

**Mant-dGppNHp**

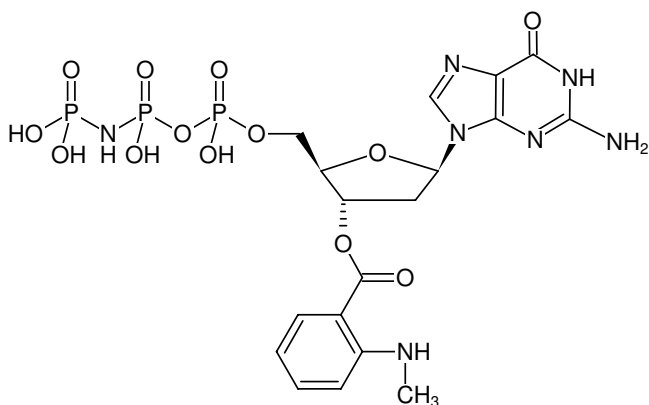
(Mant-dGMPPNP)

3'-O-(N-Methyl-anthraniloyl)-2'-deoxyguanosine-5'- [(β,γ)-imido]triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-208S	10 µl (10 mM)
NU-208L	5 x 10 µl (10 mM)

Selected References:

Scheidig *et al.* (1995) X-ray crystal-structure analysis of the catalytic domain of the oncogene product p21 (H-Ras) complexed with caged GTP and Mant dGppNHp. *J. Mol. Biol.* **253** (1):132.



Structural formula of Mant-dGppNHp

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:** C₁₈H₂₄N₇O₁₃P₃ (free acid)**Molecular Weight:** 639.35 g/mol (free acid)**Exact Mass:** 639.06 g/mol (free acid)**CAS#:** 170032-90-3**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 252/355 nm, ε 22.6/5.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 355 nm, λ_{em} 448 nm