

Whatman Mini-UniPrep G2 Syringeless Filters 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.

Safety Precaution



IMPORTANT

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. Mini-UniPrep™ G2 starter packs contain 100 Mini-UniPrep G2 syringeless filters and a Hand Compressor. If a Hand Compressor or Multi 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 (Product code. 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 devices at one time.

¹ Hand Compressor (catalogue no. MUPG2HCPWC1) and Multi Compressor (catalogue no. MUPG2MCPWC8) is available to order separately. See [cytiva.com](https://www.cytiva.com) for further details.

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

29020242 AC V:4 02/2021

