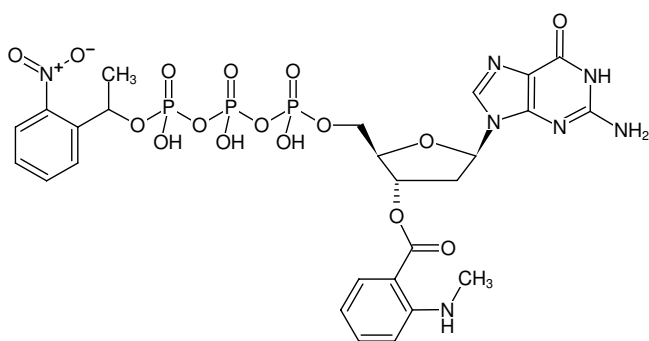


**NPE-caged-Mant-dGTP**

3'-O-(N-Methyl-anthraniloyl)-2'-deoxyguanosine-5'-triphosphate,  
P<sup>3</sup>-(1-(2-nitro-phenyl)-ethyl)-ester, Triethylammonium salt

Cat. No.	Amount
NU-303S	10 µl (10 mM)
NU-303L	5 x 10 µl (10 mM)



Structural formula of NPE-caged-Mant-dGTP

**For research use only!**

**Shipping:** shipped on blue ice

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** store dark

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

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>26</sub>H<sub>30</sub>N<sub>7</sub>O<sub>16</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 789.48 g/mol (free acid)

**Exact Mass:** 789.10 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** colorless to slightly yellow solution in water

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 253/355 nm, ε 25.4/5.7 L mmol<sup>-1</sup> cm<sup>-1</sup>  
(Tris-HCl pH 7.5), λ<sub>exc</sub> 355 nm, λ<sub>em</sub> 448 nm