

Cell Activation Cocktail

The Cell Activation Cocktail is a pre-mixed reagent that includes PMA (phorbol 12-myristate-13-acetate) and ionomycin. The mixture does not include Brefeldin A. This reagent is used for T cell activation and increases cell surface expression of certain cell surface markers (e.g. CD69).

Materials Provided

Catalog Number:	92-1334
Quantity Supplied:	1 vial(s), 100 µl per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	N/A
Source:	Synthetic
Purity:	
Endotoxin Level:	N/A
Formulation:	Each vial of this cocktail contains a solution comprised of phorbol-12-myristate 13-acetate
Storage:	-70°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	500 X
Reconstitution Volume:	N/A
Reconstitution Solvent:	N/A

Handling Instructions:

Dilute cocktail 1:5 in 100% DMSO, then 1:10 in PBS + 0.1% BSA. Should be sufficient to activate 50 mL of cells of various types.

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com