

Follicle dermal papilla cell media

Instruction manual

Product	Size	Catalog number
Follicle Dermal Papilla Cell Growth Medium (Ready-to-use) – Includes Basal Medium and SupplementMix	500 ml	C-26501
Follicle Dermal Papilla Cell Growth Medium Kit – Includes Basal Medium and SupplementPack	500 ml	C-26502
Follicle Dermal Papilla Cell Basal Medium	500 ml	C-26500
Follicle Dermal Papilla Cell Basal Medium, phenol red-free	500 ml	C-26505
Follicle Dermal Papilla Cell Growth Medium SupplementMix	for 500 ml	C-39625
Follicle Dermal Papilla Cell Growth Medium SupplementPack	for 500 ml	C-39620

Recommended for

- Human Follicle Dermal Papilla Cells

Product description

Our Follicle Dermal Papilla Cell Growth Medium is a low-serum medium that has been developed for the *in vitro* cultivation of human follicle dermal papilla cells.

The Follicle Dermal Papilla 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.

Supplementantation details

The Follicle Dermal Papilla Cell Growth Medium contains all the growth factors and supplements necessary for the optimal growth of human follicle dermal papilla cells, e.g. basic Fibroblast Growth Factor and Insulin (for

details see the table on the next page). The Follicle Dermal Papilla 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 pipetting up and down. Then, transfer the entire contents 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 Follicle Dermal Papilla Cell Growth Medium are subjected to comprehensive quality control tests using primary human follicle dermal papilla cells. Each lot is checked for growth promoting activity, adherence rate and typical morphology of the tested follicle dermal papilla 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)	Follicle Dermal Papilla Cell Growth Medium
Fetal Calf Serum	0.004 ml/ml
Bovine Pituitary Extract	0.004 ml/ml (52 µg/ml)
Basic Fibroblast Growth Factor (recombinant human)	1 ng/ml
Insulin (recombinant human)	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 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