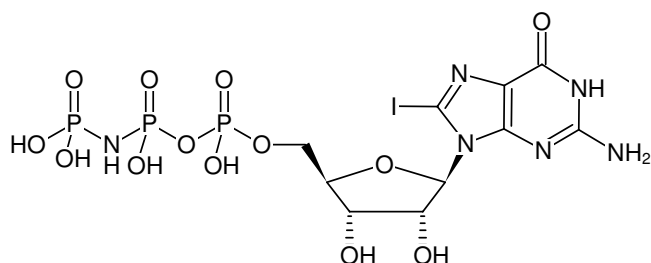


**8-Iodo-GppNHp**

(8I-GppNHp, 8I-GMPPNP)

8-Iodo-guanosine-5'-[(β,γ)-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-126	200 µl (10 mM)



Structural formula of 8-Iodo-GppNHp

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: 6 months after date of delivery**Molecular Formula:** C₁₀H₁₆N₆O₁₃P₃I (free acid)**Molecular Weight:** 648.09 g/mol (free acid)**Exact Mass:** 647.90 g/mol (free acid)**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} 261 nm, ε 17.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Schmohl *et al.* (2012) Functional analysis of Rho GTPase activation and inhibition in a bead-based miniaturized format. *Methods. Mol. Biol.* **827** : 271.Labesse *et al.* (2011) Structural and functional characterization of the Mycobacterium tuberculosis uridine monophosphate kinase: insights into the allosteric regulation. *Nucleic Acids Res.* **39** (8):3458.Ash *et al.* (2011) A suite of Switch I and Switch II mutant structures from the G-protein domain of FeoB *Acta Crystallogr. D Biol. Crystallogr.* **67** (11):973.