

# CULTREX<sup>®</sup> Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

## Mouse Laminin I, PathClear<sup>®</sup>

Catalog #: 3400-010-02

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  $\mu\text{g/ml}$  gentamycin sulfate.  
**Storage:** Store at  $-20^{\circ}\text{C}$  in a manual defrost freezer or  $-80^{\circ}\text{C}$ .  
**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  $\mu\text{g/ml}$ .

#### Sterility Testing:

- No bacterial or fungal growth detected after incubation at  $37^{\circ}\text{C}$  for 14 days following USP XXIV, Chapter 71 sterility testing.
- PathClear<sup>®</sup>: Negative by PCR test for: mycoplasma, 17 bacterial and virus strains typically included in mouse antibody production (MAP) testing, plus 13 additional murine infectious agents including LDEV, for a total of 31 organisms and viruses.
- Endotoxin concentrations < 20 EU/ml by LAL assay.

### Coating procedure:

The recommended working concentration is 0.05-10  $\mu\text{g/cm}^2$  of growth surface (0.05 -10  $\mu\text{g/ml}$ ) 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 $\mu\text{l}$
96 wells	100 $\mu\text{l}$

**C.** Incubate the plates at  $37^{\circ}\text{C}$  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.

© 2009 Trevigen, Inc. All Rights Reserved. Trevigen, Cultrex, CultreCoat and PathClear are registered trademarks, and CellSpere and 3-D Culture Matrix are trademarks of Trevigen, Inc. E7/30/09v1

## 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

### Related Products:

Catalog#	Description	Size
3455-024-K	Cultrex <sup>®</sup> 24 Well BME Cell Invasion Assay	24 inserts
3480-024-K	CultreCoat <sup>®</sup> 24 Well BME-Coated Cell Invasion Assay	24 inserts
3456-024-K	Cultrex <sup>®</sup> 24 Well Laminin I Cell Invasion Assay	24 inserts
3457-024-K	Cultrex <sup>®</sup> 24 Well Collagen I Cell Invasion Assay	24 inserts
3458-024-K	Cultrex <sup>®</sup> 24 Well Collagen IV Cell Invasion Assay	24 inserts
3455-096-K	Cultrex <sup>®</sup> 96 well BME Cell Invasion Assay	96 samples
3465-024-K	Cultrex <sup>®</sup> 24 well Migration Cell Assay	96 samples
3465-096-K	Cultrex <sup>®</sup> 96 well Migration Cell 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
3490-096-K	CultreCoat <sup>®</sup> BME 96 Well Cell Adhesion Assay	96 samples
3496-096-K	CultreCoat <sup>®</sup> 96 Well Adhesion Protein Array	96 samples

### Accessories:

Catalog#	Description	Size
3400-010-01	Cultrex <sup>®</sup> Mouse Laminin I	1 mg
3440-100-01	Cultrex <sup>®</sup> Rat Collagen I	100 mg
3410-010-01	Cultrex <sup>®</sup> Mouse Collagen IV	1 mg
3420-001-01	Cultrex <sup>®</sup> Human Fibronectin, PathClear <sup>®</sup>	1 mg
3416-001-01	Cultrex <sup>®</sup> Bovine Fibronectin, NZHD*	1 mg
3421-001-01	Cultrex <sup>®</sup> Human Vitronectin, PathClear <sup>®</sup>	50 $\mu\text{g}$
3417-001-01	Cultrex <sup>®</sup> Bovine Vitronectin, NZHD	50 $\mu\text{g}$
3438-100-01	Cultrex <sup>®</sup> Poly-L-Lysine	100 ml
3439-100-01	Cultrex <sup>®</sup> Poly-D-Lysine	100 ml
3445-048-01	Cultrex <sup>®</sup> 3-D Culture Matrix <sup>™</sup> BME	15 ml
3446-005-01	Cultrex <sup>®</sup> 3-D Culture Matrix <sup>™</sup> Laminin I	5 ml
3447-020-01	Cultrex <sup>®</sup> 3-D Culture Matrix <sup>™</sup> Collagen I	100 mg
3430-005-01	Cultrex <sup>®</sup> BME with Phenol Red	5 ml
3432-005-01	Cultrex <sup>®</sup> BME without Phenol Red	5 ml
3431-005-01	Cultrex <sup>®</sup> BME with Phenol Red; Reduced Growth Factor	5 ml
3433-005-01	Cultrex <sup>®</sup> BME; no Phenol Red; Reduced Growth Factor	5 ml
3415-001-02	Cultrex <sup>®</sup> Human BME, PathClear <sup>®</sup>	1 ml
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, Reduced Growth Factor PathClear <sup>®</sup>	5 ml
3432-005-02	Cultrex <sup>®</sup> BME, PathClear <sup>®</sup>	5 ml
3433-005-02	Cultrex <sup>®</sup> BME; no Phenol Red; Reduced Growth Factor PathClear <sup>®</sup>	5 ml
3437-100-K	Cultrex <sup>®</sup> Cell Staining Kit	100 ml
3450-048-05	CellSpere <sup>™</sup>	15 ml

\*New Zealand Herd Derived

### Reference:

Kleinman H.K. 1999. The Laminins. Proc. Indian Acad. Sci. (Chem.Sci.), 111:179-184.



**Mouse Laminin I,  
PathClear®**

Catalog #: 3400-010-02

Storage: ≤ -20°C

(Manual Defrost Freezer)

**TREVIGEN®**  
**1-800-873-8443**