

# CULTREX<sup>®</sup> Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

## Mouse Laminin I

Catalog #: 3400-010-01

Size: 1 mg

**Description:** This highly purified preparation of mouse laminin I increases cell adhesion, migration, growth, and differentiation. It is composed of  $\alpha 1\beta 1\gamma 1$  chains with a total Mr of 800,000 and is used for coating culture dishes.

**Source:** Murine Engelbreth-Holm-Swarm (EHS) tumor.

**Storage Buffer:** Dulbecco's Modified Eagle's Medium with 10  $\mu$ g/ml gentamycin sulfate.

**Storage:** Store at -20°C or -80°C in a manual defrost freezer.

**Purity:** Purity >90% by SDS-PAGE.

### Specifications:

#### Functional Assays:

- Supports the attachment (at least 50%) of MG63 human osteosarcoma cells to a laminin-coated plate at a coating solution concentration of 20  $\mu$ g/ml.

#### Sterility Testing:

- No bacterial or fungal growth detected after incubation at 37°C for 14 days following USP XXIV, Chapter 71 sterility testing.
- No mycoplasma contamination detected by PCR
- Endotoxin concentrations < 20 EU/ml by LAL assay.

### Coating procedure:

The recommended working concentration is 0.05-10  $\mu$ g/cm<sup>2</sup> of growth surface (0.05 -20  $\mu$ g/ml) depending on cell type.

**A.** Thaw stock solution on ice for several hours. In a laminar flow hood, dilute appropriately with cold serum-free, cell culture medium. Mix and transfer to the wells of tissue culture plates. Spread the solution to completely cover the bottom of the wells.

**B.** The following table is a guide for the suggested volumes required per well:

Plate type/	Volume Laminin per Well
6 wells (or 35 mm dish)	2 ml
24 wells	1 ml
48 wells	200 $\mu$ l
96 wells	100 $\mu$ l

**C.** Incubate the plates at 37°C from 4 hours to overnight. In the laminar flow hood, remove excess liquid from the wells of the tissue culture plate.

**D.** Rinse the wells once with cell culture medium and then add your cells.

### Related Products:

Catalog#	Description	Size
3455-024-K	Cultrex <sup>®</sup> 24 Well BME Cell Invasion Assay	24 inserts
3460-024-K	CultreCoat <sup>®</sup> 24 well BME-coated Cell Invasion Assay	24 inserts
3465-024-K	Cultrex <sup>®</sup> 24 Well Cell Migration Assay	24 inserts
3465-096-K	Cultrex <sup>®</sup> 96 Well Cell Migration Assay	96 samples
3455-096-K	Cultrex <sup>®</sup> BME Cell Invasion Assay	96 samples
3456-096-K	Cultrex <sup>®</sup> Laminin I Cell Invasion Assay	96 samples
3457-096-K	Cultrex <sup>®</sup> Collagen I Cell Invasion Assay	96 samples
3458-096-K	Cultrex <sup>®</sup> Collagen IV Cell Invasion Assay	96 samples

### Accessories:

Catalog#	Description	Size
3446-005-01	Cultrex <sup>®</sup> 3-D Culture Matrix <sup>™</sup> Laminin I	5 ml
3440-100-01	Cultrex <sup>®</sup> Rat Collagen I	100 mg
3442-050-01	Cultrex <sup>®</sup> Bovine Collagen I	50 mg
3447-020-01	Cultrex <sup>®</sup> 3-D Culture Matrix <sup>™</sup> Collagen I	100 mg
3410-010-01	Cultrex <sup>®</sup> Mouse Collagen IV	1 mg
3430-005-02	Cultrex <sup>®</sup> BME with Phenol Red, PathClear <sup>®</sup>	5 ml
3431-005-02	Cultrex <sup>®</sup> BME with Phenol Red, Growth Factor Reduced PathClear <sup>®</sup>	5 ml
3432-005-02	Cultrex <sup>®</sup> BME, PathClear <sup>®</sup>	5 ml
3433-005-02	Cultrex <sup>®</sup> BME Growth Factor Reduced, PathClear <sup>®</sup>	5 ml
3445-048-01	Cultrex <sup>®</sup> 3-D Culture Matrix <sup>™</sup> BME	15 ml
3430-005-01	Cultrex <sup>®</sup> BME with Phenol Red	5 ml
3432-005-01	Cultrex <sup>®</sup> BME, no Phenol Red	5 ml

### Reference:

Kleinman H.K. 1999. The Laminins. Proc. Indian Acad. Sci. (Chem.Sci.), 111:179-1841.

© 2008 Trevigen, Inc. All Rights Reserved. 3-D Culture Matrix is a trademark, and Trevigen, Cultrex, CultreCoat and PathClear are registered trademarks of Trevigen, Inc. E5/9/08v1

## TREVIGEN<sup>®</sup>

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: info@trevigen.com • www.trevigen.com



**Mouse Laminin I**  
Cat#: 3400-010-01  
Storage:  $\leq$  -20 °C  
(Manual Defrost)  
1-800-873-8443