

Whatman Mini-UniPrep G2 Syringeless Filters

Multi Compressor Product Information sheet

Introduction

Important

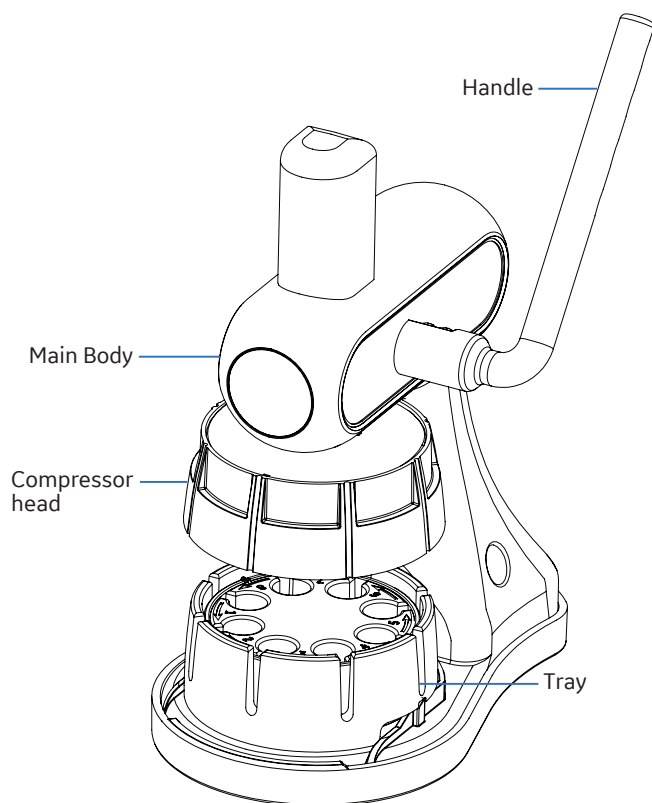
Read these instructions carefully before using the products.

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.

Safety

For use and handling of the products in a safe way, refer to the Safety Precaution section or Safety Data Sheets if applicable.



Safety Precaution

ALWAYS USE A MINI-UNIPREP G2 COMPRESSOR. Do not attempt to compress the device by fingers; there is a risk of breakage of the glass chamber which could result in personal injury if a compressor is not used. If a Multi Compressor or Hand Compressor is not available, do not use the Mini-UniPrep™ G2 Syringeless Filters until one has been obtained¹. Do not use any other device to compress the Mini-UniPrep G2 Syringeless Filters including the six-position compressor (catalogue no. CR0000006) designed to be used with the original Mini-UniPrep filters.

Do not overfill the Mini-UniPrep G2 Syringeless Filters or compress the device at an angle before the airtight seal is made. Liquid could escape from the air vent holes, resulting in the loss of sample or personal injury.

The Multi Compressor allows for the compression of up to 8 Mini-UniPrep G2 Syringeless Filters filter devices at one time.

See cytiva.com for further details.

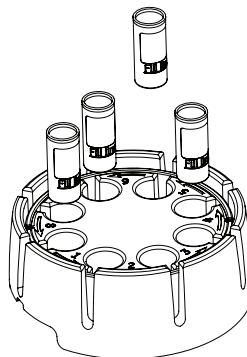
In case of sample spillage within the compressor, or should the glass chamber break during compression, user to follow their own standard operating procedures for the disposal of any such spilt samples and sharps.

The Multi Compressor is a maintenance free unit that does not contain any servicable parts.

Instructions for Use

Before you start, remember to select the Mini-UniPrep G2 Syringeless Filters that is compatible with your sample (see [Appendix, on page 3](#)).

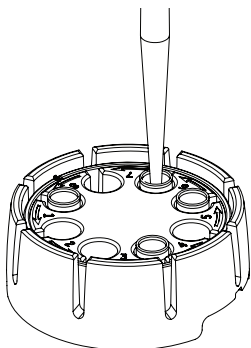
Step	Action
1	Insert up to 8 glass chambers into the Multi Compressor tray.



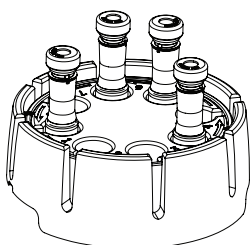
¹ A Hand Compressor (product code MUPG2HCPWC1) is available to order separately.

Step Action

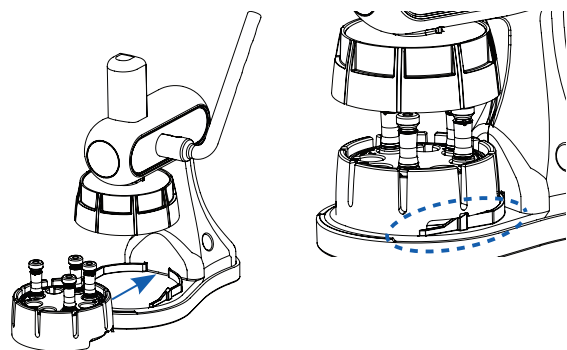
- 2 Dispense the sample into the glass chamber taking care not to overfill. The Mini-UniPrep G2 glass chamber has a maximum capacity of 500 µl indicated by the "Fill Line" on the glass chamber. Minimum sample volume that may be added to the glass chamber is 220 µl in order to collect 50 µl in the glass insert.



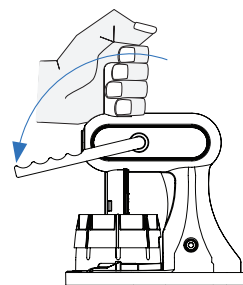
- 3 Place plunger(s) into the neck of the glass chamber(s). Do not attempt to depress further than the neck of the glass chamber



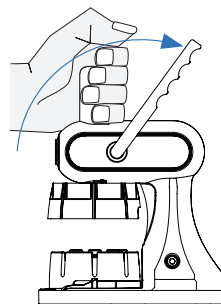
- 4 Place the tray into position on the Multi Compressor using the guide ribs for placement. Ensure tray is securely in place.



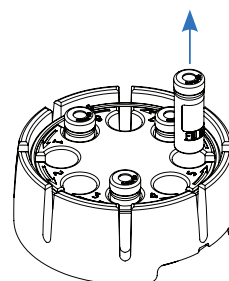
- 5 Steadying the Multi Compressor with the left hand as shown pull the handle towards the user to fully depress the plunger(s) into the chamber(s).

**Step Action**

- 6 Steadying the Multi Compressor with the left hand as shown, rotate the handle back to original position enabling the tray to be removed.



- 7 Remove Mini-UniPrep G2 device or devices.



The Mini-UniPrep G2 prepared sample is now ready for analysis. Process in either of two ways:

- Place the Mini-UniPrep G2 into an autosampler that takes standard 12 mm × 32 mm profile vials for automated sample injection into your instrument. Before your run, ensure that the autosampler needle height is set to a minimum height of 5 mm, to avoid contact with the bottom of the inner glass insert.
- Pierce the septum and draw the filtered sample into a syringe for manual injection into your instrument.

For optimum performance, the assembled and filled Mini-UniPrep G2 device should be kept in an upright position and not be subjected to excessive agitation to prevent the sample being spilled out of the protective inner glass storage reservoir.

Ordering Information

Mini-UniPrep G2 Syringeless Filters are available in different variants such as different membrane types, amber coloured version for light sensitive samples and slit septum caps.

Please visit cytiva.com for more information.

Hand Compressor (product code MUPG2HCPWC1), Multi Compressor (product code MUPG2MCPWC8) and Multi Compressor Tray (MUPG2MCWT8) may also be ordered via this website.

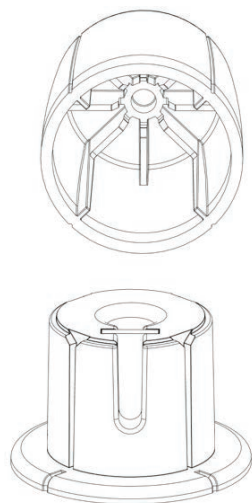


Fig 1. Hand Compressor
MUPG2HCPWC1

Appendix

Table 1. Chemical Compatibility of Membrane

Solvent	GMF	NYL	PP	DpPP	PES	PTFE	PVDF	RC
Acetic Acid 5%	R	R	R	R	R	R	R	R
Acetic Acid, Glacial	R	LR	R	R	R	R	R	NR
Acetone	R	R	R	R	NR	R	NR	R
Acetonitrile	R	R	R	R	NR	R	R	R
Ammonia, 6N	LR	R	R	R	R	R	LR	LR
Amyl Acetate	R	R	R	R	LR	R	LR	R
Amyl Alcohol	R	R	R	R	NR	R	R	R
Benzene ¹	R	LR	NR	NR	R	R	R	R
Benzyl Alcohol ¹	R	LR	R	R	NR	R	R	R
Boric Acid	R	LR	R	R	-	R	R	R
Butyl Alcohol	R	R	R	R	R	R	R	R
Butyl Chloride ¹	R	NR	NR	NR	-	R	R	-
Carbon Tetra- chloride ¹	R	LR	NR	NR	NR	R	R	R
Chloroform ¹	R	NR	LR	LR	NR	R	R	R
Chlorobenzene ¹	R	NR	LR	-	NR	R	R	R
Citric Acid	R	LR	R	-	R	R	R	R
Cresol ¹	R	NR	NR	NR	NR	R	NR	R
Cyclohexanone	R	NR	R	R	NR	R	R	R
Cyclohexane ¹	R	NR	NR	NR	NR	R	R	R
Diethyl Acetamide	R	R	R	R	-	R	NR	R
Dimethyl Forma- mide	R	R	R	R	NR	R	NR	LR
Dioxane	R	R	R	R	LR	R	LR	R
DMSO	R	R	R	R	NR	R	LR	LR
Ethanol	R	R	R	R	R	R	R	R
Ethers ¹	R	R	NR	NR	R	R	LR	R
Ethyl Acetate	R	R	R	R	NR	R	NR	R
Ethylene Glycol	R	R	R	R	R	R	R	R
Formaldehyde ¹	R	R	LR	LR	R	R	R	LR
Freon TF ¹	R	NR	NR	NR	R	R	R	-
Formic Acid	R	NR	R	R	R	R	R	LR
Hexane	R	R	R	R	R	R	R	R
Hydrochloric Acid, Conc ¹	R	NR	LR	LR	R	R	R	NR
Hydrofluoric Acid ¹	NR	NR	LR	LR	-	R	R	NR
Isobutyl Alcohol	R	R	R	R	-	R	R	R
Isopropyl Alcohol	R	R	R	R	-	R	R	R

Solvent	GMF	NYL	PP	DpPP	PES	PTFE	PVDF	RC
Methanol	R	R	R	R	R	R	R	R
Methyl Ethyl Ketone	R	R	R	R	NR	R	NR	R
Methylene Chloride ¹	R	NR	LR	LR	NR	R	R	R
Nitric Acid, Conc ¹	R	NR	NR	NR	NR	R	R	NR
Nitric Acid, 6N ¹	R	NR	LR	LR	LR	R	R	LR
Nitrobenzene ¹	R	LR	R	R	NR	R	R	R
Pentane ¹	R	R	NR	NR	R	R	R	R
Perchloro Ethylene ¹	R	LR	NR	NR	NR	R	R	R
Phenol, 0.5%	R	NR	R	R	NR	R	R	R
Pyridine	R	LR	R	R	NR	R	NR	R
Sodium Hydroxide, 6N	NR	LR	R	R	R	R	NR	NR
Sulfuric Acid, Conc ¹	R	NR	NR	NR	NR	R	NR	NR
Tetrahydrofuran	R	R	LR	LR	NR	R	R	R
Toluene ¹	R	LR	LR	LR	NR	R	R	R
Trichloroethane ¹	R	LR	LR	LR	NR	R	R	R
Trichloroethylene ¹	R	NR	LR	LR	NR	R	R	R
Water	R	R	R	R	R	R	R	R
Xylene ¹	R	LR	LR	LR	LR	R	R	R

¹ Short term resistance of housing.

Note: R = Resistant; LR = Limited Resistance; NR = Not recommended.

The table data is to be used as a guide only. Testing prior to application is recommended.

Material Abbreviations

GMF - Glass Microfiber

NYL - Nylon

PP - Polypropylene

DpPP - Polypropylene Depth Filter

PES - Polyethersulfone

PTFE - Polytetrafluoroethylene

PVDF - Polyvinylidene Difluoride

RC - Regenerated Cellulose

Membrane Guide

Type	Applications
Glass Microfiber (GMF)	High loading capacity, binder-free, glass microfiber depth filter medium.
Nylon (NYL)	Membrane for aqueous and organic samples within a pH range of 3–10.
Polyethersulfone (PES)	For aqueous-based samples. Good flow rates.
Polypropylene (Dp PP)	Aqueous and organic samples; high loading capacity depth filter.
Polypropylene (PP)	For solvent-based samples, low water breakthrough values.
Polytetrafluoroethylene (PTFE)	For aggressive samples. Great chemical compatibility.

Type	Applications
Polyvinylidene difluoride (PVDF) (Hydrophilic)	Low nonspecific protein binding membrane for samples in aqueous solutions or organic solvents.
Regenerated Cellulose (RC)	Suitable for use with either aqueous or organic solvents. Very low non-specific protein binding membrane.

cytiva.com

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

Mini-UniPrep and Whatman™ are trademarks 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

29036949 AC V:7 02/2021

