



367PR

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

A Geno Technology, Inc. (USA) brand name

Phosphatase Arrest™ III

(Cat. # 786-452)



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

INTRODUCTION

PhosphataseArrest™ III is a phosphatase inhibitor cocktail consisting of 3 phosphatase inhibitors that target alkaline and serine/threonine phosphatases. PhosphataseArrest™ III is a 100X concentrated, ready-to-use solution. If samples have high phosphatase activity then the PhosphataseArrest™ III can be used at 2-3X concentrations to ensure complete inhibition. PhosphataseArrest™ III is ideal for inhibition in tissue extractions and cell lysis experiments and is compatible with most common protein assays and IEF/2D studies.

ITEM(S) SUPPLIED (Cat. # 786-452)

Description	Size
PhosphataseArrest™ III [100X]	1.0ml

STORAGE CONDITIONS

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

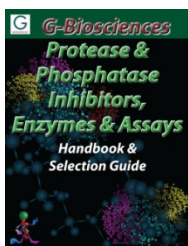
PROTOCOL

1. Vortex the PhosphataseArrest™ III vial briefly before removing the solution for use.
2. Add 10µl/ml PhosphataseArrest™ III directly into an appropriate volume of extraction buffer or protein extract to give 1X final concentration of PhosphataseArrest III.
3. For higher phosphatase inhibition, add 20-30µl/ml PhosphataseArrest III™ to give a 2-3X final concentration.

NOTE: A Protease inhibitor cocktail [FOCUS-ProteaseArrest™ (Cat# 786-108F)] 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.

This page is intentionally left blank



www.GBiosciences.com