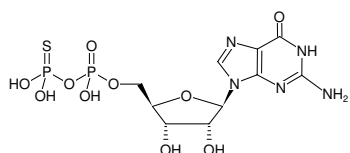


**GDPβS**

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

Cat. No.	Amount
NU-427-5	5 mg
NU-427-25	25 mg



Structural formula of GDPβS

For general laboratory use.**Shipping:** shipped on dry ice**Storage Conditions:** store at -20 °C**Shelf Life:** 6 months after date of delivery**Molecular Formula:** C₁₀H₁₅N₅O₁₀P₂S (free acid)**Molecular Weight:** 459.26 g/mol (free acid)**Exact Mass:** 459.00 g/mol (free acid)**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 252 nm, ε 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0. This can be achieved by dissolving the nucleotide in a buffer of your choice (50 - 100 mM, pH 7 - 10). Dissolve and adjust concentration photometrically.

Selected References:

Chang *et al.* (2005) Nitric Oxide-dependent Allosteric Inhibitory Role of a Second Nucleotide Binding Site in Soluble Guanylyl Cyclase. *J. Biol. Chem.* **280** (12):11513.

Balttus *et al.* (2002) Spontaneous release of GDP from Gi proteins and inhibition of adenylyl cyclase in cardiac sarcolemmal membranes. *Naunyn Schmiedebergs Arch. Pharmacol.* **365** (1):50.

Burgos *et al.* (2000) Guanine nucleotides block agonist-driven 45Ca²⁺ influx in chick embryo retinal explants. *Neuroreport.* **11** (10):2303.

Peterson *et al.* (1999) GTP (gammaS) and GDP (betaS) as electron donors: new wine in old bottles. *Life Sci.* **65** (11):1135.