



397PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# ProteCEASE™ -100, Plus EDTA

## Protease Inhibitor Cocktail for Large Volume

(Cat. # 786-336)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

*ProteCEASE™* -100, Plus EDTA protease inhibitor cocktail contains optimized concentrations of various protease inhibitors, which provide excellent inhibition of protease activities during protein purification. *ProteCEASE™* is a superior general protease inhibitor cocktail that is suitable for purification from mammalian, plant, bacteria and yeast samples. The cocktail contains both irreversible and reversible protease inhibitors to inhibit serine, cysteine and other proteases. EDTA is packed separately for inhibiting metalloproteases. At 1X concentration, *ProteCEASE™* inhibits over 90% of protease activities (e.g. Mouse Pancreas Extract 0.5mg/ml protein).

*ProteCEASE™* -100, Plus EDTA is designed for large volumes of extraction/ lysis buffers and is supplied as 10 vials per kit, with each vial suitable for 100ml buffer or extract solution. EDTA [0.5M, 100X] is also supplied with this kit as an optional metalloprotease inhibitor.

## ITEM(S) SUPPLIED (Cat. # 786-336)

Description	Size
<i>ProteCEASE™</i> -100	10 Vials
<i>ProteCEASE™</i> Buffer	2 x 2.5ml
0.5M [100X] EDTA Solution	2 x 5ml

## STORAGE CONDITIONS

The kit is shipped at ambient temperature. Upon arrival, store *ProteCEASE™* at 4°C. If stored and used as recommended, the product is stable for 1 year.

## INSTRUCTION FOR USE

*ProteCEASE™* Buffer will freeze at 4°C. Allow all components to warm to room temperature before using. Using a clean and dry pipette tip, add 0.5ml of *ProteCEASE™* Buffer to a vial of *ProteCEASE™* to reconstitute. Vortex briefly until completely dissolved. Close the *ProteCEASE™* Buffer cap tightly. Once reconstituted, the *ProteCEASE™* solution is stable for up to 1 week at 4°C.

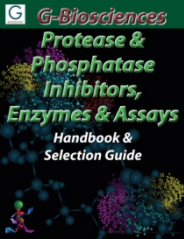
Add 0.5ml reconstituted *ProteCEASE™* -100 directly to 100ml extraction/ lysis buffer and mix it immediately for optimal protease inhibition. Once *ProteCEASE™* is added in the extraction/lysis buffer, use it within 20-30 minutes.

**Optional Protocols:**

- I. *If higher potency of protease inhibition is required, the amount of ProteCEASE™ may be doubled or sample volume reduced by half.*
- II. *If metalloproteases inhibition is needed, add 1ml of supplied 0.5M [100X] EDTA solution per 100 ml of extraction/ lysis buffer.*

**RELATED PRODUCTS**

Download our Protease & Phosphatase Inhibitors, Enzymes & Assays Handbook.



<http://info.gbiosciences.com/protease-phosphatase-inhibitors-enzymes-assay-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

Last saved: 3/5/2013 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)