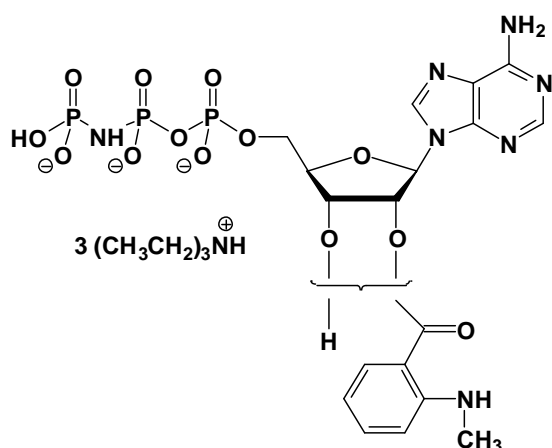


mant-AppNHp (mant-AMPPNP)

2'/3'-O-(N-Methyl-anthraniloyl)-Adenosine-5'-[(β,γ)-imido]
triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-214S	10 Units
NU-214L	50 Units



Cat. No.: NU-214

Molecular Formula: C₁₈H₂₁N₇O₁₃P₃ (Anion)

Molecular Weight: 636.32 (Anion)

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

Spectroscopic properties:

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

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 6 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution

mant-AppNHp (mant-AMPPNP)

2'/3'-O-(N-Methyl-anthraniloyl)-Adenosine-5'-[(β,γ)-imido]
triphosphate, Triethylammonium salt

Selected References:

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.

Jezewska *et al.* (1996) Interactions of Escherichia coli primary replicative helicase DnaB protein with nucleotide cofactors. *Biophys. J.* **71**:2075.

Moore *et al.* (1994) Kinetic mechanism of adenine nucleotide binding to and hydrolysis by the Escherichia coli Rep monomer. 1. Use of fluorescent nucleotide analogues. *Biochemistry* **33**:14550.

Rosenfeld *et al.* (1994) Structural and kinetic studies of the 10 S \rightleftharpoons 6 S transition in smooth muscle myosin. *J. Biol. Chem.* **269**:30187.