

ddNTP bundle

4x 10 mM (ddATP, ddCTP, ddGTP, ddTTP)

Cat. No.	Amount
NU-1019S	4x 200 μ l (4x 2 μ mol)
NU-1019L	4x 1 ml (4x 10 μ mol)

For *in vitro* use only

Quality guaranteed for 12 months

Store at -20°C, short term (up to one week) exposure to ambient temperature possible

Quality control

The dideoxynucleotide solutions have a purity of >99% and are functionally tested by chain termination sequencing reactions

Description

ddNTP Bundle contains four separate solutions of ultra-pure ddATP, ddCTP, ddGTP, and ddTTP supplied as clear aqueous solutions (pH 8.4-8.6).

ddATP

2',3'-Dideoxy-adenosine-5'-triphosphate, sodium salt

Molecular formula: C₁₀H₁₃N₅O₁₁P₃ (Anion)

Molecular weight: 472.16 (Anion)

ddCTP

2',3'-Dideoxy-cytidine-5'-triphosphate, sodium salt

Molecular formula: C₉H₁₃N₃O₁₂P₃ (Anion)

Molecular weight: 448.13 (Anion)

ddGTP

2',3'-Dideoxy-guanosine-5'-triphosphate, sodium salt

Molecular formula: C₁₀H₁₃N₅O₁₂P₃ (Anion)

Molecular weight: 488.16 (Anion)

ddTTP

2',3'-Dideoxy-thymidine-5'-triphosphate, sodium salt

Molecular formula: C₁₀H₁₄N₂O₁₃P₃ (Anion)

Molecular weight: 463.14 (Anion)

Selected References:

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