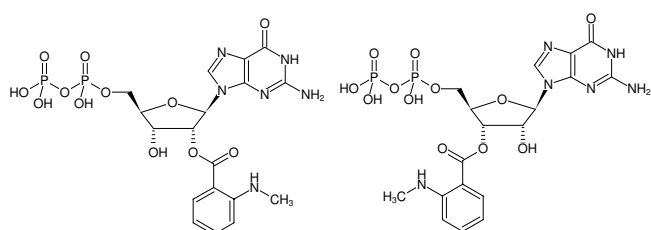




Mant-GDP

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

| Cat. No. | Amount |
|----------|--------------------|
| NU-204S | 150 µl (10 mM) |
| NU-204L | 5 x 150 µl (10 mM) |



Structural formula of Mant-GDP

For research use only!

Shipping: shipped on gel packs

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: C₁₈H₂₂N₆O₁₂P₂ (free acid)

Molecular Weight: 576.35 g/mol (free acid)

Exact Mass: 576.08 g/mol (free acid)

CAS#: 148821-02-7

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

Applications:

Conformational dynamic: DnaB/C-proteins^[1], Csk^[2]

Fluorescence stop-flow kinetics: release factor^[3], eIF5B^[4], EF-1B^[5], EF-Tu^[6]

FRET: AC^[7]

Specific Ligands:

Translation factor: eIF5B^[4], EF1B2^[5], EF-Tu^[6]

Release factor^[3]

Selected References:

[1] Roychowdhury *et al.* (2009) Interactions of Escherichia coli DnaB-DnaC protein complex with nucleotide cofactors. 1. Allosteric conformational transitions of the complex. *Biochemistry* **48**:6712.

[2] Shaffer *et al.* (2001) Nucleotide release and associated conformational changes regulate function in the COOH-terminal Src kinase, Csk. *Biochemistry* **40**:11149.

[3] Pisareva *et al.* (2006) Kinetic analysis of interaction of eukaryotic release factor 3 with guanine nucleotides. *J. Biol. Chem.* **281**:40224.

[4] Pisareva *et al.* (2007) Kinetic analysis of the interaction of guanine nucleotides with eukaryotic translation initiation factor eIF5B. *Biochemistry* **46**:2622.

[5] Pittman *et al.* (2006) Mg²⁺ and key lysine modulate exchange activity of eukaryotic translation elongation factor eIF1B. *J. Biol. Chem.* **281**:19457.

[6] Wieden *et al.* (2002) Mechanism of elongation factor (EF)-Ts-catalyzed nucleotide exchange in EF-Tu. Contribution of contacts at the guanine base. *J. Biol. Chem.* **277**:6032.

[7] Goettle *et al.* (2007) Molecular analysis of the interaction of Bordetella pertussis adenyl cyclase with fluorescent nucleotides. *Molecular Pharmacology* **72**:526.

Wehner *et al.* (2012) The guanine cap of human guanylate-binding protein 1 is responsible for dimerization and self-activation of GTP hydrolysis. *FEBS J.* **279** (2):203.

Oreb *et al.* (2011) Substrate binding disrupts dimerization and induces nucleotide exchange of the chloroplast GTPase Toc33. *Biochem J.* **436** (2):313.

Koshiba *et al.* (2011) Structure-function analysis of the yeast mitochondrial Rho GTPase, Gem1p: implications for mitochondrial inheritance. *J. Biol. Chem.* **286** (1):354.

Yang *et al.* (2010) A guanine nucleotide exchange factor is a component of the meiotic spindle pole body in Schizosaccharomyces pombe. *Mol. Biol. Cell.* **21** (7):1272.

**Mant-GDP**

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

Dahl *et al.* (2008) The Importance of P-loop and Domain Movements in EF-Tu for Guanine Nucleotide Exchange. *J. Biol. Chem.* **281** (30):21139.

Ugolev *et al.* (2008) Dissociation of Rac1 (GDP).RhoGDI Complexes by the Cooperative Action of Anionic Liposomes Containing Phosphatidylinositol 3,4,5-Trisphosphate, Rac Guanine Nucleotide Exchange Factor, and GTP. *J. Biol. Chem.* **283** (32):22257.

Ugolev *et al.* (2008) Liposomes Comprising Anionic but Not Neutral Phospholipids Cause Dissociation of Rac (1 or 2).RhoGDI Complexes and Support Amphiphile-independent NADPH Oxidase Activation by Such Complexes. *J. Biol. Chem.* **281** (28):19204.

Bos *et al.* (2007) Structure of the cyclic-AMP-responsive exchange factor Epac2 in its auto-inhibited state. *Nature* **439**:625.

Noel *et al.* (2007) Turning off the Ras switch with the flick of a finger. *Nature Structural & Molecular Biology* **4**:677.

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

Pick *et al.* (2006) Liposomes Comprising Anionic but not neutral Phospholipids cause dissociation of Rac (1 or 2)RhoGDI Complexes and support Amphiphile-independent NADPH oxidase Activation by Such Complexes. *J. Biol. Chem.* **281** (28):19204.

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.

Thanbichler *et al.* (2000) Kinetics of the interaction of translation factor SelB from *Escherichia coli* with guanosine nucleotides and selenocysteine insertion sequence RNA. *J. Biol. Chem.* **275**:20458.

Ahmadian *et al.* (1997) Individual rate constants for the interaction of Ras proteins with GTPase-activating proteins determined by fluorescence spectroscopy. *Biochemistry* **36**:4535.

Remmers *et al.* (1996) Partial G protein activation by fluorescent guanine nucleotide analogs. Evidence for a triphosphatebound but inactive state. *J. Biol. Chem.* **271**:4791.

Rensland *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.

Richter *et al.* (1995) Interferon-induced MxA protein. GTP binding and GTP hydrolysis properties. *J. Biol. Chem.* **270**:13512.

Neal *et al.* (1990) Hydrolysis of GTP by p21NRAS, the NRAS protooncogene product, is accompanied by a conformational change in the wild-type protein: use of a single fluorescent probe at the catalytic site. *Proc. Natl. Acad. Sci. USA* **87**:3562.

Related Products: