

Instruction manual

Product	Size	Catalog number
Chondrocyte Growth Medium (Ready-to-use) – Includes Basal Medium and SupplementMix	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

Our 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.

Supplementantation details

Our Chondrocyte Growth Medium contains all the components necessary for the optimal growth of human chondrocytes. The Chondrocyte 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 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 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)	Chondrocyte Growth Medium
Fetal Calf Serum	0.1 ml/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