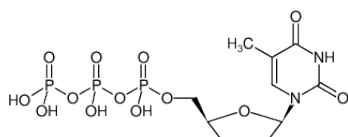




## ddTTP - high concentration

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

Cat. No.	Amount
NU-294S	10 µl (100 mM)
NU-294L	5 x 10 µl (100 mM)



Structural formula of ddTTP - high concentration

### 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>17</sub>N<sub>2</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 466.17 g/mol (free acid)

**Exact Mass:** 465.99 g/mol (free acid)

**Purity:** ≥ 98 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 100 mM - 110 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 267 nm, ε 9.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

### Selected References:

Koji *et al.* (2008) In situ detection of methylated DNA by histo endonuclease-linked detection of methylated DNA sites: a new principle of analysis of DNA methylation. *Histochem Cell Biol* **130**:917.

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