

PMA-Free Human Melanocyte Growth Supplement (HMGS-2)

Catalog Number S0165

Pub. No. MAN0001586 Rev. 3.0

Product description

PMA-Free Human Melanocyte Growth Supplement (HMGS-2) is a sterile, concentrated (100X) solution intended for use as one component in a complete culture environment for the growth of normal human epidermal melanocytes. HMGS-2 is an ionically-balanced supplement containing bovine pituitary extract (BPE), fetal bovine serum, recombinant human insulin-like growth factor-I, bovine transferrin, basic fibroblast growth factor, hydrocortisone, heparin, and endothelin-1. Each 5 mL bottle of HMGS-2 is the correct amount of supplement for a 500 mL bottle of basal Medium 254 (Cat. No. [M254500](#)).

HMGS-2 is designed for use with basal Medium 254 for the routine culture of normal human epidermal melanocytes. Additional applications may include primary isolation of melanocytes from skin.

Contents and storage

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

Product	Cat. No.	Amount	Storage
PMA-Free Human Melanocyte Growth Supplement (HMGS-2)	S0165	5 mL	-20°C; Store in a manual defrost freezer ^[1,2]

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

^[2] After long-term storage at -20°C, HMGS-2 may contain a small amount of precipitate. This precipitate is formed from cold-insoluble material in the BPE component of the HMGS-2 and will not affect the performance of the product.

Thaw the PMA-Free Human Melanocyte Growth Supplement

Place the HMGS-2 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 HMGS-2 to Medium 254, 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 www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

thermofisher.com/support | thermofisher.com/askaquestion



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. MAN0001586

Revision	Date	Description
3.0	14 December 2021	<ul style="list-style-type: none"> Replaced "bovine insulin" with "recombinant human insulin-like growth factor-I". 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.

For Research Use Only. Not for use in diagnostic procedures.