



029PR-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

# Bacterial ProteaseArrest™

## Protease Inhibitor Cocktail for Bacteria

(Cat. # 786-330, 786-432)



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

## INTRODUCTION

Bacterial ProteaseArrest™ protease inhibitor cocktail contains optimized concentrations of various bacterial protease inhibitors, which provide excellent inhibition of protease activities during protein purification from bacteria. The cocktail contains both reversible and irreversible protease inhibitors to inhibit 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.

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. Therefore Bacterial ProteaseArrest™ is supplied with an optional EDTA solution.

## ITEM(S) SUPPLIED

Description	Cat. # 786-330	Cat. # 786-432
Bacterial ProteaseArrest™ [100X]	1ml	5ml
0.5M EDTA [100X]	2ml	5ml

## STORAGE CONDITIONS

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

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

## PROTOCOL

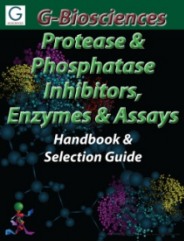
1. Allow the Bacterial ProteaseArrest™ to reach room temperature, then vortex or tap the vial before use.
2. Add 10µl 100X Bacterial ProteaseArrest™ to each 1ml of extraction buffer or extract for a 1X final concentration.
3. Mix the solution thoroughly.

**OPTIONAL:** For higher potency of protease inhibition, 20-30µl/ml Bacterial 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)