SFT Female

Opta® SFT Male

Presto HB

1.5" Sanitary Clamp

Plug for TPE tubing

All Maxicaps® MR with gamma stable filter are delivered sterile. Non gamma stable filters are delivered in the sanitary delivery condition, that means autoclaved filter capsules, irradiated lids and tub sets finally assembled in a controlled environment by trained operators for lowest possible bioburden. Sani-Tech® STHT®-R is a registered trademark of Saint-Gobain Performance Plastics Corporation. AseptiQuik® is a registered trademark of Colder Products Company.

Maxicaps® MR Virosart® Media and HF

Large-Scale Single-Use
Filtration Device:
Maxicaps® MR Virosart®

Ready-to-use, pre-
assembled, pre-sterilized

Venting

Inlet
Tubing

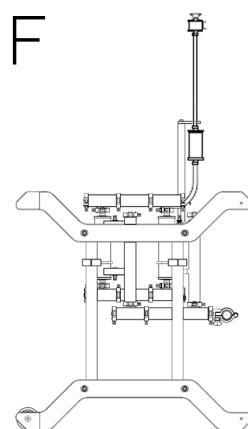
Inlet
Conne-
ctors

Virosart®
Media

Virosart®
HF

Outlet
Tubing

Outlet
Conne-
ctors



Overview Virusfilter Options

Filter	Pore Size	Material	Delivery Condition	Number of Process Modules	Filtration Area (Min.)	Filtration Area (Max.)
Virosart® Media	20 nm	Polyethersulfone Surface modified	Gamma Irradiated	3 or 6	3 m² 32.3 ft²	6 m² 64.6 ft²
Virosart® HF	20 nm	Polyethersulfone Surface modified	Gamma Irradiated	2 – 6	4.8 m² 51.7 ft²	14.4 m² 155 ft²

Connections and Tubing

Filter	Inlet Tubing	Inlet Connectors	Outlet Tubing	Outlet Connectors
Virosart® Media	Reinforced silicon tubing ¾" ID × 1 ½" OD (750 mm)	2 × AseptiQuik® G on Tri-Clamp 1 ½"	Reinforced silicon tubing ¾" ID × 1 ½" OD (750 mm)	2 × AseptiQuik® G on Tri-Clamp 1 ½"
Virosart® HF	Reinforced silicon tubing ¾" ID × 1 ½" OD (750 mm)	2 × AseptiQuik® G on Tri-Clamp 1 ½"	Reinforced silicon tubing ¾" ID × 1 ½" OD (750 mm)	2 × AseptiQuik® G on Tri-Clamp 1 ½" 1 × AseptiQuik® G on Tri-Clamp ¾", Tri-Clamp 1 ½", and flow restrictor

Venting	Tubing	Additional Reservoir	Air Filter
Venting Line Basic Maxicaps® MR Virosart®	Sani-Tech® STHT-R® ¾" × 1 ½" OD (500 mm) Sani-Tech® STHT-R® ½" ID × ⅞" OD (400 mm)	Yes – Translucent	Sartopore® Air Midicaps® 0.2 µm Size 7





Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Straße 11
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 **For more information, visit**
[sartorius.com](https://www.sartorius.com)