

**Mant-XDP**

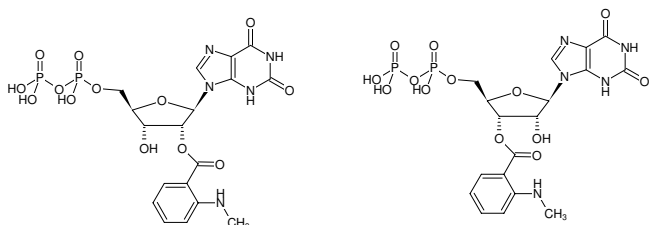
2'/3'-O-(N-Methyl-anthraniloyl)-xanthosine-5'-diphosphate, Triethylammonium salt

Cat. No.	Amount
NU-210S	10 µl (10 mM)
NU-210L	5 x 10 µl (10 mM)

Selected References:

Gille *et al.* (2004) Differential Inhibition of Adenylyl Cyclase Isoforms and Soluble Guanylyl Cyclase by Purine and Pyrimidine Nucleotides. *J. Biol. Chem.* **279**:19955.

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.



Structural formula of Mant-XDP

For research use only!**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₁₈H₂₁N₅O₁₃P₂ (free acid)**Molecular Weight:** 577.33 g/mol (free acid)**Exact Mass:** 577.06 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{\max} 254/355 nm, ϵ 19.0/5.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 355 nm, λ_{em} 448 nm