

**GELATIN**

**Catalog # 07903**

**500 mL per bottle**

**PRODUCT DESCRIPTION:**

This product contains 0.1% gelatin in water

This product is recommended for the coating of culture dishes or flasks on which embryonic stem (ES) cells or primary embryonic fibroblasts are to be grown. This product has been pre-screened for optimal performance when used with other reagents of the ES-Cult™ product line.

**STABILITY / STORAGE:**

Store at room temperature.

Stable for at least 2 years.

Contents guaranteed sterile if seal is not tampered with.

**DIRECTIONS FOR USE:**

To coat culture vessels, aseptically dispense sufficient gelatin into the dish or flask to completely cover the bottom of the vessel. Recommended volumes are:

3 mL per 25cm<sup>2</sup> flask or 60 mm tissue culture dish

7-8 mL per 100 mm tissue culture dish.

Leave gelatin solution in contact with the cultureware for at least 20 minutes at room temperature. Aspirate off all of the excess gelatin solution and leave cultureware sitting open in the biosafety cabinet until dry. These gelatinized dishes or flasks can be stored for two weeks at room temperature or at 4°C until required.

For further information, refer to Technical Manual titled “*In Vitro* Hematopoietic Differentiation of Embryonic Stem (ES) Cells”.

**THIS REAGENT IS FOR LABORATORY USE ONLY.  
IT IS NOT TO BE ADMINISTERED TO HUMANS.**