

Instruction Manual

| Product | Size | Catalog Number |
|--|-----------------------------|----------------|
| Normal Human Epidermal Melanocytes (NHEM) juvenile foreskin | 500,000 cryopreserved cells | C-12400 |
| | 500,000 proliferating cells | C-12450 |
| Normal Human Epidermal Melanocytes (NHEM) juvenile foreskin, cultured in M2 Medium | 500,000 cryopreserved cells | C-12402 |
| | 500,000 proliferating cells | C-12452 |
| Normal Human Epidermal Melanocytes (NHEM) adult donor, cultured in M2 Medium | 500,000 cryopreserved cells | C-12403 |
| | 500,000 proliferating cells | C-12453 |

Product Description

Epidermal melanocytes represent 5-10% of the cells in the epidermis. Located in the *stratum basale* with extensions to the suprabasal layers, they are specialized in the production of melanin. This protein is responsible for the pigmentation of the skin, hair and eyes and exerts a protective effect on adjacent cells by shielding them from the harmful effects of UV radiation.

PromoCell offers a range of Normal Human Epidermal Melanocytes (NHEM) produced at PromoCell's cell culture facility from donors with lightly, moderately and darkly pigmented skin (detailed pigmentation-related donor information available on request).

The cells are isolated from juvenile foreskin or adult skin from different locations, e.g. face, breast, abdomen and thighs, using either the serum free, PMA (Phorbol Myristate Acetate) containing Melanocyte Growth Medium or the unique serum-free and PMA-free Melanocyte Growth Medium M2. Since PMA is a tumor-promoting mitogen, which can interfere with experimental approaches, we recommend using Melanocytes isolated in Melanocyte

Growth Medium M2 and also using this medium for cultivation.

Shortly after isolation, all PromoCell Normal Human Epidermal Melanocytes are cryopreserved at passage 2 (P2) using PromoCell's proprietary, serum-free freezing medium, Cryo-SFM. Each cryo vial contains more than 500,000 viable cells after thawing.

Proliferating cell cultures are made from 500,000 cryopreserved cells that have been thawed and cultured for three days at PromoCell.

Quality Control

Rigid quality control tests are performed for each lot of PromoCell Normal Human Epidermal Melanocytes.

They are tested for cell morphology, adherence rate and cell viability. Furthermore, immunohistochemical tests for the cell-type specific marker, Mel-5, are carried out for each lot (see page 5). Growth performance is tested through multiple passages up to 15 population doublings (PD) under culture conditions without antibiotics and antimycotics.

In addition, all cells have been tested for the absence of HIV-1, HIV-2, HBV, HCV, HTLV-1, HTLV-2 and microbial

contaminants (fungi, bacteria, and mycoplasma).

A detailed certificate of analysis (CoA) for each lot can be downloaded at: www.promocell.com/coa

Intended Use

PromoCell Normal Human Epidermal Melanocytes are for *in vitro* research use only and not for diagnostic or therapeutic procedures.

Warning

Although tested negative for HIV-1, HIV-2, HBV, HCV, HTLV-1 and HTLV-2, the cells – like all products of human origin – should be handled as potentially infectious. No test procedure can completely guarantee the absence of infectious agents.

Follow appropriate safety precautions!

After delivery, start immediately with the protocol for cryopreserved cells (see page 2) or the protocol for proliferating cells (see page 3).

*Start immediately after delivery.
Use aseptic techniques and a laminar flow bench.*

Protocol for Cryopreserved Cells

Straight after arrival, store the cryopreserved cells in liquid nitrogen, or seed them immediately.

Note: Storage at -80°C is not sufficient for cell preservation and causes irreversible cell damage.

1. Prepare the medium

Calculate the needed culture surface area according to the plating density (see page 5) and the lot-specific cell numbers stated on the certificate of analysis. Fill the appropriate volume of PromoCell Growth Medium (at least 9 ml per vial of cells) in cell culture vessels. Place the vessels in an incubator (37°C , 5% CO_2) for 30 minutes.



2. Thaw the cells

Remove the cryovial from the liquid nitrogen container and immediately place it on dry ice – even for short transportation. Under a laminar flow bench, briefly twist the cap a quarter turn to relieve pressure, then re-tighten. Immerse the vial into a water bath (37°C) just up to the screw cap for 2 minutes. Ensure that no water enters the thread of the screw cap.



3. Disinfect the vial and seed the cells

Thoroughly rinse the cryovial with 70% ethanol under a laminar flow bench. Then, aspirate the excess ethanol from the thread area of the screw cap. Open the vial and transfer the cells to a cell culture vessel containing the prewarmed medium from step 1.



4. Incubate the cells

Place the vessel in an incubator (37°C , 5% CO_2) for cell attachment. Replace the medium after 16-24 hours. The cells should be subcultured according to the subcultivation protocol (see page 4) once they have reached 70–90% confluency.



*Start immediately after delivery.
Use aseptic techniques and a laminar flow bench.*

Protocol for Proliferating Cells

1. Incubate the cells

Unpack the culture vessel, do not open the lid, and immediately place it in an incubator (37°C, 5% CO₂) for 3 hours to allow the cells to recover from the transportation.



2. Replace the transport medium

Carefully open the vessel, rinse the inner side of the lid with 70% ethanol, and let air dry. Aspirate the transport medium from the vessel. Add 10 ml of the appropriate PromoCell Cell Growth Medium.



3. Check and incubate the cells

Check the cell density. Open the lid half a turn and place the vessel in an incubator (37°C, 5% CO₂). The cells should be subcultured according to the subcultivation protocol (see page 4) once they have reached 70–90% confluency.



Use aseptic techniques and a laminar flow bench.

Subcultivation Protocol

1. Prepare the reagents and wash the cells

Place the PromoCell DetachKit at room temperature for at least 30 minutes to adjust the temperature of the reagents. Carefully aspirate the medium from the culture vessel. Add 100 μ l Hepes BSS Solution per cm^2 of vessel surface to wash the cells and agitate the vessel carefully for 15 seconds.



2. Detach the cells

Carefully aspirate the Hepes BSS from the culture vessel. Add 100 μ l Trypsin/EDTA Solution per cm^2 of vessel surface. Note: We recommend detaching the cells at room temperature. Close the vessel and examine the cells under a microscope. When the cells start to detach, gently tap the side of the vessel to loosen the remaining cells.



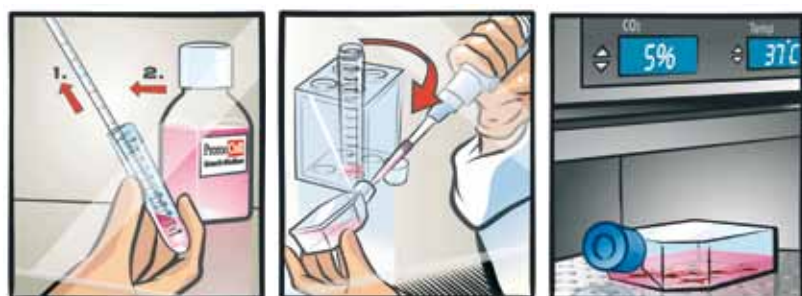
3. Neutralize the trypsin and harvest the cells

Add 100 μ l Trypsin Neutralization Solution per cm^2 of vessel surface and gently agitate. Carefully aspirate the cell suspension and transfer it to a centrifugation tube. Spin down the cells for 3 minutes at 220 x g.



4. Incubate the cells

Discard the supernatant (step 1), add 1 ml of the appropriate PromoCell Cell Growth Medium (step 2), and resuspend the cells by carefully pipetting up and down. Plate the cells according to the recommended seeding density in new cell culture vessels containing PromoCell Cell Growth Medium pre-warmed to 37°C. Place the vessels in an incubator (37°C, 5% CO_2).



Specifications

| Product | Recommended Culture Media* | Plating Density | Passage after Thawing | Marker | Population Doublings |
|---|----------------------------|--|-----------------------|--------------------|----------------------|
| Normal Human Epidermal Melanocytes (NHEM), juvenile foreskin | C-24010 | 5,000 - 10,000 cells per cm ² | P2 | Mel-5 ⁺ | > 15 |
| Normal Human Epidermal Melanocytes (NHEM), juvenile foreskin, cultured in M2 Medium | C-24300 | 5,000 - 10,000 cells per cm ² | P2 | Mel-5 ⁺ | > 15 |
| Normal Human Epidermal Melanocytes (NHEM), adult donor, cultured in M2 Medium | C-24300 | 5,000 - 10,000 cells per cm ² | P2 | Mel-5 ⁺ | > 15 |

*The catalog numbers in this table are for media in ready-to-use packaging.

Related Products

| Product | Size | Catalog Number |
|---|----------------------------|----------------|
| Melanocyte Growth Medium (Ready-to-use) | 500 ml | C-24010 |
| Melanocyte Growth Medium Kit | 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 |
| Melanocyte Growth Medium M2 (Ready-to-use) | 500 ml | C-24300 |
| Melanocyte Growth Medium M2 (Ready-to-use), phenol red-free | 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 |
| DetachKit | 30 ml | C-41200 |
| | 125 ml | C-41210 |
| | 250 ml | C-41220 |
| Cryo-SFM | 30 ml | C-29910 |
| | 125 ml | C-29912 |
| NHEM.f Pellet | 1 million cells per pellet | C-14040 |
| NHEM.f M2 Pellet | 1 million cells per pellet | C-14042 |
| NHEM M2 adult Pellet | 1 million cells per pellet | C-14043 |

PromoCell GmbH

Sickingenstr. 63/65
69126 Heidelberg
Germany

Email: 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 2015