



365PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# PhosphataseArrest™ II

(Cat. # 786-451)



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

## INTRODUCTION

PhosphataseArrest™ II is a phosphatase inhibitor cocktail consisting of 5 phosphatase inhibitors that target acid, alkaline and tyrosine phosphatases. PhosphataseArrest™ II is a 100X concentrated, ready-to-use solution. If samples have high phosphatase activity then the PhosphataseArrest™ II can be used at 2-3X concentrations to ensure complete inhibition. PhosphataseArrest™ II 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 PhosphataseArrest™ II cocktail, it is not suitable for direct IEF/2D studies.

## ITEM(S) SUPPLIED (Cat.# 786-451)

Description	Size
PhosphataseArrest™ II [100X]	1ml

## STORAGE CONDITIONS

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

## PREPARATION BEFORE USE

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

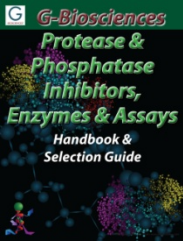
## PROTOCOL

1. Add 10µl/ml PhosphataseArrest™ II 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](http://www.GBiosciences.com) or contact us.

Last saved: 8/13/2012 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)