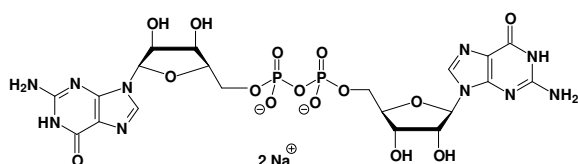


GP₂G

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

Cat. No.	Amount
NU-531S	20 Units
NU-531L	100 Units



Cat. No.: NU-531

Molecular Formula: C₂₁H₂₇N₁₀O₁₅P₂ (Anion)

Molecular Weight: 721.45 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

* 1 unit = 1 µl of a 10 mM solution

For research use only!