

Whatman Autovial

Syringeless Filters

Instructions for use

Introduction

Important

Read these instructions carefully before using the products.

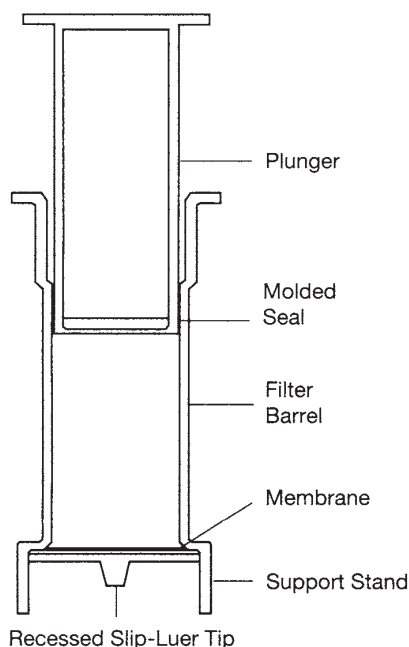
Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

Description

Autovial™ Syringeless Filters are preassembled filtration devices for removing particulates from samples. They replace syringe-coupled filtration devices with a single, disposable unit.

Each Autovial consists of two parts: A filter-barrel and a plunger. The proven design features an integral filter, built-in air purge and support stand which protects the recessed slip-luer tip.



Step	Action
1	Select the Autovial Syringeless Filters that is compatible with your sample. (See Chemical Compatibility of Membrane, on page 2). If your solvent is not listed, please contact Cytiva. Stand the Autovial filter-barrel on the bench top and pour in the sample. Each Autovial has 12 mL capacity.
2	Select an Autovial plunger and place it into the filter-barrel until the tip is securely in place. There will be a gap of air between the sample and the plunger. It is not necessary to push the plunger more than one half inch into the Autovial filter-barrel.
3	Place the tip of the Autovial into the mouth of an auto-sampler vial or container. Compress the Autovial plunger. Filtration will begin immediately. Continue to compress the plunger until it reaches the bottom of the Autovial filter-barrel to purge the membrane.

For Direct Instrument Injection

Place a needle on the Autovial slip-luer outlet for direct injection into the instrument.

Membrane Guide

Type	Applications
Cellulose Acetate (CA)	Low nonspecific protein binding membrane for samples in aqueous solutions.
Glass Microfiber (GMF)	Low nonspecific protein binding membrane for samples in aqueous or organic solutions within a pH range of 3-10.
Mixed Esters	0.65 Membrane for use in food industry applications.
Nylon (NYL)	Membrane for aqueous and organic samples within a pH range of 3-10.
Polyethersulfone (PES)	Low nonspecific protein binding membrane for samples in aqueous solutions.
polypropylene (PP)	For aggressive samples and high debris.
Polytetrafluoroethylene (PTFE)	PTFE membrane for aggressive samples or with >50% organic solvent.

Type	Applications
Polyvinylidene Fluoride (PVDF)	Low nonspecific protein binding membrane for samples in aqueous and/or organic solvents.

Chemical Compatibility of Membrane

Solvent	CA	GMF	NYL	PP	PES	PTF E	PVD F
Acetic Acid 5% +	LR	R	R	R	R	R	R
Acetic Acid, Glacial	NR	R	LR	R	R	R	R
Acetone	NR	R	R	R	NR	R	NR
Acetonitrile	NR	R	R	R	NR	R	R
Ammonia, 6N	-	LR	R	R	R	R	LR
Amyl Acetate	NR	R	R	R	NR	R	LR
Amyl Alcohol	R	R	R	R	R	R	R
Benzine ¹	R	R	LR	LR	NR	R	R
Benzyl Alcohol ¹	LR	R	LR	R	NR	R	R
Boric Acid	R	R	LR	R	R	R	R
Butyl Alcohol	R	R	R	R	R	R	R
Butyl Chloride ¹	-	R	NR	NR	-	R	R
Carbon Tetrachloride ¹	NR	R	LR	LR	NR	R	R
Chlorobenzene	-	R	NR	LR	NR	R	R
Chloroform ¹	NR	R	NR	LR	NR	R	R
Citric Acid	-	R	NR	-	-	R	R
Cresol	NR	R	NR	R	NR	R	NR
Cyclohexane	R	R	R	R	R	R	R
Cyclohexanone	NR	R	NR	R	NR	R	R
Diethyl Acetamide	NR	R	R	R	NR	R	NR
Dimethyl Formamide	NR	R	R	R	NR	R	NR
Dioxane	NR	R	R	R	NR	R	LR
DMSO	NR	R	R	R	NR	R	LR
Ethanol	R	R	R	R	R	R	NR
Ethers	LR	R	R	R	R	R	LR
Ethyl Acetate	NR	R	R	R	NR	R	LR
Ethylene Glycol	R	R	R	R	R	R	R
Formaldehyde	LR	R	R	R	R	R	R
Formic Acid	LR	R	NR	R	LR	R	R
Freon TF	R	R	R	R	R	R	R
Hexane	R	R	R	R	R	R	R
Hydrochloric Acid (Conc)	LR	R	NR	LR	R	R	R
Hydrofluoric Acid	NR	NR	NR	LR	-	R	R
Isobutyl Alcohol	R	R	R	R	R	R	R
Isopropyl Acetate	NR	R	R	R	NR	R	R
Methanol	R	R	R	R	R	R	R
Methyl Ethyl Ketone	LR	R	R	R	NR	R	LR
Methylene Chloride ¹	R	R	R	R	R	R	R
Nitric Acid (Conc)	LR	R	NR	NR	NR	R	NR
Nitric Acid, 6N	LR	R	NR	LR	LR	R	LR

Solvent	CA	GMF	NYL	PP	PES	PTF E	PVD F
Nitrobenzene ¹	NR	R	LR	R	LR	R	R
Pentane	R	R	R	LR	R	R	R
Perchloro Ethylene	R	R	R	R	NR	R	R
Phenol (0.5%)	R	R	R	R	LR	R	R
Pyridine	NR	R	LR	R	NR	R	R
Sodium Hydroxide, 6N	NR	NR	LR	R	R	R	NR
Sulfuric Acid (Conc)	NR	R	NR	R	NR	R	R
Tetrahydrofuran	NR	R	R	LR	NR	R	R
Toluene ¹	LR	R	LR	LR	NR	R	R
Trichloroethane ¹	NR	R	LR	R	NR	R	R
Trichloroethylene ¹	R	R	NR	R	NR	R	R
Water	R	R	R	R	R	R	R
Xylene ¹	R	R	LR	LR	NR	R	R

¹ Short term resistance of housing

(R = Resistant; LR = Limited Resistance; NR = Non Resistant; - = No data).

Ordering Information - Autovial Syringeless Filters

Autovial 12 - No Prefilter

PES Filtration Media

Product code AV125NPUSU

Autovial 12 Syringeless Filter, no prefilter, PES, 0.45µm 50/pk

PVDF Filtration Media

Product code AV125NPUAQU

Autovial 12 Syringeless Filter, no prefilter, PVDF, 0.45µm 50/pk

Autovial 12 - With Glass Prefilter

CA Filtration Media

Product code AV125UCA

Autovial 12 Syringeless Filter, glass prefilter, CA, 0.45 µm 50/pk

Nylon Filtration Media

Product code AV125SNAO

Autovial 12 Syringeless Filter, glass prefilter, Sterile, Nylon, 0.2µm 40/pk

Product code AV125ENAO

Autovial 12 Syringeless Filter, glass prefilter, Nylon, 0.2 µm 50/pk

Product code AV125UNAO

Autovial 12 Syringeless Filter, glass prefilter, Nylon, 0.45 µm 50/pk

Product code AV525UNAO

Autovial 12 Syringeless Filter, glass prefilter, Nylon, 0.45 µm 1000/pk

PVDF Filtration Media

Product code AV125SAQU

Autovial 12 Syringeless Filter, glass prefilter, Sterile, PVDF, 0.2 µm 40/pk

Product code AV125EAQU

Autovial 12 Syringeless Filter, glass prefilter, PVDF, 0.2 µm 50/pk

Product code AV125UAQU

Autovial 12 Syringeless Filter, glass prefilter, PVDF, 0.45 µm 50/pk

Ordering Information - Autovial Syringeless Filters

Product code AV525UAQU

Autovial 12 Syringeless Filter, glass prefilter, PVDF, 0.45 µm 1000/pk
PTFE Filtration Media

Product code AV125SORG

Autovial 12 Syringeless Filter, glass prefilter, Sterile, PTFE, 0.2 µm
40/pk

Product code AV125EORG

Autovial 12 Syringeless Filter, glass prefilter, PTFE, 0.2µm 50/pk

Product code AV125UORG

Autovial 12 Syringeless Filter, glass prefilter, PTFE, 0.45 µm 50/pk

Product code AV525UORG

Autovial 12 Syringeless Filter, glass prefilter, PTFE, 0.45 µm 1000/pk
GMF Filtration Media

Product code AV125UGMF

Autovial 12 Syringeless Filter, glass prefilter, GMF, 0.45 µm 50/pk
GF/B Filtration Media

Product code AV525BGMF

Autovial 12 Syringeless Filter, glass prefilter, GF/B, 1.0 µm 1000/pk

Autovial 12 - With Polypropylene Prefilter

PP Filtration Media

Product code AV125EPP

Autovial 12 Syringeless Filter, Polypropylene Prefilter, PP, 0.2 µm
50/pk

Product code AV125UPP

Autovial 12 Syringeless Filter, Polypropylene Prefilter, PP, 0.45 µm
50/pk

cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

Autovial and, Whatman™ are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

All other third-party trademarks are the property of their respective owners.

© 2020–2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit cytiva.com/contact

91985 AB V:4 02/2021

