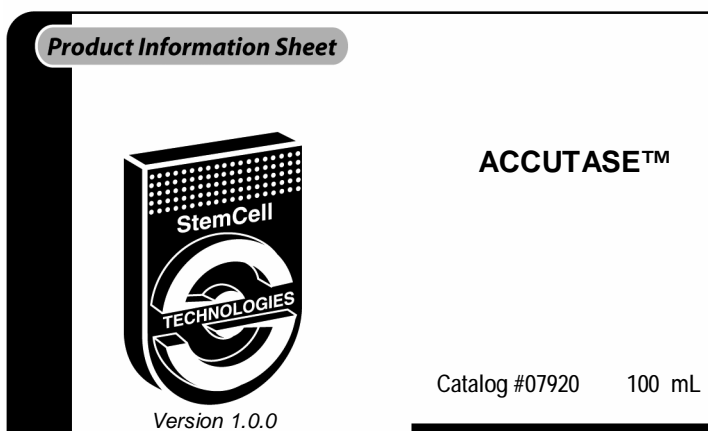


## PRODUCT DESCRIPTION:

A cell detachment solution of proteolytic and collagenolytic enzymes. Useful for the routine detachment of cells from standard tissue culture plasticware and adhesion coated plasticware. ACCUTASE™ does not contain mammalian or bacterial derived products.

Each lot of ACCUTASE™ is tested for Sterility (by USP membrane filtration method), enzymatic activity (tested with synthetic chromagenic tetrapeptides) and cell detachment from tissue culture plastic.

A Certificate of Analysis is available upon request.



## APPLICATIONS:

ACCUTASE™ has been shown to be efficiently detach or dissociate fibroblasts, neurospheres, keratinocytes, vascular endothelial cells, hepatocytes, vascular smooth muscle cells, hepatocyte progenitors, bone marrow stem cells, macrophages, 293 cells, L929 cells, immortalized mouse testicular germ cells, 3T3, Vero, COS, HeLa, NT2, MG63, M24 and A375 metastatic melanomas, gliomas U251 and D54 and T1080 fibrosarcoma cells.

## FORMAT:

1X ACCUTASE™ enzymes in Dulbecco's PBS (0.2 g/L KCl, 0.2 g/L KH<sub>2</sub>PO<sub>4</sub>, 8 g/L NaCl, and 1.15 g/L Na<sub>2</sub>HPO<sub>4</sub>) containing 0.5 mM EDTA•4Na and 3 mg/L Phenol Red.

## STABILITY AND STORAGE:

Stable when stored at -20°C. Refer to lot expiration date on bottle. Recommended storage temperature upon receipt is -20°C. After thawing, ACCUTASE™ may be stored for up to 2 months at 2 - 8°C. May be stored in smaller aliquots to avoid repeated freeze thaw cycles.

**DO NOT STORE AT ROOM TEMPERATURE (15 – 25°C).**

## PROCEDURE:

1. Thaw ACCUTASE™ in the refrigerator overnight, in a vessel of cool water or at room temperature (15 – 25°C). Do not thaw at 37°C.
2. Wash plate, flask or beads with sterile PBS.
3. Add ACCUTASE™ to culture dish or flask using aseptic procedures at 10 mL per 75 cm<sup>2</sup> surface area.
4. Return culture to 37°C incubator and allow cells to detach for about 5 - 10 minutes.
5. Count cells and passage as usual; no additional washes or enzyme inhibitors are required.

ACCUTASE™ is a registered trademark of Innovative Cell Technologies, Inc. San Diego, CA

**THIS REAGENT IS FOR RESEARCH ONLY.  
IT IS NOT TO BE ADMINISTERED TO HUMANS.**

## StemCell Technologies

In North America  
Tel: 1.604.877.0713  
Fax: 1.604.877.0704  
Toll Free Tel: 1.800.667.0322  
Toll Free Fax: 1.800.567.2899  
e-mail: info@stemcell.com  
www.stemcell.com

In the United Kingdom  
Tel: +44.(0).20.7691.3561  
Fax: +33.(0).4.76.18.99.63  
Toll Free within United Kingdom:  
Tel: 0800.731.27.14  
Fax: 0800.731.27.13  
e-mail: info@stemcellgb.com

In Europe  
Tel: +33.(0).4.76.04.75.30  
Fax: +33.(0).4.76.18.99.63  
e-mail: info@stemcellfrance.com

Revised:  
September 2007