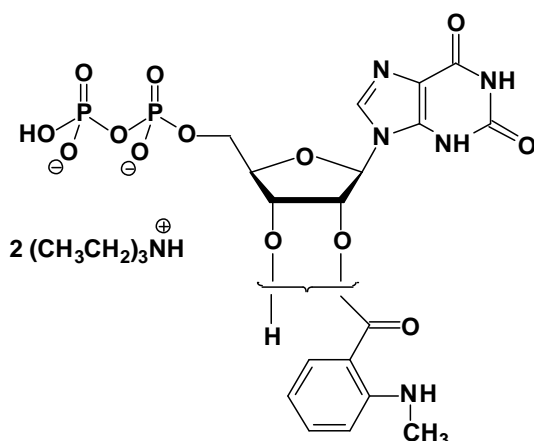


mant-XDP

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

Cat. No.	Amount
NU-210S	10 Units
NU-210L	50 Units



Selected References:

Gille (2003) Differential Inhibition of Adenylyl Cyclase Isoforms and Soluble Guanylyl Cyclase by Purine and Pyrimidine Nucleotides. *J. Biol. Chem.* **279**:19955.

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.* **369**:141.

Cat. No.: NU-210

Molecular Formula: C₁₈H₁₉N₅O₁₃P₂ (Anion)

Molecular Weight: 575.32 (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 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution