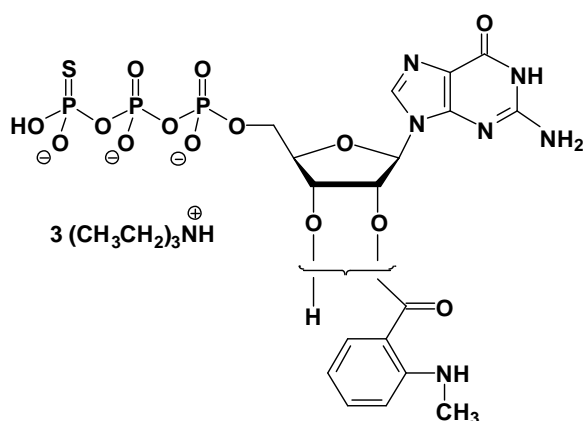


mant-GTP γ S

2'/3'-O-(N-Methyl-anthraniloyl)-guanosine-5'-(γ -thio)-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-209S	10 Units
NU-209L	50 Units



Cat. No.: NU-209

Molecular Formula: C₁₈H₂₀N₆O₁₄P₃S (Anion)

Molecular Weight: 669.36 (Anion)

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

Spectroscopic properties:

λ_{\max} 252/355 nm; ϵ 22600/5700; λ_{exc} 355 nm;
 λ_{em} 448 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 6 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution

Selected References:

Herzig *et al.* (2007) Functional Adenylyl Cyclase Inhibition in Murine Cardiomyocytes by 2'(3')-O-(N-Methylanthraniloyl)-Guanosine 5'-[gamma-Thio]triphosphate. *JPET* **321**:608.

Gottle *et al.* (2007) Molecular analysis of the interaction of Bordetella pertussis adenylyl cyclase with fluorescent nucleotides. *Molecular Pharmacology Fast Forward*, June 6, 2007; DOI: 10.1124/mol.107.034413 (published online).

Celda *et al.* (2007) Structural insights into GTPase domain of *Escherichia coli* MnmE protein. *Proteins* **66** (3):726.

Armengod *et al.* (2005) Effects of Mutagenesis in Switch I Region and Conserved Arginines of *Escherichia coli* MnmE Protein, A GTPase Involved in tRNA Modification. *J. Biol. Chem.* **280** (35):30660

Gille *et al.* (2003) 2'(3')-O-(N-methylanthraniloyl)-substituted GTP analogs: a novel class of potent competitive adenylyl cyclase inhibitors. *J. Biol. Chem.* **278**:12672.

Remmers *et al.* (1999) Interdomain interactions regulate GDP release from heterotrimeric G proteins. *Biochemistry* **38** (42):13795.

Remmers *et al.* (1998) Detection and quantitation of heterotrimeric G proteins by fluorescence resonance energy transfer. *Anal. Biochem.* **257** (1):89.