



395PR

**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, EDTA Free

Protease Inhibitor Cocktail for Large Volume

(Cat. # 786-328)



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

## INTRODUCTION

ProteCEASE™ -100, EDTA Free 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 is EDTA free and contains both irreversible and reversible protease inhibitors to inhibit serine, cysteine and other proteases. At 1X concentration, ProteCEASE™ inhibits over 90% of protease activities (e.g. Mouse Pancreas Extract 0.5mg/ml protein).

ProteCEASE™ -100, EDTA Free 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.

## ITEM(S) SUPPLIED

Description	For 1000ml
ProteCEASE™ -100	10 Vials
ProteCEASE™ Buffer	2 x 2.5ml

## 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.

## PROTOCOL

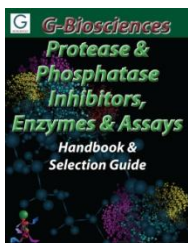
1. ProteCEASE™ Buffer will freeze at 4°C. Allow all components to warm to room temperature before using.
2. Using a clean and dry pipette tip, add 0.5ml of ProteCEASE™ Buffer to a vial of ProteCEASE™ -100 to reconstitute.
3. Vortex briefly until completely dissolved. Close the ProteCEASE™ Buffer cap tightly. Once reconstituted, the ProteCEASE™ solution is stable for up to 1 week at 4°C.
4. Add 0.5ml reconstituted ProteCEASE™ -100 directly to 100ml extraction/lysis buffer and mix immediately for optimal protease inhibition.
5. Once ProteCEASE™ is added in the extraction/lysis buffer; use it within 20-30 minutes.

**OPTIONAL:** *If higher potency of protease inhibition is required, the amount of ProteCEASE™ may be doubled or sample volume reduced by half.*

**NOTE:** *ProteCEASE™ -100, EDTe Free is an EDTA free preparation. If metalloproteases inhibition is needed, add appropriate volume of 0.5M EDTA; e.g. 1ml of 0.5M EDTA per 100 ml of sample or extraction/ lysis buffer or order our ProteCEASE™ -100, Plus EDTA kit, Cat # 786-336.*

## RELATED PRODUCTS

Download our Protease & Phosphatase Inhibitors, Enzyme & 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: 5/5/2016 IA



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