

**PRODUCT SPECIFICATION SHEET****POLY-D-LYSINE HYDROBROMIDE, HIGH MOLECULAR WEIGHT**

Poly-D-Lysine is a synthetic molecule used as a thin coating to enhance the attachment of cells to plastic and glass surfaces. It has been used to culture a wide variety of cell types including human amniotic fluid cells and chick embryo cells.

CATALOG NUMBER: 354210 LOT NUMBER: \_\_\_\_\_

SOURCE: Synthetic

QUANTITY & PHYSICAL FORM: 20 milligrams per vial, lyophilized from aqueous solution.

RECONSTITUTION AND USE: Equilibrate vial to room temperature. Resuspend in sterile distilled water and swirl to dissolve. If entire amount of material is not to be used immediately, transfer aliquots to sterile plastic tubes and store at -20°C. It is recommended that solubilized product is used within 3 months. **DO NOT STORE IN FROST-FREE FREEZER. AVOID REPEATED FREEZE THAWS.**

Poly-D-Lysine is generally used in the concentration range of 2.5 to 5.0 micrograms per square cm of growth surface.

Please see attached for coating directions.

MOLECULAR WEIGHT: 500,000 - 550,000 daltons

PURITY: \_\_\_\_\_

QUALITY CONTROL: Poly-D-Lysine is lyophilized from a membrane filtered (0.2 micron) solution. Tested and found negative for the presence of bacteria, fungi and mycoplasma.

STABILITY: Stable for a minimum of 3 months from day of shipment when stored at 2-8°C.

REFERENCES:

1. McKeehan, W.L., et.al., In Vitro, **13**:399 (1977).
2. McKeehan, W.L., Methods for Preparation of Media, Supplements, and  
150 Fifth Avenue, NY,NY 10011, p 209-213 (1984).

**BD Biosciences**  
Clontech  
Discovery Labware  
Immunocytometry Systems  
Pharmingen

BD Biosciences Discovery Labware

Two Oak Park, Bedford, MA 01730, tel: 800.343.2035, fax: 800.743.6200, www.bdbiosciences.com



Use these recommendations as guidelines to determine the optimal coating conditions for your culture system.

#### Coating procedure

1. Dilute poly-d-lysine to desired concentration using serum-free culture medium. The final solution should be sufficiently dilute to evenly coat the growth surface.  
  
e.g.:           If the final coating concentration will be 5 ug/ cm<sup>2</sup>, dilute the material to 50 ug/ml and add 1 ml/35 mm dish, 3 ml/60 mm dish, etc.
2. Add appropriate amount of diluted poly-d-lysine to culture surface.
3. Incubate at room temperature for 1 hour.
4. Aspirate remaining material.
5. Rinse plates carefully - avoid scraping bottom surface. Unbound poly-d-lysine may be inhibitory to cell growth.
6. Plates are ready for use. They may also be stored at 4°C damp or air dried if sterility is maintained.

**FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.**