

GW9508

GW 9508 is a free fatty acid receptor (FFAR) agonist that potentiates glucose and KCl-stimulated insulin secretion in MIN6 cells. It is selective for FFAR1 (GPR40) and FFAR4 (GPR120). pEC50 values are 7.32 and 5.46 for FFAR1 and FFAR4, respectively. Developed by GlaxoSmithKline.

Materials Provided

Catalog Number:	92-1039
Quantity Supplied:	1 vial(s), 1 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:	347.4 g/mol
Source:	Synthetic
Purity:	≥ 98%
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:	100 mM
Reconstitution Volume:	0.029 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