

R- α -Methylhistamine

(R)(-)- α -Methylhistamine is a potent histamine receptor agonist, particularly H3 (81, 185, and 270 times more potent than histamine for H1, H2, and H3 receptors, respectively). Also binds H4 with $K_i=23\text{nM}$. Inhibits histamine synthesis and release.

Materials Provided

Catalog Number:	92-1170
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:	198.09 g/mol
Source:	Synthetic
Purity:	$\geq 98\%$
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	23-25°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	50 mM
Reconstitution Volume:	0.505 mL
Reconstitution Solvent:	H ₂ O

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