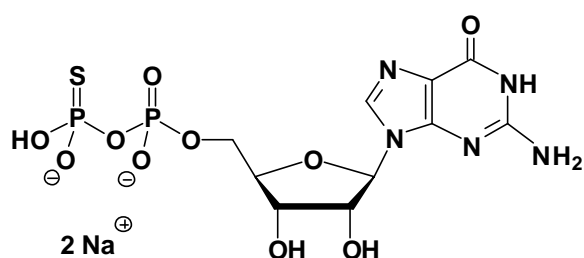


GDP β S

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

Cat. No.	Amount
NU-427S	50 Units
NU-427L	250 Units



Cat. No.: NU-427

Molecular Formula: C₁₀H₁₃N₅O₁₀P₂S (Anion)

Molecular Weight: 457.24 (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 μ l of a 10 mM solution

Selected References:

Baltus *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.