

MatTek Permeation Fixture (EPI-100-FIX) User Note

- a) The fixture is a 2 piece device with 4 screws connecting the top to the bottom. Remove the screws and place the Millicell in the bottom piece.
- b) Slide the tube (annulus) of the top piece into the Millicell.
- c) Return the screws to close the device.
- d) Tighten snugly by hand, and then further tighten an additional half turn with pliers or an Allen wrench. **The screws should not be over-tightened.** Any tightening which will deform the plastic device is too tight and will most likely compromise the seal.
- e) Place the fixture in a 6-well plate and add the donor solution into the annulus of the top piece, with the receiver solution in the 6-well plate.

Note: Care should be taken to ensure no air bubbles are trapped beneath the fixture. By submerging and tilting the device the air bubbles should release.

Dimensions and Volumes: The diameter of the inner annulus is 0.57 cm and the area is 0.256 cm². The receiver solution should be 5.0 ml and the donor solution can be up to 0.5 ml.

Re-use of MatTek Permeation Fixtures: The MatTek Permeations Fixtures can be sterilized by soaking in 70% ethanol for 30 minutes. However, re-sterilization by autoclaving should NOT be done since autoclaving will result in deformation of the device.

