



# 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	29
Sartobran® P Family	31
Sartoclean® Family	32
Sartolon Family	34
Sartopure® Family	35
Sartofluor® Family	39
Midisart®	40
Virosart®	41
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.

■ recommended  
 ■ alternatively recommended

## Biotech

Unit Operation		Particle Reduction		Prefilters   Bioburden Reduction				Mycoplasma Retentive Filters	Sterile Filters							
		Sartopure®		Sartoclear®		Sartoguard			Sartoclean®		Sartopore® 2		Sartopore® Platinum		Sartobran® P	Sartolon
		PP3	GF Plus	GF	GF	PES	GF		GF	CA	0.2 µm	XLG	XLI	0.2 µm	HB	0.2 µm
Application		Material ▶	GF		PES	PES + GF	CA + GF	CA		PES	PES	mod. PES	mod. PES	CA		
Media Filtration	Serum-containing		■		■	■			■		■					
	Serum-free				■				■		■					
	Protein-free								■		■					
	Chemically defined								■		■	■				
Buffer & Auxillaries 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							■				■		■		

# Application | Filter Matrix

■ recommended  
 ■ alternatively recommended

## Pharma

Material ▶	Particle Reduction			Prefilters   Bioburden Reduction				Mycoplasma Retentive Filters	Sterile Filters							
	Sartopure®		Sartoclear®	Sartoguard		Sartoclean®			Sartopore® 2			Sartopore® Platinum		Sartobran® P	Sartolon	Sartofluor®
	PP3	GF Plus		PES	GF	GF	CA		0.2 µm	XLG	XLI	0.2 µm	HB	0.2 µm	PA	LG
	PP	GF		PES	PES + GF	CA + GF	CA	PES	PES	PES	mod. PES	mod. PES	CA	PA	PTFE	
<b>Classical Pharma</b>																
SVP	■	■		■							■					
LVP	■			■												
API	Antibiotics (solvents)	■												■		■
	Water-based	■			■									■		
Ophthalmics	■	■		■						■	■		■			
<b>Blood &amp; Plasma</b>																
Albumin & Globulines	Intermediate Process Filtration		■	■		■										
	Filtration after Virus Inactivation	■	■			■										
	Form & Fill				■											
Clotting Factors	Intermediate Process Filtration						■						■			■
	Form & Fill												■			■
<b>Others</b>																
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](http://www.sartorius.com/validation-service)

# Application | Filter Matrix

■ recommended  
 ■ alternatively recommended

## Air & Gas Filtration

		Midisart®		Sartopore® Air	Sartopure®	Sartofluor®	
		Sartopore® Air	2000	PES	GA <sup>1</sup>	GA	HR
		PES	PTFE	PES	Hydrophob. GF	PTFE	PTFE
<b>Compressed Gas</b>							
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					■	

<sup>1</sup> Non-sterilizing grade

## Virus Retention

		Prefilters		Virus Retentive Filters			
		Virosart® Max	Sartopore® 2 XLM	Virosart® HF	Virosart® CPV	Virosart® Media	Virosart® HC
		Polyamide	PES	mod. PES	PES	PES	mod. PES
<b>Biotech</b>							
mAb, rec. Proteins	Virus Retention DSP	■		■	■		
	Virus Retention for Media Filtration		■			■	
Viral Vaccines, Viral Vectors and Cell Therapy	Virus Retention for Media Filtration		■			■	
<b>Blood &amp; Plasma</b>							
Plasma Derivates	Virus Retention DSP	■					■

# Our Portfolio

## Sartopore® Platinum

Sterilizing Grade 0.45 | 0.2 µm

Filter Material		Filter Designs	Sterilization Methods
Prefilter Membrane:	modified Polyethersulfone	Cartridges	In-line steamable or autoclavable
Final Membrane:	modified Polyethersulfone	T-Style Maxicaps®	Autoclavable or γ-irradiatable (hybrid version) <span style="border: 1px solid black; padding: 0 2px;">S</span>
		Maxicaps®	Autoclavable or γ-irradiatable (hybrid version) <span style="border: 1px solid black; padding: 0 2px;">S</span>
		Midicaps®	Autoclavable <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable
		Capsules Size 4	Autoclavable <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable

## Sartopore® 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) <span style="border: 1px solid black; padding: 0 2px;">S</span>
		Maxicaps®	Autoclavable or γ-irradiatable (hybrid version) <span style="border: 1px solid black; padding: 0 2px;">S</span>
		Midicaps®	Autoclavable <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable
		Capsules Size 4	Autoclavable <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable

## 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 <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable
		Capsules Size 4	Autoclavable <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable
		Capsules Size 5	Autoclavable <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable

## 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 <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable
		Capsules Size 4	Autoclavable <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable

## 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 <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable
		Capsules Size 4	Autoclavable <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable

## 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 <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable
		Capsules Size 4	Autoclavable <span style="border: 1px solid black; padding: 0 2px;">S</span> γ-irradiatable

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 <span style="float: right;">[S]</span>
			γ-irradiatable
		Capsules Size 4	Autoclavable <span style="float: right;">[S]</span>
			γ-irradiatable

## Sartoguard GF

Bioburden Reduction | 0.8 | 0.1 µm (nom.)  
 Particle Retention 1.2 | 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 Size 4	Autoclavable

[S] delivered in sterile condition (autoclaved)

## 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 <span style="float: right;">[S]</span>
		Maxicaps®	Autoclavable <span style="float: right;">[S]</span>
		Midicaps®	Autoclavable <span style="float: right;">[S]</span>
		Capsules Size 4	Autoclavable <span style="float: right;">[S]</span>
		Capsules Size 5	Autoclavable <span style="float: right;">[S]</span>

## Sartoclean® CA

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

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 <span style="float: right;">[S]</span>
		Maxicaps®	Autoclavable <span style="float: right;">[S]</span>
		Midicaps®	Autoclavable <span style="float: right;">[S]</span>

## Sartoclean® GF

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

Filter Material		Filter Designs	Sterilization Methods
Filter Active Fleece:	Glass Fiber	Mini Cartridges	In-line steamable or autoclavable
Prefilter Membrane:	Cellulose Acetate	Cartridges	In-line steamable or autoclavable
Final Membrane:	Cellulose Acetate	T-Style Maxicaps®	Autoclavable <span style="float: right;">[S]</span>
		Maxicaps®	Autoclavable <span style="float: right;">[S]</span>
		Midicaps®	Autoclavable <span style="float: right;">[S]</span>

## Sartolon

Sterilizing Grade 0.45 | 0.2 µm

Filter Material		Filter Designs	Sterilization Methods
Prefilter Membrane:	Polyamide (Nylon)	Mini Cartridges	In-line steamable or autoclavable
Final Membrane:	Polyamide (Nylon)	Cartridges	In-line steamable or autoclavable
		Maxicaps®	Autoclavable <span style="float: right;">[S]</span>
		Midicaps®	Autoclavable <span style="float: right;">[S]</span>

## Sartopure® PP3

Particle Retention <sup>1</sup> (for all)	100 µm	3 µm
	50 µm	1.2 µm
	20 µm	0.65 µm
	8 µm	0.45 µm
	5 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 Size 4	Autoclavable
	Capsules Size 5	Autoclavable

## Sartopure® GF Plus

Particle Retention <sup>1</sup> (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

<sup>1</sup> nom.

## Sartopore® Air

Sterilizing Grade	0.2 µm
-------------------	--------

Filter Material	Filter Designs	Sterilization Methods
Membrane: hydrophobic Polyethersulfone	Maxicaps®	γ-irradiatable
	Midicaps®	γ-irradiatable
	Capsules Size 4	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 <span style="float: right;">[S]</span>
	Midicaps®	Autoclavable <span style="float: right;">[S]</span>
	Capsules Size 4	Autoclavable <span style="float: right;">[S]</span>
	Capsules Size 5	Autoclavable <span style="float: right;">[S]</span>

## 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-irradiated

## 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-irradiated

## 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 <span style="float: right;">[S]</span>
		Midicaps®	Autoclavable <span style="float: right;">[S]</span>
		Capsules Size 4	Autoclavable <span style="float: right;">[S]</span>

[S] delivered in sterile condition (autoclaved)





# Flexibility in Design

**Capsules**

**Connector Type O**  
 ¼" multiple stepped hose barb

**Connector Type S**  
 ¾" Tri-Clamp  
 25 mm







**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

**T-Style Maxicaps®**

**Connector Type S**  
 1½" Tri-Clamp  
 50.5 mm

**Connector Type O**  
 ½" single stepped hose barb







**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

**Hollow Fiber Lab Modules**

**Connector Inlet Type M**  
 Male Luer Lock



**Connector Outlet Type L**  
 Female Luer Lock



**Hollow Fiber Mid Scale Modules**

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


**Connector Outlet Type O**  
 ½" single stepped hose barb

**Hollow Fiber Process Modules**

**Connector Inlet Type S**  
 Sanitary 1½" Tri-Clamp  
 50.5 mm


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



**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



## Small Filters for Special Cases


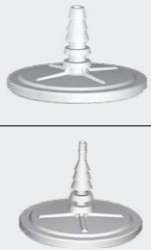


**Midisart®**

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

**Small Stepped Hose Barb**  
 4 - 12 mm

**Thread**  
 ⅜" NPT

**Tri-Clamp**  
 ¾" Tri-Clamp  
 25 mm









■ Multi-use    ■ Single-use    \* On request

**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

**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

- 83 T-Style Maxicaps®
- 53 Midicaps®
- 13 Capsule Size 4 & 5
- 73 Maxicaps®
- Hollow Fiber Modules

## Cartridge Adapter Type

- 25 | 27 | 28

## Mini Cartridge Adapter Type

- 15 | 18
- 14 (only Sartofluor® Junior)

## Pore Size Prefilter

- A No prefilter
- P Polypropylene layer
- T Single layer PTFE membrane
- E 3 µm
- F 1.2 µm
- G 0.8 µm
- D 0.65 µm
- H 0.45 µm
- I 0.35 µm
- M 0.2 µm
- V Virosart® CPV/HC
- N Virosart® Max

## Sterilization

- (T-Style) Maxicaps® | Midicaps® | Capsules
- G - γ-irradiatable | hybrid\*
- autoclavable
- Hollow Fiber Modules**
- C - prepared for γ-irradiation
- CG γ-irradiated

## Special Additions

- P Pharmaceutical grade
- GA Gas Filtration
- LG Liquid Filtration
- IT Vent valve for connecting Sartocheck® (only Maxicaps®)
- HB High Bubble Point

549	73	07	H	1	G-	SS	IT	
-----	----	----	---	---	----	----	----	--

## Filter Type

- 505 Sartopure® PP3
- 510 Sartolon
- 517 Sartofluor® HR
- 518 Sartofluor®
- 519 Sartopure® Air
- 523 Sartobran® P
- 539 Virosart® HC
- 544 Sartopure® 2
- 545 Virosart® CPV
- 547 Sartoguard PES
- 548 Sartoguard GF
- 549 Sartopure® Platinum
- 555 Sartopure® GF Plus
- 559 Sartopure® GA
- 560 Sartoclean® GF
- 562 Sartoclean® CA
- 54A Virosart® Max
- 3VI Virosart® HF
- 3V2 Virosart® Media

## Pore Size Final Filter

- 50 = 50 µm
- 20 = 20 µm
- 01 = 8 µm
- 42 = 5 µm
- 02 = 3 µm
- 03 = 1.2 µm
- 04 = 0.8 µm
- 05 = 0.65 µm
- 06 = 0.45 µm
- 07 = 0.2 µm
- 58 = 0.1 µm
- 28 = 20 nm

## Filter Size

- Cartridges | (T-Style) Maxicaps®**
- 0 5" (only cartridges)
- 1 10"
- 2 20"
- 3 30"
- 4 40" (only cartridges)
- Capsules**
- 4
- 5
- Mini Cartridges | Midicaps®**
- 7
- 8
- 9
- 0 (only Midicaps®)
- Sartoscale**
- S Sartoscale 47
- V Sartoscale 25
- Hollow Fiber Modules**
- T 1.7 cm<sup>2</sup>
- B 5 cm<sup>2</sup>
- D 200 cm<sup>2</sup>
- F 0.2 m<sup>2</sup>
- G 0.3 m<sup>2</sup>
- H 0.8 m<sup>2</sup>
- I 1 m<sup>2</sup>
- M 2.4 m<sup>2</sup>

## Connectors

- Mini | Standard Cartridges**
- Silicone O-Rings
- E EPDM O-Rings
- V Fluoroelastomer O-Rings
- (T-Style) Maxicaps® | Midicaps® | Hollow Fiber Process + mid scale Modules**
- S Tri-Clamp 50.5 mm (1½")
- O ½" single stepped hose barb
- F Tri-Clamp 25 mm (¾")
- Capsules**
- O ¼" multiple stepped hose barb
- S Tri-Clamp 25 mm (¾")
- Sartoscale 47**
- F Tri-Clamp 25mm (¾")
- Sartoscale 25**
- L Female Luer Lock
- X Male Luer Slip
- Hollow Fiber Lab Modules**
- L Female Luer Lock
- M Male Luer Slip

## Pack Size

- blank box of 1
- V box of 2
- M box of 3
- A box of 4
- B box of 5
- C box of 6
- D box of 10
- G box of 25
- H box of 40
- K box of 50

\* hybrid = autoclavable or γ-irradiatable

Note: the first code letter represents the inlet, the second the outlet.

# Sartopore® Family

## Sartopore® Platinum

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

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	<b>Pore Size</b>	<b>1.0 m<sup>2</sup>   10.8 ft<sup>2</sup></b>	<b>2.0 m<sup>2</sup>   21.5 ft<sup>2</sup></b>	<b>3.0 m<sup>2</sup>   32.3 ft<sup>2</sup></b>
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-***x	5497307H2G-***x	5497307H3G-***x

		Size 7	Size 8	Size 9	Size 0
	<b>Pore Size</b>	<b>0.065 m<sup>2</sup>   0.7 ft<sup>2</sup></b>	<b>0.13 m<sup>2</sup>   1.4 ft<sup>2</sup></b>	<b>0.26 m<sup>2</sup>   2.8 ft<sup>2</sup></b>	<b>0.52 m<sup>2</sup>   5.6 ft<sup>2</sup></b>
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
	<b>Pore Size</b>	<b>0.021 m<sup>2</sup>   0.22 ft<sup>2</sup></b>
Capsules	0.45 + 0.2	5491307H4-**-B
Gamma Capsules	0.45 + 0.2	5491307H4G-**-B

		25	47
	<b>Pore Size</b>	<b>4.5 cm<sup>2</sup></b>	<b>17.3 cm<sup>2</sup></b>
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 | D = 10 pieces  
 \*\* = IT vent valve design suitable for connection of Sartocheck® integrity tester

## Sartopore® Platinum HB

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	<b>Pore Size</b>	<b>1.0 m<sup>2</sup>   10.8 ft<sup>2</sup></b>	<b>2.0 m<sup>2</sup>   21.5 ft<sup>2</sup></b>	<b>3.0 m<sup>2</sup>   32.3 ft<sup>2</sup></b>
Cartidges	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
	<b>Pore Size</b>	<b>0.065 m<sup>2</sup>   0.7 ft<sup>2</sup></b>	<b>0.13 m<sup>2</sup>   1.4 ft<sup>2</sup></b>	<b>0.26 m<sup>2</sup>   2.8 ft<sup>2</sup></b>	<b>0.52 m<sup>2</sup>   5.6 ft<sup>2</sup></b>
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
	<b>Pore Size</b>	<b>0.021 m<sup>2</sup>   0.22 ft<sup>2</sup></b>
Capsules	0.45 + 0.2	5491307H4- -**HBB
Gamma Capsules	0.45 + 0.2	5491307H4G- -**HBB

		25	47
	<b>Pore Size</b>	<b>4.5 cm<sup>2</sup></b>	<b>17.3 cm<sup>2</sup></b>
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 | D = 10 pieces

## Sartopore® 2

		Size 0 (5")	Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.3 m <sup>2</sup>   3.2 ft <sup>2</sup>	0.6 m <sup>2</sup>   6.5 ft <sup>2</sup>	1.2 m <sup>2</sup>   12.9 ft <sup>2</sup>	1.8 m <sup>2</sup>   19.4 ft <sup>2</sup>
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 <sup>2</sup>   0.5 ft <sup>2</sup>	0.1 m <sup>2</sup>   1.1 ft <sup>2</sup>	0.2 m <sup>2</sup>   2.2 ft <sup>2</sup>	0.45 m <sup>2</sup>   4.8 ft <sup>2</sup>
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 <sup>2</sup>   0.16 ft <sup>2</sup>	0.03 m <sup>2</sup>   0.32 ft <sup>2</sup>
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 <sup>2</sup>	17.3 cm <sup>2</sup>
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 | D = 10 pieces  
 \*\* = 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 <sup>2</sup>   8.6 ft <sup>2</sup>	1.6 m <sup>2</sup>   17.2 ft <sup>2</sup>	2.4 m <sup>2</sup>   25.8 ft <sup>2</sup>
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 <sup>2</sup>   0.7 ft <sup>2</sup>	0.13 m <sup>2</sup>   1.4 ft <sup>2</sup>	0.26 m <sup>2</sup>   2.8 ft <sup>2</sup>	0.52 m <sup>2</sup>   5.6 ft <sup>2</sup>
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 <sup>2</sup>   0.22 ft <sup>2</sup>
Capsules	0.8 + 0.2	5441307G4- ** -B
Gamma Capsules	0.8 + 0.2	5441307G4G- ** -B

		25	47
	Pore Size	4.5 cm <sup>2</sup>	17.3 cm <sup>2</sup>
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 | D = 10 pieces

## Sartopore® 2 XLI

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size	0.8 m <sup>2</sup>   8.6 ft <sup>2</sup>	1.6 m <sup>2</sup>   17.2 ft <sup>2</sup>	2.4 m <sup>2</sup>   25.8 ft <sup>2</sup>
Cartridges	0.35 + 0.2	544250711	544250712	544250713
Maxicaps®	0.35 + 0.2	544730711- -** <sup>xx</sup>	544730712- -** <sup>xx</sup>	544730713- -** <sup>xx</sup>
Gamma Maxicaps®	0.35 + 0.2	544730711G- -** <sup>xx</sup>	544730712G- -** <sup>xx</sup>	544730713G- -** <sup>xx</sup>
T-Style Maxicaps® γ-irradiatable or autoclavable	0.35 + 0.2	544830711G- -**	544830712G- -**	544830713G- -**

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.065 m <sup>2</sup>   0.7 ft <sup>2</sup>	0.13 m <sup>2</sup>   1.4 ft <sup>2</sup>	0.26 m <sup>2</sup>   2.8 ft <sup>2</sup>	0.52 m <sup>2</sup>   5.6 ft <sup>2</sup>
Midicaps®	0.35 + 0.2	544530717- -** -A	544530718- -** -A	544530719- -** -A	544530710- -** -V
Gamma Midicaps®	0.35 + 0.2	544530717G- -** -A	544530718G- -** -A	544530719G- -** -A	544530710G- -** -V

		Size 4
	Pore Size	0.021 m <sup>2</sup>   0.22 ft <sup>2</sup>
Capsules	0.35 + 0.2	544130714- -** -B
Gamma Capsules	0.35 + 0.2	544130714G- -** -B

		25	47
	Pore Size	4.5 cm <sup>2</sup>	17.3 cm <sup>2</sup>
Sartoscale	0.35 + 0.2	54453071V- -LX- -C	54453071S- -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 | D = 10 pieces  
<sup>xx</sup> = 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 <sup>2</sup>   4.3 ft <sup>2</sup>	0.8 m <sup>2</sup>   8.6 ft <sup>2</sup>	1.6 m <sup>2</sup>   17.2 ft <sup>2</sup>	2.4 m <sup>2</sup>   25.8 ft <sup>2</sup>
Cartridges	0.2 + 0.1	5442558M0	5442558M1	5442558M2	5442558M3
Maxicaps®	0.2 + 0.1		5447358M1- -** <sup>xx</sup>	5447358M2- -** <sup>xx</sup>	5447358M3- -** <sup>xx</sup>
Gamma Maxicaps®	0.2 + 0.1		5447358M1G- -** <sup>xx</sup>	5447358M2G- -** <sup>xx</sup>	5447358M3G- -** <sup>xx</sup>
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 <sup>2</sup>   0.7 ft <sup>2</sup>	0.13 m <sup>2</sup>   1.4 ft <sup>2</sup>	0.26 m <sup>2</sup>   2.8 ft <sup>2</sup>	0.52 m <sup>2</sup>   5.6 ft <sup>2</sup>
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 <sup>2</sup>   0.22 ft <sup>2</sup>
Capsules	0.2 + 0.1	5441358M4- -** -B
Gamma Capsules	0.2 + 0.1	5441358M4G- -** -B

		25	47
	Pore Size	4.5 cm <sup>2</sup>	17.3 cm <sup>2</sup>
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 | D = 10 pieces  
<sup>xx</sup> = IT vent valve design suitable for connection of Sartocheck® integrity tester

## Sartopore® Air

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

		Size 7	Size 8	Size 9	Size 0
	Pore Size	0.06 m <sup>2</sup>   0.54 ft <sup>2</sup>	0.1 m <sup>2</sup>   1.1 ft <sup>2</sup>	0.17 m <sup>2</sup>   1.83 ft <sup>2</sup>	0.34 m <sup>2</sup>   3.66 ft <sup>2</sup>
Midicaps® γ-irradiatable	0.2	5195307A7G-**-A	5195307A8G-**-A	5195307A9G-**-A	5195307A0G-**-V

		Size 4
	Pore Size	0.016 m <sup>2</sup>   0.17 ft <sup>2</sup>
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 | C = 6 pieces | V = 2 pieces | M = 3 pieces | D = 10 pieces

## Sartoguard Family

### Sartoguard PES

		Size 1 (10")	Size 2 (20")	Size 3 (30")			
	Pore Size <sup>1</sup>	0.8 m <sup>2</sup>   8.6 ft <sup>2</sup>	1.6 m <sup>2</sup>   17.2 ft <sup>2</sup>	2.4 m <sup>2</sup>   25.8 ft <sup>2</sup>			
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 <sup>1</sup>	0.065 m <sup>2</sup>   0.7 ft <sup>2</sup>	0.13 m <sup>2</sup>   1.4 ft <sup>2</sup>	0.26 m <sup>2</sup>   2.8 ft <sup>2</sup>	0.52 m <sup>2</sup>   5.6 ft <sup>2</sup>
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 <sup>1</sup>	0.021 m <sup>2</sup>   0.22 ft <sup>2</sup>
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 <sup>1</sup>	4.5 cm <sup>2</sup>	17.3 cm <sup>2</sup>
Sartoscale	0.8 + 0.1	5475358GV- -LX- -C	5475358GS- -FF- -M
	1.2 + 0.2	5475307FV- -LX- -C	5475307FS- -FF- -M

<sup>1</sup>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 | D = 10 pieces

## Sartoguard GF

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

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

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

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

<sup>1</sup> 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 | D = 10 pieces

## Sartobran<sup>®</sup> P Family

### Sartobran<sup>®</sup> P

		Size 0 (5")	Size 1 (10")	Size 2 (20")	Size 3 (30")
	<b>Pore Size</b>	<b>0.3 m<sup>2</sup>   3.2 ft<sup>2</sup></b>	<b>0.6 m<sup>2</sup>   6.5 ft<sup>2</sup></b>	<b>1.2 m<sup>2</sup>   12.9 ft<sup>2</sup></b>	<b>1.8 m<sup>2</sup>   19.4 ft<sup>2</sup></b>
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 <sup>®</sup>	0.65 + 0.45		5237306D1- -** <sup>xx</sup>	5237306D2- -** <sup>xx</sup>	5237306D3- -** <sup>xx</sup>
	0.45 + 0.2		5237307H1- -** <sup>xx</sup>	5237307H2- -** <sup>xx</sup>	5237307H3- -** <sup>xx</sup>
T-Style Maxicaps <sup>®</sup>	0.65 + 0.45		5238306D1- -**	5238306D2- -**	5238306D3- -**
	0.45 + 0.2		5238307H1- -**	5238307H2- -**	5238307H3- -**

		Size 7	Size 8	Size 9	Size 0
	<b>Pore Size</b>	<b>0.05 m<sup>2</sup>   0.54 ft<sup>2</sup></b>	<b>0.1 m<sup>2</sup>   1.1 ft<sup>2</sup></b>	<b>0.2 m<sup>2</sup>   2.2 ft<sup>2</sup></b>	<b>0.45 m<sup>2</sup>   4.8 ft<sup>2</sup></b>
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 <sup>®</sup>	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
	<b>Pore Size</b>	<b>0.015 m<sup>2</sup>   0.16 ft<sup>2</sup></b>	<b>0.03 m<sup>2</sup>   0.32 ft<sup>2</sup></b>
Capsules	0.65 + 0.45		5231306D5- -OO- -B
	0.45 + 0.2	5231307H4- ----B	5231307H5- -OO- -B

		25	47
	<b>Pore Size</b>	<b>4.5 cm<sup>2</sup></b>	<b>17.3 cm<sup>2</sup></b>
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 | D = 10 pieces

<sup>xx</sup> = IT vent valve design suitable for connection of Sartocheck<sup>®</sup> integrity tester



# Sartoclean® Family

## Sartoclean® CA

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size <sup>1</sup>	0.6 m <sup>2</sup>   6.5 ft <sup>2</sup>	1.2 m <sup>2</sup>   12.8 ft <sup>2</sup>	1.8 m <sup>2</sup>   19.3 ft <sup>2</sup>
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 <sup>1</sup>	0.05 m <sup>2</sup>   0.5 ft <sup>2</sup>	0.1 m <sup>2</sup>   1.1 ft <sup>2</sup>	0.2 m <sup>2</sup>   2.2 ft <sup>2</sup>	0.45 m <sup>2</sup>   4.8 ft <sup>2</sup>
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 <sup>1</sup>	4.5 cm <sup>2</sup>	17.3 cm <sup>2</sup>
Sartoscale	3.0 + 0.8	5625304EV- -LX- -C	5625304ES- -FF- -M
	0.8 + 0.65	5625305GV- -LX- -C	5625305GS- -FF- -M

<sup>1</sup> 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 | D = 10 pieces

## Sartoclean® GF

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	Pore Size <sup>1</sup>	0.47 m <sup>2</sup>   5.1 ft <sup>2</sup>	0.94 m <sup>2</sup>   10.2 ft <sup>2</sup>	1.41 m <sup>2</sup>   15.3 ft <sup>2</sup>
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 <sup>1</sup>	0.05 m <sup>2</sup>   0.54 ft <sup>2</sup>	0.08 m <sup>2</sup>   0.85 ft <sup>2</sup>	0.16 m <sup>2</sup>   1.7 ft <sup>2</sup>	0.32 m <sup>2</sup>   3.4 ft <sup>2</sup>
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 <sup>1</sup>	4.5 cm <sup>2</sup>	17.3 cm <sup>2</sup>
Sartoscale	3.0 + 0.8	5605304EV- -LX- -C	5605304ES- -FF- -M
	0.8 + 0.65	5605305GV- -LX- -C	5605305GS- -FF- -M

<sup>1</sup> 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 | D = 10 pieces

# Sartolon Family

## Sartolon

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

Size 9		
	<b>Pore Size</b>	<b>0.2 m<sup>2</sup>   2.2 ft<sup>2</sup></b>
Mini Cartridges	0.45 + 0.2	5101507H9- - - - -B
Midicaps <sup>®</sup>	0.45 + 0.2	5105307H9- -** -A

		25	47
	<b>Pore Size</b>	<b>4.5 cm<sup>2</sup></b>	<b>17.3 cm<sup>2</sup></b>
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 | V = 2 pieces | M = 3 pieces | D = 10 pieces  
 \*\* = IT vent valve design suitable for connection of Sartocheck<sup>®</sup> integrity tester

# Sartopure<sup>®</sup> Family

## Sartopure<sup>®</sup> PP3 0.45 μm

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

		Size 7	Size 8	Size 9	Size 0
	<b>Retention Rate<sup>1</sup></b>	<b>0.04 m<sup>2</sup>   0.43 ft<sup>2</sup></b>	<b>0.08 m<sup>2</sup>   0.86 ft<sup>2</sup></b>	<b>0.13 m<sup>2</sup>   1.4 ft<sup>2</sup></b>	<b>0.26 m<sup>2</sup>   2.8 ft<sup>2</sup></b>
Midicaps <sup>®</sup>	0.45	5055306P7- -** -A	5055306P8- -** -A	5055306P9- -** -A	5055306P0- -** -V

## Sartopure<sup>®</sup> PP3 0.65 – 3 μm

		Size 1 (10")	Size 2 (20")	Size 3 (30")	Size 4 (40")
	<b>Retention Rate<sup>1</sup></b>	<b>0.45 m<sup>2</sup>   4.8 ft<sup>2</sup></b>	<b>0.9 m<sup>2</sup>   9.7 ft<sup>2</sup></b>	<b>1.35 m<sup>2</sup>   14.5 ft<sup>2</sup></b>	<b>1.8 m<sup>2</sup>   19.4 ft<sup>2</sup></b>
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 <sup>®</sup>	3	5058302P1- -**	5058302P2- -**	5058302P3- -**	
	1.2	5058303P1- -**	5058303P2- -**	5058303P3- -**	
	0.65	5058305P1- -**	5058305P2- -**	5058305P3- -**	
Maxicaps <sup>®</sup>	3	5057302P1- -**	5057302P2- -**	5057302P3- -**	
	1.2	5057303P1- -**	5057303P2- -**	5057303P3- -**	
	0.65	5057305P1- -**	5057305P2- -**	5057305P3- -**	

		Size 7	Size 8	Size 9	Size 0
	<b>Retention Rate<sup>1</sup></b>	<b>0.05 m<sup>2</sup>   0.54 ft<sup>2</sup></b>	<b>0.09 m<sup>2</sup>   0.97 ft<sup>2</sup></b>	<b>0.15 m<sup>2</sup>   1.61 ft<sup>2</sup></b>	<b>0.30 m<sup>2</sup>   3.23 ft<sup>2</sup></b>
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 <sup>®</sup>	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

<sup>1</sup> 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 | D = 10 pieces

## Sartopure® PP3 5 – 100 µm

		Size 1 (10")	Size 2 (20")	Size 3 (30")	Size 4 (40")
	Retention Rate <sup>1</sup>	0.65 m <sup>2</sup>   7 ft <sup>2</sup>	1.3 m <sup>2</sup>   14 ft <sup>2</sup>	1.95 m <sup>2</sup>   21 ft <sup>2</sup>	2.6 m <sup>2</sup>   28 ft <sup>2</sup>
Cartridges	100	505**98P1	505**98P2	505**98P3	
	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®	100	5058398P1- **	5058398P2- **	5058398P3- **	
	50	5058350P1- **	5058350P2- **	5058350P3- **	
	20	5058320P1- **	5058320P2- **	5058320P3- **	
	8	5058301P1- **	5058301P2- **	5058301P3- **	
	5	5058342P1- **	5058342P2- **	5058342P3- **	
Maxicaps®	100	5057398P1- **	5057398P2- **	5057398P3- **	
	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 <sup>1</sup>	0.065 m <sup>2</sup>   0.7 ft <sup>2</sup>	0.12 m <sup>2</sup>   1.3 ft <sup>2</sup>	0.21 m <sup>2</sup>   2.26 ft <sup>2</sup>	0.42 m <sup>2</sup>   4.52 ft <sup>2</sup>
Mini Cartridges	20			5051520P9- - - - -B	
	8			5051501P9- - - - -B	
	5	5051542P7- - - - -B	5051542P8- - - - -B	5051542P9- - - - -B	
Midicaps®	100	5055398P7- ** -A	5055398P8- ** -A	5055398P9- ** -A	5055398P0- ** -V
	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

<sup>1</sup>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 | D = 10 pieces

## Sartopure® PP3 Capsules & Sartoscale

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

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

		Size 4	Size 5
	Retention Rate <sup>1</sup>	0.018 m <sup>2</sup>   0.19 ft <sup>2</sup>	0.036 m <sup>2</sup>   0.39 ft <sup>2</sup>
Capsules	100	5051398P4- -SS- -B	
	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 <sup>1</sup>	4.5 cm <sup>2</sup>	17.3 cm <sup>2</sup>
Sartoscale	0.45	5055306PV- -LX- -C	5055306PS- -FF- -M
	0.65	5055305PV- -LX- -C	5055305PS- -FF- -M
	1.2	5055303PV- -LX- -C	5055303PS- -FF- -M
	3	5055302PV- -LX- -C	5055302PS- -FF- -M
	5	5055342PV- -LX- -C	5055342PS- -FF- -M
	8	5055301PV- -LX- -C	5055301PS- -FF- -M
	20	5055320PV- -LX- -C	5055320PS- -FF- -M
	50	5055350PV- -LX- -C	5055350PS- -FF- -M
	100	5055398PV- -LX- -C	5055398PS- -FF- -M

<sup>1</sup>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 | D = 10 pieces

## Sartopure® GF Plus

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

		Size 7	Size 8	Size 9	Size 0
	<b>Retention Rate<sup>1</sup></b>	<b>0.04 m<sup>2</sup>   0.4 ft<sup>2</sup></b>	<b>0.08 m<sup>2</sup>   0.9 ft<sup>2</sup></b>	<b>0.12 m<sup>2</sup>   1.3 ft<sup>2</sup></b>	<b>0.25 m<sup>2</sup>   2.7 ft<sup>2</sup></b>
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
	<b>Retention Rate<sup>1</sup></b>	<b>4.5 cm<sup>2</sup></b>	<b>17.3 cm<sup>2</sup></b>
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")
	<b>Retention Rate<sup>1</sup></b>	<b>0.7 m<sup>2</sup>   7.5 ft<sup>2</sup></b>	<b>1.4 m<sup>2</sup>   15.1 ft<sup>2</sup></b>	<b>2.1 m<sup>2</sup>   22.6 ft<sup>2</sup></b>
Cartridges	0.2	559**07P1----GA	559**07P2----GA	559**07P3----GA

		Size 9
	<b>Retention Rate<sup>1</sup></b>	<b>0.2 m<sup>2</sup>   2.2 ft<sup>2</sup></b>
Mini Cartridges	0.2	5591507P9---GAB

<sup>1</sup> 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 | D = 10 pieces

# Sartofluor® Family

## Sartofluor®

		Size 0 (5")	Size 1 (10")	Size 2 (20")	Size 3 (30")
	<b>Pore Size</b>	<b>0.375 m<sup>2</sup>   4.05 ft<sup>2</sup></b>	<b>0.75 m<sup>2</sup>   8.1 ft<sup>2</sup></b>	<b>1.5 m<sup>2</sup>   16.1 ft<sup>2</sup></b>	<b>2.25 m<sup>2</sup>   24.2 ft<sup>2</sup></b>
Cartridges GA	0.2	518**07T0----GA	518**07T1----GA	518**07T2----GA	518**07T3----GA
	<b>Pore Size</b>		<b>0.5 m<sup>2</sup>   5.4 ft<sup>2</sup></b>	<b>1.0 m<sup>2</sup>   10.8 ft<sup>2</sup></b>	<b>1.5 m<sup>2</sup>   16.1 ft<sup>2</sup></b>
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
	<b>Pore Size</b>	<b>0.05 m<sup>2</sup>   0.5 ft<sup>2</sup></b>	<b>0.1 m<sup>2</sup>   1.1 ft<sup>2</sup></b>	<b>0.2 m<sup>2</sup>   2.2 ft<sup>2</sup></b>	<b>0.45 m<sup>2</sup>   4.8 ft<sup>2</sup></b>
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
	<b>Pore Size</b>	<b>0.015 m<sup>2</sup>   0.16 ft<sup>2</sup></b>	<b>0.03 m<sup>2</sup>   0.32 ft<sup>2</sup></b>
Capsules	0.2	5181307T4-**-B	5181307T5--OO--D

## Sartofluor® HR

		Size 1 (10")	Size 2 (20")	Size 3 (30")
	<b>Pore Size</b>	<b>0.56 m<sup>2</sup>   6.7 ft<sup>2</sup></b>	<b>1.12 m<sup>2</sup>   12.1 ft<sup>2</sup></b>	<b>1.68 m<sup>2</sup>   18.1 ft<sup>2</sup></b>
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 <sup>2</sup>   8.6 ft <sup>2</sup>	2.4 m <sup>2</sup>   25.8 ft <sup>2</sup>
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 <sup>2</sup>   0.22 ft <sup>2</sup>	0.2 m <sup>2</sup>   2.15 ft <sup>2</sup>
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 <sup>2</sup>	5 cm <sup>2</sup>
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 <sup>2</sup>   10.8 ft <sup>2</sup>
Process module γ-irradiatable	20 nm	3V2--28-IV-SF
Sterile (γ-irradiated)	20 nm	3V2--28-IVGSF

	Pore Size	0.3 m <sup>2</sup>   3.2 ft <sup>2</sup>
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 <sup>2</sup>
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 <sup>2</sup>   7.5 ft <sup>2</sup>	1.4 m <sup>2</sup>   15 ft <sup>2</sup>	2.1 m <sup>2</sup>   22.6 ft <sup>2</sup>
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 <sup>2</sup>   2.15 ft <sup>2</sup>
Midicaps® (Autoclavable)	20 nm	5455328V9--FF--V

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

		5 cm <sup>2</sup>
	Pore Size	
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 <sup>2</sup>   9.7 ft <sup>2</sup>	1.8 m <sup>2</sup>   19.4 ft <sup>2</sup>	2.7 m <sup>2</sup>   29 ft <sup>2</sup>
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 <sup>2</sup>   2.9 ft <sup>2</sup>
Midicaps® (Autoclavable)	20 nm	5395328V9--FF--V

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

		5 cm <sup>2</sup>
	Pore Size	
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")
	<b>Pore Size</b>	0.7 m <sup>2</sup>   7.5 ft <sup>2</sup>	1.4 m <sup>2</sup>   15 ft <sup>2</sup>	2.1 m <sup>2</sup>   22.6 ft <sup>2</sup>
Cartridge (Autoclavable   Steamable)	0.1 μm	54A2558N1	54A2558N2	54A2558N3
T-Style Maxicaps® (Autoclavable)	0.1 μm	54A8358N1--SS	54A8358N2--SS	54A8358N3--SS

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

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

		5 cm <sup>2</sup>	21 cm <sup>2</sup>
Minisart®   Sartolab Sterile (Autoclaved)	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)

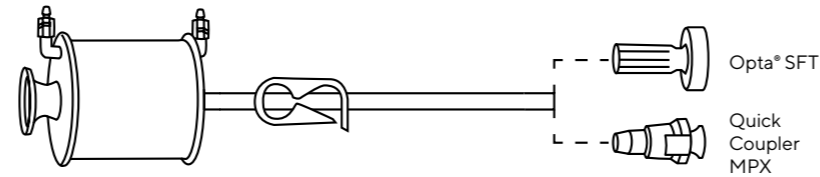
Sartopore® Platinum

Sartopore® 2, XLM and XLG

500 mm TPE tubing

Opta®

MPX Connector



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

## 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® Platinum 0.45 | 0.2 µm

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

## Sartopore® 2 XLG 0.8 | 0.2 µm

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

## Sartopore® 2 0.45 | 0.2 µm

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

## Sartopore® 2 XLM 0.2 | 0.1 µm

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

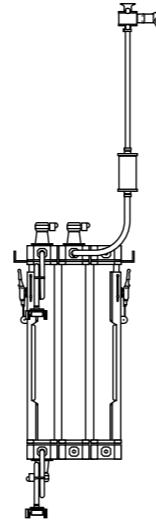
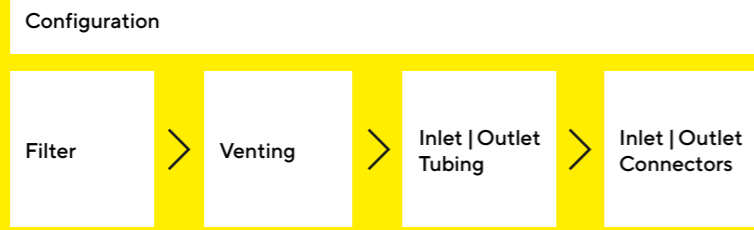
C-Flex® is a registered trademark of Saint-Gobain Performance Plastics Corporation.



# Maxicaps® MR

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

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



## 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	Polyethersulfone	Sterile	2-9	4.8 m²   51.6 ft²	21.6 m²   232.2 ft²
Sartoguard GF	0.1 µm   0.2 µm	Polyethersulfone	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	Polyethersulfone	Sterile	2-9	3.6 m²   38.8 ft²	16.2 m²   174.6 ft²
Sartopore® 2 XLG	0.2 µm	Polyethersulfone	Sterile	2-9	4.8 m²   51.6 ft²	21.6 m²   232.2 ft²
Sartopore® 2 XLI	0.2 µm	Polyethersulfone	Sterile	2-9	4.8 m²   51.6 ft²	21.6 m²   232.2 ft²
Sartopore® 2 XLM	0.1 µm	Polyethersulfone	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

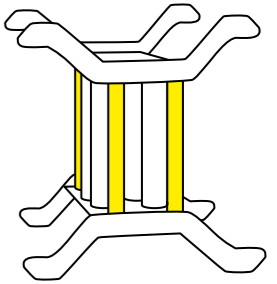
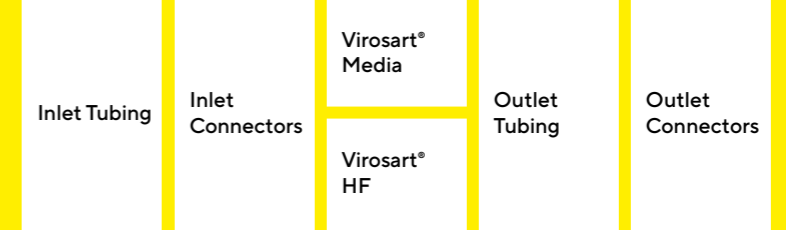
Venting	Inlet   Outlet Tubing	Inlet   Outlet Connectors
Basic	Tuflux® SIL   750 mm	AseptiQuik® G
Advanced	Sani-Tech® SHT-R®   750 mm	AseptiQuik® L
	Pharma-50   750 mm	Opta® SFT Female
	Tuflux® TPE   1,500 mm	Opta® SFT Male
	C-Flex® 374   1,500 mm	Presto HB
		1.5" Sanitary Clamp
		Plug for TPE tubing

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. C-Flex® and Sani-Tech® SHT-R are registered trademarks 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



## 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	2 x AseptiQuik® G on Tri-Clamp 1½"	Reinforced silicon tubing	2 x AseptiQuik® G on Tri-Clamp 1½"
Virosart® HF	Reinforced silicon tubing	2 x AseptiQuik® G on Tri-Clamp 1½"	Reinforced silicon tubing	2 x AseptiQuik® G on Tri-Clamp 1½" 1 x AseptiQuik® G on Tri-Clamp ¾", Tri-Clamp 1½", and flow restrictor

AseptiQuik® is a registered trademark of Colder Products Company.




**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 further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)