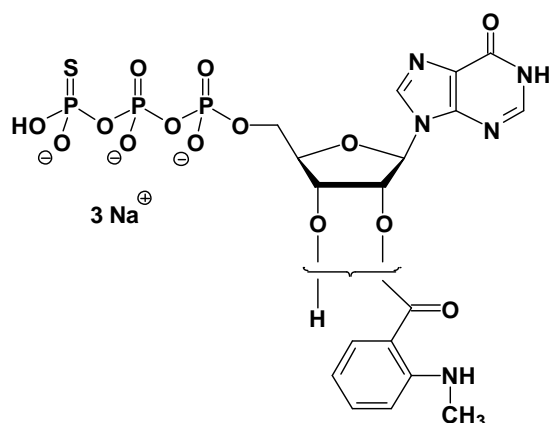


mant-ITP γ S

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

Cat. No.	Amount
NU-215S	10 Units
NU-215L	50 Units



Selected References:

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).

Gille *et al.* (2004) Differential Inhibition of Adenylyl Cyclase Isoforms and Soluble

Guanylyl Cyclase by Purine and Pyrimidine Nucleotides. *THE JOURNAL OF BIOLOGICAL CHEMISTRY* **279** (19):19955.

Cat. No.: NU-215

Molecular Formula: $C_{18}H_{19}N_5O_{14}P_3S$ (Anion)

Molecular Weight: 654.35 (Anion)

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

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!

1 unit = 1 μ l of a 10 mM solution