

PD0325901

PD0325901 is a synthetic orally active MEK1 and MEK2 inhibitor ($K_i=1.1$ and 0.79 , respectively) that research shows inhibits the growth of melanoma cell lines in-vitro and in-vivo. Also inhibits production of VEGF proangiogenic cytokines, and suppresses the phosphorylation of ERK and the RAS/RAF/MEK/ERK signaling pathway, which is frequently active in tumors.

Materials Provided

Catalog Number:	92-1184
Quantity Supplied:	1 vial(s), 5 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:	482.19 g/mol
Source:	Synthetic
Purity:	≥ 99%
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:	10 mM
Reconstitution Volume:	1.037 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