

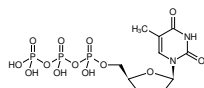
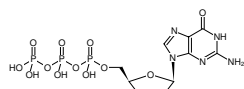
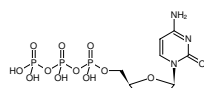
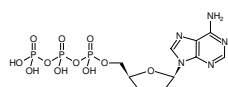


ddNTP Bundle

ddATP, ddCTP, ddGTP and ddTTP 10 mM solutions

2',3'-Dideoxyadenosine-5'-triphosphate, Lithium salt; 2',3'-Dideoxycytidine-5'-triphosphate, Lithium salt; 2',3'-Dideoxyguanosine-5'-triphosphate, Lithium salt; 2',3'-Dideoxythymidine-5'-triphosphate, Lithium salt

Cat. No.	Amount
NU-1019S	4 x 100 µl (10 mM)
NU-1019L	4 x 500 µl (10 mM)



Structural formula of ddNTP Bundle

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:

ddATP: C₁₀H₁₆N₅O₁₁P₃ (free acid)

ddCTP: C₉H₁₆N₃O₁₂P₃ (free acid)

ddGTP: C₁₀H₁₆N₅O₁₂P₃ (free acid)

ddTTP: C₁₀H₁₇N₂O₁₃P₃ (free acid)

Molecular Weight:

ddATP: 475.18 g/mol (free acid)

ddCTP: 451.16 g/mol (free acid)

ddGTP: 491.18 g/mol (free acid)

ddTTP: 466.17 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:

ddATP: λ_{max} 259 nm, ε 15.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

ddCTP: λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

ddGTP: λ_{max} 252 nm, ε 14.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

ddTTP: λ_{max} 267 nm, ε 9.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



ddNTP Bundle

ddATP, ddCTP, ddGTP and ddTTP 10 mM solutions

2',3'-Dideoxyadenosine-5'-triphosphate, Lithium salt; 2',3'-Dideoxycytidine-5'-triphosphate, Lithium salt; 2',3'-Dideoxyguanosine-5'-triphosphate, Lithium salt; 2',3'-Dideoxythymidine-5'-triphosphate, Lithium salt

Description:

ddNTP Bundle contains four separate solutions of ddATP, ddCTP, ddGTP, and ddTTP at a concentration of 10-11 mM supplied as clear aqueous solutions (pH 7.5 ±0.5).

ddATP (#NU-1015)

ddCTP (#NU-1016)

ddGTP (#NU-1017)

ddTTP (#NU-1018)

Selected References:

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