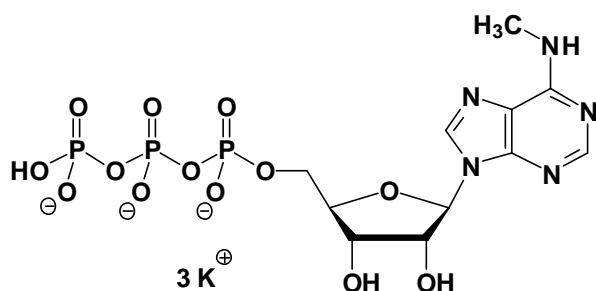


N⁶-Methyl-ATP

N⁶-Methyl-adenosine-5'-triphosphate, Potassium salt

Cat. No.	Amount
NU-1101S	150 Units
NU-1101L	750 Units



Cat. No.: NU-1101

Molecular Formula: C₁₁H₁₅N₅O₁₃P₃ (Anion)

Molecular Weight: 518.18 (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