

## Instruction Manual

Product	Size	Catalog Number
Accutase-Solution, primary human cell culture tested	100 ml	C-41310

### Product Description

Accutase is a sterile filtered ready-to-use solution specially developed for gentle and effective cell dissociation. It is a unique mixture of proteolytic and collagenolytic enzymes from an invertebrate species. Accutase is supplied in Dulbecco's PBS containing 0.5 mM EDTA.

### Purity

Accutase has been ultra-filtered (100,000 dalton cut off) before lyophilization and contains only purified enzymes. No stabilisers or other proteins have been added.

### Instruction of use

Use Accutase in the same manner that trypsin EDTA solutions are used to detach adherent cells from surfaces. Always handle cells with standard aseptic techniques if the cells are subcultured or propagated further.

1. Wash the cell substrate with a small volume of sterile PBS. Remove the wash liquid.

2. Add 8-10 ml Accutase per 75 cm<sup>2</sup> surface area to the culture vessel.

3. Return the culture vessel to an incubator at 37 °C and allow the cells to detach for 5 to 45 minutes.

4. When the cells have completely detached from the cell culture vessel carefully aspirate the cell suspension and transfer it to a centrifugation tube. Spin down the cells for 3 minutes at 220 x g. Inhibition of the enzyme reaction is not required.

5. Discard the supernatant, resuspend the cells in the appropriate medium and seed the cells in new cell culture vessels.

### Storage and Stability

Store at -20°C immediately after arrival. If stored properly, the product is stable until the expiry date stated on the label. Once thawed, Accutase is stable for 2 months when stored at 2 to 8 °C.

### Quality Control

All lots of PromoCell Accutase-Solution are subjected to comprehensive quality control tests. Each lot is routinely tested for biological function, absence of cytotoxicity, and physical parameters such as osmolality 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](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