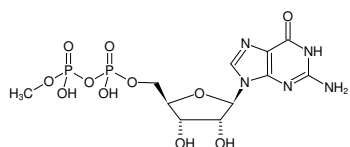


**GDP β methyl**GDP- β -methyl ester, Triethylammonium salt

Cat. No.	Amount
NU-977	200 μ l (10 mM)

Selected References:

Wilke *et al.* (2012) Activity-Based Probe for Histidine Kinase Signaling *J. Am. Chem. Soc.* **134** (22):9150.

Structural formula of GDP β methyl**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:** C₁₁H₁₇N₅O₁₁P₂ (free acid)**Molecular Weight:** 457.23 g/mol (free acid)**Exact Mass:** 457.04 g/mol (free acid)**CAS#:** 57817-66-0 (free acid)**Purity:** \geq 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{\max} 252 nm, ϵ 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)