# **Cascade Biologics**™

invitrogen cell culture



## Medium 254

Cat. no. M-254-500

500 ml

## **Product Description**

Medium 254 is a sterile, liquid tissue culture medium prepared with 200 µM calcium chloride for the culture of normal human epidermal melanocytes. This medium is intended for use as one component in a complete Medium 254 is a basal medium culture environment. containing essential and non-essential amino acids, vitamins, other organic compounds, trace minerals, and This medium does not contain inorganic salts. antibiotics, antimycotics, hormones, growth factors, or proteins. This 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. To support plating and long-term proliferation of normal human melanocytes, Medium 254 must be supplemented with Human Melanocyte Growth Supplement (HMGS, cat. no. S-002-Human PMA-Free Melanocyte Growth Supplement-2 (HMGS-2, cat. no. S-016-5). Each of these supplements contains all of the growth factors, hormones, and tissue extracts necessary for growth of melanocytes in Medium 254.

#### **Intended Use**

Medium 254 is intended for use in the routine culture of normal human epidermal melanocytes. When supplemented with HMGS or HMGS-2, Medium 254 will support the plating and proliferation of melanocytes at varying culture densities from 5 x  $10^3$  cells/cm<sup>2</sup> to 1 x  $10^5$  cells/cm<sup>2</sup>. Additional applications for use may include primary isolation of melanocytes from skin.

This product is for research use only. Not for use in animals, humans, or diagnostic procedures.

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

Medium 254 is stored at 4° C in our facility and is shipped at ambient temperature. Upon receipt, the medium should be stored at 4° C and should <u>not</u> be frozen. **Protect from light.** Several components of this tissue culture medium are light-labile, and we recommend that the medium not be exposed to light for lengthy periods of time. If the medium is warmed prior to use, do not exceed 37° C. When stored in the dark at 4° C, the product is stable until the expiration date on the label.

Follow the supplementation instructions below to prepare the medium for use.

#### **Preparation of Supplemented Medium 254**

**Note**: For information on HMGS or HMGS-2, please refer to the product sheets that accompany those products.

- Thaw one bottle of HMGS or one bottle of HMGS-2 according to the instructions provided with those products. Make sure that the cap of the bottle is tight. Gently swirl the bottle of supplement. Avoid splashing the supplement into the cap of the bottle or causing the supplement to foam.
- Wipe the outside of the containers with a disinfecting solution such as 70% ethanol or isopropanol.
- To add the HMGS or HMGS-2, transfer the entire contents of the bottle of supplement to the bottle of medium using sterile technique in a laminar flow culture hood.
- 4. 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 of Supplemented Medium 254

Once Medium 254 has been supplemented with HMGS or HMGS-2, the supplemented medium should be stored in the dark at 4° C and should <u>not</u> be frozen. When stored in the dark at 4° C, the supplemented medium is stable for 1 month.

## Cascade Biologics™

invitrogen cell culture



#### 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. ©2009 Life Technologies Corporation. All rights reserved.

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