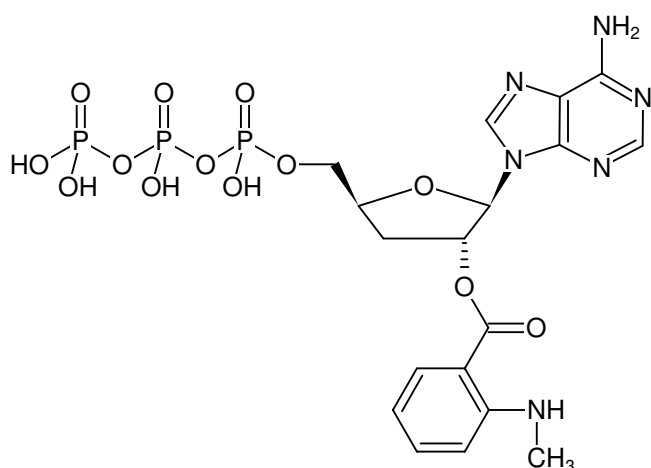


**2'-Mant-3'-dATP**

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

Cat. No.	Amount
NU-223S	20 µl (10 mM)
NU-223L	5 x 20 µl (10 mM)



Structural formula of 2'-Mant-3'-dATP

**For research use only!****Shipping:** shipped on blue ice**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:** C<sub>18</sub>H<sub>23</sub>N<sub>6</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 624.33 g/mol (free acid)**Exact Mass:** 624.05 g/mol (free acid)**CAS#:** 313378-46-0**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<sub>max</sub> 255/355 nm, ε 23.3/5.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5), λ<sub>exc</sub> 355 nm, λ<sub>em</sub> 448 nm**Selected References:**Suryanarayana *et al.* (2009) Distinct interactions of 2'- and 3'-O-(N-methyl)anthraniloyl-isomers of ATP and GTP with the adenyl cyclase toxin of *Bacillus anthracis*, edema factor. *Biochemical Pharmacology* **78** (3):224.Kainov *et al.* (2008) Structural Basis of Mechanochemical Coupling in a Hexameric Molecular Motor. *J. Biol. Chem.* **283** (6):3607.Goettle *et al.* (2007) Molecular analysis of the interaction of Bordetella pertussis adenyl cyclase with fluorescent nucleotides. *Molecular Pharmacology* **72** (3):526.Lisal *et al.* (2005) Cooperative Mechanism of RNA Packaging Motor. *J. Biol. Chem.* **280** (24):23157.Tuma *et al.* (2005) Cooperative Mechanism of RNA Packaging Motor. *J. Biol. Chem.* **280** (249):23157.