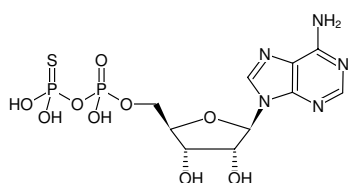


**ADPβS**

Adenosine-5'-(β-thio)-diphosphate, Lithium salt

Cat. No.	Amount
NU-433-5	5 mg
NU-433-25	25 mg



Structural formula of ADPβS

For general laboratory use.**Shipping:** shipped on dry ice**Storage Conditions:** store at -20 °C**Shelf Life:** 6 months after date of delivery**Molecular Formula:** C₁₀H₁₅N₅O₉P₂S (free acid)**Molecular Weight:** 443.26 g/mol (free acid)**Exact Mass:** 443.01 g/mol (free acid)**CAS#:** 35094-45-2 (free acid), 73536-95-5 (lithium salt)**Purity:** ≥ 85 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically.

Selected References:

Isfort *et al.* (2011) Real-time imaging reveals that P2Y2 and P2Y12 receptor agonists are not chemoattractants and macrophage chemotaxis to complement C5a is phosphatidylinositol 3-kinase (PI3K)- and p38 mitogen-activated protein kinase (MAPK)-independent. *J. Biol. Chem.* **286** (52):44776.

Chang *et al.* (2005) Nitric Oxide-dependent Allosteric Inhibitory Role of a Second Nucleotide Binding Site in Soluble Guanylyl Cyclase. *J. Biol. Chem.* **280** (12):11513.

Goody *et al.* (1971) Thiophosphate Analogs of Nucleoside Di- and Triphosphates. *J. Amer. Chem. Soc.* **93**:6252.