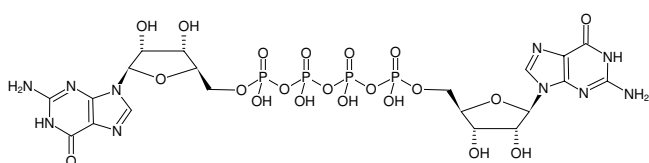


**GP₄G**P¹-(5'-Guanosyl)-P⁴-(5'-guanosyl)-tetraphosphate, Sodium salt

Cat. No.	Amount
NU-533S	20 µl (10 mM)
NU-533L	5 x 20 µl (10 mM)

Structural formula of GP₄G**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:** C₂₀H₂₈N₁₀O₂₁P₄ (free acid)**Molecular Weight:** 868.39 g/mol (free acid)**Exact Mass:** 868.04 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} 252 nm, ε 25.0 L mol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)