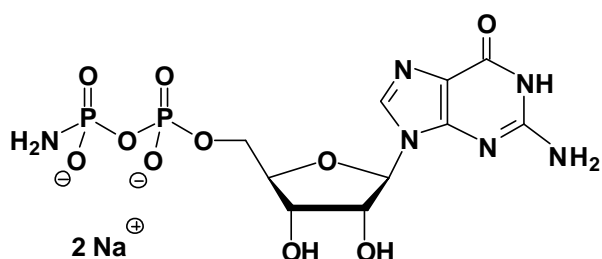


GppNH₂

Guanosine-5'-(β-amino)-diphosphate, Sodium salt

Cat. No.	Amount
NU-1136S	150 Units
NU-1136L	750 Units



Cat. No.: NU-1136

Molecular Formula: C₁₀H₁₄N₆O₁₀P₂ (Anion)

Molecular Weight: 440.20 (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 6 months after date of delivery.

For research use only!

1 unit = 1 μl of a 10 mM solution