

PromoExQ Fibroblast Growth Medium D-ACF



Instruction manual

Product	Size	Catalog number
PromoExQ Fibroblast Growth Medium D-ACF, phenol red-free	500 ml	EQ-C-23017

Recommended for

- Normal Human Dermal Fibroblasts (NHDF), juvenile donor

Product description

Our PromoExQ Fibroblast Growth Medium D-ACF, phenol red-free (prf) is a defined and animal component-free medium that has been developed for the GMP-compliant¹ cultivation of juvenile Normal Human Dermal Fibroblasts.

The optimized phenol red-free medium also supports the growth of adult Normal Human Dermal Fibroblasts. The medium consists of a 450 ml bottle of PromoExQ basal medium and one vial of PromoExQ SupplementMix. Adding the PromoExQ SupplementMix to the basal medium results in the complete growth medium.

Supplementation details

Our PromoExQ Fibroblast Growth Medium D-ACF (prf) contains all the growth factors and supplements necessary for the optimal growth of Normal Human Dermal Fibroblasts.

Note: The medium does not contain anti-

biotics 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 15–25°C. Aseptically mix the supplement solutions by carefully pipetting up and down. Then, transfer the entire content of the supplement to the basal medium. Close the bottle and swirl gently until a homogenous mixture is formed.

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 PromoExQ 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 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.

Note: 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.

Quality control

All lots of our PromoCell PromoExQ Fibroblast Growth Medium D-ACF (prf) are subjected to comprehensive quality control tests using Normal Human Dermal Fibroblasts from juvenile donors. Each lot of PromoExQ Fibroblast Growth Medium D-ACF (prf) is 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.

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.

¹ according to international guidelines

Download corresponding documents: www.promocell.com/eq-certificates

If you require special media modifications, 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

09/2025