

Whatman™ VACU-GUARD Disposable Filter

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.

Background

Description

Whatman™ VACU-GUARD is a disposable filter designed to protect vacuum systems. This filter contains a hydrophobic polytetrafluoroethylene (PTFE) membrane, which provides retention for air particulates and aerosols.

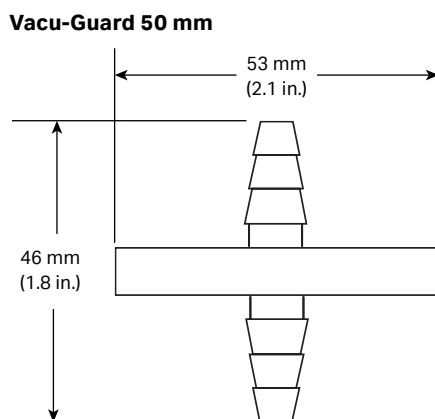
Multiple reuse is the responsibility of the operator who should protect the device from cross contamination and detect loss of integrity by appropriate testing.

Typical applications

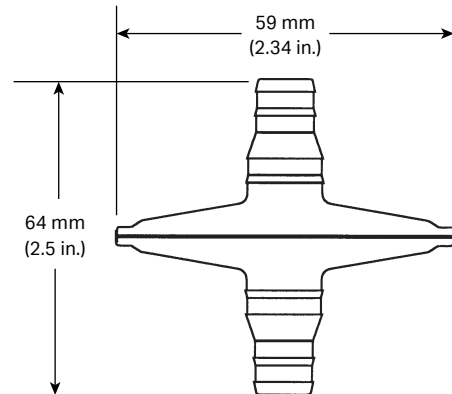
VACU-GUARD is intended to be used to prevent aerosol and fluid contamination in vacuum pumps or aspiration suction systems.

Technical information

Illustration of VACU-GUARD



Vacu-Guard 60 mm

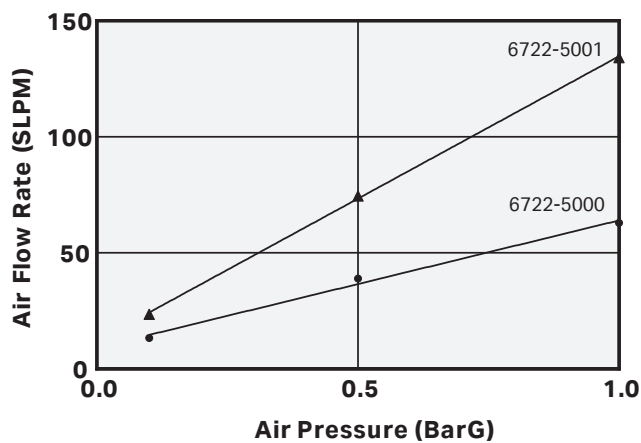


Technical data

Housing and support:	Polypropylene	
Filter media:	PTFE membrane	
Pore size:	0.45 µm	
Air particle retention:	99.9% retention of all particles ≥ 0.1 µm	
Effective filtration area:	Specification	Product code
	16 cm ²	6722-5000
	25 cm ²	6722-5001
Inlet/outlet connections:	Specification	Product code
	SB	6722-5000
	½ SB	6722-5001
	SB: 6 to 10 mm (¼ to ⅜ in.) stepped barb	
	½ SB: 10 to 12 mm (⅜ to ½ in.) stepped barb	
Dimensions (W×L):	Specification	Product code
	53 × 46 mm (2.1 × 1.8 in.)	6722-5000
	59 × 64 mm (2.34 × 2.5 in.)	6702-5001
Sealing method:	Heat-fused	
Autoclavable:	121°C (250°F) for 20 minutes at 0.1 MPa (1.0 bar, 15 psi)	
Maximum operating pressure:	0.1 MPa (1.0 bar, 15 psi)	

Operating temperature:	Ambient
Flow direction:	Inlet to outlet
Biosafety:	Materials pass USP Class VI

Typical air flow rate



Operating Instructions

Safety

When considering the specific factors of your application, refer to Technical data for correct use. Make sure not to exceed the Maximum operating pressure and follow temperature or chemical compatibility recommendations.



CAUTION

If the Maximum operating pressure is exceeded, bursting of the device can occur resulting in loss of sample or personal injury.



Give feedback on this document

Visit cytiva.com/techdocfeedback or scan the QR code.



cytiva.com

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corp. or an affiliate doing business as Cytiva.

Whatman is a trademark of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

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

© 2020–2022 Cytiva

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

90400 AC V:8 07/2022

Venting

For venting applications, connect the inlet port of the VACU-GUARD to the vessel, leaving the outlet open to the atmosphere. The connection is made by securing the tubing to the filter ports using band clamps.

Note: Change filter if there is condensation or contact with fluid preventing sufficient air flow.

In-line

To use the VACU-GUARD for in-line application, securely connect both ports of the filter into the flow stream such that the orientation flows from inlet to outlet. The connections are made by securing the tubing to the filter ports using band clamps.

Note: Change filter if there is condensation or contact with fluid preventing sufficient air flow.

Ordering information

Product Code	Product Name	Qty./Pk.
6722-5000	VACU-GUARD 50 mm	10
6722-5001	VACU-GUARD 60 mm	10

