

Troglitazone

Troglitazone is a synthetic antidiabetic and anti-inflammatory drug previously prescribed to patients with type 2 diabetes mellitus, but was withdrawn from the market in 2000 due to toxicity issues. Selective agonist of the peroxisome proliferator-activated receptor- γ (PPAR γ , EC50=555nM) with no activity for PPAR α or PPAR δ .

Materials Provided

Catalog Number:	92-1138
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:	441.54 g/mol
Source:	Synthetic
Purity:	\geq 98% by HPLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	\leq -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:	1.132 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