

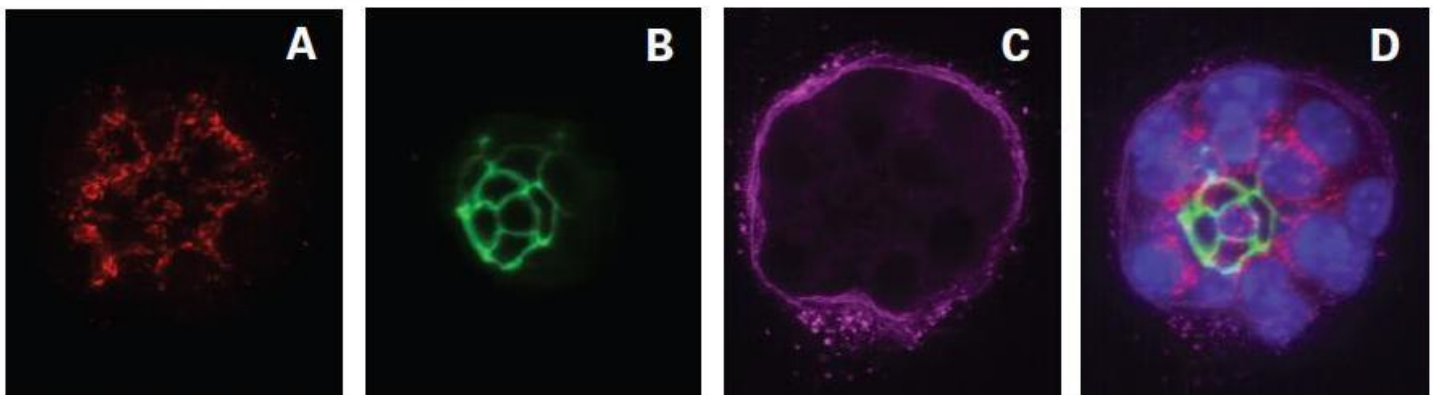


Cultrex® 3D Culture Products

All Cultrex® Brand products are designed, developed and delivered by Trevigen, Inc.

3D Culture Matrix™ Basement Membrane Matrix (BME), RGF

Trevigen's 3D Culture Matrix BME, RGF has been developed, produced and qualified specifically for use in 3D culture studies. Cultrex 3D Culture Matrix RGF BME is an extracellular matrix hydrogel that directs cells to grow in three dimensions and assemble into organotypic structures *in vitro*. Cultrex 3D Culture Matrix RGF BME is a standardized basement membrane matrix for use in 3D cultures, and a special process has been employed to reduce growth factors and provide a protein concentration of approximately 15 mg/ml.



Non transgenic primary mammary cells grown on Cultrex 3D Culture Matrix develop into a polarized acinus. Confocal microscopy (5 μ m projection) demonstrates epithelial polarity: DAPI stain, blue; GM130, red (Golgi protein, apical marker: panel A); ZO1, green (tight junction, apical; panel B); Integrin α 6, magenta (baso-lateral: panel C), overlay shown in panel D. Image courtesy of Martin Jecklinger.

Name	Catalog Number	Size
Cultrex® 3D Culture Matrix™ Reduced Growth Factor BME	3445-011-01	1 ml
	3445-005-01	5 ml
	3445-010-01	10 ml

3D Culture Matrix™ BME Coated Plate

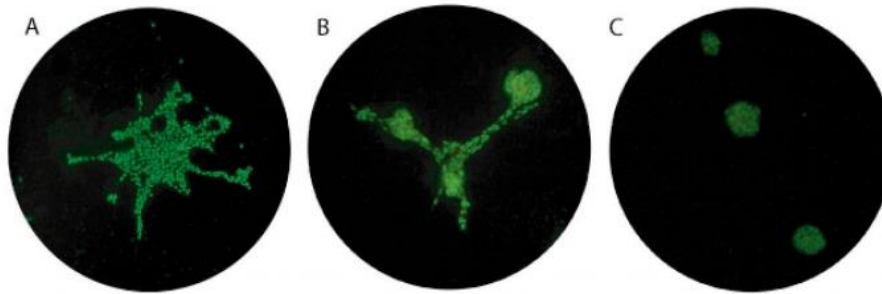
The 3D Culture Matrix BME Coated Plate is coated with a reconstituted basement membrane matrix to provide the environment that is common for epithelial or endothelial cell types. This plate provides a convenient, standardized, physiologically predictive format for evaluating the pharmacological effects of compounds in these cell culture models.

Name	Catalog Number	Size
Cultrex® 3D Culture Matrix™ BME Coated Plate	3445-096-CP	1 plate



3D Culture Matrix™ Rat Collagen I

Trevigen's 3D Culture Matrix Rat Collagen I is a purified stromal ECM protein that has been developed, produced and qualified specifically for use in 3D culture studies. 3D Culture Matrix Collagen I is an extracellular matrix hydrogel that directs cells to grow in three dimensions and assemble into organotypic structures *in vitro*. 3D Culture Matrix® Rat Collagen I may also be used to supplement customized hydrogel or medium formulations for cell culture.



Mammary epithelial cells, MCF-10A cultured on 3D Culture Matrix Rat Collagen I are induced to differentiate with the addition of 3D Culture Laminin I at : a) 9mg/ml, b) 1 mg/ml, and c) 2mg/mL.

Name	Catalog Number	Size
Cultrex® 3D Culture Matrix™ Rat Collagen I	3447-020-01	20 ml

3D Culture Matrix™ Laminin I

The 3D Culture Matrix Laminin I is a purified basement membrane protein that has been developed, produced and qualified specifically for use in 3D culture studies. Cultrex 3D Culture Matrix Laminin I is an extracellular matrix hydrogel that directs cells to grow in three dimensions and assemble into organotypic structures *in vitro*. Cultrex 3D Culture Matrix Laminin I may also be used to supplement customized hydrogel or medium formulations for cell culture.

Name	Catalog Number	Size
Cultrex® 3D Culture Matrix™ Laminin I	3446-005-01	30 mg

3D Culture Cell Harvesting Kit

The 3D Culture Cell Harvesting Kit provides an optimized and standardized solution for the isolation of intact 3D structures from Basement Membrane Extract (BME) Matrix or Laminin I hydrogels and preparation of cell lysates for subsequent biochemical analysis.

Name	Catalog Number	Size
Cultrex® 3D Culture Cell Harvesting Kit	3448-020-K	20 tests

Trevigen, Inc.

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Phone: 301-216-2800 • Fax: 301-560-4973 • Email: info@trevigen.com • <http://www.trevigen.com>