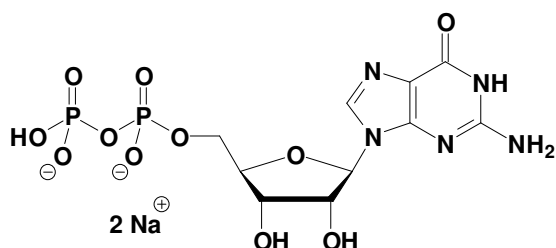


GDP

Guanosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1172S	1 g
NU-1172L	5 g



Cat. No.: NU-1172

Molecular Formula: C₁₀H₁₃N₅O₁₁P₂ (Anion)

Molecular Weight: 441.18 (Anion)

Purity: > 95%, lyophilized powder

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.

For research use only!