

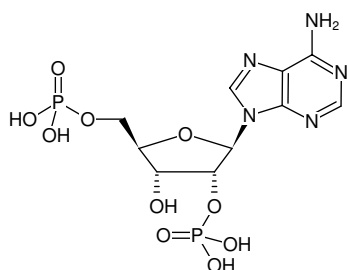


Adenosine-2',5'-bisphosphate

2',5'-pAp

Adenosine-2',5'-bisphosphate, Sodium salt

Cat. No.	Amount
NU-1113S	100 µl (10 mM)
NU-1113L	5 x 100 µl (10 mM)



Structural formula of Adenosine-2',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#: 3805-37-6

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:

Bierau *et al.* (2007) Pharmacogenetic significance of inosine triphosphatase. *Pharmacogenomics* **8** (9):1221.

Toth-Zsamboki *et al.* (2001) The P2Y₁ (1) receptor antagonist adenosine-2',5'-diphosphate non-selectively antagonizes the platelet P2X₁ (1) ion channel. *Thromb. Haemostasis* **86** (5):1338.

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.

Gu *et al.* (2000) Crystal structures of the complexes of trichosanthin with four substrate analogs and catalytic mechanism of RNA N-glycosidase. *Proteins* **39** (1):37.

Muruganandam *et al.* (1994) Isolation of nitric-oxide synthase from human platelets. *BBA-Gen. Subjects* **1200** (1):1.

Forstermann *et al.* (1991) Calmodulin-dependent endothelium-derived relaxing factor nitric-oxide synthase activity is present in the particulate and cytosolic fractions of bovine aortic endothelial-cells. *P. Natl. Acad. Sci. USA* **88** (5):1788.

Schmidt *et al.* (1991) Purification of a soluble isoform of guanylyl cyclase-activating-factor synthase. *P. Natl. Acad. Sci. USA* **88** (2):365.

Buitrago *et al.* (1986) Occurrence of Adenosine 2',5'-bisphosphate in rat-liver. *Biochim. Biophys. Acta* **881** (2):276.

Johnston *et al.* (1980) Interferon-induced (2'-5')oligoadenylate synthetase - adsorption to and assay on Adenosine 2',5'- diphosphate-sepharose. *Biochemistry-US* **19** (24):5580.

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

Hojeberg *et al.* (1976) Affinity chromatography and binding studies on immobilized Adenosine 5'-monophosphate and Adenosine 2',5'-bisphosphate of nicotinamide nucleotide transhydrogenase from *Pseudomonas-aeruginosa*. *Eur. J. Biochem.* **66** (3):467.