

Wortmannin

Wortmannin is a naturally occurring steroid metabolite that potently inhibits phosphatidylinositol 3-kinases (PI3Ks, IC₅₀=1-10nM). Also inhibits PLK1 (IC₅₀=24nM), PLK3 (IC₅₀=49nM), MTOR, DNAPK, PI4Ks, MLCK, and MAPK at higher concentrations. Helps enhance apoptotic effects by inhibiting the PI3K/Akt cascade and blocking the antiapoptotic effects of cytokines.

Materials Provided

Catalog Number:	92-1095
Quantity Supplied:	1 vial(s), 1 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:	428.4 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:	1.5 mM
Reconstitution Volume:	1.556 mL
Reconstitution Solvent:	DMSO

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