

Whatman GD/X Syringe Filters

Introduction

Product codes

6869-2502

6869-2504

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 Whatman GD/X™ Syringe Filters are designed to enable easy filtration of hard-to-filter samples up to 10mL. They are available in a wide variety of filter choices with a polypropylene housing and utilize the most advanced construction methods and design features available today. This level of engineering provides for the finest disposable filtration devices possible.

Each unit contains a unique prefiltration stack of Whatman Multigrade GMF 150 (10:1 µm) and GF/F. This advanced Prefiltration Technology combined with the favorable apparent pressure generated within the device allows you to filter even the most difficult samples with less hand force than ever before possible.

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

Operating Instructions

Safety:

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; 10mL, 140 psi; 5mL, 180 psi; 3mL, 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.

Pre-Filter Media:

The Whatman GD/X Syringe Filters contain a unique prefiltration stack of Whatman Multigrade GMF 150 (10:1 µm) and Grade GF/F prefilters. Multigrade GMF 150 is a combination of two glass microfiber filters in one. Manufactured from 100% Borosilicate glass, its construction consists of a coarse layer on top, meshed with a fine layer below. The Grade GF/F Filter will retain fine particles down to 0.7µm. Whatman Prefiltration Technology allows you to filter even the most difficult samples with less hand force than ever before possible. Compared to an unprotected membrane, the volume of sample filtered can be three to seven times greater.

Efficiency:

To maximize filtration throughput, use the largest pore size filter that will provide the required cleanliness. To extend filter life use low flow rates or pressures.

To Use With a Syringe:

Step	Action
1	Fill the syringe with the solution to be filtered.
2	Secure the filled syringe to the FII inlet of the syringe filter with a twisting motion.
3	With the outlet pointed upward, gradually apply thumb pressure to the syringe plunger to initiate flow.
4	Continue thumb pressure until all the air in the device is displaced with liquid.
5	Once liquid starts to exit syringe filter outlet, stop applying pressure, point the device downward and away from user.
6	Position syringe filter over suitable collection container or other apparatus and apply pressure again to filter sample. Change filters when flow becomes too slow or resistance becomes excessive.

Air Locks:

Seriously hamper flow rates. To eliminate, point the outlet of the filter device upward during the initiation of liquid flow.

Whatman GD/X Syringe Filters

Product code	Product Description	Qty/Box
6869-2502	25 mm GD/X 0.2 m NYL	150
6869-2504	25 mm GD/X 0.45 m NYL	150

For Laboratory Use Only.

cytiva.com

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

Whatman GD/X 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

91811 AB V:5 02/2021

