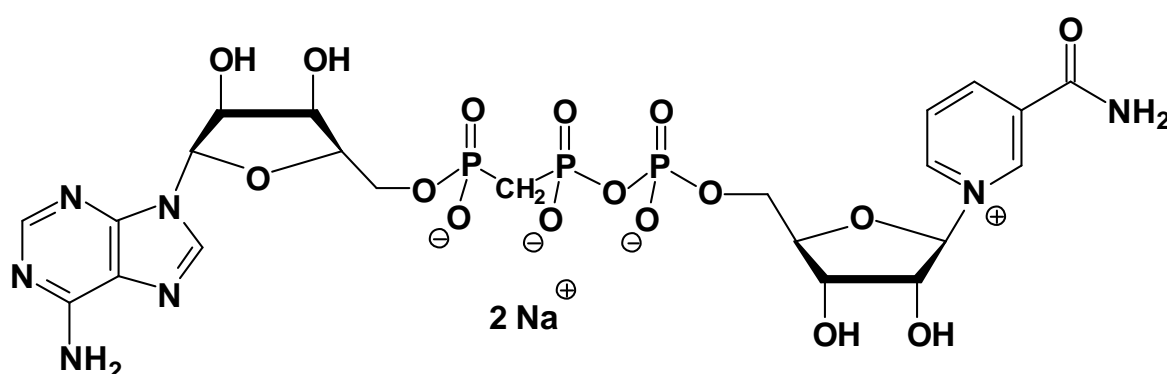


ApC_{pp}(Nic)

P¹-(5'-Adenosyl) P³-[5'-(1-β D-Ribofuranosyl-nicotinamide)]

[(α,β)-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-515	300 Units



Cat. No.: NU-515

Molecular Formula: C₂₂H₂₈N₇O₁₆P₃ (Anion)

Molecular Weight: 739.42 (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