

Whatman™ HEPA-CAP 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-CAP 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-CAP 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.

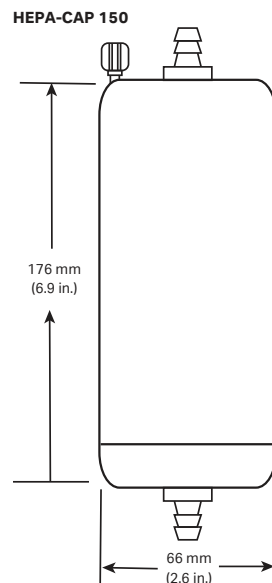
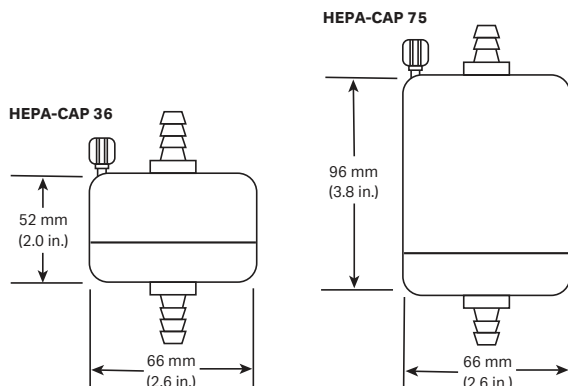
Typical applications

HEPA-CAP 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-CAP

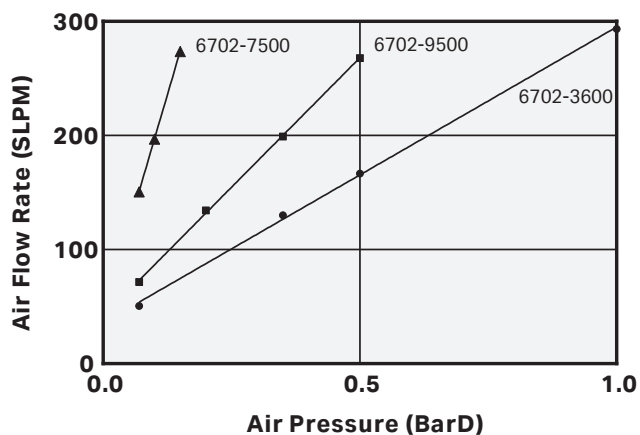


Technical data

Housing and support:	Polypropylene	
Filter media:	GMF	
Particle retention:	99.97% retention of all particles $\geq 0.3 \mu\text{m}$	
Effective filtration area:	550 cm ² – HEPA-CAP 36 1240 cm ² – HEPA-CAP 75 2440 cm ² – HEPA-CAP 150	
Inlet/outlet connections:	Specification	Product code
	SB	6702-3600
	FNTP	2609T, 2709T, 2809T
	½ SB	6702-7500, 6702-9500
	SB: 6 to 10 mm (¼ to ⅜ in.) stepped barb	
	FNTP: ⅜ in. female NPT threaded connection	
	½ SB: 10 to 12 mm (⅜ to ½ in.) stepped barb	

Dimensions with inlet/outlet connections (W×L):	Specification	Product code
	66 × 92 mm (2.6 × 3.6 in.)	6702-3600
	66 × 56 mm (2.6 × 2.2 in.)	2609T
	66 × 162 mm (2.6 × 6.4 in.)	6702-7500
	66 × 101 mm (2.6 × 4.0 in.)	2709T
	66 × 223 mm (2.6 × 8.8 in.)	6702-9500
	66 × 180 mm (2.6 × 7.1 in.)	2809T
Sealing method:	Heat-fused	
Autoclavable:	121°C (250°F) for 20 minutes at 0.1 MPa (1.0 bar, 15 psi)	
Maximum operating pressure:	0.41 MPa (4.1 bar, 60 psi)	
Operating temperature:	Ambient	
Flow direction:	Bidirectional	
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.

Venting

Step	Action
1	Securely connect the inlet port to the vessel, leaving the outlet open to the atmosphere. <ul style="list-style-type: none"> a. For filters with stepped hose barbs, secure tubing to the capsule using band clamps. b. For filters with threaded connections, hand-tighten the connection to the vessel.
2	Change filter if there is condensation or contact with fluid preventing sufficient air flow.

In-line use

Step	Action
1	Securely connect both ports of the filter into the flow stream using flow arrows to guide orientation of the filter system. <ul style="list-style-type: none"> a. For filters with stepped hose barbs, secure tubing to the capsule ports using band clamps. b. For filters with threaded connections, hand-tighten the connections of both ports.
2	Change filter if there is condensation or contact with fluid preventing sufficient air flow.

Ordering information

Product Code	Product Name	Qty./Pk.
6702-3600	HEPA-CAP 36	1
2609T	HEPA-CAP 36	5
6702-7500	HEPA-CAP 75	1
2709T	HEPA-CAP 75	5
6702-9500	HEPA-CAP 150	1
2809T	HEPA-CAP 150	5



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

90318 AC V:8 08/2022

