

Melanocyte Media M2

Instruction Manual

| Product | Size | Catalog Number |
|---|------------|----------------|
| Melanocyte Growth Medium M2 (Ready-to-use) – Includes Basal Medium and SupplementMix | 500 ml | C-24300 |
| Melanocyte Growth Medium M2 (Ready-to-use), phenol red-free – Includes phenol red-free Basal Medium and SupplementMix | 500 ml | C-24305 |
| Melanocyte Basal Medium M2 | 500 ml | C-24400 |
| Melanocyte Basal Medium M2, phenol red-free | 500 ml | C-24405 |
| Melanocyte Growth Medium M2 SupplementMix | for 500 ml | C-39420 |

Recommended for

- Normal Human Epidermal Melanocytes (NHEM), juvenile foreskin, cultured in M2 Medium
- 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 Medium M2 is a serum-free and PMA (Phorbol Myristate Acetate)-free Medium that was developed for the optimized in vitro cultivation of human melanocytes. It is also suitable for human pigment epithelial cells, neuronal cells, and melanoma cells.

The medium is optimized for primary human cells, but can also be used for equine, porcine, mouse, and rat melanocytes.

Our Melanocyte Growth Medium M2 is available as Medium (Ready-to-use) which consists of a 500 ml bottle of Basal Medium and one vial of SupplementMix. Adding the SupplementMix to the Basal Medium results in the complete Growth Medium.

Basal Medium (with or without phenol red) and the SupplementMix can also be purchased separately.

Supplementation Details

Our Melanocyte Growth Medium M2 contains all the growth factors and supplements necessary for the optimal growth of human melanocytes. Further, the Medium lacks the tumor promoting mitogen Phorbol Myristate Acetate (PMA), which may interfere with some experimental approaches. Melanocyte Growth Medium M2 does not contain antibiotics or antimycotics and is 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. Aseptically mix the SupplementMix by carefully pipetting up and down. 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 4–8°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 SupplementMix 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.

Quality Control

All lots of PromoCell Melanocyte Growth Medium M2 are subjected to comprehensive quality control tests using primary human melanocytes. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested melanocytes. 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 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 Neutralizing 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