

CULTREX[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Mouse Collagen IV

Catalog #: 3410-010-01

Size: 2 x 0.5 mg

Description: Collagen IV is a major component of basement membranes, which underlies epithelial and endothelial cells and surrounds muscle, fat and nerve cells. It can be used as a thin coating on tissue culture surfaces to promote cell attachment and proliferation and to study its effects on cell behavior.

Specifications:

Concentration: 0.5 mg/ml

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

Storage Buffer: 50 mM Hydrochloric Acid, 1.0 mM TCEP

Storage/Stability: Product is stable for 3 months from date of shipment when stored at -80 °C. **For optimal stability, store in liquid nitrogen. Keep Frozen; repeated freeze-thaws will destroy product integrity.**

Purity: >90% by SDS-PAGE.

Materials Qualification:

Functional Assay:

- ◆ Tested for ability to promote attachment and spreading of rat PC-12 pheochromocytoma cells.

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 concentration < 20 EU/ml by LAL assay.

Coating Procedure:

The recommended working concentration is 1-10 µg/cm² (1-10 µg/ml) of growth surface depending on cell type.

- A.** Thaw stock solution on ice for several hours. Position plates on ice and prechill pipette tips. Dilute the Collagen IV stock solution appropriately with cold sterile deionized water. Mix and transfer to the wells of cold tissue culture plates. Spread the solution to completely cover the bottom of the wells.

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

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

Plate Type /Volume	Collagen IV per Well
6 wells (or 35 mm dish)	1 ml
24 wells	200 µl
48 wells	50 µl
96 wells	20 µl

C. Dry down the Collagen IV solution at room temperature for 1 to 2 hours.

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

Related Products:

Catalog#	Description
3465-024-K	Cultrex® 24 well BME Cell Invasion Assay
3465-096-K	Cultrex® 96 well BME Cell Invasion Assay
3456-096-K	Cultrex® Laminin I Cell Invasion Assay
3457-096-K	Cultrex® Collagen I Cell Invasion Assay
3458-096-K	Cultrex® Collagen IV Cell Invasion Assay

Accessories:

Catalog#	Description	Size
3400-010-01	Cultrex® Mouse Laminin I	1 mg
3440-100-01	Cultrex® Rat Collagen I	100 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 µg
3417-001-01	Cultrex® Bovine Vitronectin, NZHD	50 µg
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	5 ml
3446-005-01	Cultrex® 3-D Culture Matrix™ Laminin I	5 ml
3447-020-01	Cultrex® 3-D Culture Matrix™ Collagen I	5 ml
3430-005-01	Cultrex® BME with Phenol Red	5 ml
3431-005-01	Cultrex® BME with Phenol Red, Growth Factor Reduced	5 ml
3432-005-01	Cultrex® BME, no Phenol Red	5 ml
3430-005-02	Cultrex® BME with Phenol Red, PathClear®	5 ml
3431-005-02	Cultrex® BME with Phenol Red, Growth Factor Reduced, PathClear®	5 ml
3432-005-02	Cultrex® BME PathClear®	5 ml
3433-005-02	Cultrex® BME, Growth Factor Reduced, PathClear®	5 ml
3437-100-K	Cultrex® Cell Staining Kit	100 ml
3450-048-05	CellSpense™	15 ml

*New Zealand Herd Derived

Reference:

Hudson B.G., S.T. Reeders , and Tryggvason K. 1993. Type IV collagen: structure, gene organization, and role in human diseases. Molecular basis of Goodpasture and Alport syndromes and diffuse leiomyomatosis. J. Biol. Chem. 268, 26033-36.



Collagen IV

Cat#: 3410-010-01

Storage: ≤ -80 °C

1-800-873-8443