

Airway Epithelial Cell Media

PromoCell

Instruction Manual

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

Recommended for

- Human Nasal Epithelial Cells (HNEpC)
- Human Bronchial Epithelial Cells (HBEpC)
- Human Tracheal Epithelial Cells (HTEpC)

Product Description

PromoCell Airway Epithelial Cell Growth Medium is a serum-free medium optimized for the *in vitro* cultivation of epithelial cells from large air passages. The medium is optimized for primary human cells, but can also be used for equine, porcine, murine, bat and rat airway epithelial cells.

PromoCell Airway Epithelial Cell 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

PromoCell Airway Epithelial Cell Growth Medium contains all the growth factors and supplements necessary for the optimal growth of human airway epithelial cells, e.g. Epidermal Growth Factor, Insulin, and Transferrin (for details see table on page 2). PromoCell Airway Epithelial Cell Growth 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 pi-

petting 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.

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, prewarm only an aliquot of the complete medium and keep the remaining medium refrigerated at 4–8°C.

Quality Control

All lots of PromoCell Airway Epithelial Cell Growth Medium are subjected to comprehensive quality control tests using primary human airway epithelial cells. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested airway epithelial cells. 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)

■ Bovine Pituitary Extract	0.004 ml / ml
■ Epidermal Growth Factor (recombinant human)	10 ng / ml
■ Insulin (recombinant human)	5 µg / ml
■ Hydrocortisone	0.5 µg / ml
■ Epinephrine	0.5 µg / ml
■ Triiodo-L-thyronine	6.7 ng / ml
■ Transferrin (recombinant human)	10 µg / ml
■ Retinoic Acid	0.1 ng / ml

If you require special media modifications, we offer a Custom Media Service starting at 10 bottles per order. Please ask for details.

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