

Whatman Polydisc HD Disposable Filter Device

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.

Description

The Polydisc HD disposable filter devices are designed to provide disposable filters with high flow rates and long life for fine filtration. These filters are ideal for removing particulates in the range of 5.0 µm to 10.0 µm. They provide the capability to filter in the range between gross filter papers and microporous membranes used for absolute filtration.

Polydisc HD is an all polypropylene (PP) filter device. The filter media is MAPP (Monofilament Anisotropic Polypropylene) a microfine continuous PP material formed into a variety of media by varying the filament's diameter and the layer density.

Disposable filtration devices provide great labor efficiency while insuring superior filtration when compared to hand assembled reusable filter housings.

This bulletin provides general information on the products listed.

- Total Polypropylene (PP) Construction. Housing, Filter Media and Support System
- Bonded & Sealed with Advanced Fusion Technology. No mold releases, glues, powdered metals, adhesives, epoxies or other extraneous materials are used.
- High Flow
- Long Life
- Autoclavable
- Lightweight – won't cause collapsed tubing
- Integral & disposable: reduces equipment costs, eliminates clean up time, insures installation integrity and purity
- Biosafe materials USP Class VI

Polydisc HD – 50mm PP Filters

Product code	Product	Pore Size (µm)	Qty./Box
6728-5050	Polydisc HD	5.0	10
6728-5100	Polydisc HD	10.0	10

Typical Applications for Polydisc HD Disposable Filter Devices

General Fine Filtration: Most liquids/gases requiring 10.0 µm down to 5.0 µm filtration retention

Prefiltration: extends life of RO/UF/MF media

Sample Prep: Ground Water • Biological Solutions • Effluents

Equipment: Clean air/gas

Food & Beverage: Oils & Syrup • Spirits • Beverages • Wine & Juices • Bottled water

Miscellaneous: Water to wash/rinse • Vials, Glassware & Tanks • Make-up water • General Prefiltration

OPERATING INSTRUCTIONS

Safety: Considering the special factors of your application consult the table of Technical Data to determine the correctness of use. Do not exceed the pressure, temperature or chemical compatibility recommendations.

Filter Selection: Select filter based on filter efficiency and other information presented. This information is published as a general guide. The user must evaluate a specific product to determine the appropriateness of use.

Filtration Installations and Use: Small volumes may be filtered using a syringe; larger volumes may be filtered by connecting a hose to the outlet of a pressure vessel, or "in line" using gravity vacuum or peristaltic pump for the pressure required to filter. Use hose clamps for higher pressures. To maximize recovery of filtered solution purge the housing with air at the end of filtration.

TECHNICAL DATA: Polydisc HD Disposable Filter Device

Product code	Product Name	Rated ¹ Microns	Filter Media	Connections	Flow Rates ² at 1.0 bar (14.5 psi)		Qty./ Pkg.
					AIR SLFM	WATER Liters/min.	
6728-5050	Polydisc HD	5.0	Polypropylene	6-10mm (1/4-3/8") SB	110	1.5	10
6728-5100	Polydisc HD	10.0	Polypropylene	6-10mm (1/4-3/8") SB	140	2.2	10

¹ Liquid Rating. (Efficiency in gas streams is significantly higher.)

² Typical Values

Dimensions:	53 mm (2.1 in.) x 46 mm (1.8 in.)
Weight:	11.5 grams
Filtration Area:	16 cm ² / 2.48 in. ² / 0.02 ft ²
Maximum Pressure:	4.1 bar (60 psi)
Retention:	>99% of particles > the µm rating
Flow Direction:	Flow should enter from the inlet
Housing:	Polypropylene
Connectors:	SB Stepped Hose Barbs for 6 to 10 mm (1/4" to 3/8") ID tubing (Accepts male luer)
Volume "Hold Up":	Full Housing, 1.0 ml With Air Purge, <0.1 ml
Biosafe:	All Materials Pass USP Class VI
Sterilization:	Autoclave at 121°C (132°C Max) for 20 minutes

CHEMICAL RESISTANCE SUMMARY

Classes of Substances	Guide for use ¹
	Polypropylene Stability at 20°C (68°F)
Acids, dilute	Long Term
Acids, concentrated	Not Usable
Alcohols	Long Term
Aldehydes	Long Term
Bases	Long Term
Esters	Not Usable
Hydrocarbons, aliphatic	Varies
Hydrocarbons, aromatic	Not Usable
Hydrocarbons, halogenated	Not Usable
Ketones	Not Usable
Water	Long Term

¹ Published as a general guide only. Due to time, temperature, and stress variations the user must evaluate the specific product and application to determine the appropriateness of use.

cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

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

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

© 2020–2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

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

90808 AC V:5 08/2021

