

# Medium 154

Catalog Number: M-154-500  
500 ml

## Instructions for storage and use

### Product Description

Medium 154 is a sterile, liquid tissue culture medium prepared with 200  $\mu$ M calcium chloride for the growth of normal human epidermal keratinocytes. This medium is intended for use as one component in a complete culture environment. Medium 154 is a basal medium containing essential and non-essential amino acids, vitamins, other organic compounds, trace minerals, and inorganic salts. This medium does not contain 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 keratinocytes, Medium 154 must be supplemented with Human Keratinocyte Growth Supplement (HKGS, Cat. S-001-5) or Human Keratinocyte Growth Supplement Kit (HKGS Kit, Cat. S-001-K).

### Intended Use

Medium 154 is intended for use in the routine culture of normal human epidermal keratinocytes. When supplemented with HKGS or HKGS Kit, Medium 154 will support the plating and proliferation of keratinocytes at varying culture densities from clonal (25 cells/cm<sup>2</sup>) to high density (8 x 10<sup>4</sup> cells/cm<sup>2</sup>). Additional applications for use may include primary isolation of keratinocytes from skin, and clonal growth assays. For optimal results when performing primary isolations, Medium 154 should be used in conjunction with Coating Matrix Kit (Cat. R-011-K).

***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 154 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 not 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.

Please use the supplementation instructions (page 2) to prepare the medium for use.

## Preparation of Supplemented Medium 154

**Note:** For information on HKGS or HKGS Kit, please refer to the product sheet that accompanies those products.

1. Thaw the frozen components of the HKGS Kit, or one bottle of HKGS according to the instructions provided with those products. Make sure that the caps of all of the bottles are tight. Gently swirl the bottle(s) of supplement. Avoid splashing the supplement into the cap of the bottle or causing the supplement to foam.
2. Wipe the outside of the containers with a disinfecting solution such as 70% ethanol or isopropanol.
3. To add the HKGS, transfer the entire contents of the bottle of supplement to the bottle of medium using sterile technique in a laminar flow culture hood. To add the HKGS Kit, transfer the desired amount of each component of the HKGS Kit to the bottle of medium using sterile technique in a laminar flow culture hood. **Note:** addition of less than the entire amount of any component may affect the performance of the supplemented medium. If antibiotics/antimycotics are desired, add the antibiotic/antimycotic solution included in HKGS Kit using the same technique as above.
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 154

Once Medium 154 has been supplemented with HKGS or HKGS Kit, the supplemented medium should be stored in the dark at 4° C and should not be frozen. When stored in the dark at 4° C, the supplemented medium is stable for 1 month.

## Selected References

The Medium 154 formulation is based on medium MCDB 151, with trace elements as in medium MCDB 104, and the high amino acid modifications of Pittelkow.

Cook, Pittelkow, and Shipley; J. Cell. Physiol. 146:277-289, 1991

Peehl and Ham; In Vitro 16: 526-540, 1980

McKeehan et al.; In Vitro 13: 399-416, 1977

Pittelkow and Scott; Mayo Clin. Proc. 61: 771-777, 1986

## Terms and Conditions of Sale

Cascade Biologics, Inc. (hereinafter, CBI) warrants that its products will perform according to the information provided in various publications that it distributes, and as described herein for the intended shelf life of the product when stored under the conditions prescribed by CBI. If any product does not perform according to the published information provided by us, CBI will replace the product free of charge to the original customer. The remedy of product replacement shall be the customers' sole and exclusive remedy for defective product, unless CBI is unable to deliver replacement product, in which case CBI shall reimburse the customer for the purchase price of the defective product. Customer understands that the foregoing limited warranty is in lieu of all other warranties and CBI hereby disclaims all other warranties including, but not limited to, implied warranty of merchantability or fitness for adequacy for any particular purpose or use. CBI shall not be liable to customer or to any party claiming through customer for any incidental or consequential damages, including, but not limited to any lost profits, lost savings, or lost business, whether arising out of contract, tort, or otherwise. Customer acknowledges that products are intended for research use only and are not to be used in animals, humans, or diagnostic procedures.