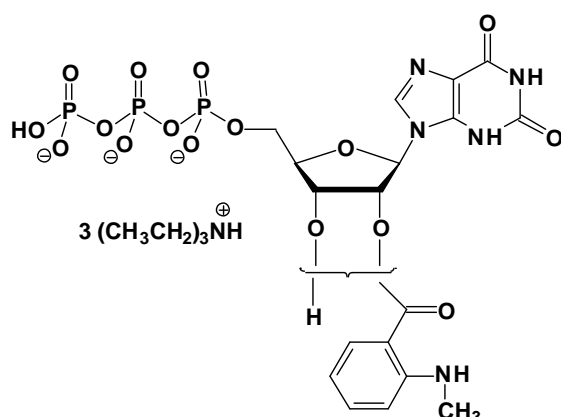


mant-XTP

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

Cat. No.	Amount
NU-213S	10 Units
NU-213L	50 Units



Selected References:

Tung-Chung Mou *et al.* (2005) Structural Basis for the Inhibition of Mammalian Membrane Adenylyl Cyclase by 2'(3')-O-(N-Methylanthraniloyl)- guanosine 5'-Triphosphate. *J. Biol. Chem.* **280** (8) :7253.

Gille *et al.* (2004) Xanthine nucleotide-specific G-protein α -subunits: a novel approach for the analysis of G-protein-mediated signal transduction *Naunyn-Schmiedeberg's Arch Pharmacol* **369** :141.

Aronsson *et al.* (2003) Unusual nucleotide-binding properties of the chloroplast protein import receptor, atToc33 *FEBS Letters* **544** :79.

Cat. No.: NU-213

Molecular Formula: C₁₈H₁₉N₅O₁₆P₃ (Anion)

Molecular Weight: 654.28 (Anion)

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

Spectroscopic properties:

λ_{\max} 254/355 nm; ϵ 19000/5800; λ_{exc} 355 nm;
 λ_{em} 448 nm

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 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution