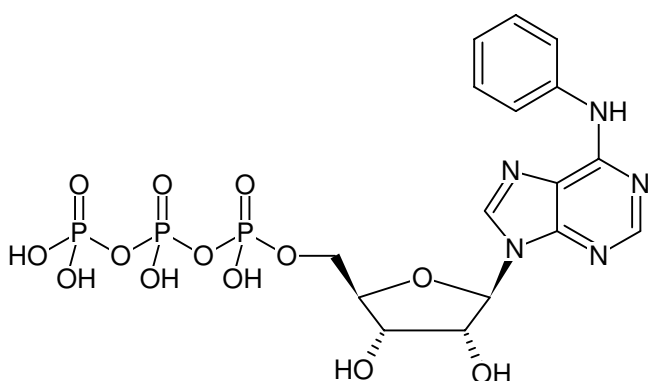


**N⁶-Phenyl-ATP**N⁶-Phenyladenosine-5'-O-triphosphate, Sodium salt

| Cat. No. | Amount |
|----------|--------------------|
| NU-879S | 250 µl (10 mM) |
| NU-879L | 5 x 250 µl (10 mM) |

Structural formula of N⁶-Phenyl-ATP**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 after date of delivery**Molecular Formula:** C₁₆H₂₀N₅O₁₃P₃ (free acid)**Molecular Weight:** 583.28 g/mol (free acid)**Exact Mass:** 583.03 g/mol (free acid)**CAS#:** 105740-47-4**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 288 nm, ε 20.8 L mmol⁻¹ cm⁻¹
(Tris-HCl pH 7.5)