Cascade Biologics™

invitrogen cell culture



Coating Matrix Kit

Animal Origin-Free

Cat. no. R-011-K

Product Description

Coating Matrix Kit is an animal origin-free kit containing materials suitable for coating up to 750 cm² of tissue culture plasticware. Coating Matrix Kit contains 0.5 ml sterile recombinant human Type-1 collagen as a Coating Matrix (Item # 50-9700) and 50 ml Dilution Medium (Item # 50-9701).

Intended Use

Coating Matrix Kit enhances the growth of human keratinocytes *in vitro*, and is intended for use in conjunction with EpiLife[®] Medium (cat. no. M-EPI-500-CA, M-EPICF-500, or M-EPICF/PRF-500) supplemented with either Supplement S7 (cat. no. S-017-5) or EDGS (cat. no. S-012-5) for the routine culture of normal human keratinocytes. Coating Matrix Kit may also be used for the primary isolation of keratinocytes in an animal-origin-free environment.

This product is for research use only. Not intended for human or animal therapeutic or diagnostic use.

Caution: If handled improperly, some components of this product may present a health hazard. Take appropriate precautions when handling this product, including the wearing of protective clothing and eyewear. Dispose of properly.

Storage and Stability

Coating Matrix Kit is stored at 2° to 8° C at our facility and is shipped at ambient temperature. When stored at 2° to 8° C, the product is stable until the expiration date on the label. Do not freeze.

Instructions for Coating Flasks

- Using sterile technique in a laminar flow culture hood, add Dilution Medium (50-9701) to each flask (5 ml per each 75 cm² flask, or 1.7 ml per each 25 cm² flask).
- 2. Add Coating Matrix (50-9700) directly to the Dilution Medium in each flask (50 μ l per each 75 cm² flask, or 17 μ l per each 25 cm² flask). Rock back and forth vigorously to ensure uniform distribution of the coating matrix over the surface of the flask.
- 3. Cap the flasks and incubate for 30 minutes at room temperature.
- Remove excess Coating Matrix/Dilution Medium from each flask. The flasks may be used immediately, or may be stored at 2° to 8° C for short periods.

EpiLife[®] is a Registered Trademark of Life Technologies.

Limited Use Label License No. 5: Invitrogen Technology

The purchase of this product conveys to the buyer the non-trnsferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies Corporation which cover this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Life Technologies is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500. Email: outlicensing@invitrogen.com.

©2009 Life Technologies Corporation. All rights reserved.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.

For research use only.