

TH588

TH588 is a potent selective inhibitor of MTH1 (NUDT1). It can also induce DNA damage and triggers an ATM-p53-mediated death response and DNA repair in U2OS cells.

Materials Provided

Catalog Number:	92-1325
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:	295.17 g/mol
Source:	Synthetic
Purity:	>99% 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:	20 mg/mL
Reconstitution Volume:	0.250 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