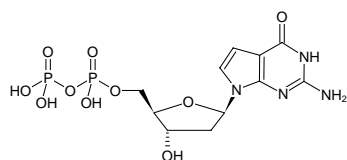


**7-Deaza-dGDP**

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

**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.

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>11</sub>H<sub>16</sub>N<sub>4</sub>O<sub>10</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 426.21 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>abs</sub> 259 nm; ε 12.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**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.