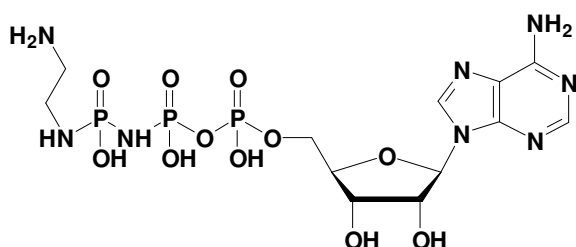


γ -[(2-Aminoethyl)-imido]-AppNHp (AMPPNP)

γ -[(2-Aminoethyl)-imido]-adenosine-5'-[(β,γ)-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-819S	30 Units
NU-819L	150 Units



Cat. No.: NU-819

Molecular Formula: $C_{12}H_{23}N_8O_{11}P_3$ (free acid)

Molecular Weight: 548.28 (free acid)

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

Spectroscopic Properties: λ_{max} 259nm;
 ϵ 15.400 $M^{-1}cm^{-1}$ (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 6 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution