



401PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

ProteCEASE™ -50, Plus EDTA

Protease Inhibitor Cocktail for Large Volume

(Cat. # 786-334, 786-335)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

ProteCEASE™ -50, 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™ -50, Plus EDTA is designed for large volumes of extraction/ lysis buffers and is supplied as 10 vials /kit (Cat# 786-334) or 20 vials/kit (Cat# 786-335), with each vial suitable for 50ml sample. EDTA [0.5M, 100X] is also supplied with this kit as an optional metalloprotease inhibitor.

ITEM(S) SUPPLIED

Description	Cat# 786-334 [For 500ml]	Cat# 786-335 [For 1000ml]
ProteCEASE™ -50	10 Vials	20 Vials
ProteCEASE™ Buffer	2 x 2.5ml	4 x 2.5ml
0.5M [100X] EDTA Solution	5ml	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.

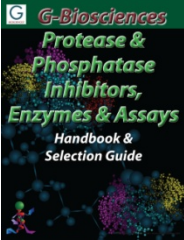
Add 0.5ml reconstituted ProteCEASE™ -50 directly to 50ml extraction/ lysis buffer and mix it immediately for optimal protease inhibition. Once ProteCEASE™ is added in the extraction/ lysis buffer, use it within 20-30 minutes.

OPTIONAL PROTOCOLS:

- If higher potency of protease inhibition is required, the amount of ProteCEASE™ may be doubled or sample volume reduced by half.
- If metalloproteases inhibition is needed, add 0.5ml of supplied 0.5M [100X] EDTA solution per 50 ml of extraction/ lysis buffer.

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 or contact us.

Last saved: 9/18/2012 CMH



www.GBiosciences.com