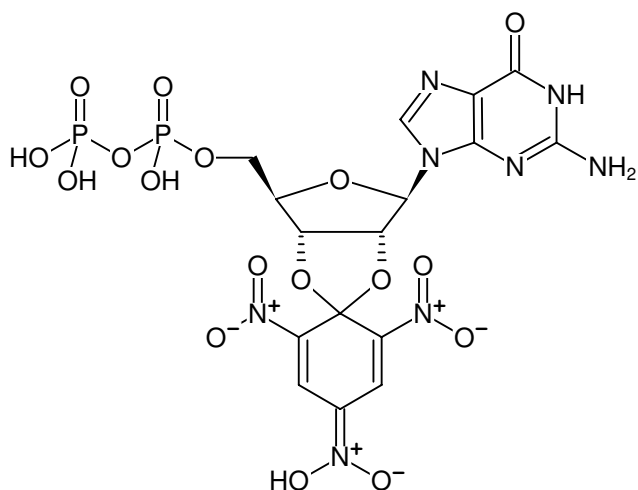


**TNP-GDP**

2',3'-O-Trinitrophenyl-guanosine-5'-diphosphate, Triethylammonium salt

Cat. No.	Amount
NU-217S	1 μ mol
NU-217L	5 x 1 μ mol



Structural formula of TNP-GDP

For research use only!**Shipping:** shipped on gel packs**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:** C₁₆H₁₆N₈O₁₇P₂**Molecular Weight:** 654.29 g/mol**Exact Mass:** 654.01 g/mol**Purity:** \geq 95 % (HPLC)**Form:** solid**Color:** orange

Spectroscopic Properties: λ_{\max} 252/408/470 nm,
 ϵ 24.1/26.5/18.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 408/470 nm,
 λ_{em} 552 nm

Selected References:

Suryanarayana *et al.* (2009) Differential Inhibition of Various Adenylyl Cyclase Isoforms and Soluble Guanylyl Cyclase by 2',3'-O- (2,4,6-Trinitrophenyl)-Substituted Nucleoside 5'-Triphosphates. *J. Pharmacol. Exp. Ther.* **330** (3):687.

Sprang *et al.* (2006) Broad Specificity of Mammalian Adenylyl Cyclase for Interaction with 2',3'-Substituted Purine- and Pyrimidine Nucleotide Inhibitors. *Mol. Pharmacol.* **70**:878.

Hiratsuka (1985) A chromophoric and fluorescent analog of GTP, 2',3'-O-(2,4,6-trinitrocyclohexadienylidene)-GTP, as a spectroscopic probe for the GTP inhibitory site of liver glutamate dehydrogenase. *J. Biol. Chem.* **260** (8):4784.