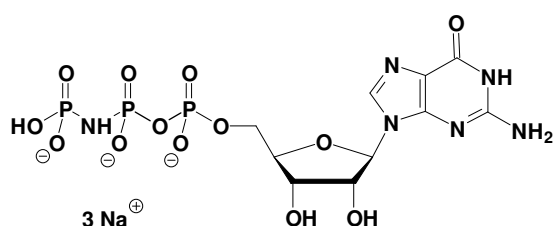


GppNHp (GMPPNP)

Guanosine-5'-[(β,γ)-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-401-10	10 mg
NU-401-50	50 mg



Cat. No.: NU-401

Molecular Formula: C₁₀H₁₄N₆O₁₃P₃ (Anion)

Molecular Weight: 519.17 (Anion)

Purity: > 95%

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!

Selected References:

Tarricone *et al.* (2001) The structural basis of Arfaptin-mediated cross-talk between Rac and Arf signalling pathways. *Nature* **411** (6834):215.

Esters *et al.* (2001) High-resolution crystal structure of *S. cerevisiae* Ypt51(Delta C15)-GppNHp, a small GTP-binding protein involved in regulation of endocytosis. *J. Mol. Biol.* **298** (1):111.

Chook *et al.* (1999) Structure of the nuclear transport complex karyopherin-beta 2-Ran center dot GppNHp. *Nature* **399** (6733):230.

Rudolph *et al.* (1999) Nucleotide binding to the G12V-mutant of Cdc42 investigated by x-ray diffraction and fluorescence spectroscopy: Two different nucleotide states in one crystal. *Protein Sci.* **8** (4):778.

Vetter *et al.* (1999) Structure of a Ran-binding domain complexed with Ran bound to a GTP analogue: implications for nuclear transport. *Nature* **398** (6722): 39.

Rittinger *et al.* (1997) Structure at 1.65 angstrom of RhoA and its GTPase-activating protein in complex with a transition-state analogue. *Nature* **389** (6652):758.

Rittinger *et al.* (1997) Crystal structure of a small G protein in complex with the GTPase-activating protein rhoGAP. *Nature* **388** (6643):693.

Hirshberg *et al.* The crystal structure of human rac1, a member of the rho-family complexed with a GTP analogue. *Nat. Struct. Biol.* **4** (2):147.

Nassar *et al.* (1995) The 2.2-Angstrom crystal-structure of the ras-binding domain of the serine threonine kinase c-Raf1 in complex with Rap1a and a GTP analog. *Nature* **375** (6532):554.

Tolkovsky (1980) 2'-deoxyadenosine functionally uncouples adenylate-cyclase from the guanyl nucleotide subunit without altering simultaneous gppnhp occupancy. *FEBS Lett.* **116** (2):165.