

PromoExQ Pericyte Growth Medium 2

Instruction manual

Product	Size	Catalog number
PromoExQ Pericyte Growth Medium 2, phenol red-free	500 ml	EQ-C-28042

Recommended for

- Human Pericytes

Our PromoExQ Pericyte Growth Medium 2, phenol red-free (prf) has been developed for the GMP-compliant¹ isolation and *in vitro* expansion of human pericytes from all tissues. The medium 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.

PromoExQ Pericyte Growth Medium 2 (prf) contains all components necessary for the optimal growth of human pericytes.

PromoExQ Pericyte Growth Medium 2 (prf) does not contain antibiotics or antimycotics and is formulated for the use in an incubator with an atmosphere of 5% CO₂.

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.

For routine subculture of pericytes the use of Accutase is recommended with the PromoExQ Pericyte Growth Medium 2 (prf).

The basal medium is shipped at room temperature and should be stored at 4 to 8°C in the dark upon arrival. The SupplementMix is shipped on dry ice and should be stored at -20 to -16°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.

All lots of PromoExQ Pericyte Growth Medium 2 (prf) are subjected to comprehensive quality control tests using primary human placenta-derived pericytes. Each lot of PromoExQ Pericyte Growth Medium 2 (prf) is checked for growth promoting activity and for the absence of microbial contaminants (fungi, bacteria, and mycoplasma) in the basal media and the supplements. In addition, endotoxin levels are determined in basal media. For details see lot specific Certificate of Analysis.

The regulated products are intended for research use or further manufacturing. They are not intended for direct administration to humans or animals. For safety precautions please see appropriate MSDS.

¹ according to international guidelines

If you require customization of our media, we offer Custom GMP Services.
Contact us at gmp@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