

Instructions for Use

Accutase

A sterile filtered ready-to-use solution specially developed for gentle and effective cell dissociation.



2729369-000



SARTORIUS

A sterile filtered ready-to-use solution specially developed for gentle and effective cell dissociation.

REF

03-073-1B (100 mL)



Store at: -20 to -10°C

Contents

1 Product Description4

2 Purity4

3 Instructions for Use4

4 Storage and Stability5

5 Quality Control5

6 Intended Use5

1 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.

2 Purity

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

3 Instructions for 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 cm² surface area to the cultureware.
3. Return the cultureware to an incubator (+37 °C) and allow cells to detach for 5–45 minutes.
4. When the cells have completely detached from the cell cultureware carefully aspirate the cell suspension and transfer it to a centrifugation tube.

5. Spin down the cells for 3 minutes at 220 xg. Inhibition of the enzyme reaction is not required.
6. Discard the supernatant, resuspend the cells in the appropriate medium and seed the cells in new cell cultureware.

4 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 2–8°C.

5 Quality Control

All lots 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).

6 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.

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone: +49 551 308 0
www.sartorius.com

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote the other gender as well.

Copyright notice:

This instruction manual, including all of its components, is protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

© Sartorius Germany

Last updated:

02 | 2021

Printed in the EU on paper bleached
without chlorine. | MB
DIR: 2729369-000-00
Revision 01