

TC-O 9311

TC-O 9311 is a synthetic small molecule and potent agonist of the G protein-coupled receptor 139 (GPR139, EC50=39nM). Lacks activity against a selected group of 90 other diverse targets.

Materials Provided

Catalog Number:	92-1232
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:	378.89 g/mol
Source:	Synthetic
Purity:	≥ 98% HPLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	2-8°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	20 mM
Reconstitution Volume:	1.320 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