

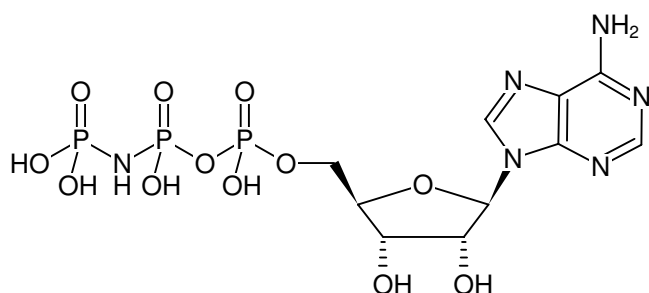
**AppNHp**

(AMPPNP)

AMPPNHP

Adenosine-5'-[( $\beta,\gamma$ )-imido]triphosphate, Tetralithium salt

Cat. No.	Amount
NU-407-10	10 mg
NU-407-50	50 mg



Structural formula of AppNHp

**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:** C<sub>10</sub>H<sub>17</sub>N<sub>6</sub>O<sub>12</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 506.19 g/mol (free acid)**Exact Mass:** 506.01 g/mol (free acid)**CAS#:** 72957-42-7**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:**  $\lambda_{\max}$  259 nm,  $\epsilon$  15.4 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Applications:**X-ray analysis<sup>[1, 2]</sup>

Hydrolyse studies<sup>[3, 4]</sup> Agonistic ligand, mainly for nucleoside receptor A<sub>1</sub>  
 Nucleosidephosphates stabilized against hydrolytic degradation can directly bind to nucleoside receptors.

**Specific Ligands:**Thymidylate kinase<sup>[1, 2]</sup>for P2Y<sub>2</sub> receptor<sup>[5]</sup>

**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.

When stored at -20 °C, product may hydrolyze, thereby forming AppNH<sub>2</sub> at a rate of up to 1 % per month!

**Selected References:**

[1] Segura-Pena *et al.* (2007) Quaternary structure change as a mechanism for the regulation of thymidine kinase 1-like enzymes. *Structure* **15**:1555.

[2] Ostermann *et al.* (2003) Structure of human thymidylate kinase in complex with prodrugs: Implications for the structure-based design of novel compounds. *Biochemistry* **42**:2568.

[3] Shimizu *et al.* (1997) Hydrolysis of AMPPNP by the motor domain of NCD, a kinesin-related protein. *Mol. Biol. Cell* **8**:1497.

[4] Suzuki *et al.* (1997) Hydrolysis of AMPPNP by the motor domain of ncd, a kinesin-related protein. *FEBS Lett.* **409** (1):29.

[5] Lazarowski *et al.* (1995) Pharmacological selectivity of cloned human P2U-purinoreceptor: potent activation by diadenosine tetrakisphosphate. *Br. J. Pharmacol.* **116** (1):1619.

Bodey *et al.* (2009) 9-Angström structure of a microtubule-bound mitotic motor. *J. Mol. Biol.* **388** (2):218.

Karow *et al.* (2009) A conformational change in the helicase core is necessary but not sufficient for RNA unwinding by the DEAD box helicase YxiN. *Nucleic Acids Research* **37** (13):4464.

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.

Menendez-Arias *et al.* (2005) Suppression of Multidrug-resistant HIV-1 Reverse Transcriptase Primer Unblocking Activity by a-Phosphate-modified Thymidine Analogues. *J. Mol. Biol.* **349** (3):451.

**AppNHp**

(AMPPNP)

AMPPNHP

Adenosine-5'-[( $\beta,\gamma$ )-imido]triphosphate, Tetralithium salt

Menz *et al.* (2001) The structure and nucleotide occupancy of bovine mitochondrial F-1-ATPase are not influenced by crystallisation at high concentrations of nucleotide. *FEBS Lett.* **494** (1):11.

Tereshko *et al.* (2001) Crystal structures of the catalytic domain of human protein kinase associated with apoptosis and tumor suppression. *Nat. Struct. Biol.* **8** (10):899.

Prodromou *et al.* (2000) The ATPase cycle of Hsp90 drives a molecular 'clamp' via transient dimerization of the N-terminal domains. *EMBO J.* **19** (16):4383.

Gil *et al.* (1999) N-Acetyl-L-glutamate kinase from Escherichia coli: cloning of the gene, purification and crystallization of the recombinant enzyme and preliminary X-ray analysis of the free and ligand-bound forms. *Acta Cryst. D* **55**:1350.

Shi *et al.* (1999) EDC cross-linking of actin and myosin S1. II. AMPPNP induces an allosteric transition from pre-hydrolysis state to posthydrolysis state. *Biophys. J.* **76** (1):A163.

Williams *et al.* (1986) Effects of purine nucleotides on the binding of [<sup>3</sup>H]cyclopentyladenosine to adenosine A1-receptors in rat brain membranes. *J. Neurochem.* **47** (1):88.