

## 2-Methylthio-ATP

2-(Methylthio)adenosine 5'-triphosphate (2-methylthio-ATP) is an ATP analog that is a more selective agonist of the P2y purinergic receptor. Adenosine 5'-triphosphate is a nucleoside triphosphate that is key to intracellular energy transfer. It is responsible for transporting chemical energy within cells for metabolism, is an end product of phosphorylation, cellular respiration, and fermentation, and is used in many cellular processes.

### Materials Provided

|                           |                           |
|---------------------------|---------------------------|
| <b>Catalog Number:</b>    | <b>92-1214</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> | 641.2 g/mol                                     |
| <b>Source:</b>           | Synthetic                                       |
| <b>Purity:</b>           | ≥ 90%   |
| <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>    | 10 mM            |
| <b>Reconstitution Volume:</b>  | 1.56 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)