

TO901317

TO901317 is a small molecule Liver X receptor (LXR) agonist, with EC₅₀ of ~50nM for LXR α and LXR β . This activation increases the efflux of cholesterol from enterocytes, inhibiting the absorption of cholesterol.

Materials Provided

Catalog Number:	92-1126
Quantity Supplied:	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

Molecular Weight:	481.3 g/mol
Source:	Synthetic
Purity:	≥ 98% by TLC
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:	2.077 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