

# Fibroblast Growth Media D-ACF

## Instruction manual

| Product   | Size   | Catalog number |
|---|--------|----------------|
| Fibroblast Growth Medium D-ACF, phenol red-free | 500 ml | C-23017        |
| Fibroblast Growth Medium D-ACF                  | 500 ml | C-23027        |

### Recommended for

- Normal Human Dermal Fibroblasts (NHDF), juvenile donor

### Product description

Our Fibroblast Growth Media D-ACF are defined and animal-component-free media. They have been developed for the *in vitro* cultivation of juvenile Normal Human Dermal Fibroblasts, but also the growth of adult Normal Human Dermal Fibroblasts is supported. Our Fibroblast Growth Medium D-ACF and Fibroblast Growth Medium D-ACF, phenol red-free are available as Ready-to-use media. The media (Ready-to-use) consist of a 450 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

Our Fibroblast Growth Media D-ACF contain all the growth factors and supplements necessary for the optimal growth of Normal Human Dermal Fibroblasts. The media do

not contain antibiotics or antimycotics and are formulated for use in an incubator with an atmosphere of 5% CO<sub>2</sub>.

### 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 each supplement to the Basal Medium. Close the bottle and swirl gently until a homogenous mixture is formed.

### 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 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.

### Quality control

All lots of PromoCell Fibroblast Growth Media D-ACF are subjected to comprehensive quality control tests using Normal Human Dermal Fibroblasts from juvenile donors. Each lot is checked for growth promoting activity and for the absence of microbial contaminants (fungi and bacteria).

### Intended use

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

**Note:** Due to the absence of serum, PromoCell media are not suitable for trypsin neutralization (e.g. when splitting the cells). Instead we recommend using our DetachKit (C-41200, C-41210, C-41220), which contains HEPES BSS, Trypsin/EDTA and Trypsin Neutralizing Solution.

---

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

[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