

Instruction Manual

Product	Size	Catalog Number
Fibroblast Growth Medium 2 (Ready-to-use)	500 ml	C-23020
Fibroblast Growth Medium 2 Kit	500 ml	C-23120
Fibroblast Basal Medium 2	500 ml	C-23220
Fibroblast Basal Medium 2, phenol red-free	500 ml	C-23225
Fibroblast Growth Medium 2 SupplementMix	for 500 ml	C-39325
Fibroblast Growth Medium 2 SupplementPack	for 500 ml	C-39320

Recommended for

- Normal Human Dermal Fibroblasts (NHDF), adult donor
- Human Pulmonary Fibroblasts (HPF)
- Human Aortic Adventitial Fibroblasts (HAoAF)
- Human Uterine Fibroblasts (HUF)

Product Description

PromoCell Fibroblast Growth Medium 2 is a low-serum medium and has been developed for the *in vitro* cultivation of adult Human Fibroblasts. The medium is optimized for primary human cells, but can also be used for bovine, porcine, and rat fibroblasts, as well as fibroblast cell lines.

PromoCell Fibroblast Growth Medium 2 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

PromoCell Fibroblast Growth Medium 2 contains all the growth factors and supplements necessary for the optimal growth of human adult fibroblasts, e.g. basic Fibroblast Growth Factor, Insulin, and Fetal Calf Serum (for details see table on page 2). PromoCell Fibroblast 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 Fibroblast Growth Medium 2 are subjected to comprehensive quality control tests using primary human adult fibroblasts. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested adult fibroblasts. 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)

▪ Fetal Calf Serum	0.02 ml / 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. 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