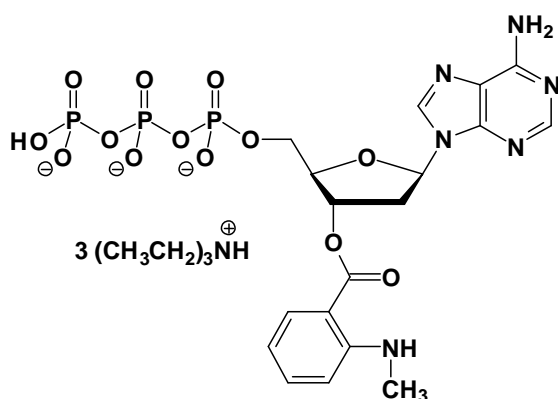


mant-dATP

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

Cat. No.	Amount
NU-203S	50 Units
NU-203L	250 Units



Cat. No.: NU-203

Molecular Formula: C₁₈H₂₀N₆O₁₃P₃ (Anion)

Molecular Weight: 621.31 (Anion)

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

Spectroscopic properties:

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

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).

Tuma *et al.* (2005) Cooperative Mechanism of RNA Packaging Motor. *J. Biol. Chem.* **280** (249): 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.

Churchich (1997) Conformational changes at the nucleotide binding of GroEL induced by binding of protein substrates - Luminescence studies. *J. Biol. Chem.* **272** (32):19645.

Bruce *et al.* (1997) Characterization of the molecular-chaperone function of the heat-shock-cognate-70-interacting protein. *Eur. J. Biochem.* **245** (3):738.

Schultz *et al.* (1997) Structural and catalytic properties of CMP kinase from *Bacillus subtilis*: A comparative analysis with the homologous enzyme from *Escherichia coli*. *Arch. Biochem. Biophys.* **340** (1):144.

Churchich (1995) Binding of a fluorescent nucleotide analog to hsc70 - the effect of peptide protein interactions on the luminescence properties of the probe. *Eur. J. Biochem.* **231** (3):736.

Bujalowski *et al.* (1994) Structural characteristics of the nucleotide-binding site of *Escherichia coli* primary replicative helicase DNAb protein - studies with ribose and base-modified fluorescent nucleotide analogs. *Biochemistry-US* **33** (15):4682.

Pineda *et al.* (1993) Interaction of ATP analogs with yeast 3-phosphoglycerate kinase - affinity labeling of the hinge region. *Eur. J. Biochem.* **212** (3):719.

Sarfati *et al.* (1991) New fluorescent and photoactivable analogs acting on nucleotide binding enzymes. *Tetrahedron Lett.* **32** (36):4699.