

## Embelin

Embelin is a natural cell-permeable benzoquinone inhibitor of the X-linked inhibitor of apoptosis (XIAP) protein (IC<sub>50</sub>=4.1μM), acting as an anti-tumor, anti-inflammatory, and analgesic. Activates caspases, blocks growth, and promotes apoptosis in tumor cells with high levels of XIAP. Also prevents XIAP- and caspase-dependent inflammation, and inhibits IKK, preventing NF-κB activation.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1202</b>
<b>Quantity Supplied:</b>	1 vial(s), 10 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

<b>Molecular Weight:</b>	294.39 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 98% by HPLC
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	2-8°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

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

<b>Stock Concentration:</b>	100 mM
<b>Reconstitution Volume:</b>	0.340 mL
<b>Reconstitution Solvent:</b>	DMSO

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)