

Whatman Autovial 5 Syringeless Filters

Product Information sheet

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.

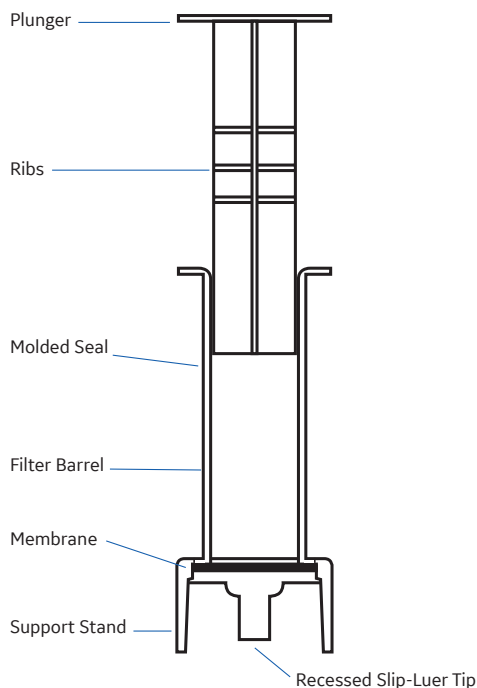
Safety

For use and handling of the products in a safe way, refer to the Safety Data Sheets.

Description

Autovial™ 5 syringeless filters are preassembled filtration devices for removing particulates from samples up to 5 mL in volume. 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.

Autovial is for laboratory use only.



Step	Action
1	Select the Autovial 5 Syringeless Filter that is compatible with your sample (See Chemical Compatibility Guide). If your solvent is not listed, please contact our technical services staff. Stand the Autovial filter-barrel on the bench top and pour in the sample. Each Autovial has 5 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 or air between the sample and the plunger. Push the plunger approximately 13 mm into the Autovial filter-barrel.
3	Place the tip of the Autovial into the opening 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 air-purge the membrane.

Membrane Guide

Type	Applications
Glass Microfiber (GMF)	High loading capacity, binder-free, glass microfiber depth filter medium. Broad chemical compatibility. Fast flow rates.
Nylon (NYL)	Membrane for aqueous and organic samples within a pH range of 3-10.
Polytetrafluoroethylene (PTFE)	PTFE membrane for samples with >50% organic solvent, and aggressive samples.
Polyvinylidene Fluoride (PVDF)	Low nonspecific protein binding membrane for samples in aqueous solutions or organic solvents.

Chemical Compatibility Guide

Solvent	GMF	NYL	PTFE	PVDF
Acetic Acid 5% +	R	R	R	R
Acetic Acid, Glacial	R	LR	R	R
Acetone	R	R	R	NR
Acetonitrile	R	R	R	R
Ammonia, 6N	LR	R	R	LR
Amyl Acetate	R	R	R	LR
Amyl Alcohol	R	R	R	R
Benzene ¹	R	LR	R	R
Benzyl Alcohol ¹	R	LR	R	R
Boric Acid	R	LR	R	R
Butyl Alcohol	R	R	R	R
Butyl Chloride ¹	R	NR	R	R
Carbon Tetrachloride ¹	R	LR	R	R
Chloroform ¹	R	NR	R	R
Cyclohexanone	R	NR	R	R
Chlorobenzene	R	*	R	R
Citric Acid	R	R	R	R
Cresol	R	NR	R	NR
Cyclohexane	R	R	R	R
Diethyl Acetamide	R	R	R	NR
Dimethyl Formamide	R	R	R	NRN
Dioxane	R	R	R	LR
DMSO	R	R	R	LR
Ethanol	R	R	R	R
Ethers	R	R	R	LR
Ethyl Acetate	R	R	R	LR
Ethylene Glycol	R	R	R	R
Formaldehyde	R	R	R	R
Freon TF	R	R	R	R
Formic Acid	R	NR	R	R
Hydrochloric Acid (Conc)	R	NR	R	R
Hydrofluoric Acid	NR	NR	R	R
Hexane	R	R	R	R
Isobutyl Alcohol	R	R	R	R
Isopropyl Acetate	R	R	R	R
Methanol	R	R	R	R
Methyl Ethyl Ketone	NR	R	R	R
Methylene Chloride ¹	NR	LR	R	LR
Nitric Acid (Conc)	R	NR	R	NR
Nitric Acid, bN	R	NR	R	LR
Nitrobenzene ¹	R	LR	R	R
Pentane	R	R	R	R
Perchloro Ethylene	R	R	R	R

Solvent	GMF	NYL	PTFE	PVDF
Pyridine	R	LR	R	R
Phenol (0.5%)	R	R	R	R
Sodium Hydroxide, 6N	NR	LR	R	NR
Sulfuric Acid (Conc)	R	NR	R	NR
Tetrahydrofuran	R	R	R	R
Toluene ¹	R	LR	R	R
Trichloroethane*	R	LR	R	R
Trichloroethylene ¹	R	NR	R	R
Water	R	R	R	R
Xylene ¹	R	LR	R	R

¹ Short term resistance of housing

Table abbreviations

R = Resistant; LR = Limited Resistance; NR = Non Resistant; + = Insufficient Data

Ordering Information

Product code	Description	Size	Qty/Box
AV115NPEORG	PTFE	0.2 µm	50
AV115NPUORG	PTFE	0.45 µm	50
AV115NPUAQU	PVDF	0.45 µm	50
AV115NPUNYL	NYL	0.45 µm	50
AV115UGMF	GMF	0.45 µm	50

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

29249519 AB V:4 01/2021

