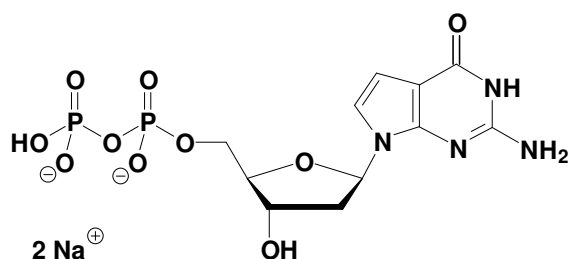


## 7-Deaza-dGDP

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

Cat. No.	Amount
NU-1178S	100 Units
NU-1178L	500 Units



**Cat. No.:** NU-1178

**Molecular Formula:** C<sub>11</sub>H<sub>14</sub>N<sub>4</sub>O<sub>10</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 424.20 (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.