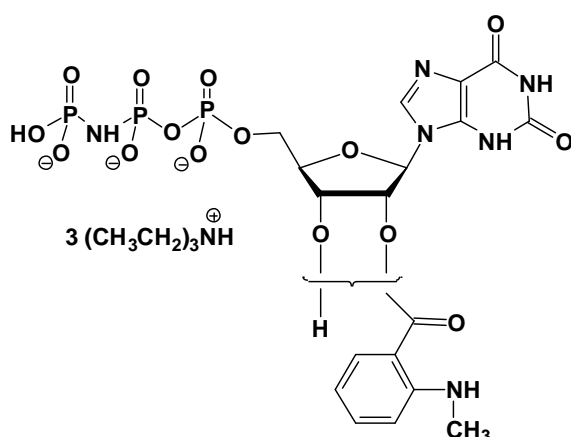


mant-XppNHp (mant-XMPPNP)

2'/3'-O-(N-Methyl-anthraniloyl)-xanthosine-5'-[(β,γ)-imido]triphosphate,
Triethylammonium salt

Cat. No.	Amount
NU-211S	5 Units
NU-211L	25 Units



Selected References:

Seifert (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* **269**:141.

Gille (2004) Differential Inhibition of Adenylyl Cyclase Isoforms and Soluble Guanylyl Cyclase by Purine and Pyrimidine Nucleotides. *THE JOURNAL OF BIOLOGICAL CHEMISTRY* **279**:19955.

Marlovits (2002) The membrane protein FeoB contains an intramolecular G protein essential for Fe(II) uptake in bacteria. *PNAS* **99**:16243.

Cat. No.: NU-211

Molecular Formula: C₁₈H₂₀N₆O₁₅P₃ (Anion)

Molecular Weight: 653.30 (Anion)

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

Spectroscopic properties:

λ_{\max} 254/355 nm; ϵ 19000/5800; λ_{exc} 355nm;
 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 6 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution