



dNTP Mix dUTP - Solution

Premix of 10 mM dATP, dCTP, dGTP and 20 mM dUTP

2'-Deoxyadenosine-5'-triphosphate, Sodium salt; 2'-Deoxycytidine-5'-triphosphate, Sodium salt; 2'-Deoxyguanosine-5'-triphosphate, Sodium salt; 2'-Deoxyuridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1020S	1 ml
NU-1020L	5x 1 ml

Description:

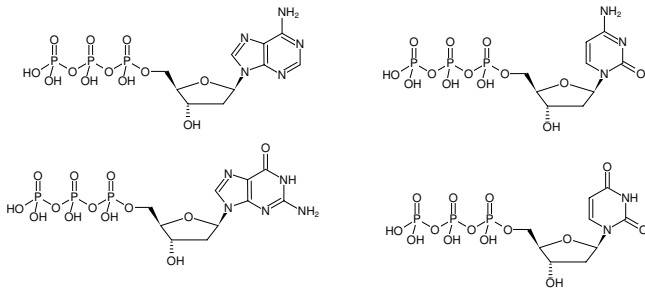
dNTP Mix incl. dUTP is a mixture of 10 mM ultrapure dATP, dCTP, and dGTP, and 20 mM dUTP supplied as clear aqueous solution (pH 8.5). dUTP can be used in place of dTTP in PCR and RT-PCR protocols to prevent carry-over contaminations from previous amplifications.

Selected References:

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.



Structural formula of dNTP Mix dUTP - Solution

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Shelf Life: 12 months

Molecular Formula:

dATP: C₁₀H₁₆N₅O₁₂P₃ (free acid)

dCTP: C₉H₁₆N₃O₁₃P₃ (free acid)

dGTP: C₁₀H₁₆N₅O₁₃P₃ (free acid)

dUTP: C₉H₁₅N₂O₁₄P₃ (free acid)

Molecular Weight:

dATP: 491.18 g/mol (free acid)

dCTP: 467.15 g/mol (free acid)

dGTP: 507.18 g/mol (free acid)

dUTP: 468.14 g/mol (free acid)

Purity: ≥ 99 % (HPLC)

Form: clear aqueous solution

pH: 8.5 ± 0.2 (22 °C)