

## IBMX

IBMX, also referred to as 3-isobutyl-1-methylxanthine, is a competitive phosphodiesterase inhibitor which raises cAMP levels, activates PKA, inhibits TNF-alpha, synthesizes leukotrienes, and reduces a number of inflammation and innate immune mechanisms. IBMX also is known among the class of compounds to act as a nonselective adenosine receptor antagonist.

### Materials Provided

<b>Catalog Number:</b>	<b>92-0007</b>
<b>Quantity Supplied:</b>	1 vial(s), 55.55 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>	222.25 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 99% by HPLC
<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>	500 mM
<b>Reconstitution Volume:</b>	0.5 mL
<b>Reconstitution Solvent:</b>	DMSO

**Handling Instructions:**

It is recommended that IBMX be used at a final concentration of 500  $\mu$ M

***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)