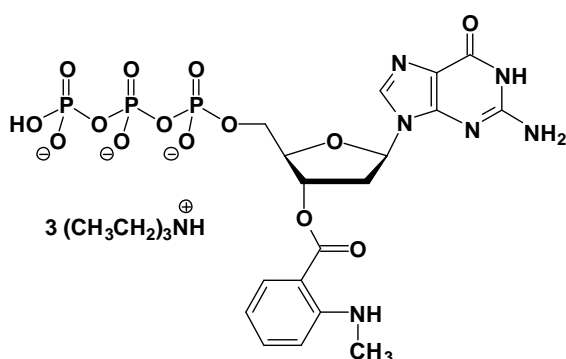


mant-dGTP

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

Cat. No.	Amount
NU-212S	10 Units
NU-212L	50 Units



Selected References:

Rodnina *et al.* (1995) Codon-dependent conformational change of elongation-factor tu preceding GTP hydrolysis on the ribosome *EMBO J.* **14 (11)**:2613.

Hazlett *et al.* (1993) Solution dynamics of p21(Ras) proteins bound with fluorescent nucleotides - a time-resolved fluorescence study. *Biochemistry-US* **32 (49)**:13575.

Cat. No.: NU-212

Molecular Formula: C₁₈H₂₀N₆O₁₄P₃ (Anion)

Molecular Weight: 637.30 (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