

A68930

A68930 is a synthetic molecule developed by Abbott that is selective for D1-like dopamine receptor over D2-like (EC50=2.1 and 3910nM, respectively). Produces antidepressant and anorectic effects in animals.

Materials Provided

Catalog Number:	92-1043
Quantity Supplied:	1 vial(s), 10 mg per vial
Reconstitution Buffer:	2 mL

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:	307.78 g/mol
Source:	Synthetic
Purity:	≥ 99%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	Desiccate at +4°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	25 mM
Reconstitution Volume:	1.3 mL
Reconstitution Solvent:	H2O (updated 2/26/16)

Handling Instructions:

Desiccate at +4°C

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