

Forskolin

Forskolin is a labdane diterpene produced by the Indian Coleus plant, *Coleus forskohlii*. It is commonly used to raise levels of cyclic AMP (cAMP) associated with the study of cell physiology mechanisms, and it activates the enzyme adenylyl cyclase to increase intracellular levels of this second messenger, which is necessary for the proper biological response of cells to hormones and other extracellular signals.

Materials Provided

Catalog Number:	92-0005
Quantity Supplied:	2 vial(s), 0.25 mg 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:	410.5 g/mol
Source:	Synthetic
Purity:	≥ 98% by HPLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	5 mM
Reconstitution Volume:	0.120 mL
Reconstitution Solvent:	DMSO

Handling Instructions:

Protect from light and moisture

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