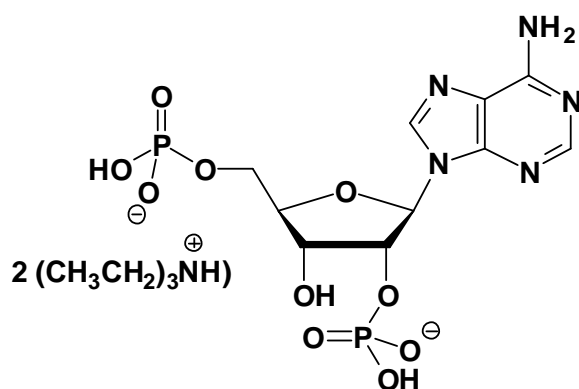


## Adenosine-2',5'-bisphosphate

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

Cat. No.	Amount
NU-1113S	100 Units
NU-1113L	500 Units



**Cat. No.:** NU-1113

**Molecular Formula:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>10</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 425.18 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

### For research use only!

1 unit = 1 µl of a 10 mM solution

### Selected References:

Toth-Zsomboki *et al.* (2001) The P2Y(1) receptor antagonist adenosine-2',5'-diphosphate non-selectively antagonizes the platelet P2X(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.