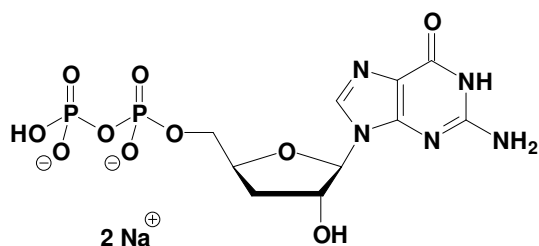


3'-dGDP

3'-Deoxy-Guanosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1144S	20 Units
NU-1144L	100 Units



Cat. No.: NU-1144

Molecular Formula: C₁₀H₁₃N₅O₁₀P₂ (Anion)

Molecular Weight: 425.18 (Anion)

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

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.

* 1 unit = 1 µl of a 10 mM solution

For research use only!

Selected References:

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.