

Preadipocyte / Adipocyte Media

PromoCell

Instruction Manual

■ Preadipocyte Growth Medium

Product	Size	Catalog Number
Preadipocyte Growth Medium (Ready-to-use)	500 ml	C-27410
Preadipocyte Growth Medium Kit	500 ml	C-27417
Preadipocyte Basal Medium	500 ml	C-27411
Preadipocyte Basal Medium, phenol red-free	500 ml	C-27415
Preadipocyte Growth Medium SupplementMix	for 500 ml	C-39425
Preadipocyte Growth Medium SupplementPack	for 500 ml	C-39427

■ Preadipocyte Differentiation Medium

Product	Size	Catalog Number
Preadipocyte Differentiation Medium (Ready-to-use)	500 ml	C-27436
Preadipocyte Differentiation Medium Kit	500 ml	C-27437
Preadipocyte Differentiation Medium SupplementMix	for 500 ml	C-39436
Preadipocyte Differentiation Medium SupplementPack	for 500 ml	C-39437

■ Adipocyte Nutrition Medium

Product	Size	Catalog Number
Adipocyte Nutrition Medium (Ready-to-use)	500 ml	C-27438
Adipocyte Nutrition Medium Kit	500 ml	C-27439
Adipocyte Basal Medium	500 ml	C-27431
Adipocyte Basal Medium, phenol red-free	500 ml	C-27435
Adipocyte Nutrition Medium SupplementMix	for 500 ml	C-39438
Adipocyte Nutrition Medium SupplementPack	for 500 ml	C-39439

Recommended for

- Human White Preadipocytes (HWP)
- Human Adipocytes

Product Description

PromoCell Preadipocyte Growth Medium is a low-serum medium developed for the undifferentiated expansion of human preadipocytes. The Preadipocyte Differentiation Medium is a serum-free medium, which induces the differen-

tiation of preadipocytes into mature adipocytes. The mature adipocytes can be cultivated using the Adipocyte Nutrition Medium. The media are optimized for primary human cells, but can also be used for murine preadipocytes/adipocytes.

PromoCell Preadipocyte/Adipocyte Media are available as Medium (Ready-to-use) or as Medium Kit.

The Medium (Ready-to-use) consists of a 500ml bottle of Basal Medium and

one vial of SupplementMix. The Medium Kit consists of a 500ml 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.

Final supplement concentrations (after addition to the medium)

	Preadipocyte Growth Medium	Preadipocyte Differentiation Medium	Adipocyte Nutrition Medium
■ Fetal Calf Serum	0.05 ml / ml	—	0.03 ml / ml
■ Endothelial Cell Growth Supplement	0.004 ml / ml	—	—
■ Epidermal Growth Factor (recombinant human)	10 ng / ml	—	—
■ Hydrocortisone	1 µg / ml	—	—
■ d-Biotin	—	8 µg / ml	8 µg / ml
■ Insulin (recombinant human)	—	0.5 µg / ml	0.5 µg / ml
■ Dexamethasone	—	400 ng / ml	400 ng / ml
■ IBMX	—	44 µg / ml	—
■ L-Thyroxine	—	9 ng / ml	—
■ Ciglitazone	—	3 µg / ml	—
■ Heparin	90 µg / ml	—	—

Supplementation Details

PromoCell Preadipocyte Growth Medium contains all the growth factors and supplements necessary for the optimal growth of undifferentiated human preadipocytes. PromoCell Preadipocyte Differentiation Medium is a serum-free medium containing all supplements for effectively inducing the differentiation of preadipocytes into mature adipocytes, e.g. Ciglitazone. PromoCell Adipocyte Nutrition Medium is a low-serum medium containing all the growth factors and supplements necessary for the maintenance of mature adipocytes (for details see the table above). The Media do not contain antibiotics or antimycotics and are 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°C–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 ECGS. This does not affect the activity. Optionally, the precipitate can be removed by centrifugation under sterile conditions.

Instructions for Use of PromoCell Preadipocyte Differentiation Medi- um / Adipocyte Nutrition Medium

For differentiating human preadipocytes, culture the cells in Preadipocyte Growth Medium to total confluency stage. Replace the Growth Medium with Preadipocyte Differentiation Medium for 72 hours. Then change to PromoCell Adipocyte Nutrition Medium to complete the differentiation process. Feed the cells every 2–3 days with fresh Adipocyte Nutrition Medium. The differentiation process to mature adipocytes is completed after 12–14 days.

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 Preadipocyte/Adipocyte Media are subjected to comprehensive quality control tests using primary human white preadipocytes. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested preadipocytes. In addition, the Preadipocyte Differentiation Medium/Adipocyte Nutrition Medium are tested for the capacity to induce the terminal differentiation of preadipocytes into mature adipocytes. 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.

If you require special media modifications, we offer a Custom Media Service starting at 10 bottles per order. Please ask for details.

PromoCell GmbH

Sickingenstr. 63/65
69126 Heidelberg
Germany

Email: 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