

CDDO Methyl Ester

CDDO Methyl Ester (bardoxolone methyl) is a semi-synthetic triterpenoid developed from the oleanolic acid scaffold that induces apoptosis and differentiation in tumor cells. Agonist of the KEAP1-Nrf2 pathway, thus inducing expression of antioxidant proteins and protection against oxidative damage from injury and inflammation. Also an antagonist of the NF- κ B pro-inflammatory transcription factor. Clinical trials are currently underway for type 2 diabetes treatment, due to its potent antidiabetic activity, but was confirmed to enhance cardiac disease during a phase 2 trial in October 2012.

Materials Provided

Catalog Number:	92-1147
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:	505.69 g/mol
Source:	Synthetic
Purity:	N/A
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	2-8°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	100 mM
Reconstitution Volume:	0.099 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