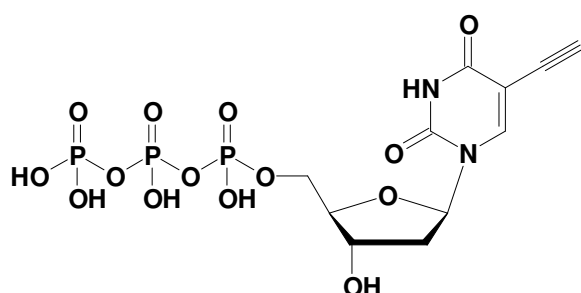


5-Ethynyl-dUTP (5-EdUTP), 100 mM

5-Ethynyl-2'-deoxyuridine 5'-triphosphate, Sodium salt

Cat. No.	Amount
CLK-T07-S	5 μ l (0,5 μ mol)
CLK-T07-M	25 μ l (2,5 μ mol)
CLK-T07-L	50 μ l (5 μ mol)



Cat. No.: CLK-T07

Molecular formula: $C_{11}H_{15}N_2O_{14}P_3$ (free acid)

Molecular weight: 492.16 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic Properties: λ_{max} 288nm;
 ϵ 12,000 $M^{-1}cm^{-1}$ Tris HCl 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!

Applications:

Reaction with profluorogenic 3-azido-7-hydroxycoumarin^[1]

Measurement of cells in S-phase^[1]

Selected References:

[1] Kai *et al.* (2010) Fluorogenic "click" reaction for labeling and detection of DNA in proliferating cells. *Biotechniques* **49(1)**:525.

Chehrehasa *et al.* (2009) EdU, a new thymidine analogue for labelling proliferating cells in the nervous system. *J. Neurosci. Methods* **177**:122.

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.