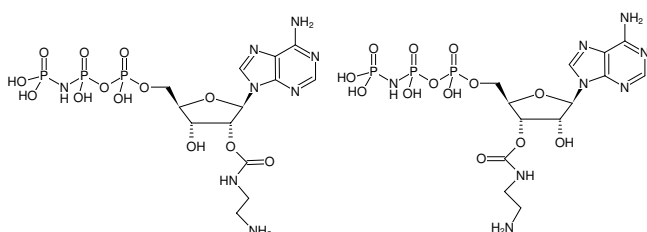


**EDA-AppNHp**

(EDA-AMPPNP)

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

Cat. No.	Amount
NU-810	500 μ l (10 mM)



Structural formula of EDA-AppNHp

For research use only!**Shipping:** shipped on dry ice**Storage Conditions:** store at -20 °C**Shelf Life:** 6 months after date of delivery**Molecular Formula:** C₁₃H₂₃N₈O₁₃P₃ (free acid)**Molecular Weight:** 592.29 g/mol (free acid)**Exact Mass:** 592.06 g/mol (free acid)**Purity:** \geq 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{max} 259 nm, ϵ 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Applications:**X-ray analysis^[1, 2]**Specific Ligands:**Kinesin^[3]Hsp70^[4]

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. When stored at -20 °C, product may hydrolyze, thereby forming EDA-AppNH₂ at a rate of up to 1 % per month!

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.