

Renal Epithelial Cell Medium 2

Instruction manual

Product	Size	Catalog number
Renal Epithelial Cell Growth Medium 2 (Ready-to-use) – includes Basal Medium and SupplementMix	500 ml	C-26030
Renal Epithelial Cell Growth Medium 2 Kit – includes Basal Medium and SupplementPack	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 SupplementMix	for 500 ml	C-39606
Renal Epithelial Cell Growth Medium 2 SupplementPack	for 500 ml	C-39605

Recommended for

- Human Renal Epithelial Cells (HREpC)
- Human Renal Cortical Epithelial Cells (HRCEpC)

Product description

Our Renal Epithelial Cell Growth Medium 2 is an optimized low-serum medium that additionally contains Epinephrine. The combination of low serum and selected growth factors ensures successful growth *in vitro* making the Renal Epithelial Cell Growth Medium 2 the first choice for our Renal Epithelial Cells.

The Renal Epithelial Cell Medium 2 is optimized for primary human cells, but can also be used for porcine and mouse renal epithelial cells.

Our Renal Epithelial Cell Media are 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 SupplementMix and SupplementPacks can also be purchased separately.

Supplementation details

The Renal Epithelial Cell Growth Medium 2 is a low-serum medium that additionally contains Epinephrine (for details see the table above). The Renal 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 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.

Storage and stability

Store the basal medium at 4 to 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 to 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 to 8°C.

Quality control

All lots of PromoCell Renal Epithelial Cell Medium are subjected to comprehensive quality control tests using primary human renal epithelial cells. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested renal 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 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)	Renal Epithelial Cell Medium 2
Fetal Calf Serum	0.05 ml/ml
L-Epidermal Growth Factor (recombinant human)	10 ng/ml
Insulin (recombinant human)	5 µg/ml
Epinephrine	0.5 µg/ml
Hydrocortisone	900 ng/ml
Transferrin (recombinant human)	5 µg/ml
Triiodo-L-thyronine	4 pg/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 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
www.promocell.com

© PromoCell GmbH