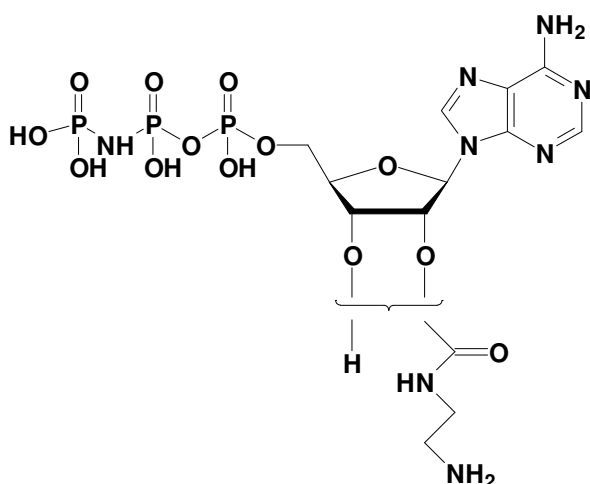


EDA-AppNHp (EDA-AMPPNP)

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, Sodium salt

Cat. No.	Amount
NU-810S	20 Units
NU-810L	100 Units



Cat. No.: NU-810

Molecular Formula: C₁₃H₂₃N₈O₁₃P₃ (free acid)

Molecular Weight: 592.29 (free acid)

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 6 months after date of delivery.

For research use only!

* 1 unit = 1 μ l of a 10 mM solution

Applications:

X-ray analysis^[1, 2]

Specific Ligands:

Kinesin^[3]

Hsp70^[4]

Selected References:

- [1] Terakado *et al.* (2010) Deleting two C-terminal α -helices is effective to crystallize the bacterial ABC transporter Escherichia coli MsbA complexed with AMP-PNP. *Acta Cryst. D* **D66**:319.
- [2] Pakhomova *et al.* (2008) Crystal structure of fosfomycin resistance kinase FomA for streptomyces wedmorensis. *J. Biol. Chem.* **283**:28518.
- [3] Sugata *et al.* (2009) Nucleotide-induced flexibility change in neck linkers of dimeric kinesin as detected by distance measurements using spin-labeling EPR. *J. Mol. Biol.* **386**:626.
- [4] Shida *et al.* (2010) Direct inter-subdomain interactions switch between the closed and open forms of the Hsp70 nucleotide binding domain in the nucleotide-free state. *Acta Cryst. D* **D66**:223.
- Mertens *et al.* (2012) Stepwise motion of a microcantilever driven by the hydrolysis of viral ATPases. *Nanotechnology* **23** (1):015501.
- Lansky *et al.* (2011) Studying kinesin's enzymatic cycle using a single-motor confocal motility assay, employing Förster resonance energy transfer. *Methods Mol. Biol.* **778**:19.
- Matamoras *et al.* (2005) Suppression of Multidrug-resistant HIV-1 Reverse Transcriptase Primer Unblocking Activity by α -Phosphate-modified Thymidine Analogues. *J. Mol. Biol.* **349**:451.