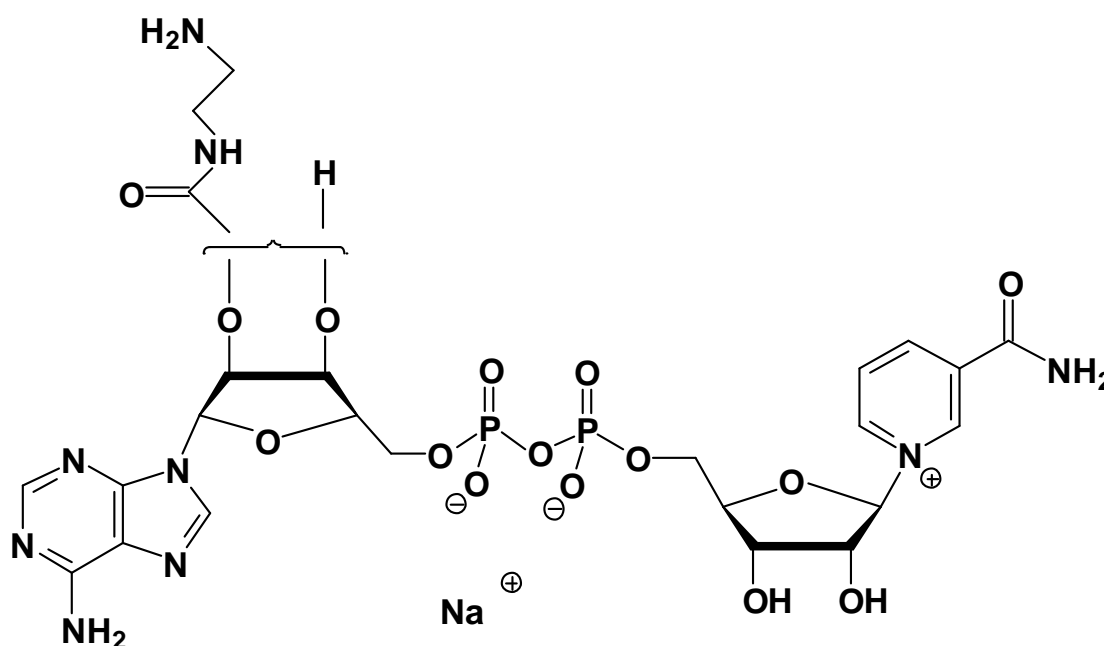


EDA-AP₂(Nic)

P¹-5'-[2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosyl]
P²-[5'-(1-β D-Ribofuranosyl-nicotinamide)] diphosphate,

Cat. No.	Amount
NU-525	300 Units



Cat. No.: NU-525

Molecular Formula: C₂₄H₃₂N₉O₁₅P₂ (Anion)

Molecular Weight: 748.51 (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