

# EZBlock™ Universal Protease and Phosphatase Inhibitor Cocktail, EDTA-Free

02/15

**CATALOG NO:** K284-1 1 ml  
K284-set 5 x 1 ml

**FORM:** 10X Liquid (in H<sub>2</sub>O)

**STORAGE:** -20 °C. Hygroscopic

**DESCRIPTION:** A broad-spectrum cocktail 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. This formulation is EDTA-free.

**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)
Aprotinin (Cat#4690)	6512	Serine proteases and esterases	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	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

## 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 (Cat. No. K283-1, 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)

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.**