

GW7647

GW 7647 is a synthetic small molecule agonist of the peroxisome proliferator-activated receptor-alpha (PPAR α) transcription factor, which is involved in lipid homeostasis regulation. Binds PPAR α , PPAR γ , and PPAR δ with EC50 of 1-10nM, ~1 μ M, and 1-10 μ M respectively. Exhibits anti-inflammatory properties, reduces NO production in macrophages, and has lipid-lowering effects.

Materials Provided

Catalog Number:	92-1134
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:	502.75 g/mol
Source:	Synthetic
Purity:	≥ 98% by HPLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	23-25°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:	0.995 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