

## User Guide

# Swinny Stainless Steel 13 mm Filter Holder

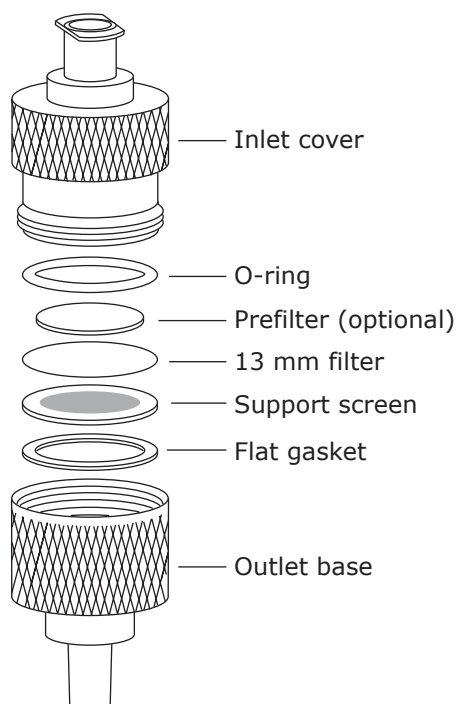
**XX3001200**

## Introduction

The Swinny Stainless Steel 13 mm Filter Holder is a syringe-operated filter holder designed to sterilize or clarify small volumes of liquids by pressure filtration. It can be used to sterilize tissue culture media or clarify HPLC samples, solvents, etc. The filter holder can be autoclaved with the filter in place for clarification/sterilization applications that do not require removing the filter intact for further use.

## Filter Loading

1. Unscrew the inlet cover from the outlet base.
2. Remove the O-ring, support screen, and flat gasket.
3. Inspect the gasket, O-ring, and sealing surfaces for damage; imperfections can cause leaks during filtration. Replace damaged parts.
4. Place the flat gasket in the base and place the support screen on top of it.
5. Using smooth-tip forceps, place a 13 mm diameter filter on the support screen, then place the O-ring on top of the filter. If a depth-type prefilter (10 mm diameter) is required to extend the life of the final filter, center the prefilter on top of the final filter before adding the O-ring.
6. Screw the cover down into the base and finger tighten firmly so that the filter is securely sealed.



## Operation

⚠ **Caution:** Do not exceed the maximum inlet and differential pressure of 6.9 bar (100 psi).

When using a syringe smaller than 10 mL, the syringe should have a Luer-Lok™ connection (not Luer-slip) to avoid leakage at the higher pressures generated by small syringes.

1. Draw the unfiltered sample into the syringe.
2. Attach the Luer-Lok™ inlet of the filter holder to the syringe. If required, attach a hypodermic needle to the filter holder outlet for dispensing the filtered liquid.
3. Press the syringe plunger to filter the sample.

## Autoclaving (Optional)

1. Make sure that all inner surfaces of the filter holder are completely dry, then assemble the holder with a filter in place. Tighten partially.
2. Wrap the filter holder with lint-free autoclave paper, and autoclave for 15 minutes at 121 °C and 1 bar (15 psi).
3. After the unit has cooled, fully tighten it, being careful not to touch the outlet tube.
4. Use aseptic technique when making outlet connections.

## Maintenance

1. Disassemble the holder and inspect all components for signs of damage. Grit embedded in PTFE surfaces can sometimes be dislodged without damaging the surface, but it is preferable to replace the O-ring or gasket.
2. Wash all components thoroughly with a sponge, hot water, and a nonabrasive cleanser. Scrub threaded parts, recesses, and orifices with a stiff bristle brush and use a pipe cleaner to remove contaminants from inlet and outlet orifices.

⚠ **Caution:** To avoid component damage, never use steel wool or abrasive materials on any part of the filter holder.

3. Rinse all parts thoroughly with hot running water, then rinse several times with cold, laboratory-grade water.
4. Allow the components to air dry while disassembled.

⚠ **Caution:** Do not wipe dry with paper or cloth, as this leaves fibers and lint on the surfaces and also generates electrostatic forces that can attract more dirt. Filtered compressed air can be used to facilitate drying.

## Specifications

<b>Maximum pressure</b>	6.9 bar (100 psi) inlet and differential
<b>Filter diameter</b>	13 mm diameter
<b>Prefilter diameter</b>	10 mm diameter (thick depth prefilter)
<b>Effective filter area</b>	Approximately 0.8 cm <sup>2</sup> (0.13 in <sup>2</sup> )
<b>Weight</b>	23 g (0.8 oz)
<b>Connections</b>	Female Luer-Lok™ inlet, male Luer-slip outlet
<b>Dimensions</b>	
Diameter	16.0 mm (0.63 in.)
Length	36.6 mm (1.44 in.)
<b>Materials of construction</b>	
Inlet cover	316 stainless steel
Outlet base	316 stainless steel
Support screen	316 stainless steel
Flat gasket	Polytetrafluoroethylene (PTFE)
O-ring	PTFE

## Statement Regarding Compliance with the Pressure Equipment Directive, 2014/68/EU

We certify that this product complies with the European Pressure Equipment Directive, 2014/68/EU of 15 May 2014. This product is classified under Article 4 § 3 of the Pressure Equipment Directive. It has been designed and manufactured in accordance with sound engineering practices to ensure safe use. The product is accompanied by user instructions and bears markings to permit identification us as the manufacturer or authorized representative of this product within the European Community. In compliance with Article 4 § 3 of the Pressure Equipment Directive, this product does not bear the CE mark.

## Product Information

Purchase products on-line at [SigmaAldrich.com](https://SigmaAldrich.com).

Description	Catalogue Number
Swinny Stainless Steel 13 mm Filter Holder  Includes inlet cover (1), outlet base (1), O-ring (1), support screen (1), flat gasket (1)	XX3001200
Swinny 13 mm Filter Holder Replacement Parts Kit  Includes O-ring (4), flat gasket (4), filter holder support screen (4)	XX30012RK
Support screen (1/pk)	XX3001210
Flat gasket, PTFE (10/pk)	XX3001202
Filter forceps, stainless, blunt end (3/pk)	XX6200006P
Plastic syringe, 20 mL capacity (12/pk)	XX1102012
Plastic syringe, 50 mL capacity (5/pk)	XX1105005

## 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 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.  
© 2015-2024 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.

