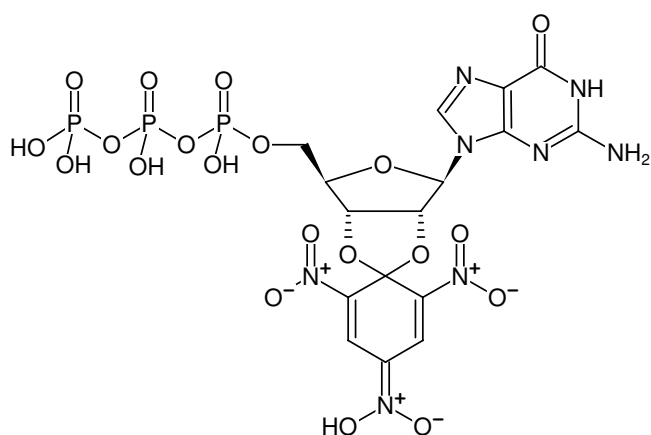


**TNP-GTP**

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

Cat. No.	Amount
NU-220S	1 μ mol
NU-220L	5 x 1 μ mol



Structural formula of TNP-GTP

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:** C₁₆H₁₇N₈O₂₀P₃**Molecular Weight:** 734.27 g/mol**Exact Mass:** 733.98 g/mol**CAS#:** 612837-26-0**Purity:** \geq 95 % (HPLC)**Form:** solid**Color:** orange

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

Selected References:

Spangler *et al.* (2011) Interaction of the diguanylate cyclase YdeH of *Escherichia coli* with 2', (3')-substituted purine and pyrimidine nucleotides. *J. Pharmacol. Exp. Ther.* **336** (1):234.

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.

Goettle *et al.* (2007) Molecular analysis of the interaction of *Bordetella pertussis* adenylyl cyclase with fluorescent nucleotides. *Molecular Pharmacology* **72** (3):526.

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.

Golczak *et al.* (2002) GTP-induced membrane binding and ion channel activity of annexin VI: is annexin VI a GTP biosensor? *Biophys. J.* **82**:2737.

Chordia *et al.* (1998) Distances between the paclitaxel, colchicine, and exchangeable GTP binding sites on tubulin. *Biochemistry* **37**:6636.

Randak *et al.* (1996) A recombinant polypeptide model of the second predicted nucleotide binding fold of the cystic fibrosis transmembrane conductance regulator is a GTP-binding protein. *FEBS Lett.* **398**:97.

Baubichon-Cortay *et al.* (1994) Overexpression and purification of the carboxyl-terminal nucleotide-binding domain from mouse P-glycoprotein. Strategic location of a tryptophan residue. *J. Biol. Chem.* **269**:22983.

Hoyer *et al.* (1986) Synthesis of 2',3'-O-(2,4,6-trinitrocyclohexadienylidene)guanosine 5'-triphosphate and a study of its inhibitory properties with adenylyl cyclase. *Arch. Biochem. Biophys.* **245**:369.

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**:4784.