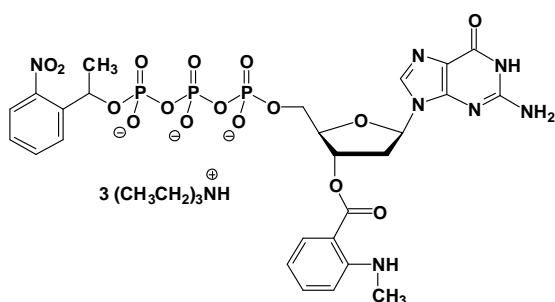


NPE-caged-mant-dGTP

3'-O-(N-Methyl-anthraniloyl)-2'-deoxy-guanosine-5'-triphosphate,
P³-(1-(2-nitrophenyl)-ethyl)-ester,

Cat. No.	Amount
NU-303S	10 Units
NU-303L	50 Units



Cat. No.: NU-303

Molecular Formula: C₂₆H₂₇N₇O₁₆P₃ (Anion)

Molecular Weight: 786.45 (Anion)

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

Spectroscopic properties:

λ_{\max} 253/355 nm; ϵ 25400/5700;

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. Store in a dark place and protect from light!

For research use only!

1 unit = 1 μ l of a 10 mM solution