

EZCell™ Protein Transport Inhibitor Cocktail (500X)

CATALOG NO: K932-500 500 µl

FORM: Liquid

STORAGE: -20 °C

DESCRIPTION:

The EZCell™ Protein Transport Inhibitor Cocktail (500X) is a cocktail composed of of Brefeldin A and Monensin, formulated in DMSO. Useful in assays that require prevention of protein transport to the extracellular space is required. Addition of this cocktail to cells in culture inhibits the intracellular protein secretory/transport pathway resulting in the accumulation of secreted proteins in the lumen of the endoplasmic reticulum and in the Golgi apparatus. This accumulation can then be detected by intracellular staining and flow cytometric analysis or Western blotting.

CONCENTRATION:

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

Product	Cat. No.	M.W.	Concentration (500X)
Brefeldin A	1560	280.36	5.3 mM
Monensin Methyl Ester	1808	684.9	1 mM

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.