

Whatman Puradisc 25 GF/F

Disposable Filter Device

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.

Description

The Puradisc 25 GF/F disposable filtration device was designed to provide a disposable filter with a high flow rate and long life, for fine filtration. This filter is ideal for removing fine particulates down to 0.7 μm . Unlike membrane filters with a comparable retention value, this product has a rapid flow rate and a high loading capacity.

Puradisc 25 GF/F has a polypropylene (PP) housing with Glass Microfiber filter medium. This filter medium is very effective with fine particulates such as precipitated proteins. It can also be used as a prefilter in conjunction with smaller pore size filtration devices.

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

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

- Whatman™ Brand GF/F Filter Media
- Extremely High Flow Rate
- Long Life - High Loading Capacity
- Chemical & Physical Stability
- Binder Free Glass Microfiber
- Particulate Retention Down to 0.7 μm
- Autoclavable
- Inlet: Female Luer Lock (FLL)
- Outlet: Male Slip Luer (ML)
- No adhesives used in construction

Puradisc 25 GF/F - 25 mm Filters

Catalog Number	Product	Pore Size	Qty./Pkg.
6825-2517	Puradisc 25 GF/F	0.7	50
6753-2510	Puradisc 25 GF/F	0.7	200

Typical Applications for Puradisc 25 GF/F Filter Devices

- TCLP Samples
- Prefiltration: Extends Life of Other Filter Media or Membranes
- Clarification of 'Difficult' Biochemical Solutions and Fluids
- Fine Filtration: Liquids requiring particulate Removal Down to 0.7 μm For Gases Retention Efficiency is Significantly Greater.
- Filtration of Precipitated Proteins
- Liquid Scintillation Counting
- Radioimmunoassay
- Water and Wastewater Analysis
- Sample Preparation

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: 20 mL, 80 psi; 10 mL, 140 psi; 5 mL, 180 psi; 3 mL, 200 psi; 1 mL, 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.

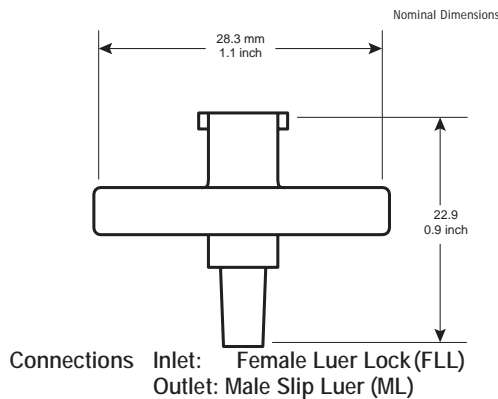
GF/F Media Considerations: The glass microfiber filter medium is produced entirely from the finest grades of borosilicate glass microfiber and contains no binders either as manufacturing aids or as wet strengthening additives. Glass microfiber filters are generally resistant to weakening or disruption of the fibrous matrix by inorganic or organic liquids.

Autoclaving: Autoclave at 121°C (131°C max.) for 20 minutes.

To use with a syringe: (3 mL, 5 mL, 10 mL)

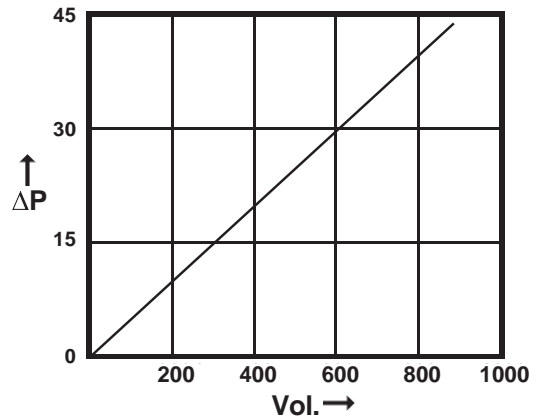
Step	Action
1	Fill the syringe with the solution to be filtered.
2	Secure the filled syringe to the FLL on the inlet with a twisting motion.
3	With outlet pointed upward, gradually apply thumb pressure to the syringe plunger to initiate flow.
4	Change filters when flow becomes too slow or resistance becomes excessive.

Technical Data: Puradisc 25 GF/F Disposable Filter Devices



Dimensions:	28.3 mm (1.1 in) x 22.9 mm (0.9 in)
Weight:	2.7 grams
Filtration Area:	4.2 cm ²
Maximum Pressure:	5.2 bar (75 psi)
Housing:	Polypropylene
Hold Up Volume:	Full Housing 0.16 mL With air purge <0.1 mL
Filter Media:	Whatman GF/F Glass Microfiber
Flow Direction:	Flow should enter from the inlet
Connectors:	Inlet – Female Luer Lock (FLL) Outlet – Male Slip Luer (ML)
Autoclaving:	Autoclave at 121°C (131°C max) at 15 psi for 20 minutes.

Water Flow Rates



Liquid Flow Rate (mL/minute)

Catalog No. Qty./Pk	Product	Media	Nominal Pore Size Microns	Water Flow Rate ¹ at 1.0 bar (14.5 psi)	
6825-2517	Puradisc 25 GF/F	Glass Microfiber	0.7	270 mL/min	50
6825-2527	Puradisc 25 GF/F	Glass Microfiber	0.7	270 mL/min	200

¹ Typical values

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

90920 AC V:5 03/2021

