

**Mant-cGMP**

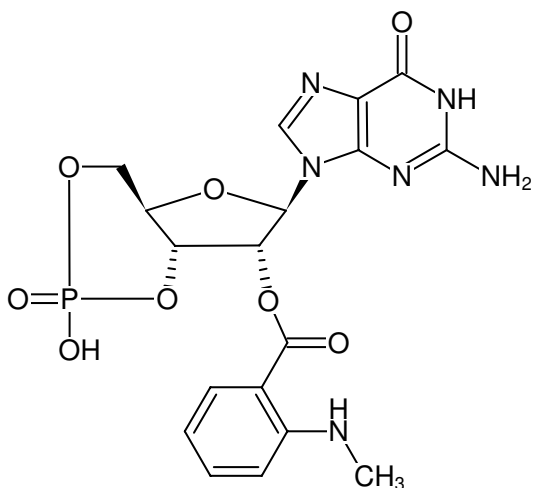
2'-O-(N-Methyl-anthraniloyl)-guanosine-3',5'-cyclic monophosphate, Sodium salt

Cat. No.	Amount
NU-240S	10 $\mu\text{mol}$
NU-240L	5 x 10 $\mu\text{mol}$

**Selected References:**

Newton *et al.* (2010) A real-time fluorescent assay of the purified nitric oxide receptor, guanylyl cyclase. *Anal. Biochem.* **402** (2):129.

Johnson *et al.* (1987) A continuous fluorescence assay for cyclic nucleotide phosphodiesterase hydrolysis of cyclic GMP. *Anal. Biochem.* **162** (1):291.



Structural formula of Mant-cGMP

**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>19</sub>N<sub>6</sub>O<sub>8</sub>P (free acid)**Molecular Weight:** 478.35 g/mol (free acid)**Exact Mass:** 478.10 g/mol (free acid)**CAS#:** 83707-15-7**Purity:**  $\geq$  98 % (HPLC)**Form:** solid**Color:** colorless to slightly yellow**Spectroscopic Properties:**  $\lambda_{\text{max}}$  252/355 nm,  $\epsilon$  22.6/5.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5),  $\lambda_{\text{exc}}$  355 nm,  $\lambda_{\text{em}}$  448 nm