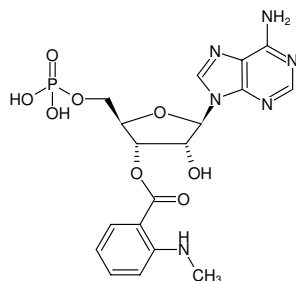
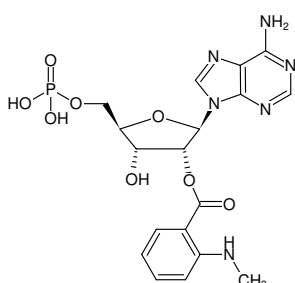


**Mant-AMP**

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

Cat. No.	Amount
NU-236S	150 µl (10 mM)
NU-236L	5 x 150 µl (10 mM)



Structural formula of Mant-AMP

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:** C₁₈H₂₁N₆O₈P (free acid)**Molecular Weight:** 480.37 g/mol (free acid)**Exact Mass:** 480.12 g/mol (free acid)**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:**Stopped-flow kinetic studies^[1, 2]Analysis of conformational changes^[1]**Specific Ligands:**Pyrophosphatase mtCBS^[1]Helicase DnaB^[2]**Selected References:**

[1] Jaemsen *et al.* (2010) Nucleotide- and substrate-induced conformational transitions in the CBS domain-containing pyrophosphatase of *Moorella thermoacetica*. *Biochemistry* **49**:1005.

[2] Bujalowski *et al.* (2000) Kinetic mechanism of nucleotide cofactor binding to *Escherichia coli* replicative helicase DnaB protein. Stopped-flow kinetic studies using fluorescent, ribose-, and base-modified nucleotide analogues. *Biochemistry* **39**:2106.

Wang *et al.* (2011) Charge isomers of myelin basic protein: structure and interactions with membranes, nucleotide analogues, and calmodulin. *PLoS One*. **6** (5):e19915.

Lísal *et al.* (2008) Cooperative Mechanism of RNA Packaging Motor. *The Journal of biological chemistry* **280** (24):23157.

Hiratsuka *et al.* (1983) New Ribose-modified Fluorescent Analogs of Adenine and Guanine Nucleotides Available as Substrates for Various Enzymes. *Biochim Biophys Acta* **742**:496.