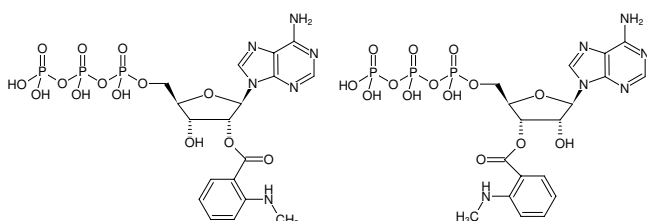


**Mant-ATP**

2'/3'-O-(N-Methyl-anthraniloyl)-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-202S	150 µl (10 mM)
NU-202L	5 x 150 µl (10 mM)



Structural formula of Mant-ATP

For research use only!**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:** C₁₈H₂₃N₆O₁₄P₃ (free acid)**Molecular Weight:** 640.33 g/mol (free acid)**Exact Mass:** 640.05 g/mol (free acid)**CAS#:** 151481-86-6**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} 255/355 nm, ε 23.3/5.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 355 nm, λ_{em} 448 nm**Applications:**Conformational dynamic: DnaB/C-protein^[1], Csk^[2]Inhibition of AC-isoforms^[3]Fluorescence stop-flow kinetics: PKA^[4]FRET: AC^[5], myosin V^[6]**Specific Ligands:**DnaC-protein^[7]Factor Rho^[8]**Selected References:**

- [1] Roychowdhury *et al.* (2009) Interactions of Escherichia coli DnaB-DnaC protein complex with nucleotide cofactors. 1. Allosteric conformational transitions of the complex. *Biochemistry* **48**:6712.
- [2] Shaffer *et al.* (2001) Nucleotide release and associated conformational changes regulate function in the COOH-terminal Src kinase, Csk. *Biochemistry* **40**:11149.
- [3] Gille *et al.* (2004) Differential inhibition of adenylyl cyclase isoforms and soluble guanylyl cyclase by purine and pyrimidine nucleotides. *J. Biol. Chem.* **279**:19955.
- [4] Ni *et al.* (2000) Insights into nucleotide binding in protein kinase A using fluorescent adenosine derivatives. *Protein Science* **9**:1818.
- [5] Goettle *et al.* (2007) Molecular analysis of the interaction of Bordetella pertussis adenylyl cyclase with fluorescent nucleotides. *Molecular Pharmacology* **72** (3):526.
- [6] Sun *et al.* (2006) Dynamics of the upper 50-kDa domain of myosin V examined with fluorescence resonance energy transfer. *J. Biol. Chem.* **281**:5711.
- [7] Galletto *et al.* (2005) The nucleotide-binding site of the Escherichia coli DnaC protein: Molecular topography of DnaC protein-nucleotide cofactor complex. *Cell Biochem. And Biophys.* **43**:331.
- [8] Jeong *et al.* (2004) Nucleotide binding induces conformational changes in Escherichia coli transcription termination factor Rho. *J. Biol. Chem.* **279**:1837.
- Seifert *et al.* (2012) Inhibitors of membranous adenylyl cyclases. *Trends Pharmacol. Sci.* **33** (2):64.
- Wang *et al.* (2011) Charge isomers of myelin basic protein: structure and interactions with membranes, nucleotide analogues, and calmodulin. *PLoS One.* **6** (5):e19915.
- Pinto *et al.* (2011) Structure-activity relationships for the interactions of 2'- and 3'- (O)- (N-methyl)anthraniloyl-substituted purine and pyrimidine nucleotides with mammalian adenylyl cyclases. *Biochem. Pharmacol.* **82** (4):358.



Mant-ATP

2'/3'-O-(N-Methyl-anthraniloyl)-adenosine-5'-triphosphate, Triethylammonium salt

Biophys. Acta **742**:496.

Hübner *et al.* (2011) Effect of MANT-nucleotides on L-type calcium currents in murine cardiomyocytes. *J. Naunyn Schmiedebergs Arch. Pharmacol.* **383** (4):573.

Spangler *et al.* (2011) Interaction of the diguanylate cyclase YdeH of *Escherichia coli* with 2', (3')-substituted purine and pyrimidine nucleotides. *J. Pharmacol. Exp. Ther.* **336** (1):234.

Erdorf *et al.* (2011) Pharmacological characterization of adenylyl cyclase isoforms in rabbit kidney membranes. *J. Naunyn Schmiedebergs Arch. Pharmacol.* **383** (4):357.

Goettle *et al.* (2010) Cytidylyl and uridylyl cyclase activity of bacillus anthracis edema factor and *Bordetella pertussis* CyaA. *Biochemistry* **49** (26):5494.

Chen *et al.* (2009) Structural Basis of Mechanochemical Coupling in a Hexameric Molecular Motor. *J. Biol. Chem.* **283** (6):3607.

Kainov *et al.* (2008) ADP but Not Pi Dissociation Contributes to Rate Limitation for *Escherichia coli* Rho. *J. Biol. Chem.* **284** (49):33773.

Kambach *et al.* (2007) Human OLA1 Defines an ATPase subfamily in the Obg Family of GTP-binding proteins. *J. Biol. Chem.* **282** (27):19928.

Scott *et al.* (2006) Chain-Terminating Dinucleoside Tetraphosphates are substrates for DNA Polymerization by Human Immunodeficiency Virus Type 1 Reverse Transcriptase with Increased Activity against Thymidine Analogue-Resistant Mutants. *AAC* **50** (11):3607.

Lísal *et al.* (2005) Cooperative Mechanism of RNA Packaging Motor. *J. Biol. Chem.* **280** (24):23157.

Tung-Chung Mou *et al.* (2005) Structural Basis for the Inhibition of Mammalian Membrane Adenylyl Cyclase by 2' (3')-O- (N-Methylanthraniloyl)-guanosine 5'-Triphosphate. *J. Biol. Chem.* **280** (8):7253.

Fujita *et al.* (1999) Fluorescence changes of a label attached near the myosin active site on nucleotide binding in rat skeletal muscle fibres. *J. Physiol.* **515**:869.

Thoenges *et al.* (1999) Tight binding of bulky fluorescent derivatives of adenosine to the low affinity E2ATP site leads to inhibition of Na⁺/K⁺-ATPase. Analysis of structural requirements of fluorescent ATP derivatives with a Koshland-Nemethy-Filmer model of two interacting ATP sites. *J. Biol. Chem.* **274**:1971.

Cheng *et al.* (1998) Interaction of mant-adenosine nucleotides and magnesium with kinesin. *Biochemistry* **37**:5288.

Mocz *et al.* (1998) Probing the nucleotide binding sites of axonemal dynein with the fluorescent nucleotide analogue 2' (3')-O- (-N-Methylanthraniloyl)-adenosine 5'-triphosphate. *Biochemistry* **37**:9862.

Jameson *et al.* (1997) Fluorescent nucleotide analogs: synthesis and applications. *Methods Enzymol.* **278**:363.

Hiratsuka (1983) New ribose-modified fluorescent analogs of adenine and guanine nucleotides available as substrates for various enzymes. *Biochim.*