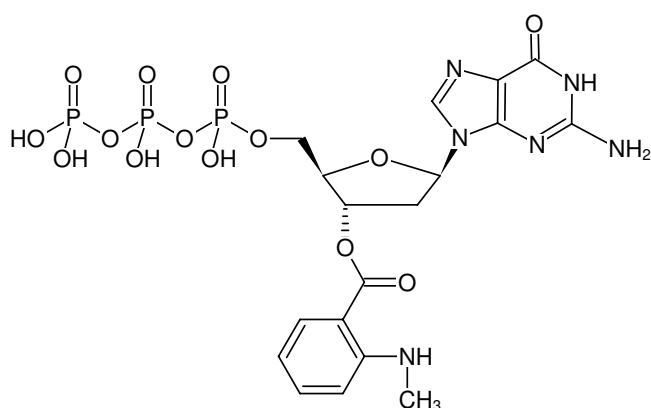


**Mant-dGTP**

3'-O-(N-Methyl-anthraniloyl)-2'-deoxyguanosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-212S	50 µl (10 mM)
NU-212L	5 x 50 µl (10 mM)



Structural formula of Mant-dGTP

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: 12 months after date of delivery**Molecular Formula:** C₁₈H₂₃N₆O₁₄P₃ (free acid)**Molecular Weight:** 640.33 g/mol (free acid)**Exact Mass:** 640.05 g/mol (free acid)**CAS#:** 124615-99-2**Purity:** ≥ 95 % (HPLC)**Form:** colorless to slightly yellow solution in water**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**Applications:**Resonance energy transfer studies (FRET)^[1]Enzyme kinetic studies^[2]**Specific Ligands:**Bacillus anthrax toxin edema factor^[1]EF-Tu^[3]**Selected References:**

[1] Suryanarayana *et al.* (2009) Distinct interactions of 2'- and 3'-O-(N-methyl)anthraniloyl-isomers of ATP and GTP with the adenylyl cyclase toxin of *Bacillus anthracis*, edema factor. *Biochemical Pharmacology* **78** (3):224.

[2] Kettlun *et al.* (2005) Potato tuber isoapyrases: Substrate specificity, affinity labelling, and proteolytic susceptibility. *Phytochemistry* **66**:975.

[3] Rodnina *et al.* (1995) Codon-dependent conformational change of elongation-factor tu preceding GTP hydrolysis on the ribosome. *EMBO J.* **14** (11):2613.

Hazlett *et al.* (1993) Solution dynamics of p21 (Ras) proteins bound with fluorescent nucleotides - a time-resolved fluorescence study. *Biochemistry-US* **32** (49):13575.