

# TREVIGEN® Product Data

*For Research Use Only. Not For Use In Diagnostic Procedures*

## Cultrex® Poly-L-Ornithine Solution

**Catalog #:** 3436-100-01

**Size:** 100 mL

**Description:** Poly-L-Ornithine, a highly positively charged amino acid chain, is generally used as a coating reagent to promote cell adhesion in culture. Poly-L-Ornithine along or in combination with Fibronectin or Laminin-I is frequently used to enhance attachment and differentiation of various types of neuronal and neural stem cells.

This product is provided ready-to-use, as 0.01% solution in sterile water, and contains polymers in the 30,000 - 70,000 Da range.

**Concentration:** 0.01% w/v (0.1 mg/ml) in sterile water.

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

### Specifications:\*

- Functional Assay: Tested for ability to promote attachment of rat PC-12 cells.
- Sterility Testing: No bacterial or fungal growth detected after incubation at 37°C for 14 days.
- Endotoxin concentration  $\leq$  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 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.

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

<u>Plate Type</u>	<u>Volume Poly-L-Ornithine/Well</u>
10 cm dish (or 75 cm <sup>2</sup> flask)	10 mL
6 cm dish (or 25 cm <sup>2</sup> flask)	3-5 mL
6 wells (or 35 mm dish)	1 mL
12 well	500 $\mu$ L

- B. Pipette the appropriate amount of Poly-L-Ornithine solution in each well. Swirl the plate to ensure that it is fully coated.
- C. Incubate the plate for 1-2 hours at 37°C.
- D. Remove excess reagent. Rinse the wells three times with cell culture medium.
- E. Plate cells on Poly-L-Ornithine-coated dish.

# 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](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

**Related Products:**

Catalog#	Description	Size
3400-010-01	Cultrex® Mouse Laminin I	1 mg
3400-010-02	Cultrex® Mouse Laminin I, PathClear®	1 mg
3401-010-02	Cultrex® Antibiotic-Free Mouse Laminin I, PathClear®	1 mg
3446-005-01	Cultrex® 3-D Culture Matrix™ Laminin I	5 ml
3400-010-03	Cultrex® Stem Cell Qualified Laminin I, PathClear®	1 mg
3434-005-02	Cultrex® Stem Cell Qualified RGF BME, PathClear®	5 ml
3420-001-03	Cultrex® Stem Cell Qualified Human Fibronectin, PathClear®	1 mg
3421-001-03	Cultrex® Stem Cell Qualified Human Vitronectin, PathClear®	200 µg
3432-005-01	Cultrex® Basement Membrane Extract, PathClear®	5 ml
3433-005-01	Cultrex® Reduced Growth Factor BME, PathClear®	5 ml
3532-005-02	Cultrex® Basement Membrane Extract, Type 2, PathClear®	5 ml
3533-005-02	Cultrex® Reduced Growth Factor BME, Type 2, PathClear®	5 ml
3632-005-02	Cultrex® Basement Membrane Extract, Type 3, PathClear®	5 ml
3445-005-01	Cultrex® 3-D Culture Matrix™ BME, PathClear®	5 ml
3447-020-01	Cultrex® 3-D Culture Matrix™ Collagen I	100 mg
3438-100-01	Cultrex® Poly-L-Lysine	100 ml
3439-100-01	Cultrex® Poly-D-Lysine	100 ml



**Cultrex® Poly-L-Ornithine**

Catalog# 3436-100-01

Storage: 2-8°C

**TREVIGEN®**  
1-800-873-8443