



363PR

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

A Geno Technology, Inc. (USA) brand name

PhosphataseArrest™ I

For Complete Inhibition of Phosphatase Activities

(Cat. # 786-450, 786-647, 786-782, 786-783 & 786-784)



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

INTRODUCTION

PhosphataseArrest™ I is a broad spectrum phosphatase inhibitor cocktail consisting of 5 phosphatase inhibitors that target all the phosphatase categories: serine/threonine (Ser/Thr) specific, tyrosine specific and dual specificity phosphatases.

PhosphataseArrest™ I is a 100X concentrated, ready-to-use solution. If samples have high phosphatase activity then the PhosphataseArrest™ I can be used at 2-3X concentrations to ensure complete inhibition. PhosphataseArrest™ I cocktail is ideal for inhibition in tissue extractions and cell lysis experiments and is compatible with most common protein assays. Due to the salt concentrations of the cocktail, PhosphataseArrest™ I is not suitable for direct IEF/2D studies.

ITEM(S) SUPPLIED

Description	Cat# 786-450	Cat# 786-647	Cat# 786-782	Cat# 786-783	Cat# 786-784
PhosphataseArrest™ I [100X]	1ml	24 x 100 µl	2ml	5ml	10ml

STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store PhosphataseArrest™ I at 4°C. If stored properly, it is stable for 1 year.

PREPARATION BEFORE USE

Vortex the PhosphataseArrest™ I vial briefly before removing the solution. If crystals develop after storage at 4°C, bring the solution to room temperature and vortex.

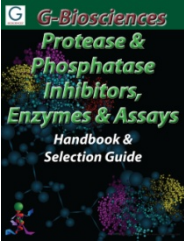
PROTOCOL

1. Add 10µl/ml PhosphataseArrest™ I directly in to an appropriate volume of extraction buffer or protein extract to a 1X final concentration.
2. For higher potency of phosphatase inhibition, add 20-30µl/ml PhosphataseArrest™ I to give a 2-3X final concentration.

NOTE: A protease inhibitor cocktail [ProteaseArrest™ (Cat. # 786-108)] and a Nuclease cocktail [FOCUS Nuclease™ (Cat. # 786-039F)] may be added in the same reaction mixture.

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: 6/24/2014 IA



www.GBiosciences.com