

User Guide

Millex®-HA Non-Sterile 25 mm Syringe Filter

with MF-Millipore™ Membrane

 **SLHA025NB, SLHA025NK**

- For research use only
- Single use only

Introduction

The Millex®-HA Non-sterile 25 millimeter (mm) syringe filter with male Luer-slip outlet is recommended for filtering 10–100 milliliter (mL) volumes to remove particles prior to instrumentation analysis. The single-use, disposable filter removes particles larger than the membrane's rated pore size.

The Millex®-HA syringe filter consists of a 0.45 µm mixed cellulose esters (MCE) membrane sealed in a polyvinylidene chloride (PVC) housing and is intended for filtration of aqueous solutions.

How to Use Millex® 25 mm Syringe Filter

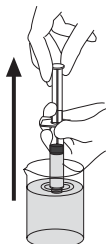
WARNINGS

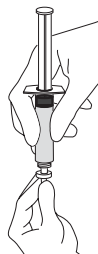
- Do not use the Millex® 25 mm syringe filter for direct patient care applications; it is designed for research use only.
- Do not use with syringes smaller than 10 mL because pressures in excess of the maximum pressure rating may be reached, potentially causing damage to the syringe filter and/or personal injury.
- Sudden loss of pressure could indicate failure of filter.
- Make sure to wet the filter membrane thoroughly before injecting the solution; improperly wetted filters can become airlocked.
- Single use only; do not re-use.
- Do not use this filter as an in-line filter.
- Discard appropriately after single use. See "Disposal" section.

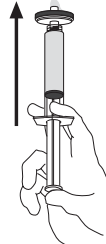
CAUTIONS

- Do not use the 25 mm syringe filter at temperatures above 45 °C (113 °F).
- Do not use the same Millex® syringe filter to filter solutions in both directions.
- Do not use the syringe filter to filter emulsions or suspensions.
- Perform a binding study before use if there is a concern about loss of analyte (proteins, nucleic acids, active pharmaceuticals) due to binding.

Instructions for Use

- 

1 Fill syringe with solution to be filtered.
- 

2 Attach syringe to the Millex® filter.
- 

3 Hold syringe with filter pointing up and "top off" by pushing a few drops through. ⚠ Excess fluid may be hazardous and should be disposed of with care.
- 

4 Push the syringe plunger to deliver filtered solution.

Optional: To purge the syringe filter and maximize sample throughput, remove the Millex® filter from the syringe, draw air into the syringe, reattach the Millex® filter, and push the plunger to force some of the air through the filter.

Specifications

Materials

Membrane	Hydrophilic MF-Millipore™ mixed cellulose esters (MCE)
Pore size	0.45 µm
Housing	Polyvinyl chloride (PVC)

Dimensions

Inlet to outlet	25 mm (0.98 in.)
Diameter	29 mm (1.14 in.)
Filtration area	4 cm ² (0.62 in ²)
Temperature limit	45 °C (113 °F)
Pressure Limit at 21 °C	75 psi (5.2 bar) inlet and differential
Filtration volume	10-100 mL
Hold-up volume	≤ 0.1 mL after air purge
Connections	Female Luer-Lok™ inlet; male Luer-slip outlet
Water flow rate at 30 psi (2.1 bar), 21 °C	≥ 180 mL/min

Chemical Compatibility

Millex-HA® syringe filters are only compatible with aqueous solutions and gases. Compatibility information was developed from technical publications, materials suppliers, and laboratory tests, and is believed to be accurate and reliable. However, because of variability in temperature, concentrations, exposure time, and other factors beyond our control that may affect the use of the filter, no warranty is provided or implied with respect to such information.

Product Ordering

Purchase products online at SigmaAldrich.com/Products.

Description	Pore Size	Membrane	250/pk	1000/pk
Millex®-HA	0.45 µm	MF-Millipore™ MCE	SLHA025NB	SLHA025NK

Disposal

Follow precautions for disposal of items contaminated with 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.

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at SigmaAldrich.com/Terms.

Contact Information

For the location of the office nearest you, go to 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.