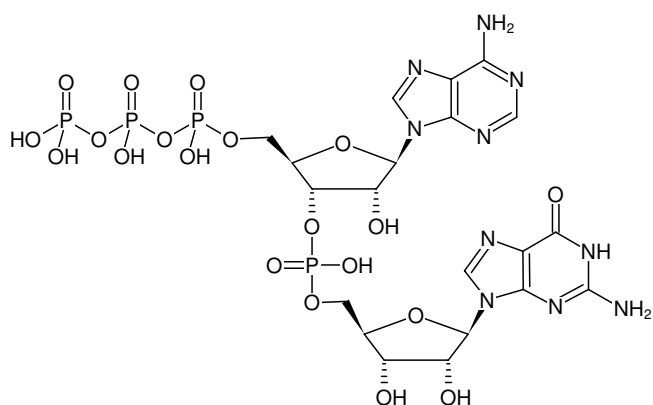




pppApG

Sodium salt

Cat. No.	Amount
NU-980	10 µl (100 mM)



Structural formula of pppApG

For research use only!**Shipping:** shipped on blue ice**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:** C₂₀H₂₈N₁₀O₂₀P₄ (free acid)**Molecular Weight:** 852.39 g/mol (free acid)**Exact Mass:** 852.04 g/mol (free acid)**Purity:** ≥ 95 % (RP-HPLC)**Form:** clear aqueous solution**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 260 nm, ε 25.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Whelan *et al.* (2004) Affinity assays using fluorescence anisotropy with capillary electrophoresis separation. *Anal. Chem.* **76** (24):7380.Korlach *et al.* (2004) Spontaneous nucleotide exchange in low molecular weight GTPases by fluorescently labeled gamma-phosphate-linked GTP analogs. *Proc. Natl. Acad. Sci. USA.* **101** (9):2800.