

EZBlock™ Protease Inhibitor Cocktail, EDTA-Free

02/15

CATALOG NO: K272-1 1 ml
 K272-5 5 ml
 K272-set 5 x 1 ml

FORM: Liquid

STORAGE: -20 °C. Hygroscopic

DESCRIPTION: A cocktail of six broad specificity protease inhibitors (see table below) that can provide maximal protection to prevent degradation of proteins in mammalian tissue extracts as well as both mammalian and bacterial cell extracts by inhibition of various serine, cysteine, aspartic proteases and aminopeptidases. This cocktail is EDTA-free making it broadly applicable for any downstream application that is affected by the presence of metal chelators.

USE: 1 ml cocktail is recommended for 20 g tissue. A dilution of 1:200 is a good starting point but the working concentration should be optimized for each individual application.

HANDLING: Do not take internally. Wear gloves and mask when handling the product! Avoid contact by all modes of exposure.

Protease Inhibitor	M.W.	Target Protease	Concentration in the vial
AEBSF, HCl (Cat#1644)	239.7	Serine proteases	100 mM
Aprotinin (Cat#4690)	6512	Broad spectrum, serine proteases	80 µM
Bestatin (Cat#1733)	308.4	Aminopeptidases	5 mM
E-64 (Cat#1739)	357.4	Cysteine Proteases	1.5 mM
Leupeptin (Cat#1648)	475.6	Serine/Cysteine proteases	2 mM
Pepstatin A (Cat#1732)	685.9	HIV proteases, Aspartic proteases, Acid proteases	1 mM

RELATED PRODUCTS:

AEBSF, HCl (Cat. No. 1644-200, 1G)

Aprotinin (Cat. No. 4690-5, 100, 1000)

Bestatin (Cat. No. 1733-25, -100)

Calyculin A (Cat. No. 1562-025)

E-64 (Cat. No. 1739-5, 25)

EZBlock™ Protease Inhibitor Cocktail II (Cat. No. K277-set)

EZBlock™ Protease Inhibitor Cocktail III (Cat. No. K278-set)

EZBlock™ Protease Inhibitor Cocktail IV (Cat. No. K279-1, set)

EZBlock™ Protease Inhibitor Cocktail V, EDTA-Free (Cat. No. K290-1, set)

EZBlock™ Protease Inhibitor Cocktail VI, General Use (Cat. No. K291-1, set)

EZBlock™ Protease Inhibitor Cocktail VII (Cat. No. K292-1, set)

EZBlock™ Protease Inhibitor Cocktail VIII (Cat. No. K293-1, set)

EZBlock™ Protease Inhibitor Cocktail IX (Cat. No. K294-1, set)

EZBlock™ Universal Protease and Phosphatase Inhibitor Cocktail (Cat. No. K283-1, set)

EZBlock™ Universal Protease and Phosphatase Inhibitor Cocktail, EDTA-Free (Cat. No. K284-1, set)

EZBlock™ Protease Inhibitor Cocktail V, EDTA-Free (Cat. No. K290-1, set)

EZLys™ Bacterial Protein Extraction Reagent (Cat. No. 8001-100, 500)

EZLys™ Lysozyme, Human (Cat. No. 8005-1G, 5G)

Leupeptin, Hemisulfate (Cat. No. 1648-25, 50, 100)

Pepstatin A (Cat. No. 1732-25, 100)

Protease Inhibitor Cocktail (Cat. No. K271-500)

PMSF (Cat. No. 1548-5)

FOR RESEARCH USE ONLY! Not to be used on humans.