

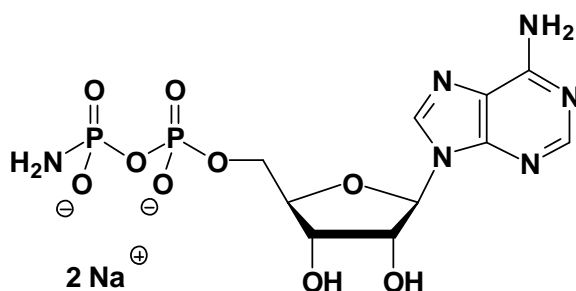
AppNH₂ (AMPPN)

Adenosine-5'-(β-amido)diphosphate, sodium salt

Cat. No.	Amount
NU-1126S	150 Units
NU-1126L	750 Units

Selected References:

Stournaras *et al.* (1988) Thiol group reactivity and polymerization of actin in the presence of ATP analogs. *Biochem Biophys Res Commun.* **155(2)**:962.



Cat. No.: NU-1126

Molecular Formula: C₁₀H₁₄N₆O₉P₂ (Anion)

Molecular Weight: 424.20 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{max} 259nm;
ε 15400

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 6 months after date of delivery.

For research use only!

1 unit = 1 μl of a 10 mM solution