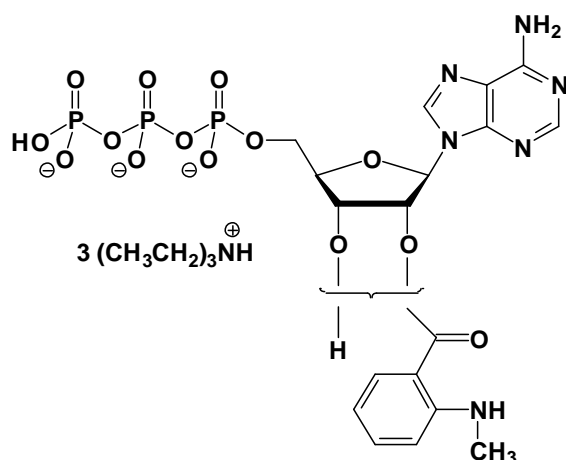


mant-ATP

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

Cat. No.	Amount
NU-202S	150 Units
NU-202L	750 Units



Cat. No.: NU-202

Molecular Formula: C₁₈H₂₀N₆O₁₄P₃ (Anion)

Molecular Weight: 637.30 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{\max} 255/355 nm; ϵ 23300/5800; λ_{exc} 355nm;
 λ_{em} 448 nm

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 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution

Selected References:

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

Gottle *et al.* (2007) Molecular analysis of the interaction of Bordetella pertussis adenylyl cyclase with fluorescent nucleotides. *Molecular Pharmacology Fast Forward*, June 6, 2007; DOI: 10.1124/mol.107.034413 (published online).

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.

Sprang *et al.* (2006) Broad and Specificity of Mammalian Adenylyl Cyclase for Interaction with 2',3'-Substituted Purine- and Pyrimidine Nucleotide Inhibitors. *Mol. Pharmacol.* **70**:878.

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.

Booth *et al.* (2000) Analysis of the properties of the N-terminal nucleotide-binding domain of human P-glycoprotein. *Biochemistry* **39**:5518.

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.

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.

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. Biophys. Acta* **742**:496.