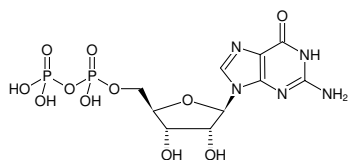


**GDP - Solid**

Guanosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1172-1G	1 g
NU-1172-5G	5 g



Structural formula of GDP - Solid

For research use only!**Shipping:** shipped on blue ice**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:** C₁₀H₁₅N₅O₁₁P₂ (free acid)**Molecular Weight:** 443.20 g/mol (free acid)**Exact Mass:** 443.02 g/mol (free acid)**CAS#:** 43139-22-6**Purity:** ≥ 90 % (HPLC)**Form:** lyophilised**Spectroscopic Properties:** λ_{max} 252 nm, ε 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Specific Ligands:**Engineered GPCR-Gα-complexes^[1]**Selected References:**[1] Jastrzebska *et al.* (2010) Complex between photoactivated rhodopsin and transducin: progress and questions. *Biochem. J.* **428**:1.Kowluru (2010) Small G-proteins in islet beta-cell function. *Endocrine reviews* **31**:52.Yang *et al.* (2010) A guaninine nucleotide exchange factor is a component of the meiotic spindle pole body in *Schizosaccharomyces pombe*. *Mol. Biol. Cell.* **21** (7):1272.