

User Guide

OptiScale®-25 Capsule Filter

Introduction

The OptiScale®-25 capsule filter is a single use filter designed for process development and scaling. It is not designed or validated as a sterilizing grade filter.

Specifications

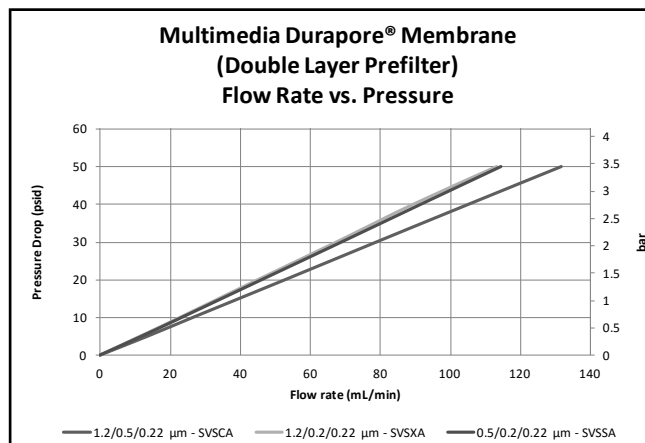
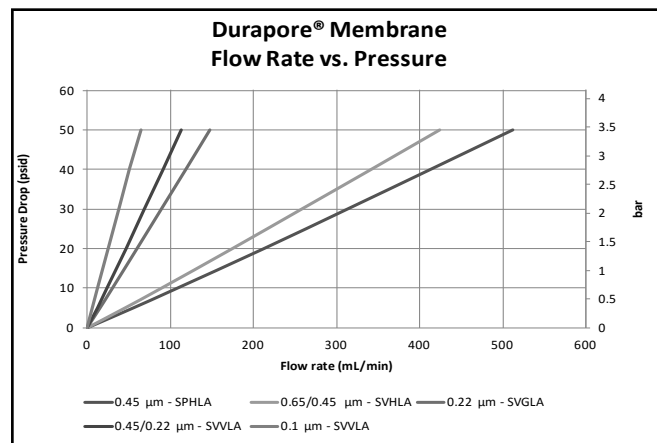
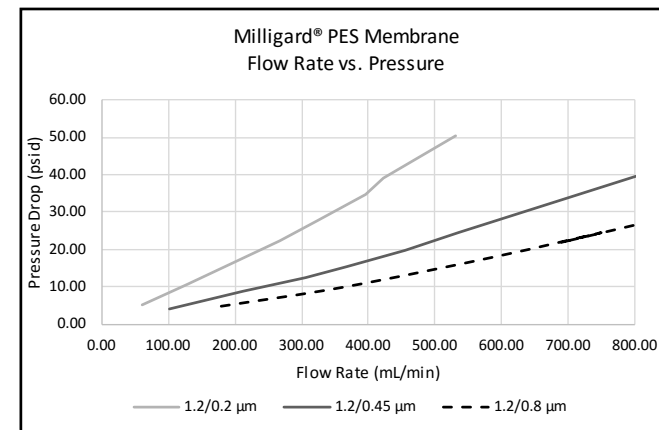
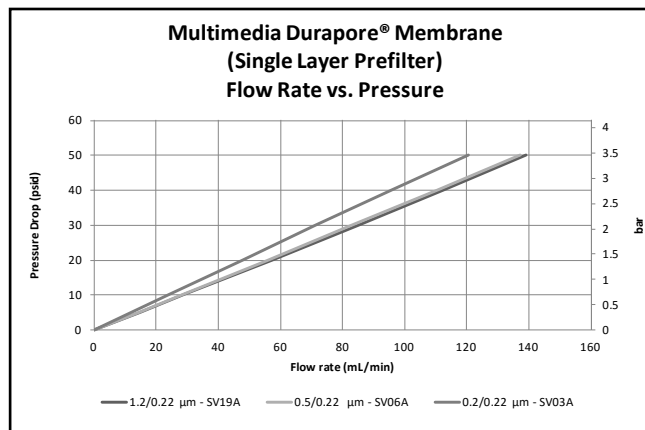
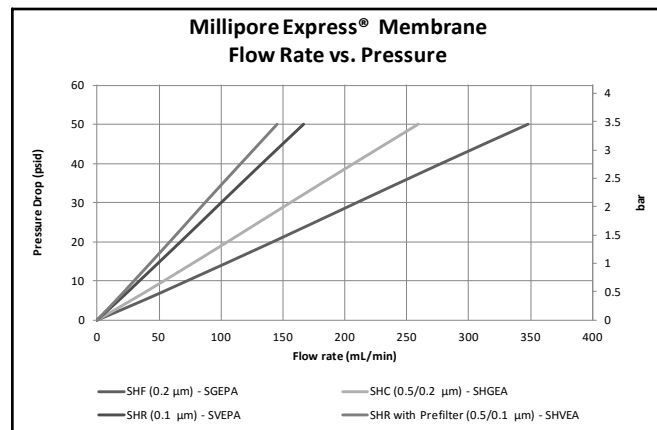
Please refer to the Certificate of Quality for complete specifications.

Maximum forward operating pressure	4.1 bar (60 psi) at ambient temperature. Do not apply reverse pressure to the filter.
Storage	Store in original packaging at room temperature.
Sterilization OptiScale®-25 capsule filters are single use devices.	Filter may be autoclaved one time for 60 minutes at 123 °C. Do not clean in place. Do not steam sterilize. Do not expose to gamma radiation.
Connection Fittings	Inlet: Female Luer-Lok™ Outlet: Male Luer slip Vent: Female Luer-Lok™ with cap
Effective Surface Area	3.50 cm ²
Labeling	Each filter is labeled with membrane type and pore size. Box labels contain product and lot data.



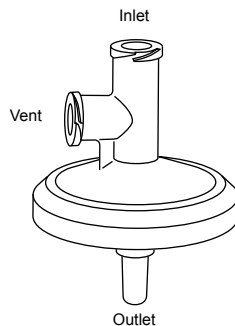
Flow Rate vs. Pressure

Refer to typical water flow rates below to verify full wetting of the membrane prior to use. Flow rates shown are at 23°C.



Installation

1. Remove filter from packaging.
2. Connect the filter inlet to the inlet tubing.
3. Connect the filter outlet to the outlet tubing.
4. Install the filter in the vertical position with the inlet on top and outlet on bottom to avoid air lock.
5. Autoclave if required.



Wetting

Notes

- Avoid introducing air into the filter.
 - The filter must be completely wetted. Any residual air in the filter may impair fluid flow and reduce the capacity of the filter.
 - Install the filter in the vertical position with the inlet on top and outlet on bottom to avoid air lock.
1. If a Luer-Lok™ valve is installed on the filter outlet, close the valve to prevent filtrate flow.
 2. Slowly introduce the process fluid through the capsule inlet port.
 3. Open the vent until no air escapes through the vent and process fluid is visible.
 4. Close the vent.

Standard Warranty

The applicable warranty for the products listed in this publication may be found at www.emdmillipore.com/terms (within the "Terms and Conditions of Sale" applicable to your purchase transaction).

Technical Assistance

For more information, contact the office nearest you or visit the Technical Service page at www.emdmillipore.com/techservice. Worldwide contact information is available at www.emdmillipore.com/offices.

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

For more information and documentation please contact:
Merck KGaA, Darmstadt, Germany
Corporation with General Partners
Frankfurter Str. 250
64293 Darmstadt, Germany
Phone: + 49 6151-72 0

For requests from USA and Canada please contact:
EMD Millipore Corporation
A subsidiary of Merck KGaA, Darmstadt, Germany
290 Concord Road
Billerica, MA 01821, USA
Phone: 1-800-645-5476

www.emdmillipore.com

The Vibrant M, Milligard, Millipore Express, Durapore and OptiScale are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2010-2017 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

00113893PU, Rev. 4 11/2017

