

Features

- Triple Cell Culture Model
- Optional incorporation of macrophages
- Highly reproducible
- Easily-handled cell culture inserts
- Cost effective alternative to animal and clinical testing

Key Applications

- Inhalation Toxicity
- Inhaled Drug Delivery
- Pulmonary Fibrosis
- Viral Infection

The EpiAlveolar Model

Mattek's EpiAlveolar System consists of normal, human-derived alveolar epithelial cells co-cultured with donor-matched fibroblasts and human microvascular endothelial cells on specially prepared, optically clear cell culture inserts. Human-derived macrophages can also be incorporated into the three-dimensional, highly differentiated model.

A



B

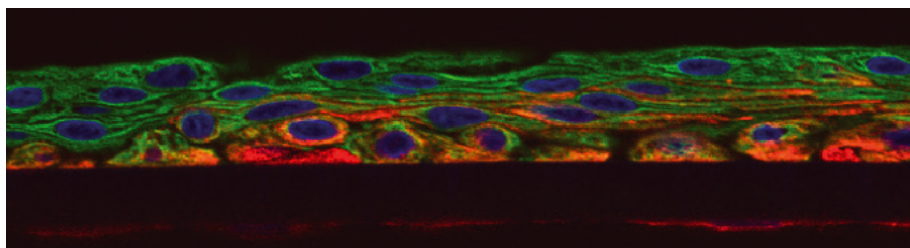
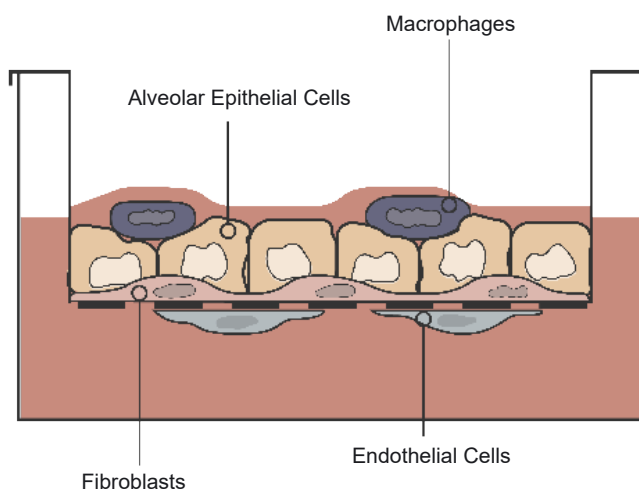
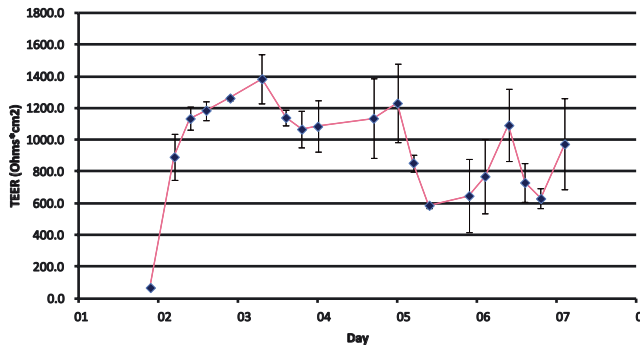


Figure 1. Cross-sections of EpiAlveolar tissue stained with: a) H&E and b) Cytokeratin 19 (green, epithelial cells), Vimentin (red, fibroblasts), DAPI (blue, nuclei.)

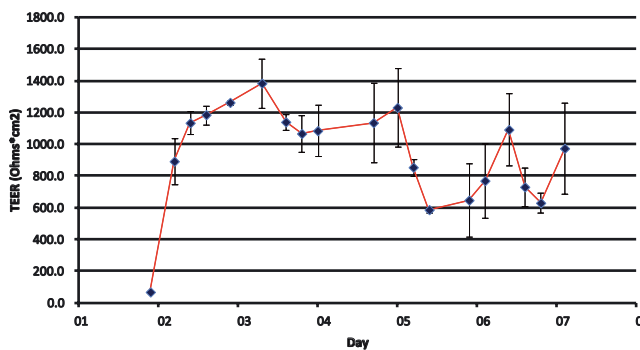
Schematic diagram of EpiAlveolar tetra-culture system containing primary endothelial cells, fibroblasts, epithelial cells and optional macrophages.



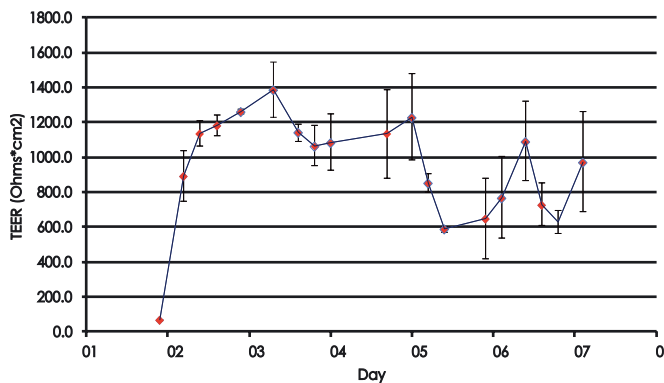
EpiAlveolar Barrier Function



Alv-100-FT-PE12 Transepithelial Electrical Resistance



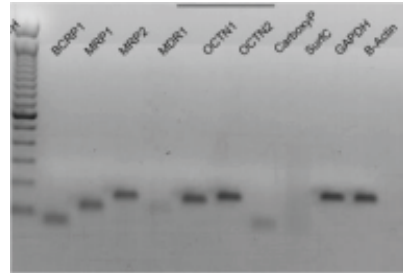
EpiAlveolar Barrier Function



EpiAlveolar maintains its barrier function over extended culture periods: Transepithelial electrical resistance (TEER) >300 Ohms*cm² for up to 60 days. Transepithelial electrical resistance (TEER) >300 Ohms*cm² for up to 60 days.

Inhalation Studies

EpiAlveolar expresses several ABC transporters, including BCRP, MRP1 and MRP2, as well as organic cation uptake transporters, making it a useful tool for inhaled drug delivery and inhalation toxicity studies. EpiAlveolar has also been used to study fibrosis and deep lung effects of nano-particles.



Pulmonary Fibrosis

Studies have shown that treatment of EpiAlveolar with TGF-β induces the production of the pro-fibrotic markers Collagen 1a1, α-Smooth Muscle Actin and Fibronectin, making it a useful model for the study of pulmonary fibrosis.

