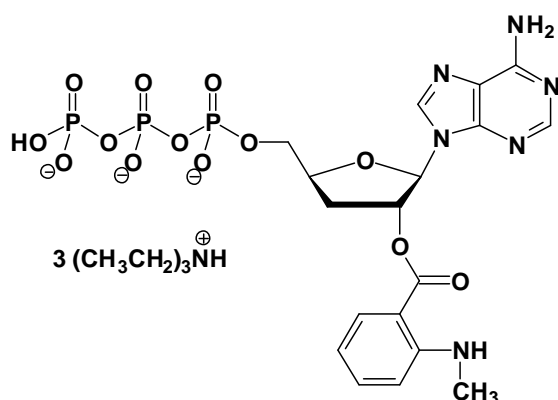


2'-Mant-3'-dATP

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

Cat. No.	Amount
NU-223S	20 Units
NU-223L	100 Units



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

Cat. No.: NU-223

Molecular Formula: C₁₈H₂₃N₆O₁₃P₃ (Anion)

Molecular Weight: 621.31 (Anion)

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

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 µl of a 10 mM solution