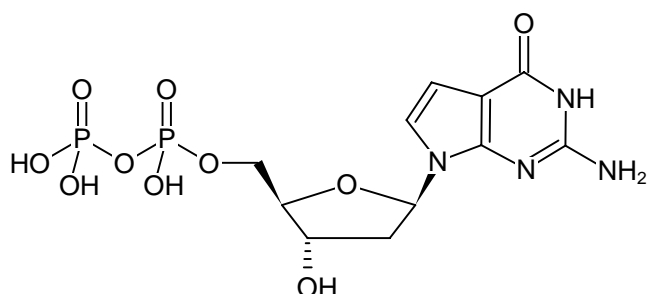


**7-Deaza-dGDP**

7-Deaza-2'-deoxyguanosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1178S	100 µl (10 mM)
NU-1178L	5 x 100 µl (10 mM)



Structural formula of 7-Deaza-dGDP

Selected References:

Mizusawa *et al.* (1986) Improvement of the dideoxy chain termination method of DNA sequencing by use of deoxy-7-deazaguanosine triphosphate in place of dGTP. *Nucleic Acids Research* **14** (3):1319.

Barr *et al.* (1986) 7-Deaza-2'-Deoxyguanosine-5'-Triphosphate: Enhanced Resolution in M13 Dideoxy Sequencing. *BioTechniques* **4**:428.

For research use only!**Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months**Molecular Formula:** C₁₁H₁₆N₄O₁₀P₂ (free acid)**Molecular Weight:** 426.21 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution, pH 7.5 ± 0.5**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{abs} 259 nm; ε 12.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)