

# Human Corneal Growth Supplement (HCGS)

Catalog Numbers S0095

Pub. No. MAN0001613 Rev. 3.0



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

#### **Product description**

Human Corneal Growth Supplement (HCGS) is a sterile, concentrated (100X) solution intended for use as one component in a complete culture environment for the growth of normal corneal epithelial cells. HCGS is an ionically-balanced supplement containing bovine pituitary extract (BPE), recombinant human insulin-like growth factor-I, hydrocortisone, bovine transferrin, and mouse epidermal growth factor. Each 5 mL bottle of HCGS is the correct amount of supplement for a 500 mL bottle of EpiLife<sup>™</sup> Medium.

HCGS is designed for use with EpiLife<sup>™</sup> Medium for the routine serum-free culture of human corneal epithelial cells.

### Contents and storage

Catalog numbers that appear as links open the web pages for those products.

Product	Cat. No.	Amount	Storage
Human Corneal Growth Supplement (HCGS)	S0095	5 mL	-20°C; Store in a manual defrost freezer <sup>[1,2]</sup>

<sup>[1]</sup> When stored at -20°C, the product is stable until the expiration date on the label.

## Thaw the Human Corneal Growth Supplement

Place the HCGS in a 37°C water bath or overnight at 4°C.

Note: If thawed in a water bath, do not leave the product at 37°C after the product has thawed.

For instructions on adding HCGS to EpiLife<sup>™</sup> Medium, see the instructions that accompany the basal medium.

### Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at <a href="https://www.thermofisher.com/us/en/home/global/terms-and-conditions.html">www.thermofisher.com/us/en/home/global/terms-and-conditions.html</a>. If you have any questions, please contact Life Technologies at <a href="https://www.thermofisher.com/support">www.thermofisher.com/support</a>.



Life Technologies Corporation | 3175 Staley Road | Grand Island, New York 14072 USA

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0001613

Revision	Date	Description	
3.0	17 November 2021	Replaced "bovine insulin" with "recombinant human insulin-like growth factor-I" in "Product description" on page 1.	
		Updated to the current document template, with associated updates to trademarks, logos, licensing, and warranty.	

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.



<sup>[2]</sup> 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.