

Sealing Foil F-RSXT-10

Gas permeable sealing foil with evaporation reduction and perforated silicone layer for the BioLector XT Microbioreactor

TECHNICAL PARAMETERS

TYPE

Two-layer sandwich silicone mat

3 mm venting holes to reduce cross section for evaporation
4 mm positioning holes for simple application

Top layer: silicone mat thickness 2 mm, including venting holes and pre-cut slits for automated or manual sampling/dosing

Bottom layer: non-gas permeable bonding sheet (glue layer) including venting holes, already glued to top layer

Separate layer: gas permeable non-woven membrane (part no. F-GP-10), to be ordered separately

Self resealing pre-cut slits for penetration by syringe (manual) or robotic tips (automated)

Suitable for foil application with the Foil Applicator XT

For non-robotic experiments (per run): 1x F-RSXT-10 + 1x F-GP-10¹

For robotic experiments (per run): 2x F-RSXT-10 + 1x F-GP-10¹

DIMENSION

Length: 121.9 mm

Width: 80.4 mm

STERILIZATION

20 kGy (β-radiation)

Sterile packed in single bags

OTHER INFORMATION

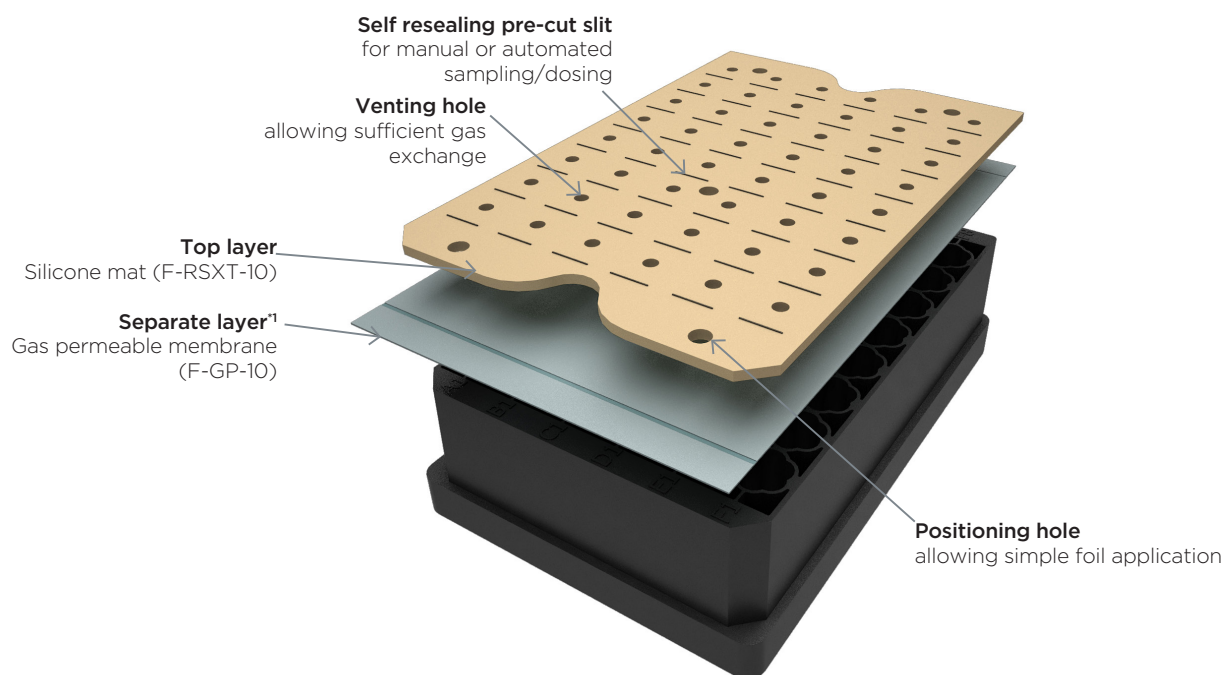
Autoclavability: no

For single use only

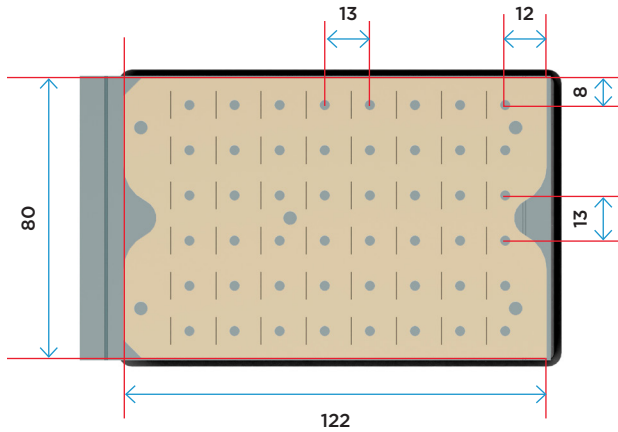
Unit size: 1x 10 pcs. in sterilized bag

Medical grade adhesive

Temperature range: -20 - 80°C



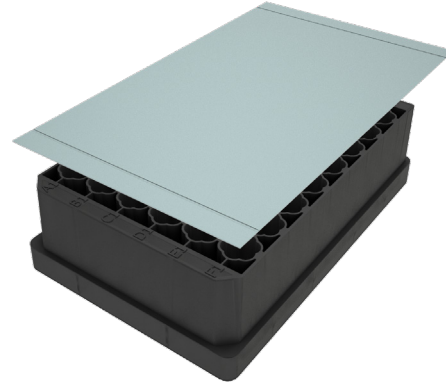
TOP VIEW DIMENSIONS FOR AUTOMATION [mm]



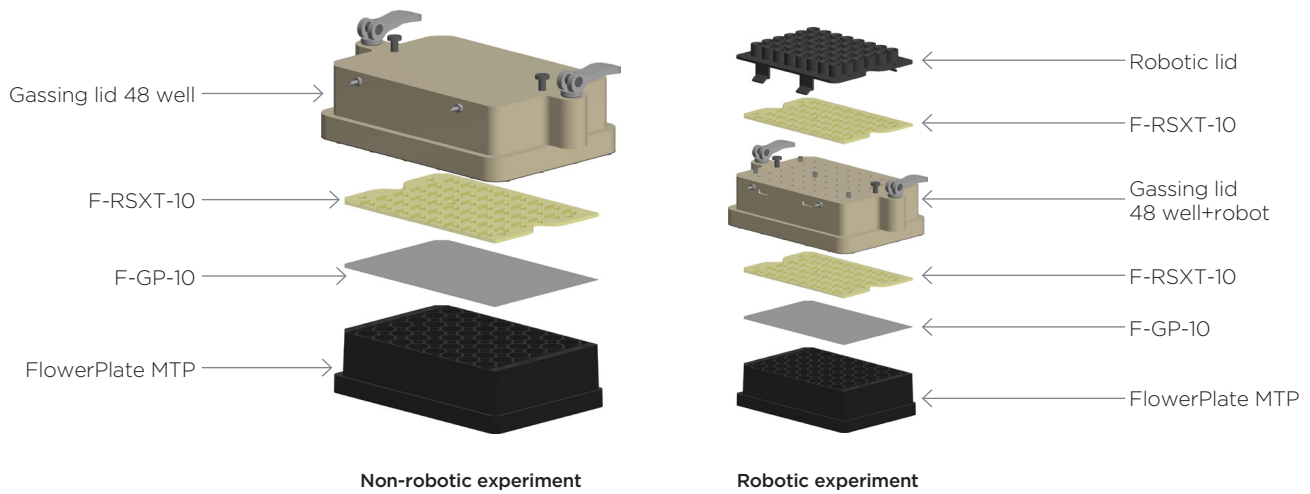
ALSO AVAILABLE SPECIALIZED SEALING FOILS

PART NO. F-GP-10

Non-woven, gas permeable membrane which serves as sterile barrier



ASSEMBLY OF MICROTITER PLATE, FOILS AND GASSING HEAD



*To be ordered separately

m2p and the m2p logo are the trademarks or registered trademarks of m2p-labs GmbH in the US and other countries. m2p-labs is a Beckman Coulter company. Information contained in this document or drawing is proprietary to m2p-labs GmbH. This document may not be reproduced in the whole or in parts for any reason without written permission from m2p-labs GmbH. All rights of design, invention and copyright are reserved.

© 2021 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at www.m2p-labs.com, beckman.com 21.04.1437.PCC

