

## Instruction Manual

Product	Size	Catalog Number
Pericyte Growth Medium 2 (Ready-to-use) – Includes Basal Medium and SupplementMix	500 ml	C-28041

### Recommended for

- Human Pericytes

### Product Description

Our 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 500 ml bottle of Basal Medium and one vial of SupplementMix.

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

### Supplementation Details

The Pericyte Growth Medium contains all components necessary for the optimal growth of human pericytes. Our 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, myco- plasma).

### Intended Use

The product is for *in vitro* use only and not for diagnostic or therapeutic procedures. For safety precautions please see appropriate MSDS.

---

If you require special media modifications, we offer a custom media service starting at 10 bottles per order.  
Contact us at [info@promocell.com](mailto: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](mailto:info@promocell.com)  
[www.promocell.com](http://www.promocell.com)

© PromoCell GmbH

05/2022