

Instruction Manual

| Product | Size | Catalog Number |
|---|------------|----------------|
| Melanocyte Growth Medium (Ready-to-use) – Includes Basal Medium and SupplementMix | 500 ml | C-24010 |
| Melanocyte Growth Medium Kit – Includes Basal Medium and SupplementPack | 500 ml | C-24110 |
| Melanocyte Basal Medium | 500 ml | C-24210 |
| Melanocyte Basal Medium, phenol red-free | 500 ml | C-24215 |
| Melanocyte Growth Medium SupplementMix | for 500 ml | C-39415 |
| Melanocyte Growth Medium SupplementPack | for 500 ml | C-39410 |

Recommended for

- Normal Human Epidermal Melanocytes (NHEM), juvenile foreskin

Product Description

Our Melanocyte Growth Medium is a PMA (Phorbol Myristate Acetate)-containing Medium and has been developed for the *in vitro* cultivation of juvenile melanocytes. Since PMA is a tumor promoting mitogen, it may interfere with some experimental approaches. In these cases, we recommend using the serum-free and PMA-free Melanocyte Growth Medium M3 (C-24310). The Melanocyte Growth Medium is optimized for primary human cells, but can also be used for equine, porcine, murine, and rat melanocytes. The Melanocyte Growth Medium is available as Medium (Ready-to-use) or as Medium Kit.

The Medium (Ready-to-use) consists of a 500 ml bottle of Basal Medium and one vial of SupplementMix. The Medium Kit consists of a 500 ml bottle of Basal Medium and the SupplementPack (a set of individual vials with pre-measured supplements) allowing the user full control over the media formulation. Adding the SupplementMix or the SupplementPack to the Basal Medium results in the complete Growth Medium. Basal Medium (with or without phenol red) as well as the SupplementMix and SupplementPack can also be purchased separately.

Supplementation Details

Our Melanocyte Growth Medium contains all

the growth factors and supplements necessary for the optimal growth of human melanocytes, e.g. Bovine Pituitary Extract, basic Fibroblast Growth Factor, and Insulin (for details see table on page 2). While the supplements are serum-free, the Basal Medium is manufactured with 0.2% serum. The Medium 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 or SupplementPack at 15–25°C. Aseptically mix the supplement solutions by carefully pipetting up and down. Then, transfer the entire content of each supplement to the Basal Medium. Close the bottle and swirl gently until a homogenous mixture is formed.

Note: Light flocculation may be seen upon thawing the supplements containing BPE. This does not affect the activity. Optionally, the precipitate can be removed by centrifugation under sterile conditions.

Storage and Stability

Store the Basal Medium at 4–8°C in the dark and the SupplementMix or SupplementPack 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 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 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.

| Final supplement concentrations (after addition to the medium) | Melanocyte Growth Medium |
|--|--------------------------|
| Bovine Pituitary Extract | 0.004 ml/ml |
| Basic Fibroblast Growth Factor (recombinant human) | 1 ng/ml |
| Insulin (recombinant human) | 5 µg/ml |
| Hydrocortisone | 0.5 µg/ml |
| Phorbol Myristate Acetate | 10 ng/ml |

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

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

info@promocell.com
www.promocell.com

© PromoCell GmbH

07/2022