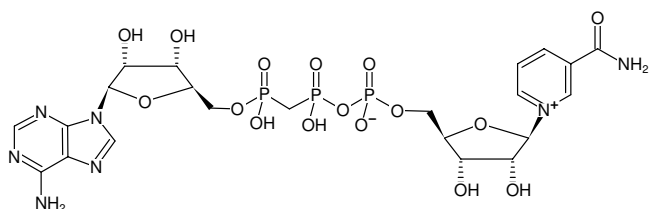


**ApCpp(Nic)**

P¹-(5'-Adenosyl) P³-[5'-(1-β D-ribofuranosyl-nicotinamide)] [(α,β)-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-515	300 µl (10 mM)



Structural formula of ApCpp(Nic)

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: C₂₂H₃₀N₇O₁₆P₃ (free acid)

Molecular Weight: 741.43 g/mol (free acid)

Exact Mass: 741.10 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 16.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)