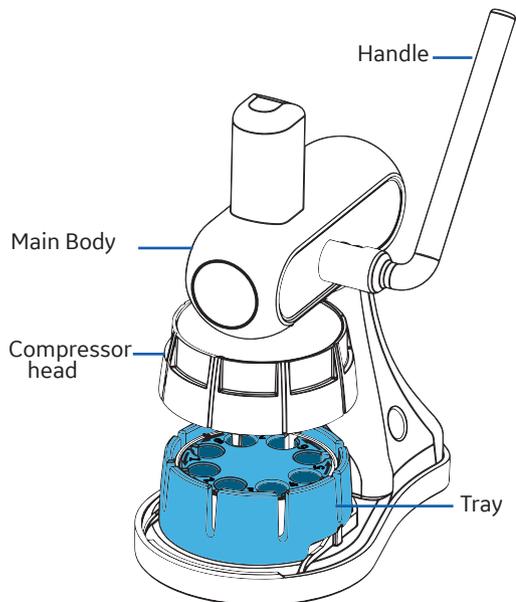


Whatman Mini-UniPrep Syringeless Filters

Multi Compressor Product Information sheet

Introduction

About



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

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 Data Sheets.

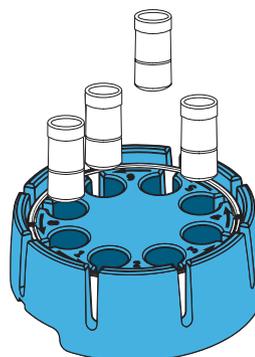
Instructions for Use

Before you start, remember to use the blue tray designed for use with the plastic Mini-UniPrep™ and select the Mini-UniPrep filter that is compatible with your sample. Chemical compatibility information can be found at: cytiva.com

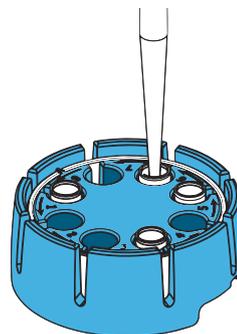
Note: Please do not use a white Multi Compressor tray to plunge the Mini-UniPrep plastic vials as this product will not be fully compressed. Always use blue Multi Compressor tray.

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

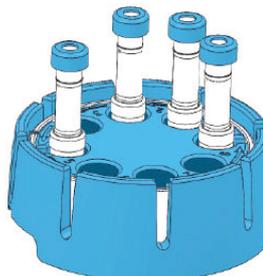
1	Insert up to 8 Mini-UniPrep plastic chambers into the Multi Compressor tray.
---	--



2	Dispense the sample into the chamber up to the fill line taking care not to overfill. The Mini-UniPrep chamber has a capacity of 400 µL indicated by the fill line. Minimum sample volume that may be added to the chamber is 135 µL in order to collect 50 µL in the plunger.
---	--

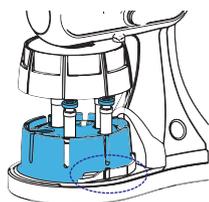
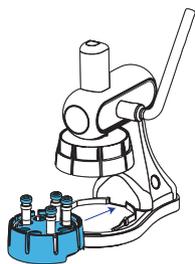


3	Place plunger(s) into the neck of the chamber(s).
---	---

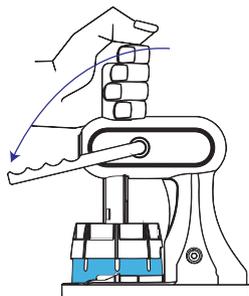


Step Action

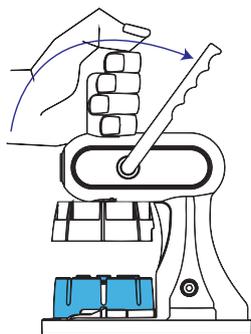
- 4 Place the tray into position on the Multi Compressor using the guide ribs for placement. Ensure tray is securely in place.



- 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).

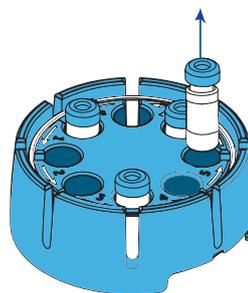


- 6 Steadying the Multi Compressor with the left hand as shown, rotate the handle back to original position enabling the tray to be removed.



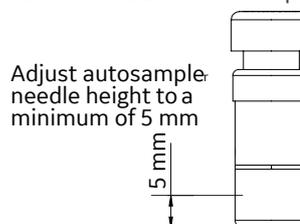
- 7 Remove Mini-UniPrep device or devices.

Step Action



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

- Place the Mini-UniPrep into an autosampler that takes standard 12 mm x 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 5mm, to avoid contact with the bottom of the plunger.



- 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 device should be kept in an upright position.

Ordering Information

Mini-UniPrep 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.

Multi Compressor for plastic Mini-UniPrep (Product code MUPMCPBC8) and Multi Compressor Tray for plastic Mini-UniPrep (MUPMCBT8) may also be ordered via this website.

Membrane Guide

Type	Applications
Glass Microfiber (GMF)	High loading capacity, binderfree, 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.
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/](https://www.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](https://www.cytiva.com/contact)

29344968 AB V:5 02/2021

