

## Histamine

Histamine is a neurotransmitter that regulates inflammatory physiological functions in the gut. Stimulates gastric secretion, bronchial smooth muscle constriction, and vasodilation. Increases the permeability of capillaries for white blood cells and other proteins to engage foreign pathogens. Potent agonist of the G protein coupled histamine receptors H1-4. Released by mast cells and basophils, and is derived by enzymatic carboxylation of the amino acid histidine.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1169</b>
<b>Quantity Supplied:</b>	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

<b>Molecular Weight:</b>	111.15 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 97%
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	≤ -20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

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

<b>Stock Concentration:</b>	25 mM
<b>Reconstitution Volume:</b>	3.6 mL
<b>Reconstitution Solvent:</b>	H <sub>2</sub> O

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)