

Trichostatin A

Trichostatin A (TSA) is a naturally occurring antibiotic that inhibits cell cycle growth of eukaryotes. Potent and selective inhibitor of histone deacetylases (HDAC, IC₅₀=70nM for HDAC1), thus inhibiting removal of acetyl groups from histones and resulting in hyperacetylation. Leads to chromatin relaxation and gene expression modulation.

Materials Provided

Catalog Number:	92-1246
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:	302.4 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:	1.653 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