

# Mammary Epithelial Cell Media

## Instruction Manual

Product	Size	Catalog Number
Mammary Epithelial Cell Growth Medium (Ready-to-use) – Includes Basal Medium and SupplementMix	500 ml	C-21010
Mammary Epithelial Cell Growth Medium Kit – Includes Basal Medium and SupplementPack	500 ml	C-21110
Mammary Epithelial Cell Basal Medium	500 ml	C-21210
Mammary Epithelial Cell Basal Medium, phenol red-free	500 ml	C-21215
Mammary Epithelial Cell Growth Medium SupplementMix	for 500 ml	C-39115
Mammary Epithelial Cell Growth Medium SupplementPack	for 500 ml	C-39110

### Recommended for

- Human Mammary Epithelial Cells

### Product Description

Our Mammary Epithelial Cell Growth Medium is a serum-free medium developed for the *in vitro* cultivation of epithelial cells from the mammary gland. This medium has been validated using primary human mammary epithelial cells tested for adherence rate and typical morphology and growth performance. The medium is optimized for primary human cells, but can also be used for rat and murine mammary epithelial cells as well as mammary carcinoma cell lines.

The Mammary 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

Our Mammary Epithelial Cell Growth Medium

contains all the growth factors and supplements necessary for the optimal growth of human mammary epithelial cells, e.g. Epidermal Growth Factor and Insulin (for details see the table on page 2). The Mammary 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, pre-warm only an aliquot of the complete medium and keep the remaining medium refrigerated at 4–8°C.

### Quality Control

All lots of PromoCell Mammary Epithelial Cell Media are subjected to comprehensive quality control tests. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested 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)	Mammary Epithelial Cell Growth 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

---

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**  
Sickingenstr. 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