

# Whatman™ HEPA-VENT 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™ HEPA-VENT is a disposable, in-line filter that will retain 99.97% of all particles  $\geq 0.3 \mu\text{m}$ . The filter media used in HEPA-VENT is used throughout scientific, research, and industrial environments to provide clean, particulate-free air and gasses.

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

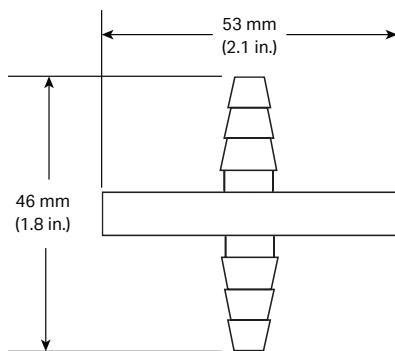
#### Typical applications

HEPA-VENT with high flow rate glass microfiber (GMF) is treated to provide mild hydrophobicity and is suitable for:

- Venting vessels used for mixing, filling, fermenting, storage, or transport by allowing particulate free air to move freely in a bidirectional manner.
- In-line filtration of a gas or air stream such as found on instruments, incubators, and culture vessels.

### Technical information

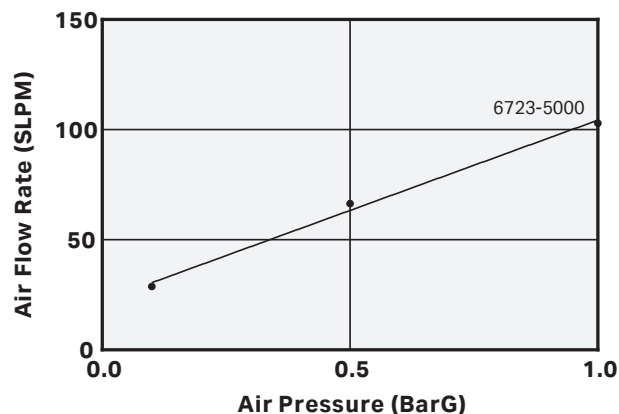
#### Illustration of HEPA-VENT



### Technical data

<b>Housing and support:</b>	Polypropylene
<b>Filter media:</b>	GMF
<b>Particle retention:</b>	99.97% retention of all particles $\geq 0.3 \mu\text{m}$
<b>Effective filtration area:</b>	16 cm <sup>2</sup>
<b>Inlet/outlet connections:</b>	6 to 10 mm (1/4 to 3/8 in.) stepped barb
<b>Dimensions (W×L):</b>	53 × 46 mm (2.1 × 1.8 in.)
<b>Sealing method:</b>	Heat-fused
<b>Autoclavable:</b>	121°C (250°F) for 20 minutes at 0.1 MPa (1.0 bar, 15 psi)
<b>Maximum operating pressure:</b>	0.41 MPa (4.1 bar, 60 psi)
<b>Operating temperature:</b>	Ambient
<b>Flow direction:</b>	Bidirectional
<b>Biosafety:</b>	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.

### Venting

For venting applications, connect the inlet port of the HEPA-VENT 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 HEPA-VENT 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.
6723-5000	HEPA-VENT	10
2216	HEPA-VENT	50



#### Give feedback on this document

Visit [cytiva.com/techdocfeedback](https://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](https://cytiva.com/contact)

90315 AC V:8 07/2022

