

**Mant-dADP**

2'-deoxy-3'-mant-ADP

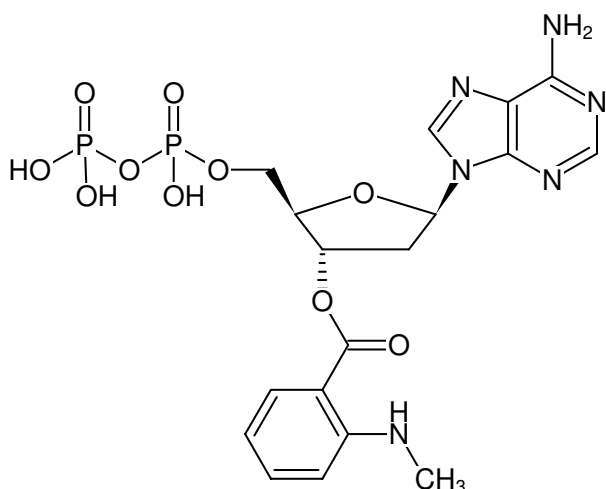
3'-O-(N-Methyl-anthraniloyl)-2'-deoxyadenosine-5'-diphosphate, Triethylammonium salt

Cat. No.	Amount
NU-238S	30 µl (10 mM)
NU-238L	5 x 30 µl (10 mM)

Selected References:

Hackney *et al.* (2016) Nucleotide-free kinesin motor domains reversibly convert to an inactive conformation with characteristics of a molten globule. *Arch. Biochem. Biophys.* **608**:42.

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* **33** (15):4682.



Structural formula of Mant-dADP

For general laboratory use.**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:** C₁₈H₂₂N₆O₁₀P₂ (free acid)**Molecular Weight:** 544.35 g/mol (free acid)**Exact Mass:** 544.09 g/mol (free acid)**CAS#:** 91828-83-0**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_{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