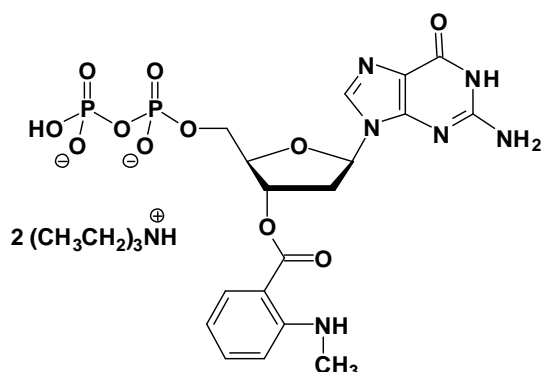


mant-dGDP

3'-(N-Methyl-anthraniloyl)-2'-deoxy-guanosine-5'-diphosphate, Triethylammonium salt

Cat. No.	Amount
NU-205S	30 Units
NU-205L	150 Units



Selected References:

Nomanbhoy *et al.* (1996) Investigation of the GTP-binding GTPase cycle of Cdc42hs using extrinsic reporter group fluorescence. *Biochemistry* **35** (14):4602.

Leonard *et al.* (1994) Investigation of the GTP-binding GTPase cycle of cdc42hs using fluorescence spectroscopy. *Biochemistry* **33** (40):12323.

Cat. No.: NU-205

Molecular Formula: C₁₈H₂₀N₆O₁₁P₂ (Anion)

Molecular Weight: 558.33 (Anion)

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

Spectroscopic properties:

λ_{\max} 252/355 nm; ϵ 22600/5700; 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