

ATP

Adenosine 5'-triphosphate (ATP) is a multifunctional nucleotide and a key component of energy storage and metabolism used in many cellular processes. Serves as a substrate for kinases and coenzymes for a wide array of enzymatic reactions in-vivo. Potent p2 purinoceptor agonist.

Materials Provided

Catalog Number:	92-1176
Quantity Supplied:	1 vial(s), 10 mg (re-packaged) 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:	551.14 g/mol
Source:	Synthetic
Purity:	≥ 99%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	100 mM
Reconstitution Volume:	181 µL with 10 mg vial
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