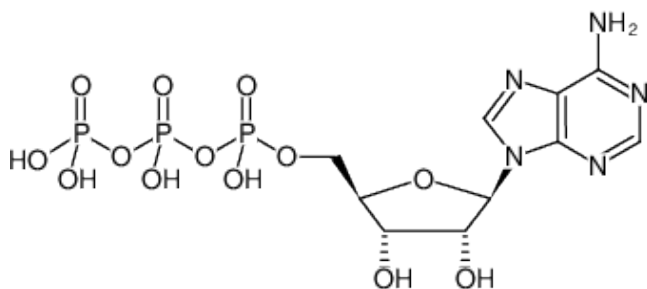


**ATP - Solid**

Adenosine 5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1049-1G	1 g
NU-1049-10G	10 g
NU-1049-100G	100 g

Quality Control Specifications: in vitro transcription (T7 RNA polymerase): visible RNA fragments after 5 min incubation, Proteases: not detectable



Structural formula of ATP - Solid

For general laboratory use.**Shipping:** Shipped on gel packs**Storage Conditions:** Store at -20 °C

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

Shelf Life: 12 months**Molecular Formula:** C₁₀H₁₆N₅O₁₃P₃ (free acid)**Molecular Weight:** 507.18 g/mol (free acid)**Exact Mass:** 507.00 g/mol (free acid)**CAS#:** 987-65-5**Purity:** ≥ 98 % (HPLC)**Form:** Powder**Color:** White**Solubility:** Soluble in water. Clear solution.**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)