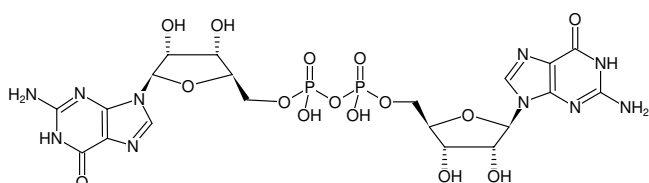


**GP₂G**P¹-(5'-Guanosyl)-P²-(5'-guanosyl)-diphosphate, Sodium salt

Cat. No.	Amount
NU-531S	20 µl (10 mM)
NU-531L	5 x 20 µl (10 mM)

Structural formula of GP₂G**Selected References:**

Stern *et al.* (2012) Studies of Mg²⁺/Ca²⁺ complexes of naturally occurring dinucleotides: potentiometric titrations, NMR, and molecular dynamics. *J. Biol. Inorg. Chem.* **17** (6):861.

Sillero *et al.* (2002) Dinucleoside polyphosphates stimulate the primer independent synthesis of poly(A) catalyzed by yeast poly(A) polymerase. *Eur. J. Biochem.* **269** (21):5323.

Jankowski *et al.* (2001) Dinucleotides as growth-promoting extracellular mediators. Presence of dinucleoside diphosphates Ap2A, Ap2G, and Gp2G in releasable granules of platelets. *J. Biol. Chem.* **276** (12):8904.

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:** C₂₀H₂₆N₁₀O₁₅P₂ (free acid)**Molecular Weight:** 708.43 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 mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)