



259PR-01

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

A Geno Technology, Inc. (USA) brand name

# Mammalian ProteaseArrest™

Protease Inhibitor Cocktail  
Specific for Mammalian Cells

(Cat. # 786-331, 786-433)



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

## INTRODUCTION

Mammalian ProteaseArrest™ protease inhibitor cocktail contains optimized concentrations of various mammalian protease inhibitors, which provide excellent inhibition of protease activities during protein purification from mammalian cells and tissues. The cocktail contains both reversible and irreversible protease inhibitors to inhibit serine, cysteine and specific inhibitors for mammalian proteases, such as aminopeptidases. A 1X concentration of Mammalian ProteaseArrest™ inhibits over 90% of protease activities.

Optional EDTA is provided to inhibit metalloproteases. Some proteins require divalent cations ( $\text{Ca}^{2+}$ ,  $\text{Mg}^{2+}$  or  $\text{Mn}^{2+}$ ) for their biological activity and the presence of EDTA may be detrimental to the protein's activity. Furthermore, EDTA inhibits the purification of proteins by immobilized metal affinity chromatography; EDTA, therefore Mammalian ProteaseArrest™ is supplied with an optional EDTA solution.

## ITEM(S) SUPPLIED

Description	Cat. # 786-331	Cat. # 786-433
Mammalian ProteaseArrest™ [100X]	1ml	5ml
0.5M EDTA [100X]	2ml	5ml

## STORAGE CONDITIONS

The kit is shipped at ambient temperature. Store refrigerated at 4°C upon arrival, stable for 1 year.

**NOTE:** Once diluted, Mammalian ProteaseArrest™ is stable for 1-2 weeks at 4°C or for 4-6 weeks at -20° C.

## PROTOCOL

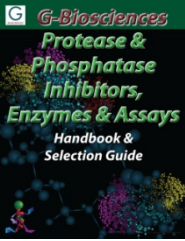
1. Allow the Mammalian ProteaseArrest™ to reach room temperature, then vortex or tap the vial before use.
2. Add 10µl 100X Mammalian ProteaseArrest™ to each 1ml of extraction buffer or extract for a 1X final concentration. Mix the solution thoroughly.

*Optional: For higher potency of protease inhibition, 20-30µl/ml Mammalian ProteaseArrest™ may be used.*

**NOTE:** For inhibition of metalloproteases, add 10µl of 100X EDTA per ml of extraction buffer or extract for a 1X final concentration. Not required if EDTA is already present.

## 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: 7/26/2012 CMH



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