

2-(methylthio)-ADP

2-(Methylthio)adenosine 5'-diphosphate (2-methylthio-ADP) induces platelet aggregation and inhibits cAMP accumulation in platelets in-vitro. It's 3-5 times more active than ADP as an aggregating agent, and 150-200 times more active as a cAMP accumulation inhibitor. Potent agonist of the P2Y1, P2Y12, and P2Y13 purinoceptors (pEC50=8.29 and 9.05nM for P2Y1 and P2Y12, and EC50=19nM for P2Y13).

Materials Provided

Catalog Number:	92-1177
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:	539.24 g/mol
Source:	Synthetic
Purity:	≥ 89%
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:	50 mM
Reconstitution Volume:	0.185 mL
Reconstitution Solvent:	H2O

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