

PromoExQ MSC Growth Media



Instruction manual

Product	Size	Catalog number
PromoExQ MSC Growth Medium 2, phenol red-free	500 ml	EQ-C-28017
PromoExQ MSC Growth Medium XF, phenol red-free	500 ml	EQ-C-28018

Recommended for

- Human Mesenchymal Stem Cells from Bone Marrow (hMSC-BM)
- Human Mesenchymal Stem Cells from Umbilical Cord Matrix (hMSC-UC)
- Human Mesenchymal Stem Cells from Adipose Tissue (hMSC-AT)

Product description

Our PromoExQ MSC Growth Media are developed for the GMP-compliant¹ *in vitro* expansion of multipotent mesenchymal stem cells (MSC) from bone marrow, the umbilical cord matrix (Wharton's Jelly), and adipose tissue. The PromoExQ MSC Growth Medium 2 (prf) (EQ-C-28017) is an optimized low-serum formulation for isolation and routine culture of human MSC. The PromoExQ MSC Growth Medium XF (prf) (EQ-C-28018) is a serum-free and xeno-free medium for the routine culture of MSC in a xeno-free environment. The PromoExQ MSC Growth Media consist of a bottle of basal medium and one vial of SupplementMix. Adding the SupplementMix to the basal medium results in the complete growth medium.

Supplementation details

Our PromoExQ MSC Growth Media contain all the growth factors and supplements necessary

for the optimal expansion of human mesenchymal stem cells.

Note: The PromoExQ MSC Growth Media do not contain antibiotics or antimycotics and are formulated for use in an incubator with an atmosphere of 5% CO₂.

Culture vessel surface coating

Our PromoExQ MSC Growth Medium XF (prf) (EQ-C-28018) does not contain attachment- and spreading factors. Therefore, culture vessels to be used with PromoExQ MSC Growth Medium XF (prf) must be precoated either with 1 µg/cm² human fibronectin or 0.5 µg/cm² human vitronectin according to the instructions of the manufacturer. When using our PromoExQ MSC Growth Medium 2 (prf), no surface coating is required.

Preparation of the supplemented medium for use

Thaw the supplements at 15–25°C. Aseptically mix the supplement solutions by carefully pipetting up and down. 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.

MSC subculture

For routine subculture of hMSC the use of Accutase is recommended with all PromoExQ MSC Growth Media.

Storage and stability

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 supplements to the basal medium, the complete growth medium is stable for 3 weeks at 4 to 8°C. Do not freeze the complete medium.

For use, pre-warm only an aliquot of the complete medium at 15 to 25°C and keep the remaining medium refrigerated at 4 to 8°C.

Note: For the SupplementMix of the PromoExQ MSC Growth Medium XF (prf) (EQ-C-28018) a concurrent stability study is being conducted in parallel with the commercial distribution to support the indicated shelf life. Although full-term data is not yet available, real-time data supports interim release, with ongoing monitoring ensuring continued compliance with stability requirements and product quality.

¹ according to international guidelines

Quality control

All lots of PromoCell PromoExQ MSC Growth Media are subjected to comprehensive quality control tests using primary human mesenchymal stem cells. Each lot of PromoExQ MSC Growth

Media are checked for growth promoting activity and for the absence of the 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.

Intended use

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.

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