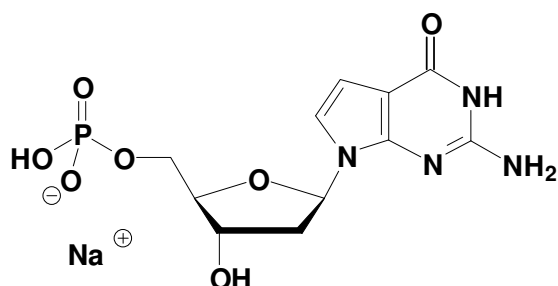


7-Deaza-dGMP

7-Deaza-2'-Deoxy-Guanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1177S	100 Units
NU-1177L	500 Units



Cat. No.: NU-1177

Molecular Formula: C₁₁H₁₄N₄O₇P (Anion)

Molecular Weight: 345.22 (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:

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.