

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Cultrex® Poly-L-Lysine

Catalog #: 3438-100-01
3438-200-01

Size: 100 mL
2 x 100 mL

Description: Poly-L-Lysine, a highly positively charged amino acid chain, is commonly used as a coating agent to promote cell adhesion in culture. This solution is provided ready to use at 0.01% and contains polymers in the 70,000 - 150,000 Da range.

Concentration: 0.01% in phosphate-buffered saline (PBS), sterile-filtered.

Storage Conditions: Product is stable for at least 6 months from the date of receipt when stored at 2 – 8 °C. Keep sterile.

Applications: Substrate for cell culture adhesion. An area of 25 cm² can be coated with 0.5 mL of a 0.1 mg/mL Poly-L-Lysine solution. Optimal conditions for attachment must be determined for each cell line and application. Slides may be dipped in the solution and air dried before applying sample. Keep sterile.

Specifications:*

- Functional Assay: Tested for ability to promote attachment 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.
- Endotoxin concentration ≤ 20 EU/mL by LAL assay.

* Mycoplasma testing: not required for synthetic product.

Coating Procedure:

The recommended working concentration is 0.1 mg/mL (as provided) but may need optimization depending on cell type.

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

<u>Plate Type</u>	<u>Volume Poly-L-Lysine/Well</u>
6 wells (or 35 mm dish)	1 mL
24 wells	200 µL
48 wells	50 µL
96 wells	20 µL

B. Pipette the appropriate amount of Poly-L-Lysine solution in each well. Swirl the plate to ensure coverage. Remove excess reagent and dry wells for 2 hours at room temperature in the biological hood to ensure sterility. **OR**

Pipette the appropriate amount of Poly-L-Lysine solution in each well. Incubate the plate for 1 - 2 hours at 37° C. Remove excess reagent.

C. Rinse the wells twice with cold sterile water, PBS, or cell culture medium. Add cells.

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
3439-100-01	Cultrex® Poly-D-Lysine	100 ml
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
3400-010-01	Cultrex® Mouse Laminin I	1 mg
3440-100-01	Cultrex® Rat Collagen I	1 ml
3442-050-01	Cultrex® Bovine Collagen I	50 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 µg
3417-001-01	Cultrex® Bovine Vitronectin, NZHD	50 µg
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



Cultrex® Poly-L-Lysine

Catalog #: 3438-100-01
3438-200-01

Storage: 2-8 °C

TREVIGEN®
1-800-873-8443