

Airway Epithelial Cell Growth Medium 2



Instruction manual

Product	Size	Catalog number
Airway Epithelial Cell Growth Medium 2	500 ml	C-21040

Recommended for

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

Product description

Our Airway Epithelial Cell Growth Medium 2 is a serum- and bovine pituitary extract (BPE)-free medium optimized for the *in vitro* cultivation of epithelial cells from large airway passages.

Our Airway Epithelial Cell Growth Medium 2 is available as a ready-to-use medium consisting of a 500 ml bottle of basal medium and one vial of SupplementMix. Adding the SupplementMix to the basal medium results in the complete growth medium.

Supplementation details

Our Airway Epithelial Cell Growth Medium 2 contains all the necessary growth factors and supplements for optimal growth of human airway epithelial cells. The Airway Epithelial Cell Growth Medium 2 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 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.

Storage and stability

Immediately after arrival, store the basal medium at 4–8°C in the dark and the SupplementMix at -20°C. 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 2 are subjected to comprehensive quality control tests using primary human airway epithelial cells (HBEPc). Each lot of basal medium and supplement is checked for growth-promoting activity and 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 the appropriate MSDS.

Note: Due to the absence of serum, this PromoCell medium is 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.

For thawing and expansion of cryopreserved epithelial cells using our Airway Epithelial Cell Growth Medium 2, please refer to the corresponding instruction manual for handling human airway epithelial cells. The instruction manual can be downloaded from our website at www.promocell.com.

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

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

09/2025