

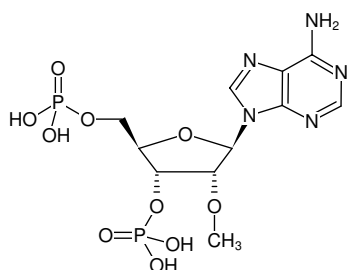


2'OMe-Adenosine-3',5'-bisphosphate

(2'OMe-pAp)

2'-O-Methyl-adenosine-3',5'-bisphosphate, Sodium salt

Cat. No.	Amount
NU-1114S	100 µl (10 mM)
NU-1114L	5 x 100 µl (10 mM)



Structural formula of 2'OMe-Adenosine-3',5'-bisphosphate

For research use only!

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: C₁₁H₁₇N₅O₁₀P₂ (free acid)

Molecular Weight: 441.23 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} 259 nm, ε 15.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)