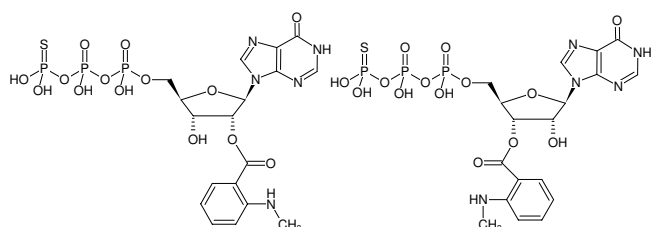




Mant-ITPyS

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

Cat. No.	Amount
NU-981	150 µl (10 mM)



Structural formula of Mant-ITPyS

For research use only!**Shipping:** shipped on dry ice**Storage Conditions:** store at -20 °C**Shelf Life:** 6 months after date of delivery**Molecular Formula:** C₁₈H₂₂N₅O₁₄P₃S (free acid)**Molecular Weight:** 657.37 g/mol (free acid)**Purity:** ≥ 90 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 279/355 nm, ε 20.5/5.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 355 nm, λ_{em} 448 nm**Selected References:**Seifert *et al.* (2012) Inhibitors of membranous adenylyl cyclases. *Trends Pharmacol. Sci.* **33** (2):64.Hübner *et al.* (2011) Effect of MANT-nucleotides on L-type calcium currents in murine cardiomyocytes. *J. Naunyn Schmiedebergs Arch. Pharmacol.* **383** (4):573.Erdorf *et al.* (2011) Pharmacological characterization of adenylyl cyclase isoforms in rabbit kidney membranes. *J. Naunyn Schmiedebergs Arch. Pharmacol.* **383** (4):357.Goettle *et al.* (2009) Characterization of Mouse Heart Adenylyl Cyclase. *J. Pharmacol. Exp. Ther.* **329** (3):1156.Goettle *et al.* (2007) Molecular analysis of the interaction of Bordetella pertussis adenylyl cyclase with fluorescent nucleotides. *Molecular Pharmacology* **72** (3):526.Gille *et al.* (2004) Differential Inhibition of Adenylyl Cyclase Isoforms and Soluble Guanylyl Cyclase by Purine and Pyrimidine Nucleotides. *J. Biol. Chem.* **279** (19):19955.