

EZBlock™ Phosphatase Inhibitor Cocktail III

02/15

CATALOG NO: K276-1 1 ml
K276-set 5 x 1 ml

FORM: Liquid

STORAGE: -20 °C. Hygroscopic

DESCRIPTION: A cocktail of three phosphatase inhibitors (see table below) that can inhibit Ser/Thr and alkaline phosphatases. This cocktail is suitable for preserving the phosphorylation state of proteins in cell lysates as well as tissue extracts from different species.

USE: Recommended dilution: 1:49. 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
(-)- <i>p</i> -Bromotetramisole Oxalate	373.2	Alkaline Phosphatases	2.5 mM
Cantharidin	196.2	Ser/Thr Phosphatases(PP2A)	500 μM
Calyculin A	1009.2	Ser/Thr Phosphatases(PP1 and PP2A)	1 μM

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 IV (Cat. No. K282-1, 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)

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.