

TREVIGEN® Product Data

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

Cultrex™ Poly-L-Lysine

Catalog #: 3438-100-01

Size: 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 kDa range.

Concentration: 0.01% in Phosphate Buffered Saline, sterile-filtered.

Storage Conditions: Store at 4°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.
- No mycoplasma contamination detected by PCR.
- Endotoxin concentration ≤ 20 EU/ml by LAL assay.

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 per 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 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 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 then add your cells.



TREVIGEN®

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-216-2801 • e-mail: info@trevigen.com • www.trevigen.com