

# Small airway epithelial cell media

## Instruction manual

Product	Size	Catalog number
Small Airway Epithelial Cell Growth Medium (Ready-to-use) – Includes Basal Medium and SupplementMix	500 ml	C-21070
Small Airway Epithelial Cell Growth Medium Kit – Includes Basal Medium and SupplementPack	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

### Recommended for

- Human Small Airway Epithelial Cells (HSAEpC)

### Product description

Our Small Airway Epithelial Cell Growth Medium is a serum-free medium developed for the *in vitro* cultivation of epithelial cells from the distal respiratory system. The medium is optimized for primary human cells, but can also be used for porcine small airway epithelial cells.

The Small 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.

### Storage and stability

Our Small Airway Epithelial Cell Growth Medium contains all the growth factors and

supplements necessary for the optimal growth of human small airway epithelial cells, e.g. L-Epidermal Growth Factor (recombinant human), Insulin, and Transferrin (for details see table on page 2). The Small Airway Epithelial Cell Growth Medium does not contain antibiotics or antimycotics and is formulated for use in an incubator with an atmosphere of 5% CO<sub>2</sub>.

### 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, prewarm only an aliquot of the complete medium at 15–25°C and keep the remaining medium refrigerated at 4–8°C.

### Quality control

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

### 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 to use 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
L-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
Bovine Serum Albumin-Fatty Acid Free (BSA-FAF)	2.5 mg/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](mailto:info@promocell.com) to find out more.

---

**PromoCell GmbH**  
 Sickingerstr. 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](mailto:info@promocell.com)  
[www.promocell.com](http://www.promocell.com)

© PromoCell GmbH

03/2024