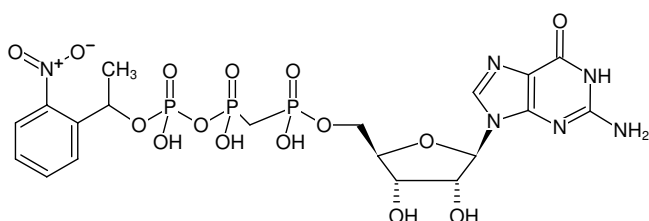


**NPE-caged-GpCpp**

(NPE-caged-GMPCPP)

Guanosine-5'-[(α,β)-methylene]triphosphate, P³-(1-(2-nitrophenyl)-ethyl)-ester, Triethylammonium salt

Cat. No.	Amount
NU-306S	10 μ l (10 mM)
NU-306L	5 x 10 μ l (10 mM)



Structural formula of NPE-caged-GpCpp

For research use only!**Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₁₉H₂₅N₆O₁₅P₃ (free acid)**Molecular Weight:** 670.35 g/mol (free acid)**Exact Mass:** 670.06 g/mol (free acid)**Purity:** \geq 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{\max} 254 nm, ϵ 16.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Derrick *et al.* (2012) The Nucleotide-Binding State of Microtubules Modulates Kinesin Processivity and Tau's Ability to Inhibit Kinesin Mediated Transport. *J. Biol. Chem.* **286** (50): 42873.Katsuki *et al.* (2011) Preparation of Dual-Color Polarity-Marked Fluorescent Microtubule Seeds *Methods. Mol. Biol.* **777**:117.Telley *et al.* (2011) Reconstitution and Quantification of Dynamic Microtubule End Tracking In Vitro Using TIRF Microscopy *Methods. Mol. Biol.* **777**:127.Kalisch *et al.* (2011) Force Generation by Dynamic Microtubules In Vitro *Methods. Mol. Biol.* **777**:147.Salvarelli *et al.* (2011) Independence between GTPase active sites in the Escherichia coli cell division protein FtsZ. *FEBS Lett.* **585** (24):3880.Khrapunovich-Baine *et al.* (2011) Hallmarks of molecular action of microtubule stabilizing agents: effects of epothilone B, ixabepilone, peloruside A, and laulimalide on microtubule conformation. *J. Biol. Chem.* **286** (13):11765.Antonio *et al.* (2011) Bacterial Tubulin Distinct Loop Sequences and Primitive Assembly Properties Support Its Origin from a Eukaryotic Tubulin Ancestor. *J. Biol. Chem.* **286** (22):19789.Yusuke Oguchi *et al.* (2011) The bidirectional depolymerizer MCAK generates force by disassembling both microtubule ends *NATURE CELL BIOLOGY*. **13** (7):846.Maurer SP *et al.* (2011) GTPyS microtubules mimic the growing microtubule end structure recognized by end-binding proteins (EBs) *Proc. Natl. Acad. Sci. U. S. A.* **108** (10):3988.Ryuzo Kawamura *et al.* (2011) Formation of motile assembly of microtubules driven by kinesins *Smart Mater. Struct.* **20** (2011): 124007.