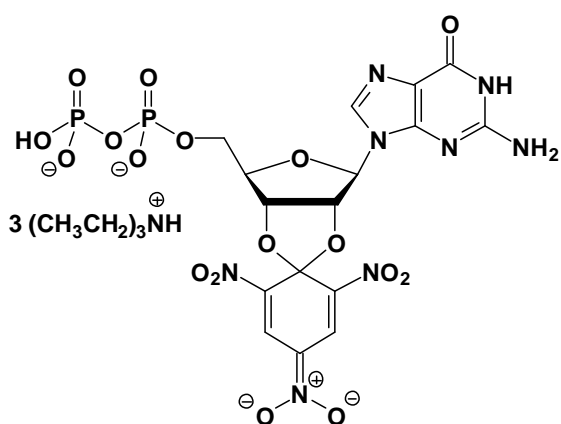


## TNP-GDP

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

Cat. No.	Amount
NU-217S	100 Units
NU-217L	500 Units



### Selected References:

Sprang *et al.* (2006) Broad and 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.

**Cat. No.:** NU-217

**Molecular Formula:** C<sub>16</sub>H<sub>15</sub>N<sub>8</sub>O<sub>17</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 653.28 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

### Spectroscopic properties:

$\lambda_{\max}$  252/408/470 nm;  $\epsilon$  24100/26500/18300;

$\lambda_{\text{exc}}$  408/470nm;  $\lambda_{\text{em}}$  552 nm

### 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 12 months after date of delivery.

### For research use only!

1 unit = 1  $\mu$ l of a 10 mM solution