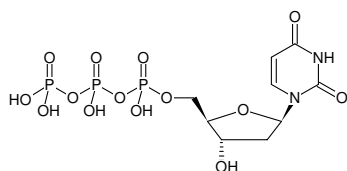


**dUTP - Solution**

100 mM Sodium salt solution

Cat. No.	Amount
NU-1008L	1 ml (100 mM)
NU-1008-10ML	10 ml (100 mM)



Structural formula of dUTP - Solution

**For in vitro 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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**Shelf Life:** 12 months**Molecular Formula:** C<sub>9</sub>H<sub>15</sub>N<sub>2</sub>O<sub>14</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 468.14 g/mol (free acid)**CAS#:** 102814-08-4**Purity:** ≥ 99 % (HPLC)**Form:** clear aqueous solution**Concentration:** 100 mM - 110 mM**pH:** 8.5 ± 0.2 (22 °C)**Spectroscopic Properties:** λ<sub>max</sub> 262 nm, ε 9.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.0)**Description:**

dTTP, PCR-grade is supplied as ultrapure aqueous solution (pH 8.5) and suitable for all molecular biology applications including PCR/qPCR, reverse transcription, DNA labeling and DNA sequencing.

**Selected References:**

Tóth *et al.* (2007) Kinetic Mechanism of Human dUTPase, an Essential Nucleotide Pyrophosphatase Enzyme. *J. Biol. Chem.* **282** (46):33572.

Erllich *et al.* (1988) Primer-directed enzymatic amplification of DNA with a thermostable DNA polymerase. *Science* **29** (239):487.

Holland *et al.* (1991) Detection of specific polymerase chain reaction product by utilizing the 5'→3' exonuclease activity of *Thermus aquaticus* DNA polymerase. *Proc. Natl. Acad. Sci. USA* **88** (16):7276.

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