



Cultrex[®]

CultreCoat[®] Vascular Permeability Assay Kits

For quantitating the degree of vascular monolayer permeability

Trevigen's Cultrex[®] 96 Well Vascular Permeability Assays accelerate the screening process for pathways and compounds that influence vascular permeability. These assays offer a flexible, standardized, high-throughput format for quantitating the degree to which genes or compounds can influence the maintenance of endothelial cell-to-cell adhesion.

Ordering Information:

Description	Size	Catalog #
CultreCoat [®] 24 Well Vascular Permeability Assay	24 Samples	3475-024-K
CultreCoat [®] 96 Well Vascular Permeability Assay	96 Samples	3475-096-K

CultreCoat[®] Vascular Permeability Assay Features:



- Pre-coated inserts save time
- Pre-coated inserts improve reproducibility of results
- Complete kit format
- Coating process assures uniform layers of BME plate to plate, and lot to lot, to minimize variability

Cultrex[®] Cell Invasion Assay Features:

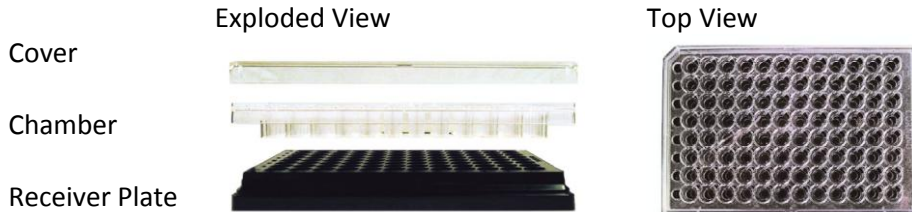
- Collagen I coated membranes.
- Kit components include vascular permeability chamber and FITC-Dextran.
- The assay has been functionally tested with HUVEC and SVEC4-10 cell lines.
- Each lot of CultreCoate Vascular Permeability Assay kits has been tested to ensure that HUVEC, primary endothelial cells, attaches, proliferates, and forms cell-cell adhesions sufficient to block 95% of FITC-Dextran permeability.

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS



CultreCoat[®] Vascular Permeability Assay Kits

Vascular Permeability Plate (96 well plate pictured)



- The receiver plate is compatible with 96 well fluorescent readers, eliminating the requirement for transfer, saving time and reducing error.
- The receiver plate was designed specifically for fluorescence detection and its black surface eliminates crosstalk between wells and minimizes background.
- The chamber is designed so that the assay may be performed without disassembly until the final analysis. This format helps prevent contamination from handling and may be adapted for high throughput screening (robotic).

Related Products:

Description	Size	Catalog#
CultreCoat [®] Cell Invasion Optimization Assay Sampler	96 wells	3484-096-K
CultreCoat [®] Cell Invasion Optimization Assay Sampler	24 wells	3484-024-K
CultreCoat [®] Adhesion Protein Array Kit	96 wells	3496-096-K
CultreCoat [®] BME 96 Well Cell Adhesion Assay	96 wells	3490-096-K
CultreCoat [®] BME 24 Well Cell Adhesion Assay	24 wells	3480-024-K
CultreCoat [®] Laminin I 96 Well Cell Adhesion Assay	96 wells	3491-096-K
CultreCoat [®] Collagen I 96 Well Cell Adhesion Assay	96 wells	3492-096-K
CultreCoat [®] Collagen IV 96 Well Cell Adhesion Assay	96 wells	3493-096-K
CultreCoat [®] Fibronectin 96 Well Cell Adhesion Assay	96 wells	3494-096-K
CultreCoat [®] Vitronectin 96 Well Cell Adhesion Assay	96 wells	3495-096-K
Cultrex [®] BME with Phenol Red	5 ml	3430-005-01
Cultrex [®] BME PathClear [®] with Phenol Red	5 ml	3430-005-02
Cultrex [®] BME with Phenol Red, Reduced Growth Factor	5 ml	3431-005-01
Cultrex [®] BME with Phenol Red, PathClear [®] Reduced Growth Factor	5 ml	3431-005-02
Cultrex [®] BME, No Phenol Red	5 ml	3432-005-01
Cultrex [®] BME, PathClear [®] No Phenol Red	5 ml	3432-005-02
Cultrex [®] BME, No Phenol Red, Reduced Growth Factor	5 ml	3433-005-01
Cultrex [®] BME, PathClear [®] No Phenol Red, Reduced Growth Factor	5 ml	3433-005-02

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS