



377PR-01

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

A Geno Technology, Inc. (USA) brand name

Plant ProteaseArrest™

Protease Inhibitor Cocktail Specific for Plant

(Cat. # 786-332, 786-434)



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

INTRODUCTION

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

ITEM(S) SUPPLIED

Description	Cat. # 786-332	Cat. # 786-434
Plant 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, Plant ProteaseArrest™ is stable for 1-2 weeks at 4°C or for 4-6 weeks at -20°C.

PROTOCOL

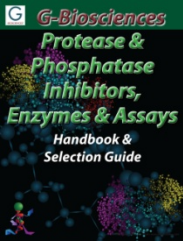
1. Allow the Plant ProteaseArrest™ to reach room temperature, then vortex or tap the vial before use.
2. Add 10µl 100X Plant ProteaseArrest™ to each 1ml of extraction buffer or extract for a 1X final concentration. Mix the solution thoroughly.

OPTIONAL: For higher potency of protease inhibition, 20-30µl/ml Plant 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, 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: 8/2/2012 CMH



www.GBiosciences.com