

EZBlock™ Universal Protease and Phosphatase Inhibitor Cocktail

Components	K283-1	K283-set	Cap Color	Part Number
	1 ml	1 set		
Cocktail Solution	1 ml	5 x 1 ml	Red	K283-1-1
0.5 M EDTA Solution	1 ml	5 x 1 ml	Clear	K283-1-2

FORM: 10X Liquid (in H₂O)

STORAGE: -20 °C. Hygroscopic

DESCRIPTION: A broad-spectrum cocktail (component part# K283-1-1) of four protease inhibitors and four phosphatase inhibitors (see table below). This cocktail is suitable for use with cell lysates as well as tissue extracts from different species. EDTA, Disodium (0.5 M solution) is provided in a separate vial (component part# K283-1-2)

USE: Recommended dilution: 1:9. A 1X solution 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.

Inhibitor	M.W.	Target Phosphatase/Phosphatase	Concentration in the vial (10X)	Component Part#
Aprotinin (Cat#4690)	6512	Serine proteases and esterases	80 µM	K283-1-1
Bestatin (Cat#1733)	308.4	Aminopeptidases	5 mM	
E-64 (Cat#1739)	357.4	Cysteine proteases	1.5 mM	
Leupeptin (Cat#1648)	475.6	Cysteine/Trypsin-like proteases	2 mM	
β-Glycerophosphate	306.1	Ser/Thr Phosphatases	10 mM	
Sodium Fluoride	42.0	Acid Phosphatases and Ser/Thr phosphatases	50 mM	
Sodium Orthovanadate	183.9	Protein Tyrosine Phosphatase, Alkaline Phosphatase	1 mM	
Sodium Pyrophosphate	446.1	Ser/Thr phosphatases	10 mM	
EDTA, Disodium	372.2	Metalloproteases	0.5 M	K283-1-2

RELATED PRODUCTS:

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

BCA Protein Quantitation Kit (Cat. No. K812-1000)

Bradford Protein Quantitation Kit (Cat. No. K810-1000)

Calyculin A (Cat. No. 1562-025)

EZBlock™ Phosphatase Inhibitor Cocktail I (50X) (Cat. No. K273-1, set)

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

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

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 EDTA-Free (Cat. No. K272-1, 5, 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, EDTA-free (Cat. No. K284-1, set)

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

Nafamostat Mesylate (Cat. No. 1760-10, -50)

Okadaic Acid (Cat. No. 1543-025)

Okadaic Acid, Ammonium Salt (Cat. No. 1766-025)

Okadaic Acid, Potassium Salt (Cat. No. 1765-025)

Okadaic Acid, Sodium Salt (Cat. No. 1764-025)

FOR RESEARCH USE ONLY!

rev. 06/16



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.