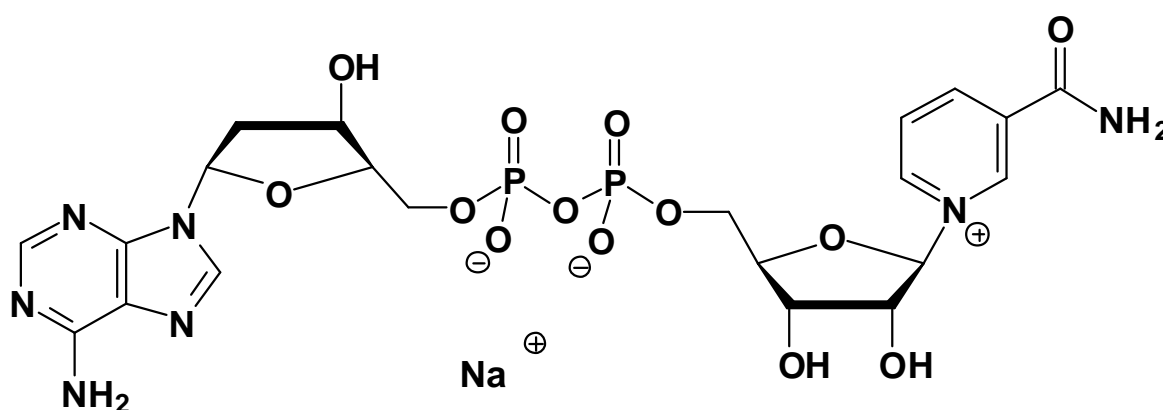


dAP₂(Nic)

P¹-[5'-(2'-Deoxy-adenosyl)] P²-[5'-(1-β D-Ribofuranosyl-nicotinamide)]
diphosphate, Sodium salt

Cat. No.	Amount
NU-517	300 Units



Cat. No.: NU-517

Molecular Formula: C₂₁H₂₆N₇O₁₃P₂ (Anion)

Molecular Weight: 646.42 (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

Selected References:

Fawcett *et al.* (1962) Preparation and properties of some nicotinamide adenine dinucleotide analogues with pentose and purine modifications. *J. of Biological Chem.* **237**(5): 1709