

Melanocyte Growth Medium M3

Instruction manual

Product	Size	Catalog number
Melanocyte Growth Medium M3 (Ready-to-use) – Includes Basal Medium and SupplementMix	500 ml	C-24310
Melanocyte Growth Medium M3, phenol red-free (Ready-to-use) – Includes Basal Medium and SupplementMix	500 ml	C-24311
Melanocyte Growth Medium M3 SupplementMix	for 500 ml	C-39430
Melanocyte Basal Medium M3, with phenol red	500 ml	C-24410

Recommended for

- Normal Human Epidermal Melanocytes (NHEM), juvenile foreskin, cultured in M3 Medium
- Normal Human Epidermal Melanocytes (NHEM), adult donor, cultured in M3 Medium

Product description

Our Melanocyte Growth Media M3 are designed for the isolation and cultivation of human melanocytes from human juvenile foreskin and human adult skin in a BPE- and serum-free environment. The media do not contain PMA (Phorbol Myristate Acetate) which could interfere with experimental settings. When using the Melanocyte Growth Media M3 the melanocytes grow well on plastic and no matrix coating is required. Isolated human juvenile and adult melanocytes show their characteristic dendritic morphology in this advanced cell culture medium. We offer different donor sites from adult skin (e.g., breast, eyelid, abdomen) isolated in Melanocyte Growth Media M3.

Our Melanocyte Growth Medium M3 (C-24310) and Melanocyte Growth Medium M3, phenol red-free (C-24311) are available as media (ready-to-use) which consist of 500 ml bottle of basal medium and one vial of SupplementMix. Adding the SupplementMix to the basal medium results in complete growth medium. Basal medium and SupplementMix of the Melanocyte Growth Medium M3 with phenol red can also be purchased separately.

Supplementation details

The Melanocyte Growth Media M3 are BPE- and serum-free and contain all growth factors and supplements necessary for the optimal growth of human melanocytes. Furthermore, the media do not contain the tumor promoting mitogen Phorbol Myristate Acetate (PMA).

The media do not contain antibiotics or antimycotics and are formulated for use in an incubator with an atmosphere of 5% CO₂.

Preparation of the supplemented medium for use

Thaw the SupplementMix at 15–25°C. Do not incubate longer than necessary. Then, transfer the entire content of the SupplementMix to the basal medium. Close the bottle and swirl gently until a homogenous mixture is formed.

Storage and stability

Store the basal medium at 2–8°C in the dark and the SupplementMix at -20°C immediately after arrival. Do not freeze the 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 2–8°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 2–8°C.

Quality and control

All lots of PromoCell Melanocyte Growth Media M3 are subjected to comprehensive quality control tests using Normal Human Epidermal Melanocytes (NHEM). Each lot is checked for growth promoting activity and for the absence of microbial contaminants (fungi and bacteria).

Intended use

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

Cell culture

Since our Melanocyte Growth Media M3 enable serum-, BPE- and PMA-free cultivation of NHEM without any cell culture plastic coating, we point out that the cells in this media require special culture conditions. To maintain the cells in a robust adherent pro-proliferative phase, we recommend passage of the cells at 70–90% confluency. According to the literature, it is known that high cell densities of NHEM can promote 3D spheroid growth.¹ Therefore, hyperconfluency should be avoided if possible.

Note: Due to the absence of serum, PromoCell Melanocyte Growth Medium M3 and PromoCell Melanocyte Growth Medium M3, phenol red-free 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.

References

1. Lin, S.-J., et al. (2005). „Formation of melanocyte spheroids on chitosan-coated surface." Biomaterials 26(12): 1413–1422

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