

# EpiLife® Basal Medium

Cat. no. UM-EPI-500

500 ml

## Product Description

EpiLife® Basal Medium is a sterile-filtered, basal liquid tissue culture medium that can be used as one component in a complete culture environment for the growth of normal human epidermal keratinocytes. EpiLife® Basal Medium is animal origin free (AOF). EpiLife® Basal Medium contains essential and non-essential amino acids, vitamins, other organic compounds, trace minerals, and inorganic salts. EpiLife® Basal Medium does not contain antibiotics, antimycotics, hormones, growth factors, or proteins. The medium is HEPES and bicarbonate buffered and is designed for use in an incubator with an atmosphere of 5% CO<sub>2</sub>/95% air.

## Intended Use

For Further Manufacturing Use.

## Storage and Stability: EpiLife® Basal Medium

Store at 2–8°C. EpiLife® Basal Medium is light-sensitive. Protect from light. Do not expose to light for lengthy periods of time. When stored in the dark at 2–8°C, the medium is stable until the expiration date shown on the label.

## Preparation of Supplemented Medium

To supplement one bottle of EpiLife® Basal Medium with the appropriate growth supplement, the following protocol is recommended:

1. Thaw the growth supplement.
2. Gently swirl the bottle of supplement. Avoid splashing the supplement into the cap of the bottle or causing the supplement to foam.
3. Obtain one bottle of EpiLife® Basal Medium.
4. Wipe the outside of the bottles with a sanitizing solution such as 70% ethanol or 70% isopropanol and move the bottles to a laminar flow hood.
5. Add 5 mL of 100X growth supplement to the bottle of medium using sterile technique.
6. Tightly cap the bottle of supplemented medium and swirl the contents to ensure a homogeneous solution. Avoid causing the medium to foam.

## Storage and Stability: Supplemented Medium

Once EpiLife® Basal Medium has been supplemented, the supplemented medium should be stored at 2–8°C. Do not freeze. When stored at 2–8°C, the supplemented medium is stable for 1 month.

*EpiLife® is a Registered Trademark of Life Technologies.*

**For research use only.**

Life Technologies Corporation • 5791 Van Allen Way • Carlsbad • CA 92008 • Tel: 800.955.6288 • [www.invitrogen.com](http://www.invitrogen.com) • E-mail: [tech\\_support@invitrogen.com](mailto:tech_support@invitrogen.com)

**Limited Use Label License No. 5: Invitrogen Technology**

The purchase of this product conveys to the buyer the non-transferable 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](mailto: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.***

Life Technologies Corporation • 5791 Van Allen Way • Carlsbad • CA 92008 • Tel: 800.955.6288 • [www.invitrogen.com](http://www.invitrogen.com) • E-mail: [tech\\_support@invitrogen.com](mailto:tech_support@invitrogen.com)