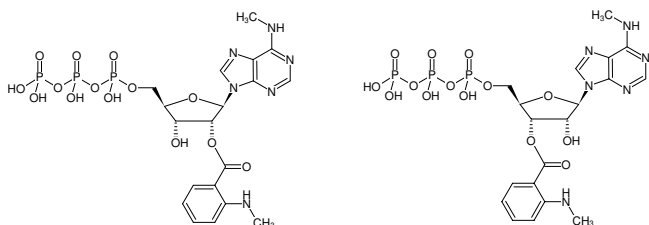




Mant-N⁶-Methyl-ATP

2'/3'-O-(N-Methyl-anthraniloyl)-N⁶-methyl-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-216S	30 µl (10 mM)
NU-216L	5 x 30 µl (10 mM)



Structural formula of Mant-N⁶-Methyl-ATP

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: C₁₉H₂₅N₆O₁₄P₃ (free acid)

Molecular Weight: 654.36 g/mol (free acid)

Exact Mass: 654.06 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