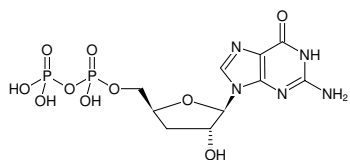


**3'-dGDP**

3'-Deoxyguanosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1144S	20 µl (10 mM)
NU-1144L	5 x 20 µl (10 mM)



Structural formula of 3'-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>10</sub>H<sub>15</sub>N<sub>5</sub>O<sub>10</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 427.20 g/mol (free acid)**Exact Mass:** 427.03 g/mol (free acid)**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 nm, ε 13.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Specific Ligands:**Affinity to RNA polymerase I and II<sup>[1]</sup>**Selected References:**[1] Shunji Izuta *et al.* (1996) 3'-Deoxyribonucleotides inhibit Eukaryotic DNA Primase1. *J. Biochem.* **119** (6):1038.Rasmussen *et al.* (1993) In vivo transcriptional pausing and cap formation on three Drosophila heat shock genes. *Proc. Natl. Acad. Sci. USA* **90**:7923.Hamel *et al.* (1976) Interactions of Guanosine Triphosphate Analogues with Elongation Factor G of Escherichia coli. *J. Biochem.* **63**:13894.