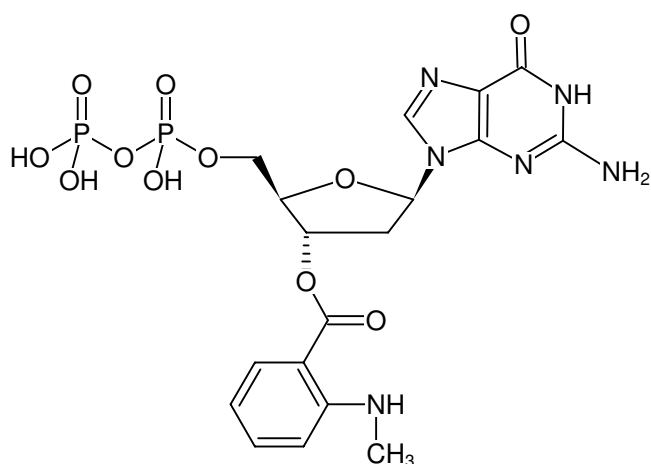


**Mant-dGDP**

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

Cat. No.	Amount
NU-205S	30 µl (10 mM)
NU-205L	5 x 30 µl (10 mM)



Structural formula of Mant-dGDP

**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:** C<sub>18</sub>H<sub>22</sub>N<sub>6</sub>O<sub>11</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 560.35 g/mol (free acid)**Exact Mass:** 560.08 g/mol (free acid)**CAS#:** 127383-33-9**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<sub>max</sub> 252/355 nm, ε 22.6/5.7 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**Applications:**Resonance energy transfer to Trp in protein cdc42hs<sup>[1]</sup>Conformational analysis of Ras-like protein<sup>[2]</sup>**Specific Ligands:**Ras-like GTP-binding protein Cdc42Hs<sup>[1, 2]</sup>**Selected References:**[1] Leonard *et al.* (1994) Investigation of the GTP-binding GTPase cycle of cdc42hs using fluorescence spectroscopy. *Biochemistry* **33** (40):12323.[2] Nomanbhoy *et al.* (1996) Investigation of the GTP-binding GTPase cycle of Cdc42Hs using extrinsic reporter group fluorescence. *Biochemistry* **35** (14):4602.Liu *et al.* (2008) Ras Is Required for the Cyclic AMP-Dependent Activation of Rap1 via Epac2. *Molecular and cellular biology* **28** (23):7109.