

Instruction Manual

Product	Size	Catalog Number
Human Nasal Epithelial Cells (HNEpC)	500,000 cryopreserved cells 500,000 proliferating cells	C-12620 C-12621
Human Tracheal Epithelial Cells (HTEpC)	500,000 cryopreserved cells 500,000 proliferating cells	C-12644 C-12645
Human Bronchial Epithelial Cells (HBEpC)	500,000 cryopreserved cells 500,000 proliferating cells	C-12640 C-12641
Human Small Airway Epithelial Cells (HSAEpC)	500,000 cryopreserved cells 500,000 proliferating cells	C-12642 C-12643
Human Renal Epithelial Cells (HREpC)	500,000 cryopreserved cells 500,000 proliferating cells	C-12665 C-12666
Human Renal Cortical Epithelial Cells (HRCEpC)	500,000 cryopreserved cells 500,000 proliferating cells	C-12660 C-12662

Product Description

The epithelium is the interface between the body and the external environment and covers all exterior surfaces and interior lumina. Depending on the tissue of origin, the functions of epithelial cells are diverse and include absorption, secretion, protection, transcellular transport, and sensation.

We offer a range of epithelial cells produced at our cell culture facility. The cells are isolated from normal human adult tissues of the airways (nasal mucosa, trachea, bronchi, and distal respiratory tract) and kidney (whole kidney and renal cortex).

Shortly after isolation, all our Human Epithelial Cells are cryopreserved at passage 2 (P2) using our proprietary, defined, animal-component free, and protein-free cryopreservation medium, Cryo-SFM. Each cryovial contains more than 500,000 viable cells after thawing. Proliferating cell cultures are made from cryopreserved cells that have been thawed and cultured for three days in our hands and shipped as growing cultures.

Quality Control

We perform rigid quality control tests for each lot of Epithelial Cells.

The cells are tested for cell morphology, adherence rate and cell viability. Furthermore, flow cytometric analyses for the cell-type specific marker cytokeratin are carried out for each lot (see page 5). Growth performance is tested through multiple passages under culture conditions without antibiotics or antimycotics. Upon request we provide HBEpC donors tested for proper epithelial cell barrier function in our Air-Liquid Interface (ALI) Medium (C-21080). ALI pre-screened HBEpC are tested by voltohmmeter for a stable trans-epithelial electrical resistance (TEER) of $>500 \Omega \cdot \text{cm}^2$ until 4 weeks of culture.

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

Our Epithelial Cells 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, cryopreserved cells should be stored in liquid nitrogen or seeded directly (see page 2). Proliferating cells must be processed immediately (see page 3).

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.

Use aseptic techniques and a laminar flow bench.

1

Prepare the medium

Calculate the required 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 retighten. Immerse the vial in a water bath (37°C) up to the height of 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 pre-warmed 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 and every two to three days thereafter. The cells should be subcultured, according to the subcultivation protocol (see page 4), once they have reached 70–90% confluency.



Protocol for Proliferating Cells

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

1

Incubate the cells

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



2

Replace the transport medium

Carefully open the vessel, rinse the inner side of the cap 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). Change the medium every two to three days. The cells should be subcultured, according to the subcultivation protocol (see page 4), once they have reached >70% confluency.



Subcultivation Protocol

Use aseptic techniques and a laminar flow bench.

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 re-suspend the cells by carefully pipetting up and down. Plate the cells according to the recommended seeding density in new cell culture vessels containing prewarmed PromoCell Growth Medium. Place the vessels in an incubator (37°C, 5% CO_2) and change the media every two or three days.



Specifications

Product	Recommended Culture Media*	Plating Density	Passage after Thawing	Marker	Population Doublings
Human Nasal Epithelial Cells (HNEpC)	C-21060	10,000 – 15,000 cells per cm ²	P2	Cytokeratin ⁺	> 15
Human Tracheal Epithelial Cells (HTEpC)	C-21060	10,000 – 15,000 cells per cm ²	P2	Cytokeratin ⁺	> 15
Human Bronchial Epithelial Cells (HBEpC)	C-21060	10,000 – 15,000 cells per cm ²	P2	Cytokeratin ⁺	> 15
Human Small Airway Epithelial Cells (HSAEpC)	C-21070	10,000 – 15,000 cells per cm ²	P2	Cytokeratin ⁺	> 15
Human Renal Epithelial Cells (HREpC)	C-26001 C-26030	10,000 – 15,000 cells per cm ²	P2	Cytokeratin ⁺	> 15
Human Renal Cortical Epithelial Cells (HRCEpC)	C-26001 C-26030	10,000 – 15,000 cells per cm ²	P2	Cytokeratin ⁺	> 15

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

Related Products

Product	Size	Catalog Number
Airway Epithelial Cell Growth Medium (Ready-to-use)	500 ml	C-21060
Airway Epithelial Cell Growth Medium Kit	500 ml	C-21160
Airway Epithelial Cell Basal Medium	500 ml	C-21260
Airway Epithelial Cell Basal Medium, phenol red-free	500 ml	C-21265
Airway Epithelial Cell Growth Medium SupplementMix	for 500 ml	C-39165
Airway Epithelial Cell Growth Medium SupplementPack	for 500 ml	C-39160
Air-Liquid Interface Medium (ALI-Airway)	500 ml	C-21080
Small Airway Epithelial Cell Growth Medium (Ready-to-use)	500 ml	C-21070
Small Airway Epithelial Cell Growth Medium Kit	500 ml	C-21170
Small Airway Epithelial Cell Basal Medium	500 ml	C-21270
Small Airway Epithelial Cell Basal Medium, phenol red-free	500 ml	C-21275
Small Airway Epithelial Cell Growth Medium SupplementMix	for 500 ml	C-39175
Small Airway Epithelial Cell Growth Medium SupplementPack	for 500 ml	C-39170
Mammary Epithelial Cell Growth Medium (Ready-to-use)	500 ml	C-21010
Mammary Epithelial Cell Growth Medium Kit	500 ml	C-21110
Mammary Epithelial Cell Basal Medium	500 ml	C-21210

Product	Size	Catalog Number
Mammary Epithelial Cell Basal Medium, phenol red-free	500 ml	C-21215
Mammary Epithelial Cell Growth Medium Supplement Mix	for 500 ml	C-39115
Mammary Epithelial Cell Growth Medium Supplement Pack	for 500 ml	C-39110
Renal Epithelial Cell Growth Medium 2 (Ready-to-use)	500 ml	C-26030
Renal Epithelial Cell Growth Medium 2 Kit	500 ml	C-26130
Renal Epithelial Cell Basal Medium 2	500 ml	C-26230
Renal Epithelial Cell Basal Medium 2, phenol red-free	500 ml	C-26235
Renal Epithelial Cell Growth Medium 2 Supplement Mix	for 500 ml	C-39606
Renal Epithelial Cell Growth Medium 2 Supplement Pack	for 500 ml	C-39605
Detach Kit	30 ml 125 ml 250 ml	C-41200 C-41210 C-41220
Cryo-SFM	30 ml 125 ml	C-29910 C-29912
HNEpC Pellet	1 million cells per pellet	C-14062
HTEpC Pellet	1 million cells per pellet	C-14064
HBEpC Pellet	1 million cells per pellet	C-14063
HSAEpC Pellet	1 million cells per pellet	C-14065
HREpC Pellet	1 million cells per pellet	C-14067
HRCEpC Pellet	1 million cells per pellet	C-14068
BPE-15	2 ml	C-30020
BPE-26	2 ml	C-30021

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