

## User Guide

# Millex®-GS/HA/AA Filter

with MF-Millipore™ Membrane

 **SLGSR33SS, SLGSR33SB, SLHAR33SS, SLHAR33SB, SLAAR33SS, SLAAR33SB**

- 33 mm
- Sterile
- Single use only
- Non-pyrogenic
- For research use only

## Introduction

This document provides compatibility information, operating steps, and specifications for the MF-Millipore™ mixed cellulose esters family of sterile Millex® filters. The Millex® filter's bidirectional support of the filter membrane enables users to filter aqueous solutions in either direction; forward (from the syringe into the container) or backward (from the container into the syringe). The Millex® filter removes microorganisms, particles, precipitates, and undissolved powders larger than the membrane's rated pore size. These single-use filters consist of a membrane filter sealed in an acrylic housing. They are non-pyrogenic and non-toxic.

## Applications

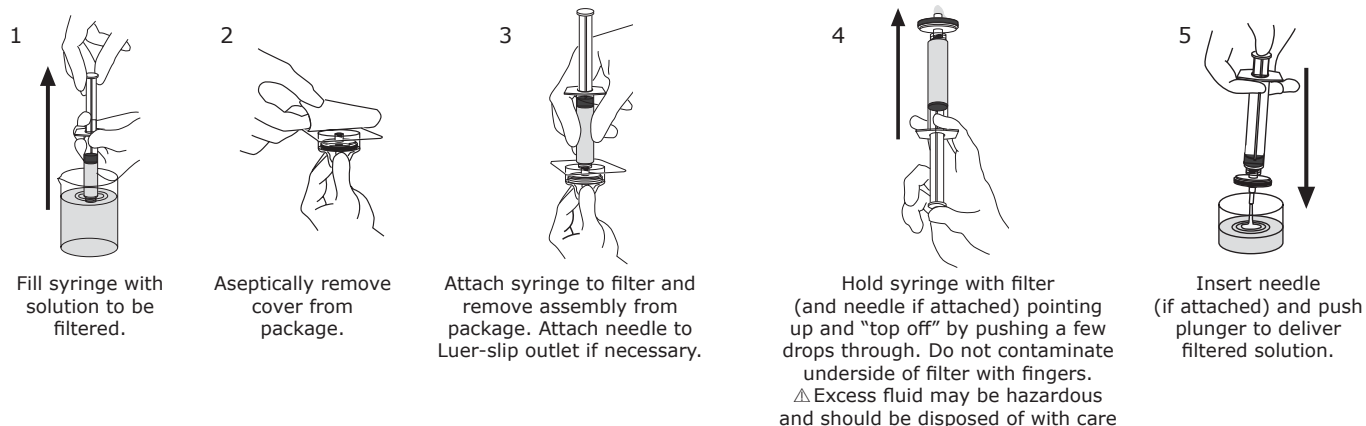
For research use only. Typical research laboratory applications include the sterile filtration (GS) and/or clarification (GS/HA/AA) of protein solutions, tissue culture media, additives, buffers, and water.

## How to Use the Millex® Sterile Filter

### WARNINGS

- To ensure sterility, do not use this product if the package is damaged.
- Do not use this product as an in-line filter; it was not designed for long-term continuous use.
- Do not use with syringes smaller than 10 mL because pressures in excess of the maximum pressure rating may be reached, potentially causing damage to the filter and/or personal injury.
- Filter solutions with a temperature limit of 45 °C (113 °F).
- Do not use the Millex® filter to filter emulsions or suspensions because it was not designed for that purpose.
- Do not use the Millex® filter to filter solutions containing 5 milligrams (mg) or less of active materials unless binding studies have been performed.
- Do not use the same Millex® filter to filter solutions in both directions.
- Single use only; do not re-use or resterilize.
- Discard appropriately after single use. See "Disposal" section.

## Instructions for Use



## Specifications

### Materials

Membrane	MF-Millipore™ mixed cellulose esters (MCE)
Pore size	Millex® -GS filter: 0.22 µm Millex® -HA filter: 0.45 µm Millex® -AA filter: 0.8 µm
Housing	Modified acrylic copolymer (MMA)

### Dimensions

Inlet to outlet	27 mm (1.06 in.)
Diameter	33 mm (1.30 in.)
Filtration area	4.52 cm <sup>2</sup> (0.70 in <sup>2</sup> )

<b>Temperature limit</b>	45 °C (113 °F) maximum
<b>Housing Pressure at 21 °C</b>	10.3 bar (150 psi) inlet maximum
<b>Filtration volume</b>	10 mL to 100 mL
<b>Hold-up volume</b>	≤0.1 mL after air purge
<b>Sterilization method</b>	Ethylene oxide gas
<b>Connections</b>	Female Luer-Lok™ inlet; male Luer-slip outlet
<b>Flow rate at 2.1 bar (30 psi), 21 °C</b>	Millex® -GS filter: ≥75 mL/min Millex® -HA filter: ≥180 mL/min Millex® -AA filter: ≥360 mL/min

## Chemical Compatibility

The Millex® filter with MF-Millipore™ membrane is compatible with most aqueous solutions. Based on information from technical publications, materials suppliers, and laboratory tests, we believe the agents listed in the following chart are safe to use with Millex® filters. However, because of the effects of variability in temperature, concentrations, duration of exposure, and other factors outside of our control, we do not provide or imply a warranty with respect to this information.

### Chemicals

Acetic acid (5%)	Guanidine hydrochloride (6 M)*	Perchloroethylene
Alconox® detergent (1%)*	Guanidine thiocyanate (5 M)*	Petroleum based oils*
Aliphatic ethers	Helium (gas)**	Petroleum ether*
Ammonium hydroxide (6 N)*	Hexane	Silicone oils*
Benzyl alcohol (1%)	Hydrochloric acid (1 N)	Sodium carbonate
Boric acid (aqueous solution)*	Hydrogen (gas)**	Sodium chloride (2 M)
Butyl alcohol*	HYPO (aqueous solution)	Sodium dodecyl sulfate*
Formic acid (50%)*	Kerosene*	Tween® 20 surfactant*
Freon® solvent (TF or PCA)*	Mineral spirits*	Urea (8 M)
Gasoline*	Nitrogen (gas)**	Water (brine)
Glycerine (glycerol)	Ozone (gas)**	Water (deionized)
(aqueous solution)	Pentane	

\* Test before using.

\*\* May leak through membrane, application dependent.

## Product Ordering

Purchase products online at [SigmaAldrich.com/Products](https://www.sigmaaldrich.com/Products).

Description	Pore Size, µm	Diameter	Membrane	50/pk	250/pk
Millex®-GS	0.22	33 mm	MF-Millipore™ MCE	SLGSR33SS	SLGSR33SB
Millex®-HA	0.45	33 mm	MF-Millipore™ MCE	SLHAR33SS	SLHAR33SB
Millex®-AA	0.8	33 mm	MF-Millipore™ MCE	SLAAR33SS	SLAAR33SB

---

## Disposal

Follow precautions for disposal of items contaminated with potentially infectious or hazardous material according to all applicable international, federal, state, and local regulations.

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

## Technical Assistance

Visit the tech service page on our web site at [SigmaAldrich.com/TechService](https://SigmaAldrich.com/TechService).

## Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/Terms](https://SigmaAldrich.com/Terms).

## Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/Offices](https://SigmaAldrich.com/Offices).

Merck, Millipore, MF-Millipore, Millex and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.  
© 2019-2025 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

The life science business of Merck operates  
as MilliporeSigma in the U.S. and Canada.

