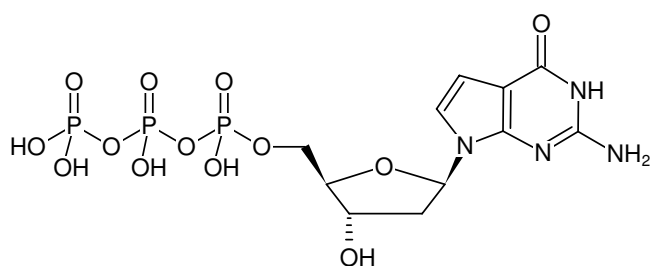


**7-Deaza-dGTP**

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

Cat. No.	Amount
NU-1179S	10 µl (100 mM)
NU-1179L	5 x 10 µl (100 mM)



Structural formula of 7-Deaza-dGTP

For general laboratory use.**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:** C₁₁H₁₇N₄O₁₃P₃ (free acid)**Molecular Weight:** 506.19 g/mol (free acid)**Exact Mass:** 506.00 g/mol (free acid)**CAS#:** 101515-08-6**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 259 nm, ε 12.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Pivonková *et al.* (2010) Direct voltammetric analysis of DNA modified with enzymatically incorporated 7-deazapurines. *Anal Chem.* **82** (16):6807.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.