

MDL 29,951

MDL 29,951 is a synthetic small molecule selective agonist of the G protein-coupled receptor 17 (GPR17) that is active in intact cells. Selective over endogenous purinergic receptors. May be used in GPR17 studies and has a role in the maturation of oligodendrocytes.

Materials Provided

Catalog Number:	92-1264
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:	302.11 g/mol
Source:	Synthetic
Purity:	≥ 98%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	RT 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:	0.827 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