

Syringe Filters – Sartorius Minisart® Selection Guide:

How to Choose the Optimal Housing and Membrane Material for Your Application.

Innovative design features, coupled with the largest surface areas and high flow rates make Minisart® syringe filters the ideal choice for all your syringe filtration needs.

- Largest effective filtration area (EFA)
- Low adsorption
- Low hold-up volume
- Gamma/EO sterilized
- High throughput
- For sterile filtration, analytical sample preparation, and clarification of media additives, buffers, chemical reagents and even gases!

1. Sample Composition

Aqueous		Aqueous Solvents		
Hydrophilic		Hydrophilic		Hydrophobic
For Use With:				
<ul style="list-style-type: none"> ▪ Buffers ▪ Protein analysis 	<ul style="list-style-type: none"> ▪ Tissue culture media 	<ul style="list-style-type: none"> ▪ Aqueous solvent mixtures solvents 	<ul style="list-style-type: none"> ▪ Solvent mixtures solvents 	<ul style="list-style-type: none"> ▪ Solvents gases acids bases
Recommended Filter Material:				
CA	PES	RC	NY	PTFE
Cellulose acetate	Polyethersulfone	Regenerated cellulose	Polyamide, nylon	Polytetrafluoroethylene



2. Pore Sizes

Sterile Filtration		Sample preparation / clarification / particle removal				Prefiltration	
For Use With:							
<ul style="list-style-type: none"> ▪ Small bacteria ▪ Mycoplasma ▪ Colloids >0.1 µm 	<ul style="list-style-type: none"> ▪ 0.2 µm – UHPLC, etc. (Columns < 3 µm particles) bacteria 	<ul style="list-style-type: none"> ▪ HPLC, etc. (Columns >3 µm particles) particles 	<ul style="list-style-type: none"> ▪ Particles ▪ Yeast cells 	<ul style="list-style-type: none"> ▪ Particles ▪ Yeast cells 	<ul style="list-style-type: none"> ▪ Particles ▪ Yeast cells ▪ Platelets 	<ul style="list-style-type: none"> ▪ Large particles ▪ Cells 	<ul style="list-style-type: none"> ▪ Glass prefilter ▪ Glass + membrane ▪ Highly particle-laden samples
Recommended Pore Size:							
0.1 µm	0.2 µm	0.45 µm	0.65 µm	0.8 µm	1.2 µm	5 µm	GF (Glass Fiber)

3. Sample Volumes

For Use With:			
<ul style="list-style-type: none"> ▪ 1 mL to 200 mL 	<ul style="list-style-type: none"> ▪ 1 mL to 100 mL 	<ul style="list-style-type: none"> ▪ 0.5 mL to 15 mL 	<ul style="list-style-type: none"> ▪ 0.05 mL to 1 mL
Recommended Diameter:			
28 mm for up to 200 mL	25 mm for up to 100 mL	15 mm for up to 15 mL	4 mm for up to 1 mL

