

EZBlock™ Protease Inhibitor Cocktail XIV

CATALOG NO: K223-1 1 vial
K223-set 5 x 1 vials

FORM: Lyophilized solid

STORAGE: -20 °C. Hygroscopic

DESCRIPTION: A cocktail of three protease inhibitors (shown in the table below) with specificity for the inhibition of metalloproteases, cysteine proteases including cathepsins and papain. This cocktail is recommended for inhibition of calcium-dependent endopeptidases, trypsin-like serine proteases, papain, and cysteine proteases. Reconstitute each vial with 1 ml D.I water to obtain a 100X concentrated stock solution.

Protease Inhibitor	Cat. No.	M.W.	Target Protease	Concentration (100X)
Antipain dihydrochloride	2834	677.62	Cathepsin A, B, and D as well as papain, trypsin and trypsin-like serine proteases	0.2 mg/ml
Leupeptin hemisulfate	1648	493.61	Serine/Cysteine proteases	0.05 mg/ml
EDTA, disodium	N/A	372.2	Metalloproteases	50 mg/ml

RELATED PRODUCTS:

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

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

Calyculin A (Cat. No. 1562-025)

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

EZBlock™ Phosphatase Inhibitor Cocktail I (Cat. No. K273-1, 1EA)

EZBlock™ Phosphatase Inhibitor Cocktail II (Cat. No. K275-1, 1EA)

EZBlock™ Phosphatase Inhibitor Cocktail III (Cat. No. K276-1, 1EA)

EZBlock™ Phosphatase Inhibitor Cocktail IV (Cat. No. K282-1, 1EA)

EZBlock™ Protease Inhibitor Cocktail EDTA-Free (Cat. No. K272-1, 5, 1EA)

EZBlock™ Protease Inhibitor Cocktail II (Cat. No. K277-1EA)

EZBlock™ Protease Inhibitor Cocktail III (Cat. No. K278-1EA)

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

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, 1EA)

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

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

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

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

PMSF (Cat. No. 1548-5)

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

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

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