

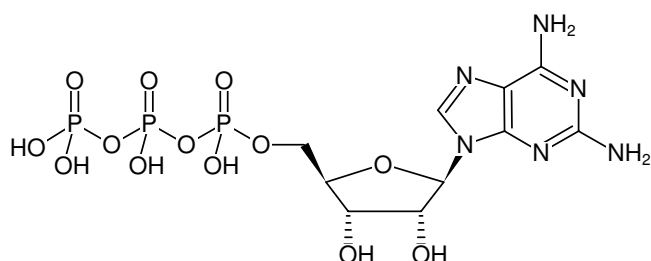
**2-Amino-ATP**

DAP-rTP

2-Aminoadenosine-5'-Triphosphate, Sodium salt

2,6-Diaminopurine-ribose-5'-Triphosphate

| Cat. No. | Amount             |
|----------|--------------------|
| NU-250S  | 10 µl (100 mM)     |
| NU-250L  | 5 x 10 µl (100 mM) |



Structural formula of 2-Amino-ATP

**Selected References:**

Lei *et al.* (2017) Structure-activity relationships of the ATP cofactor in ligase-catalysed oligonucleotide polymerisations. *Org. Biomol. Chem.* **15** (11):2349.

Ehteshami *et al.* (2013) Formation of a Quaternary Complex of HIV-1 Reverse Transcriptase with a Nucleotide-competing Inhibitor and Its ATP Enhancer. *J. Biol. Chem.* **288** (24):17336.

Höbartner *et al.* (2007) Engineering a selective small-molecule substrate binding site into a deoxyribozyme. *Angew. Chem. Int. Ed. Engl.* **468** (39):7420.

Wigle *et al.* (2007) Conformationally selective binding of nucleotide analogues to *Escherichia coli* RecA: a ligand-based analysis of the RecA ATP binding site. *Biochemistry* **45** (14):4502.

Ward *et al.* (1969) Fluorescence studies of nucleotides and polynucleotides. I. Formycin, 2-aminopurine riboside, 2,6-diaminopurine riboside, and their derivatives. *J. Biol. Chem.* **244** (5):1228.

**For research use only!****Shipping:** shipped on blue ice**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:** C<sub>10</sub>H<sub>17</sub>N<sub>6</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 522.20 g/mol (free acid)**Exact Mass:** 522.01 g/mol (free acid)**CAS#:** 18549-34-3 (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ<sub>max</sub> 279 nm, ε 9.89 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5), λ<sub>exc</sub> 280 nm, λ<sub>em</sub> 350 nm