

CULTREX® 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 μ 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 μ 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 μ g/cm² of growth surface (0.05 -20 μ 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 μ l
96 wells	100 μ 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.

© 2009 Trevigen, Inc. All Rights Reserved. 3-D Culture Matrix is a trademark, and Trevigen, Cultrex, CultreCoat and PathClear are registered trademarks of Trevigen, Inc. E7/20/09v1

TREVIGEN®

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

Accessories:

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

*New Zealand Herd Derived

Reference:

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



Mouse Laminin I

Cat#: 3400-010-01

Storage: \leq -20 °C

(Manual Defrost Freezer)

1-800-873-8443