



157PR-02

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

A Geno Technology, Inc. (USA) brand name

FOCUS ProteaseArrest™

A 2D Electrophoresis Compatible Inhibitor Cocktail
For Complete Inhibition of Protease Activities

(Cat. # 786-108F)



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

INTRODUCTION

FOCUS ProteaseArrest™ is a protease inhibitor cocktail specifically developed for sample preparation for IEF/2D-studies. No EDTA is used, which allows optimal action of nuclease activity for removing nucleic acids from the samples. FOCUS ProteaseArrest™ contains both irreversible and reversible protease inhibitors and inhibits serine, cysteine and metalloproteases etc. Due to the optimized concentration of the various inhibitors, the FOCUS ProteaseArrest™ shows excellent inhibition of protease activities and is therefore suitable for the protection of proteins extracted from animal cells / tissues, plants, yeast and bacteria etc. It is compatible with our FOCUS Nuclease™ for removing nucleic acid contamination from samples.

FOCUS ProteaseArrest™ at 1X concentration in extraction buffer at pH 7-8 inhibits over 90% of protease activities.

ITEM(S) SUPPLIED (Cat. # 786-108F)

Description	Size
FOCUS ProteaseArrest™ [100X]	1ml

STORAGE CONDITIONS

It is shipped at ambient temperature. Store FOCUS-Protease-Arrest™ at -20°C upon arrival. If stored properly, it is stable for 1 year.

PROTOCOL

1. Allow the FOCUS Protease Arrest solution to warm to room temperature and briefly vortex or tap the vial to mix it completely before removing the solution for use.
2. Add FOCUS ProteaseArrest™ 10µl/ml directly in an appropriate volume of extraction buffer or protein extract to 1X final concentration.

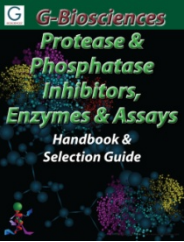
NOTE: When FOCUS ProteaseArrest™ is added in the buffer or extract, it is stable for 1-2 weeks at 4°C and 8-10 weeks at -20°C.

3. For higher potency of protease inhibition, add FOCUS ProteaseArrest™ 15µl/ml to give 1.5X final concentration.

NOTE: A Nuclease cocktail [e.g. our 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: 3/28/2014 IA



www.GBiosciences.com