

Whatman Puradisc 25 GD 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

Puradisc 25 GD disposable filter devices were designed to provide faster and more efficient prefiltration of aqueous or solvent solutions. They are made of multilayered, binder-free glass microfibre filters with a polypropylene housing using the most advanced methods and design features available today.

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

This bulletin provides general information on the products listed below. The specifications in the Technical Data section are intended to provide the basis for establishing functional use, as well as setting quality assurance test performance levels.

- Multilayered — Multiple Porosity
- High Flow Binder Free Glass Microfibre Filter Media.
- Polypropylene Housing
- Connectors

Inlet: Female Luer Lock (FLL)
Outlet: Male Luer (ML)

- Ultra-Clean with No Mold Releasing Agents or Glues Used
- Available in 2 Particle Retention Ratings, Labeled for Easy Identification
- Autoclavable

Puradisc 25 GD — 25mm Filters

Product Code	Product Name	Pore Size (µm)	Media	Qty./Pkg.
67832510	Puradisc 25 GD	1	GMF-150	100
67832520	Puradisc 25 GD	2	GMF-150	100
67922510	Puradisc 25 GD	1	GMF-150	1000

Typical Applications for Puradisc 25 GD Filter Devices

Serum Samples

Tissue Culture Media

Immunologicals
Buffer Solutions
Particle Counting Solutions
Salt Solutions
Aqueous Solutions

Biological Samples
Virus Suspensions
Air/Gas Filtration
HPLC, TLC and GC
Sample Clarification

General Fine Filtration:

Most liquids/gases requiring filtration to 1.0µm retention.

Prefiltration:

extends life of RO/UF/MF membranes.

Semiconductor/Magnetic Media:

Solvents • Acids • Bases • Etchants • Photoresists • Strippers • Air Guns/Cleaning Wands • Coatings • Oxide Dispersions • Media Chemicals • Gases

Laboratories:

Clean air/gas for instruments/environments • Analytical Reagents • Solvents • Buffers Diluents • Serum, Ascites Fluid

Sample prep:

for Ground Water • Process Solutions Effluents

Photographic:

Emulsions, Make-up, or Rinse water

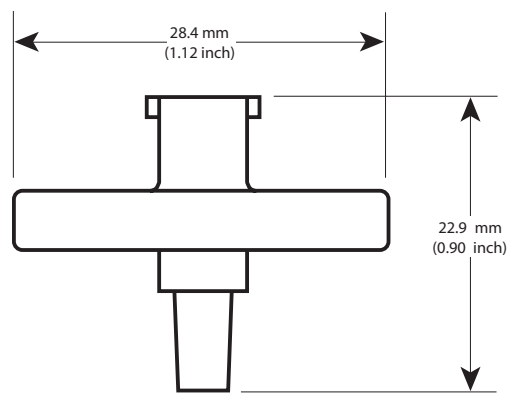
Food & Beverage Analysis:

Syrup • Oils • Spirits • Water • Beverages • Wine • Juices • Bottled Water

Pharmaceutical • Cosmetic • Personal Care:

Water to wash/rinse • Vials, glassware • General prefiltration • Filtration with requirements to 1µm nominal for clarity

TECHNICAL DATA: Puradisc 25 GD Disposable Filter Devices



Connections • Inlet: Female Luer Lock (FLL)
 Outlet: Male Luer (ML)

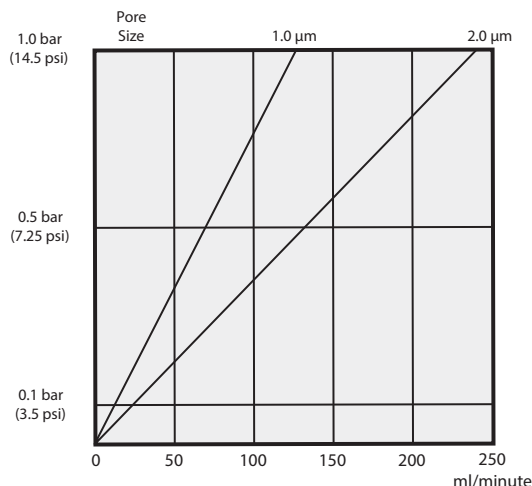
Product Code	Pore Size (µm)	Media
6783-2510	1.0	GMF 150
6783-2520	2.0	GMF 150
6792-2510	1.0	GMF 150

CHEMICAL RESISTANCE SUMMARY

Classes of Substances 20°C (68°F)	Polypropylene Guide for Use ¹
Hydrocarbons, aliphatic	One Week
Hydrocarbons, aromatic	One Week
Hydrocarbons, halogenated	One Week
Ketones	One Week
Oxidising agents, strong	One Week
Acids, dilute	Long Term
Acids concentrated	Long Term
Alcohols	Long Term
Aldehydes	One Month
Bases	Long Term
Esters	One Week

¹ For guidance only. User should determine suitability of syringe filter for particular application.

TYPICAL WATER FLOW RATES



OPERATING INSTRUCTIONS

Safety: When considering the special factors of your application, consult the Technical Data to determine correctness of use. Do not exceed the pressure, temperature or chemical compatibility recommendations. High pressures can be obtained when using syringes. The smaller the syringe the higher the pressure that can be generated. As a general guide, the following pressures can be obtained by hand with the syringes indicated: 20mL, 80 psi; 10mL, 140 psi; 5mL, 180 psi; 3mL, 200 psi; 1mL, 250 psi. Each user should determine the pressure they can generate by hand with a specific size syringe and take appropriate safety precautions not to exceed the recommended rating for the device used. If these limitations are exceeded, bursting of the device may occur resulting in loss of sample or personal injury.

Filter Media: Multigrade GMF-150 is a combination of two microfiber filters in one. Manufactured from 100% borosilicate glass microfibers, the filter is unique in its construction: A coarse layer on top, meshed with a fine layer below. This filtration technique allows for higher particulate loading capacity, faster flow rate and extended life of filter. Multigrade GMF-150, used as a prefilter, nearly doubles the volume of sample filtered compared to a single density prefilter. Compared to an unprotected membrane, the volume of sample filtered is three to seven times greater. Conventional prefilters cannot perform up to the same standard as the Multigrade GMF-150 simply because prefilters of a uniform density do not have the loading capacity afforded by Whatman's multilayered filter technology.

Efficiency: To maximize filtration throughput, use the largest pore size filter that will provide the required cleanliness. To extend filter life use low flow or pressure and prefilters.

To Use With A Syringe:

Step	Action
1	Fill the syringe with the solution to be filtered.
2	Secure the filled syringe to the FLL on the inlet with a twisting motion.
3	With outlet pointed upward, gradually apply thumb pressure to the syringe plunger to initiate flow.

Step	Action
-------------	---------------

4	Change filters when flow becomes too slow or resistance becomes excessive.
---	----------------------------------------------------------------------------

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

90958 AB V:4 02/2021

