

EZBlock™ Phosphatase Inhibitor Cocktail IV

02/15

CATALOG NO: K282-1 1 ml
K282-set 5 x 1 ml

FORM: Liquid

STORAGE: -20 °C. Hygroscopic

DESCRIPTION: A cocktail of six phosphatase inhibitors (see table below) that can inhibit both acid/alkaline phosphatases, Ser/Thr phosphatases and protein tyrosine phosphatases (PTPs). This cocktail is suitable for use with cell lysates as well as tissue extracts from different species.

USE: Recommended dilution: 1:99. 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.

Phosphatase Inhibitor	M.W.	Target Phosphatase	Concentration in the vial (100X)
β-Glycerophosphate	306.1	Ser/Thr Phosphatases	100 mM
Imidazole	68.1	Alkaline Phosphatases	200 mM
Sodium Fluoride	42.0	Acid Phosphatases and Ser/Thr phosphatases	100 mM
Sodium Molybdate	205.9	Acid Phosphatases	115 mM
Sodium Orthovanadate	183.9	Protein Tyrosine Phosphatase, Alkaline Phosphatase	100 mM
Sodium Tartrate, Dihydrate	230.1	Acid Phosphatases	400 mM

RELATED PRODUCTS:

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

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

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

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

Calyculin A (Cat. No. 1562-025)

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

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

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

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

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.