

Keratinocyte Growth Medium 3

Instruction Manual

Product	Size	Catalog Number
Keratinocyte Growth Medium 3 (Ready-to-use) – Includes Basal Medium, SupplementMix and CaCl ₂ -Solution	500 ml	C-20021

Recommended for

- Normal Human Epidermal Keratinocytes (NHEK) GM3, juvenile foreskin
- Normal Human Epidermal Keratinocytes (NHEK) GM3, adult donor

Product Description

Our Keratinocyte Growth Medium 3 is the improved animal extract free version of the BPE-containing Keratinocyte Growth Medium 2. Keratinocyte Growth Medium 3 is optimized for the isolation and expansion of human keratinocytes from juvenile or adult donors. The BPE-free medium can be used without ECM coating of the cell culture vessel and supports the growth of a fast proliferating and homogeneous cell population.

Keratinocyte Growth Medium 3 is available as ready-to-use Medium. The Medium (Ready-to-use) consists of a bottle of Basal Medium, one vial of CaCl₂ stock solution, and one vial of SupplementMix. Adding the supplement solutions to the Basal Medium results in the complete Growth Medium.

Supplementation Details

Our Keratinocyte Growth Medium 3 contains all the growth factors and supplements necessary for the optimal adherence and growth of human keratinocytes. The formulation is confidential. The Keratinocyte Growth Medium 3 does not contain antibiotics or antimycotics and is formulated for use in an incubator with an atmosphere of 5% CO₂.

Note: CaCl₂ is provided as separate stock solution. This allows users to adjust the CaCl₂ concentration individually. To ensure maximal growth performance of normal human keratinocytes, we recommend adding the entire 60 µl CaCl₂ stock solution, which results in final 0.06 mM CaCl₂ concentration.

In case you need a higher CaCl₂ concentration (e.g. for differentiation experiments), an additional CaCl₂-Solution must be added separately.

Storage and Stability

Store the 500 ml Basal Medium at 4–8°C in the dark and the SupplementMix and the CaCl₂ solution at -20°C immediately after arrival. Do not freeze the 500 ml Basal Medium. If stored properly, the products are stable until the expiry date stated on the label. After adding the supplements to the Basal Medium, the shelf life of the complete medium is 6 weeks at 4–8°C. Do not freeze the complete medium.

For use, pre-warm only an aliquot of the complete medium and keep the remaining medium refrigerated at 4–8 °C.

Note: The Keratinocyte Growth Medium 3 contains light sensitive components. Avoid longer exposition times to light e.g. store in a cabinet when pre-warming the medium.

Initiation of Keratinocyte Culture and Subcultivation Process

For detailed information about keratinocyte seeding density, detachment

and subcultivation protocol please follow our Instruction Manual *Keratinocytes in Keratinocyte Growth Medium 3* and our Application Note *Isolation of Fast Proliferating and Homogeneous Keratinocyte Populations with PromoCell's Improved Keratinocyte Growth Medium 3*. Both documents are available for download on our website www.promocell.com.

Quality Control

All lots of PromoCell Keratinocyte Growth Medium 3 are subjected to comprehensive quality control tests using primary human keratinocytes. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested keratinocytes. Approved in-house lots of media are used as a reference.

In addition, all lots of media have been tested for the absence of microbial contaminants (fungi, bacteria, mycoplasma).

Intended Use

The products are for in vitro use only and not for diagnostic or therapeutic procedures. For safety precautions please see the appropriate MSDS.

Note: Due to their low serum content or the absence of serum, PromoCell media are not suitable for trypsin neutralization (e.g. when splitting the cells.) Instead we recommend using our DetachKit (C-41200, C-41210, C-41220), which contains HEPES BSS, Trypsin/EDTA and Trypsin Neutralization Solution.

If you require special media modifications, we offer a custom media service starting at 10 bottles per order.
Contact us at info@promocell.com to find out more.

PromoCell GmbH
Sickingenstr. 63/65
69126 Heidelberg
Germany

info@promocell.com
www.promocell.com

USA/Canada

Phone: 1 - 866 - 251 - 2860 (toll free)
Fax: 1 - 866 - 827 - 9219 (toll free)

Deutschland

Telefon: 0800 - 776 66 23 (gebührenfrei)
Fax: 0800 - 100 83 06 (gebührenfrei)

France

Téléphone: 0800 - 90 93 32 (ligne verte)
Téléfax: 0800 - 90 27 36 (ligne verte)

United Kingdom

Phone: 0800 96 03 33 (toll free)
Fax: 0800 169 85 54 (toll free)

Other Countries

Phone: +49 6221 - 649 34 0
Fax: +49 6221 - 649 34 40

© PromoCell GmbH