

EDGS

EpiLife® Defined Growth Supplement

Cat. no. S-012-5

5 ml

Product Description

EpiLife® Defined Growth Supplement (EDGS) is a sterile, concentrated (100X) solution intended for use as one component in a complete culture environment for the growth of normal human epidermal keratinocytes. EDGS is an ionically-balanced supplement containing purified bovine serum albumin (BSA), purified bovine transferrin, hydrocortisone, recombinant human insulin-like growth factor type-1 (rhIGF-1), prostaglandin E₂ (PGE₂) and recombinant human epidermal growth factor (rhEGF). The bovine products in EDGS have been isolated from animals of North American or Australian origin. The isolation process for these components includes steps that inactivate viruses. All components of EDGS are greater than 95% pure. Each 5 ml bottle of EDGS is the correct amount of supplement for a 500 ml bottle of EpiLife® basal medium. When a 500 ml bottle of medium is supplemented with EDGS the final concentrations of the components in the supplemented medium are: BSA, 30 µg/ml; bovine transferrin, 5 µg/ml; hydrocortisone, 11 ng/ml; rhIGF-1, 10 ng/ml; PGE₂, 18 ng/ml; and rhEGF, 1 ng/ml.

Intended Use

EDGS is intended for use in conjunction with EpiLife® basal medium for the culture of normal human epidermal keratinocytes in a serum-free, chemically-defined environment. For optimal cell growth, EDGS should be used in conjunction with Coating Matrix Kit (Cat. R-011-K) and Defined Trypsin Inhibitor (DTI, Cat. R-007-100). **This product is for research use only. Not for use in animals, humans, or diagnostic procedures.**

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.

For research use only.

Life Technologies Corporation • 5791 Van Allen Way • Carlsbad • CA 92008 • Tel: 800.955.6288 • www.invitrogen.com • E-mail: tech_support@invitrogen.com

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

EDGS is stored at -20° C at our facility and is shipped on dry ice. Upon receipt, the product should be stored at -20° C in a freezer that is not self-defrosting. When stored at -20° C the product is stable until the expiration date shown on the label.

After long-term storage at -20° C, HCGS may contain a small amount of precipitate. This precipitate is formed from cold-insoluble material in the BPE component of the HCGS and will not affect the performance of the product.

Thawing

To thaw, place the product in a 37° C water bath or overnight at 4° C. If thawed in a water bath, do not leave the product at 37° C after the product has thawed. For instructions on adding EDGS to EpiLife® basal medium, refer to the instructions that accompany the basal medium.

EpiLife® is a Registered Trademark of Life Technologies.