



391PR

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

A Geno Technology, Inc. (USA) brand name

Protease Inhibitor Set

(Cat. # 786-207)



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

INTRODUCTION 3

ITEM(S) SUPPLIED 4

STORAGE CONDITIONS 4

PREPARATION BEFORE USE 5

Related Products 5

INTRODUCTION

The Protease Inhibitor Set contains twelve (12) ready to use individual protease inhibitors for characterization of protease activity. Each set contains 4-(2-Aminoethyl)-benzenesulfonyl-fluoride hydrochloride (AEBSF), ALLN (Calpain inhibitor), Antipain-dihydrochloride, Aprotinin, Bestatin, Chymostatin, E-64, EDTA-Na₂, Leupeptin, Pepstatin, Phosphoramidon, and PMSF in 100X concentration. The substrate specificities of various protease inhibitors are listed below. The 100X concentration of individual inhibitors in the set are carefully adjusted to inhibit >90% activity of individual protease in crude tissue extracts at 1X concentration. Various concentrations and/or combinations of protease inhibitors may be used to inhibit a broad spectrum of protease activity.

Our Protease Screening kit is supplied in conjunction with the Protease Inhibitor Set as Cat # 786-325 to determine protease activity. The Protease Screening Kit is based on a ready to use dye-labeled protein (as protease substrate). Proteases digest the substrate and release dye-labeled peptides. The absorbance of dye-labeled peptide is measured at 574nm for determination of protease activity. The kit is sufficient for 50 assays in microtiter format. The Protease Screening kit (Cat. # 786-137) may be purchased separately. Individual protease inhibitors are also offered in larger quantity for preparative scale uses.

ITEM(S) SUPPLIED (Cat. # 786-207)

Inhibitor	Specificity	Size
AEBSF [100X] 4-(2-Aminoethyl)-benzenesulfonyl-fluoride, hydrochloride	Irreversible inhibitor of serine proteases- trypsin, chymotrypsin, plasmin, plasma kallikrein, & thrombin	25µl
ALLN [100X] (N-Acetyl-Leu-Leu-Nle-CHO (Calpain Inhibitor I))	Cell-permeable, peptide aldehyde inhibitor of Calpain and other neutral cysteine proteases	25µl
Antipain-dihydrochloride [100X]	Inhibits papain, trypsin and to small extent plasmin	25µl
Aprotinin [100X]	Inhibits serine proteases- plasmin, kallikrein, trypsin, chymotrypsin	25µl
Bestatin [100X]	Inhibitor of amino-peptidases and other exopeptidases (does not inhibit carboxypeptidases)	25µl
Chymostatin [100X]	Specific inhibitor of α , β , γ , δ chymotrypsin	25µl
E-64 [100X] (N-(N-(L-3 Trans-carboxirane-2-carbonyl)-L-leucyl)-agmatine)	Inhibitor of papain and other cystein proteases like cathepsin B & L	25µl
EDTA-Na₂ [100X]	Inhibits metalloproteases	25µl
Leupeptin [100X]	Inhibits Serine & Cystein proteases e.g. Plasmin, Trypsin, Papain & Cathepsin B	25µl
Pepstatin [100X]	Inhibits Aspartic proteases, such as pepsin, renin, cathepsin D, chymosin	25µl
Phosphoramidon [100X]	Specifically inhibits thermolysin, collagenase & other metalloendoproteinases	25µl
PMSF [100X] Phenylmethylsulfonyl fluoride	Inhibits serine proteases; chymotrypsin, trypsin & thrombin, also inhibits cystein proteases such as papain (reversible by DTT)	25µl

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Upon arrival, store at 4°C. When stored & used properly, the kit components are stable for 1 year.

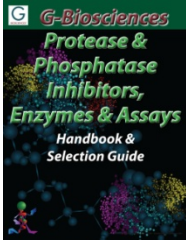
NOTE: Various concentrations and/or combinations of individual protease inhibitors may also be used for characterization of protease.

PREPARATION BEFORE USE

Allow the Protease Inhibitor solutions to warm to room temperature. Centrifuge the vials for a brief 10 seconds to collect the solution at the bottom of the vial. Vortex each vial for 15-20 seconds before use. Add appropriate volume of 100X inhibitor solution (s) in the samples to be tested to give a 1X or more (as needed) final concentration.

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: 7/26/2012 CMH

This page is intentionally left blank

This page is intentionally left blank



www.GBiosciences.com