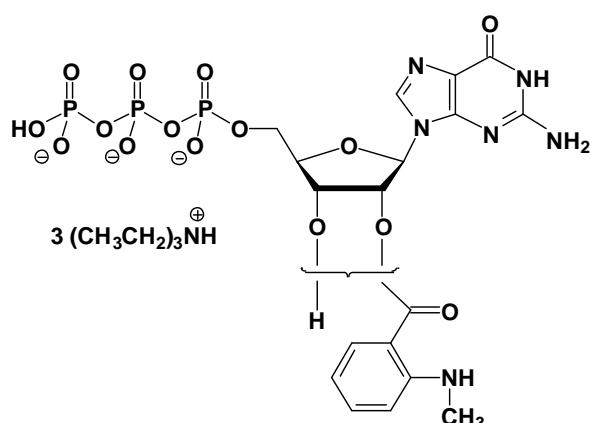


mant-GTP

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

Cat. No.	Amount
NU-206S	150 Units
NU-206L	750 Units



Cat. No.: NU-206

Molecular Formula: C₁₈H₂₀N₆O₁₅P₃ (Anion)

Molecular Weight: 653.30 (Anion)

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

Spectroscopic properties:

λ_{\max} 255/355 nm; ϵ 23300/5800; λ_{exc} 355nm;
 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 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution

Selected References:

Kambach *et al.* (2007) Human OLA1 Defines an ATPase subfamily in the Obg Family of GTP-binding proteins. *J. Biol. Chem.* **282** (27):19928.

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).

Pestova *et al.* (2006) Kinetic Analysis of Interaction of Eukaryotic Release Factor 3 with Guanine Nucleotides. *J. Biol. Chem.* **281** (52):40224

Wintermeyer *et al.* (2006) Role and timing of GTP binding and hydrolysis during EF-G-dependent RNA translocation on the ribosome. *PNAS* **103** (37):13670

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.

Tung-Chung Mou *et al.* (2005) Structural Basis for the Inhibition of Mammalian Membrane Adenylyl Cyclase by 2'(3')-O-(N-Methylantraniloyl)- guanosine 5'-Triphosphate. *J. Biol. Chem.* **280** (8):7253.

Alexandrov *et al.* (2001) Fluorescence methods for monitoring interactions of Rab proteins with nucleotides, Rab escort protein, and geranylgeranyltransferase. *Methods Enzymol.* **329**:14.

Jagath *et al.* (1998) Interaction of guanine nucleotides with the signal recognition particle from Escherichia coli. *Biochemistry* **37**:15408.
 Simon *et al.* (1996) Kinetics of interaction of Rab5 and Rab7 with nucleotides and magnesium ions. *J. Biol. Chem.* **271**:20470.

Renstrand *et al.* (1995) Substrate and product structural requirements for binding of nucleotides to H-ras p21: the mechanism of discrimination between guanosine and adenosine nucleotides. *Biochemistry* **34**:593.

Remmers *et al.* (1994) Fluorescent guanine nucleotide analogs and G protein activation. *J. Biol. Chem.* **269**:13771.

Hazlett *et al.* (1993) Solution dynamics of p21ras proteins bound with fluorescent nucleotides: a time-resolved fluorescence study. *Biochemistry* **32**:13575.

Brownbridge *et al.* (1993) Interaction of GTPase activating proteins (GAPs) with p21ras measured by a novel fluorescence anisotropy method. Essential role of Arg-903 of GAP in activation of GTP hydrolysis on p21ras. *J. Biol. Chem.* **268**:10914.

Eccleston *et al.* (1991) Fluorescence approaches to the study of the p21ras GTPase mechanism. *Biochem. Soc. Trans.* **19**:432.