

Product	Size	Catalog Number
Pericyte Growth Medium 2 (Ready-to-use)	500 ml	C-28041

## Recommended for

- Human Pericytes

## Product Description

PromoCell Pericyte Growth Medium has been developed for the *in vitro* cultivation of human pericytes from all tissues. The Medium (Ready-to-use) consists of a 500ml bottle of Basal Medium and one vial of SupplementMix.

Adding the SupplementMix to the Basal Medium results in the complete Growth Medium.

## Supplementation Details

PromoCell Pericyte Growth Medium contains all components necessary for the optimal growth of human pericytes. PromoCell Pericyte Growth Medium does not contain antibiotics or antimycotics and is formulated for the use in an incubator with an atmosphere of 5% CO<sub>2</sub>.

## Preparation of the supplemented Medium for Use

Thaw the SupplementMix in a 37°C water bath with occasional swirling. Do not

incubate longer than necessary! **In case of visible precipitates after complete thawing, mix gently until all precipitates have re-dissolved.** Then, transfer the entire content of the SupplementMix to the Basal Medium. Close the bottle and swirl gently until a homogenous mixture is formed.

## Pericyte Subculture

For routine subculture of Pericytes the use of Accutase (C-41310) is recommended with all PromoCell Pericyte Growth Media.

## Storage and Stability

Store the Basal Medium at 4–8°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 4–8°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 Pericyte Growth Medium are subjected to comprehensive quality control tests using primary human placenta-derived pericytes. Each lot is checked for growth promoting activity, adherence rate, and typical morphology of the tested pericytes. 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 product is for *in vitro* use only and not for diagnostic or therapeutic procedures. For safety precautions please see appropriate MSDS.

### 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

© PromoCell GmbH