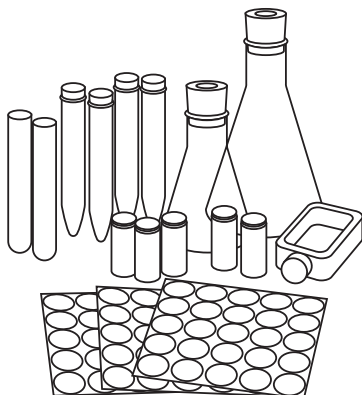


User Guide

Milliseal™ Adhesive Membranes

FWMS01800



Introduction

Milliseal™ Adhesive Membrane discs are 0.45 µm hydrophobic polytetrafluoroethylene (PTFE) membranes with all the compatibility characteristics of Fluoropore™ membranes. Milliseal™ discs are applied as a clean barrier and allow the sample to vent while preventing outside contamination. The 18 mm disc has a 4 mm band of adhesive around the edge which allows it to adhere to the top of any small volume container, including glass, stainless steel, and some plastics. The covered container can then be transported or stored in an open area without risk of outside contamination. Milliseal™ membrane discs are designed to replace metal caps, Parafilm® film, cotton plugs, stoppers, and other tube coverings, especially when venting is necessary.

How to Use Milliseal™ Membrane Discs

1. To remove the membrane disc from the sheet, place your forefinger on the top center of the disc. With membrane forceps, rub from the edge of the disc inward to lift the edge of the disc.
2. Pull the disc off the sheet with the forceps.
3. Center the disc on the top of the container, adhesive side down, then press it in place with your forefinger or thumb. Adhere securely by pressing it down with your thumb.

Ordering Information

Purchase this product on-line at SigmaAldrich.com.

| Description | Qty/Pk | Cat. No. |
|------------------------------------------|--------|-----------|
| Milliseal™ Adhesive Membrane, 18 mm disc | 375 | FWMS01800 |

Notice

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.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Technical Assistance

Visit the tech service page on our web site at SigmaAldrich.com/TechService.

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at SigmaAldrich.com/Terms.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/Offices.

Merck, Millipore, Milliseal, and Sigma-Aldrich 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. © 1999-2025 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

MERCK