



394PR-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

# Recom ProteaseArrest™

Recombinant Protein Protease Inhibitor Cocktail

(Cat. # 786-376, 786-436)



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

## INTRODUCTION

A broad range, 100X concentrated, ready-to-use protease inhibitor cocktail. Recom ProteaseArrest™ contains optimized concentrations of various bacterial protease inhibitors, which provide excellent inhibition of protease activities during recombinant protein purification from bacteria. The cocktail contains both reversible and irreversible protease inhibitors to inhibit metallo, serine and cysteine proteases. It also contains specific inhibitors for bacterial proteases, such as aspartic proteases and aminopeptidases. A 1X concentration of Bacterial ProteaseArrest™ inhibits over 90% of protease activities.

Recom ProteaseArrest™ cocktail does not use EDTA as its metalloprotease inhibitor as it would inhibit the activity of proteins that require divalent cations ( $\text{Ca}^{2+}$ ,  $\text{Mg}^{2+}$  or  $\text{Mn}^{2+}$ ) for their biological activity. In addition, EDTA would inhibit the purification of proteins using immobilized metal affinity chromatography (IMAC), for example His tagged or CBP tagged proteins. Recom ProteaseArrest™ cocktail is compatible with immobilized metal affinity chromatography.

## ITEM(S) SUPPLIED

Description	Cat. # 786-376	Cat. # 786-436
Recom ProteaseArrest™ [100X]	1ml	5ml

## STORAGE CONDITIONS

Shipped at ambient temperature. Store refrigerated at 4°C upon arrival, stable for 1 year.

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

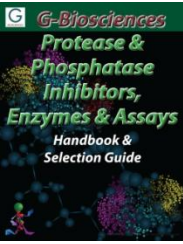
## PROTOCOL

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

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

## 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: 7/30/2014 CMH



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