

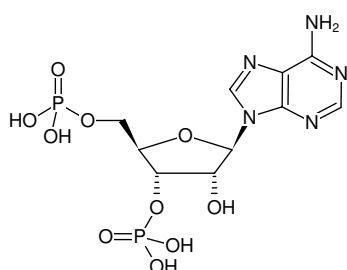


Adenosine-3',5'-bisphosphate

3',5'-pAp

Adenosine-3',5'-bisphosphate, Sodium salt

Cat. No.	Amount
NU-547S	100 µl (10 mM)
NU-547L	5 x 100 µl (10 mM)



Structural formula of Adenosine-3',5'-bisphosphate

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₀H₁₅N₅O₁₀P₂ (free acid)

Molecular Weight: 427.20 g/mol (free acid)

Exact Mass: 427.03 g/mol (free acid)

CAS#: 1053-73-2

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 259 nm, ε 15.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Jacobson *et al.* (2001) Ribose modified nucleosides and nucleotides as ligands for purine receptors. *Nucleos. Nucleot. Nucl.* **20** (4-7):333.

Jung *et al.* (2001) Platelet collagen receptor integrin alpha (2)beta (1) activation involves differential participation of ADP-receptor subtypes P2Y1 and P2Y12 but not intracellular calcium change. *Eur. J. Biochem.* **268** (12):3513.

Sheng *et al.* (2001) Bacterial expression, purification, and characterization of rat hydroxysteroid sulfotransferase STa. *Protein Expres. Purif.* **21** (1):235.

Leonidas *et al.* (2001) Mapping the ribonucleolytic active site of eosinophil-derived neurotoxin (EDN) - High resolution crystal structures of EDN complexes with adenylic nucleotide inhibitors. *J. Biol. Chem.* **276** (18):15009.

Kakuta *et al.* (1998) The sulfuryl transfer mechanism - Crystal structure of a vanadate complex of estrogen sulfotransferase and mutational analysis. *J. Biol. Chem.* **273** (42):27325.

Dichtl *et al.* (1997) Lithium toxicity in yeast is due to the inhibition of RNA processing enzymes. *EMBO J.* **16** (23):7184.

Mandon *et al.* (1994) Purification of the Golgi Adenosine 3'-phosphate 5'-phosphosulfate transporter, a homodimer within the membrane. *P. Natl. Acad. Sci. USA* **91** (22):10707.

Yang *et al.* (1994) Acyl carrier protein (ACP) import into chloroplasts - covalent modification by a stromal holoACP synthase is stimulated by exogenously added CoA and inhibited by Adenosine 3',5'-bisphosphate. *Eur. J. Biochem.* **224** (2):743.

Lee *et al.* (1992) Purification and identification of the heat-stable factor required for pregnenolone-binding protein-activity - evidence that the factor is Adenosine 3',5'-diphosphate. *J. Biol. Chem.* **267** (16):10982.

Chung *et al.* (1978) Conformation of Adenosine 2',5'-diphosphate and Adenosine 3',5'-diphosphate. *J. Carb.-Nucleos.-Nucl.* **5** (6):469.

Subbarao *et al.* (1978) Adenosine-3',5'-diphosphate and Coenzyme-A - effects on platelet-function. *Biochem. Pharmacol.* **27** (8):1193.

McDonald *et al.* (1978) Inhibition of catalytic activity of bovine pancreatic ribonuclease by Adenosine-3',5' diphosphate. *Fed. Proc.* **37** (6):1806.