

Prefilter 6000

Instructions for Use

Characteristics

The Prefilter (Code No. 18458201) filters buffers and solvents at pressures up to 4 MPa (40 bar, 569 psi).

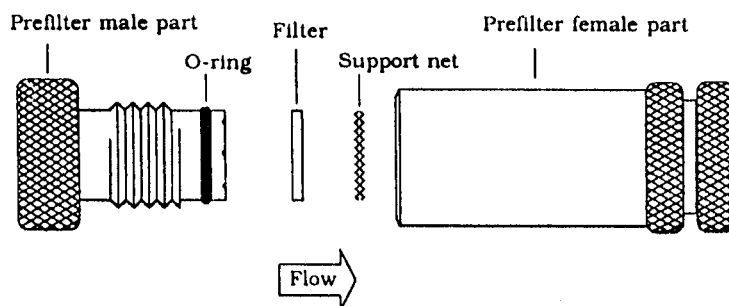


Figure 1: Disassembled prefilter.

Unpacking

Please check the contents against this list.

Designation	Number supplied
Prefilter	1
Filter kit	1

Connecting the prefilter

Install the Prefilter between the pump and the mixer.

To replace the filter

Step	Action
1	Unscrew the two sections of the Prefilter.
2	Remove the filter with a spatula.
3	Insert a new filter and press it down with a spatula.
4	Screw the two parts of the Prefilter together.

Spare parts

Designation	Number per pack	Code number
Filters + O-rings + support nets	50+5+5	18458401

Page intentionally left blank



cytiva.com

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

All other third-party trademarks are the property of their respective owners.

© 2020 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

56458300 AE V:4 10/2020