

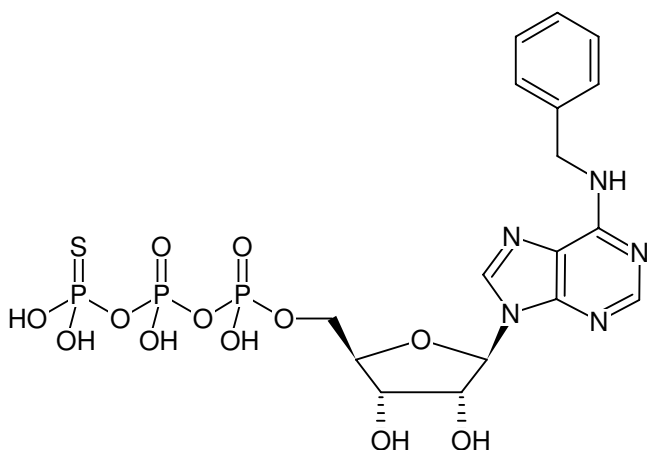
**N⁶-Benzyl-ATPyS**

6-Bn-ATPyS

N⁶-Benzyladenosine-5'-O-(3-thiotriphosphate), Sodium salt

Cat. No.	Amount
NU-241	3,4 mg (approx. 5 µmol)

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically.

Structural formula of N⁶-Benzyl-ATPyS**For research use only!****Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 6 months after date of delivery**Molecular Formula:** C₁₇H₂₂N₅O₁₂P₃S (free acid)**Molecular Weight:** 613.37 g/mol (free acid)**Exact Mass:** 613.02 g/mol (free acid)**CAS#:** 944834-42-8**Purity:** ≥ 90 % (HPLC)**Form:** solid**Spectroscopic Properties:** λ_{max} 269 nm, ε 20.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)