

Smooth Muscle Cell Media 2

Instruction manual

Product	Size	Catalog number
Smooth Muscle Cell Growth Medium 2 (Ready-to-use) – Includes Basal Medium and SupplementMix	500 ml	C-22062
Smooth Muscle Cell Growth Medium 2 Kit – Includes Basal Medium and SupplementPack	500 ml	C-22162
Smooth Muscle Cell Basal Medium 2	500 ml	C-22262
Smooth Muscle Cell Basal Medium 2, phenol red-free	500 ml	C-22267
Smooth Muscle Cell Growth Medium 2 SupplementMix	for 500 ml	C-39267
Smooth Muscle Cell Growth Medium 2 SupplementPack	for 500 ml	C-39262

Recommended for

- Human Umbilical Artery Smooth Muscle Cells (HUASMC)
- Human Coronary Artery Smooth Muscle Cells (HCASMC)
- Human Pulmonary Artery Smooth Muscle Cells (HPASMC)
- Human Aortic Smooth Muscle Cells (HAoSMC)
- Human Bronchial Smooth Muscle Cells (HBSMC)
- Human Tracheal Smooth Muscle Cells (HTSMC)
- Human Uterine Smooth Muscle Cells (HUtSMC)

Product description

Our Smooth Muscle Cell Growth Medium 2 is a low serum medium developed for the *in vitro* cultivation of smooth muscle cells from blood vessels and hollow organs. The medium is optimized for primary human cells, but can also be used for bovine, porcine, dog, murine and rat smooth muscle cells.

The Smooth Muscle 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 the SupplementMix and SupplementPack can also be purchased separately.

Supplementation details

Our Smooth Muscle Cell Growth Medium 2 contains all the growth factors and supplements necessary for the optimal growth of human smooth muscle cells. The medium contains Epidermal Growth Factor, basic Fibroblast Growth Factor, and Insulin (for details see table on page 2). The Smooth Muscle 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 4–8°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 Smooth Muscle Cell Growth Medium are subjected to comprehensive quality control tests using primary human smooth muscle cells. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested smooth muscle 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)	Size
Fetal Calf Serum	0.05 ml/ml
Epidermal Growth Factor (recombinant human)	0.5 ng/ml
Basic Fibroblast Growth Factor (recombinant human)	2 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

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

06/2024