

**TNP-GppNHp**

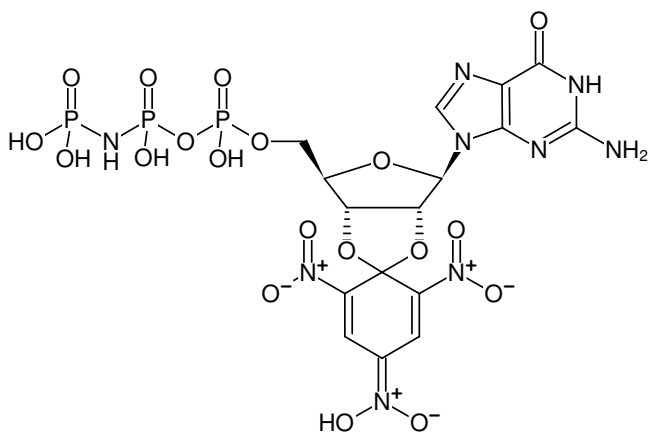
(TNP-GMPPNP)

2',3'-O-Trinitrophenyl-guanosine-5'-[(β,γ)-imido]triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-218S	0,1 μmol
NU-218L	5 x 0,1 μmol

**Selected References:**

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



Structural formula of TNP-GppNHp

**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:** 6 months after date of delivery**Molecular Formula:** C<sub>16</sub>H<sub>18</sub>N<sub>9</sub>O<sub>19</sub>P<sub>3</sub>**Molecular Weight:** 733.28 g/mol**Exact Mass:** 732.99 g/mol**Purity:** ≥ 90 % (HPLC)**Form:** orange lyophilised solid

**Spectroscopic Properties:** λ<sub>max</sub> 252/408/470 nm,  
 ε 24.1/26.5/18.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5), λ<sub>exc</sub> 408/470 nm,  
 λ<sub>em</sub> 552 nm