



465PR-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

# Yeast/Fungal ProteaseArrest™

Protease Inhibitor Cocktail Specific  
for Yeast and Fungi

(Cat. # 786-333, 786-435)



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

## INTRODUCTION

Yeast/Fungal ProteaseArrest™ protease inhibitor cocktail contains optimized concentrations of various yeast and fungal protease inhibitors, which provide excellent inhibition of protease activities during protein purification from yeast or fungi. The cocktail contains both reversible and irreversible protease inhibitors to inhibit serine, cysteine, metalloproteases and specific inhibitors for yeast or fungal proteases, such as aspartic proteases. At 1X concentration, Yeast/Fungal ProteaseArrest™ inhibits over 90% of protease activities.

## ITEM(S) SUPPLIED

Cat. #	786-333	786-435
Yeast/ Fungal ProteaseArrest™ [100X]	1ml	5ml

## STORAGE CONDITIONS

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

**Note:** *Once diluted, Yeast/Fungal ProteaseArrest™ is stable for 1-2 weeks at 4°C or for 4-6 weeks at -20°C.*

## INSTRUCTIONS FOR USE

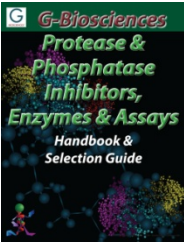
1. Allow the Yeast/Fungal ProteaseArrest™ to reach room temperature, then vortex or tap the vial before use.
2. Add 10µl 100X Yeast/Fungal ProteaseArrest™ to each 1ml of extraction buffer or extract for a 1X final concentration. Mix the solution thoroughly.

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

**Note:** Some proteins require divalent cations for their biological activity and the presence of metalloprotease inhibitor may be detrimental to the protein's activity and its purification by immobilized metal affinity chromatography.

## RELATED PRODUCTS

Download our Protease & Phosphatase: Inhibitors, Enzymes & 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)