

2-CI-IB-MECA

2-CI-IB-MECA is a potent and selective agonist of A3 adenosine receptor (ADORA3, $K_i=0.33\text{nM}$). Selective over A1 and A2 receptors by 2500- and 1400-fold, respectively. ADORA3 has been implicated in cardioprotective functions during ischemia, neuroprotection and neurodegeneration, as well as cell proliferation and cell death mediation.

Materials Provided

Catalog Number:	92-1045
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:	544.74 g/mol
Source:	Synthetic
Purity:	$\geq 98\%$
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:	10 mM
Reconstitution Volume:	1.836 mL
Reconstitution Solvent:	DMSO

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