

EZCell™ Cell Stimulation Cocktail (500X)

CATALOG NO:	K930-100	100 µl
	K930-set	(4 x 100 µl)
FORM:	Liquid	
STORAGE:	-20 °C	

DESCRIPTION: The Cell Stimulation Cocktail (500X) is a ready-to-use cocktail useful to induce activation of many cell types to produce cytokines. This cocktail can be used to induce cytokine production in *in vitro* cell culture for subsequent detection in immunoassays. Each vial contains Protein Kinase C activator Phorbol 12-Myristate 13-Acetate (PMA) (40.5 µM) and the calcium ionophore, Ionomycin (670 µM) formulated in DMSO (500X).

CONCENTRATION: 500X. Use at 1X (dilute 1:500 or to 2 µl/ml final concentration)

Product	Cat. No.	M.W.	Concentration (500X)
Phorbol 12-Myristate 13-Acetate (PMA)	1544	616.83	40.5 µM
Ionomycin	1565	709.02	670 µM

RELATED PRODUCTS:

EZCell™ Cell Stimulation Cocktail (500X) **(K930)**

EZCell™ Cell Stimulation & Protein Transport Inhibitor Cocktail (500X) **(K931)**

EZCell™ Protein Transport Inhibitor Cocktail (500X) **(K932)**

Brefeldin A **(1560)**

Ionomycin **(1565)**

Monensin Methyl Ester **(1808)**

Phorbol 12-Myristate 13-Acetate (PMA) **(1544)**

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