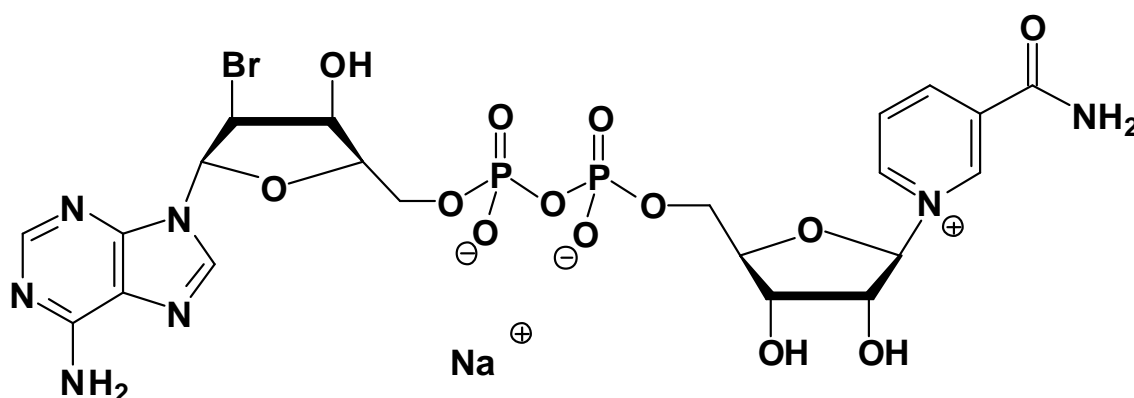


2'Br-AP₂(Nic)

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

Cat. No.	Amount
NU-524	300 Units



Cat. No.: NU-524

Molecular Formula: C₂₁H₂₅BrN₇O₁₃P₂ (Anion)

Molecular Weight: 725.32 (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