

Camptothecin

Camptothecin is a cytotoxic quinoline plant alkaloid that inhibits DNA and RNA synthesis by specifically inhibiting the DNA topoisomerase I enzyme. Shows great anti-tumor potential in the clinic by inducing DNA breaks and apoptosis, but shows in-vivo cytotoxicity. Derivatives of camptothecin have been developed and are used in chemotherapy, such as toptecan and irinotecan.

Materials Provided

Catalog Number:	92-1146
Quantity Supplied:	1 vial(s), 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:	348.36 g/mol
Source:	Synthetic
Purity:	≥ 98%
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:	5 mM
Reconstitution Volume:	14.35 mL
Reconstitution Solvent:	DMSO

Handling Instructions:

Desiccate

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