

**IppNHp**

(IMPPNP)

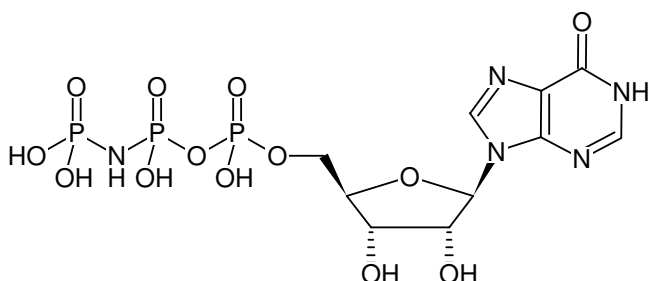
Inosine-5'-[(β,γ)-imido]triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1204-5	5 mg

Selected References:

Gille *et al.* (2003) GDP Affinity and Order State of the catalytic Site Are Critical for Function of Xanthine Nucleotide-selective G_{αs} Proteins. *J. Biol. Chem.* **278**:7822.

Schuster *et al.* (1976) Effect of inosine 5'- (beta,gamma-imido) triphosphate and other nucleotides on beef-heart mitochondrial ATPase. *J. Biol. Chem.* **251**(21):6705.



Structural formula of IppNHp

For research use only!**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: 6 months after date of delivery**Molecular Formula:** C₁₀H₁₆N₅O₁₃P₃ (free acid)**Molecular Weight:** 507.18 g/mol (free acid)**Exact Mass:** 507.00 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 249 nm, ε 12.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)