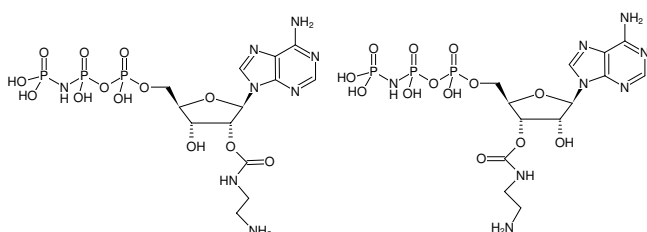


**EDA-AppNHp**

(EDA-AMPPNP)

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-[( $\beta,\gamma$ )-imido] triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-810	500 $\mu$ l (10 mM)



Structural formula of EDA-AppNHp

**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:** 6 months after date of delivery**Molecular Formula:** C<sub>13</sub>H<sub>23</sub>N<sub>8</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 592.29 g/mol (free acid)**Exact Mass:** 592.06 g/mol (free acid)**Purity:**  $\geq$  95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5  $\pm$  0.5**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>**Specific Ligands:**Kinesin<sup>[3]</sup>Hsp70<sup>[4]</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. When stored at -20 °C, product may hydrolyze, thereby forming EDA-AppNH<sub>2</sub> at a rate of up to 1 % per month!

**Selected References:**

[1] Terakado *et al.* (2010) Deleting two C-terminal  $\alpha$ -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  $\alpha$ -Phosphate-modified Thymidine Analogues. *J. Mol. Biol.* **349**:451.