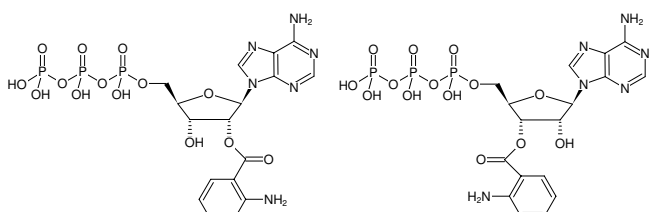


**Ant-ATP**

2'/3'-O-Anthraniloyl-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-891S	150 µl (10 mM)
NU-891L	5 x 150 µl (10 mM)



Structural formula of Ant-ATP

For general laboratory use.**Shipping:** shipped on gel packs**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:** C₁₇H₂₁N₆O₁₄P₃ (free acid)**Molecular Weight:** 626.30 g/mol (free acid)**Exact Mass:** 626.03 g/mol (free acid)**CAS#:** 289633-56-3**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 255 nm, ε 20.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 330 nm, λ_{em} 428 nm**Selected References:**

Yan *et al.* (2011) Role of protein conformational dynamics in the catalysis by 6-hydroxymethyl-7,8-dihydropterin pyrophosphokinase. *Protein Pept. Lett.* **18** (4):328.

Li *et al.* (2005) Is the critical role of loop 3 of Escherichia coli 6-hydroxymethyl-7,8-dihydropterin pyrophosphokinase in catalysis due to loop-3 residues arginine-84 and tryptophan-89? Site-directed mutagenesis, biochemical, and crystallographic studies. *Biochemistry* **44** (24):8590.

Shi *et al.* (2000) Dissecting the nucleotide binding properties of Escherichia coli 6-hydroxymethyl-7,8-dihydropterin pyrophosphokinase with fluorescent 3' (2')-o-anthraniloyladenine 5'-triphosphate. *Biochim. Biophys. Acta* **1478** (2):289.

Karasaki *et al.* (1987) Inhibition of single and double-stranded DNA-dependent ATPase of RecA protein by ATP ribose-modified analogs. *Journal of UOEH* **9** (2):141.