

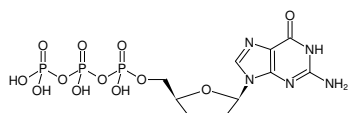
**ddGTP**

2',3'-Dideoxyguanosine-5'-triphosphate, Lithium salt

Cat. No.	Amount
NU-1017S	100 µl (10 mM)
NU-1017L	5 x 100 µl (10 mM)

**Selected References:**

Sanger *et al.* (1977) DNA sequencing with chain-terminating inhibitors. *Proc. Natl. Acad. Sci. USA* **74**:5463.



Structural formula of ddGTP

**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:** C<sub>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>12</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 491.18 g/mol (free acid)**Exact Mass:** 491.00 g/mol (free acid)**Purity:** ≥ 98 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 252 nm, ε 14.2 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)