

Cell Dissociation Solution ACF

Instruction Manual

Product	Size	Catalog Number
Cell Dissociation Solution ACF	100 ml	C-41320

Product Description

Our Cell Dissociation Solution ACF is a non-enzymatic, defined and animal-component free alternative to trypsin. The formulation is based on a proprietary blend of chelators and allows for efficient detachment of adherent cells under defined conditions in routine subculture. The gentle cell disaggregation guarantees high cell viability, even after prolonged exposure times. Unlike enzyme-based solutions the Cell Dissociation Solution ACF does not alter cell surface proteins and neutralization is not necessary.

Note: The Cell Dissociation Solution ACF is not intended for use with very strongly adherent cells, e.g., macrophages.

Instruction of use

1. Pre-warm the Cell Dissociation Solution ACF to ambient temperature.
2. Aspirate the medium from the culture vessel.
3. Wash the cell monolayer twice with Dulbecco's PBS w/o Ca²⁺/Mg²⁺ (C-40232).

4. Discard the PBS, add the Cell Dissociation Solution ACF (50–100 µl/cm²) and gently swirl the vessel to completely cover the vessel surface.

5. Incubate the cells at 37°C, and periodically check under a microscope. When the cells begin to round up tapping the side of the vessel will facilitate the detachment.

Note: The incubation time needed to dislodge the cells will vary depending on the cell type (typically 5 to 10 minutes) and culture medium used and must therefore be determined empirically by the user for each type of culture. Five minutes is a good starting point to check the status of the detachment process for the first time.

6. After detachment, disperse cells into suspension by gentle repeated pipetting.

7. Centrifuge the cells at 300 x g for 4 minutes. Remove as much of the Cell Dissociation Solution ACF as possible and resuspend the pellet in an appropriate medium.

Storage and Stability

Store at 2 to 8°C immediately after arrival. If stored properly, the product is stable until the expiry date stated on the label.

Quality Control

All lots of PromoCell Cell Dissociation Solution ACF are subjected to comprehensive quality control tests. Each lot is routinely tested for biological function, absence of cytotoxicity, and physical parameters such as osmolarity and pH level. Approved in-house lots are used as a reference.

In addition, all lots have been tested for the absence of microbial contaminants (fungi, bacteria).

Intended Use

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

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