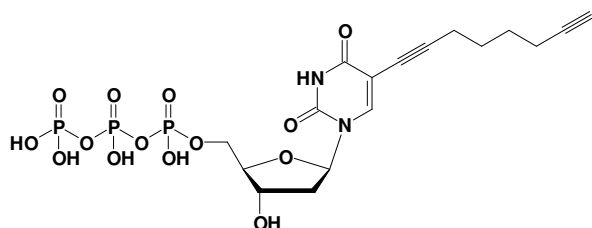


C8-Alkyne-dUTP, 100 mM

5-(Octa-1,7-diynyl)-2'-deoxyuridine 5'-triphosphate, Sodium salt

Cat. No.	Amount
CLK-T05-S	5 μ l (0,5 μ mol)
CLK-T05-M	25 μ l (2,5 μ mol)
CLK-T05-L	50 μ l (5 μ mol)



Cat. No.: CLK-T05

Molecular formula: C₁₇H₂₃N₂O₁₄P₃ (free acid)

Molecular weight: 572.29 (free acid)

Spectroscopic properties:

$\epsilon_{260\text{nm}}$ 3900 cm⁻¹ M⁻¹ (in water)

$\epsilon_{292\text{nm}}$ 11000 cm⁻¹ M⁻¹ (in water)

Purity: > 99%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

Selected References:

Carell *et al.* (2007) Synthesis of Highly Modified DNA by a Combination of PCR with Alkyne-Bearing Triphosphates and Click Chemistry. *Chem. Eur. J.* **13**:9486.

Carell *et al.* (2006) Directed DNA Metallization. *J. Am. Chem. Soc.* **128** (5):1398.