

## Instruction Manual

Product	Size	Catalog Number
Chondrocyte Growth Medium (Ready-to-use)	500 ml	C-27101
Chondrocyte Basal Medium	500 ml	C-27111
Chondrocyte Basal Medium, phenol red-free	500 ml	C-27115
Chondrocyte Growth Medium SupplementMix	for 500 ml	C-39635

### Recommended for

- Human Chondrocytes (HCH)

### Product Description

PromoCell Chondrocyte Growth Medium has been developed for the *in vitro* cultivation of human chondrocytes. The medium is optimized for primary human cells, but can also be used for primary porcine and rat chondrocytes.

The Medium (Ready-to-use) consists of a 500 ml bottle of Basal Medium and one vial of SupplementMix. Adding the SupplementMix to the Basal Medium results in the complete Growth Medium. Basal Medium (with or without phenol red) as well as the SupplementMix can also be purchased separately.

### Supplementation Details

PromoCell Chondrocyte Growth Medium contains all the components necessary for the optimal growth of hu-

man chondrocytes. PromoCell Chondrocyte Growth Medium does not contain antibiotics or antimycotics and is formulated for use in an incubator with an atmosphere of 5% CO<sub>2</sub>.

### Preparation of the supplemented Medium for Use

Thaw the SupplementMix at 15 to 25°C. Aseptically mix the SupplementMix by carefully pipetting up and down. Then, transfer the entire content of the SupplementMix 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 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 SupplementMix 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 Chondrocyte Growth Medium are subjected to comprehensive quality control tests using primary human chondrocytes. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested chondrocytes. 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).

## Final supplement concentrations (after addition to the medium)

■ Fetal Calf Serum	0.1 ml / ml
--------------------	-------------

### 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.

#### PromoCell GmbH

Sickingenstr. 63/65  
69126 Heidelberg  
Germany

Email: [info@promocell.com](mailto:info@promocell.com)  
[www.promocell.com](http://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