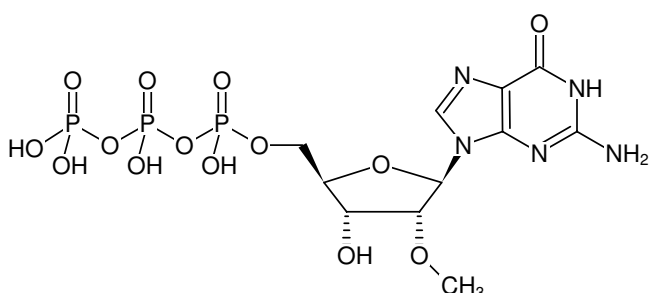


**2'OMe-GTP**

2'-O-Methylguanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1127S	50 µl (100 mM)
NU-1127L	5 x 50 µl (100 mM)



Structural formula of 2'OMe-GTP

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: 12 months after date of delivery**Molecular Formula:** C₁₁H₁₈N₅O₁₄P₃ (free acid)**Molecular Weight:** 537.21 g/mol (free acid)**Exact Mass:** 537.01 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 252 nm, ε 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

Heck *et al.* (2008) Effects of mutagenic and chain-terminating nucleotide analogs on enzymes isolated from hepatitis C virus strains of various genotypes. *Antimicrob. Agents Chemother.* **52** (6):1901.

Paeshuysse *et al.* (2007) The imidazopyrrolopyridine analogue AG110 is a novel, highly selective inhibitor of pestiviruses that targets the viral RNA-dependent RNA polymerase at a hot spot for inhibition of viral replication. *J. Virol.* **81** (20):11046.

Paeshuysse *et al.* (2006) A novel, highly selective inhibitor of pestivirus replication that targets the viral RNA-dependent RNA polymerase. *J. Virol.* **80** (1):149.

Dutarte *et al.* (2005) A relaxed discrimination of 2'-O-methyl-GTP relative to GTP between de novo and Elongative RNA synthesis by the hepatitis C RNA-dependent RNA polymerase NS5B. *J. Biol. Chem.* **280** (8):6359.

Fresno *et al.* (1980) Inhibitory effects of 'cap' analogues on globin mRNA and encephalomyocarditis RNA translation in a reticulocyte cell-free system. *Eur. J. Biochem.* **103** (1):125.