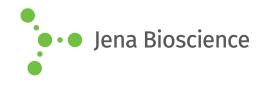
DATA SHEET





■ 5-Ethynyl-dUTP (5-EdUTP)

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

Cat. No.	Amount
CLK-T07-S	5 μl (100 mM)
CLK-T07-L	5 x 5 μl (100 mM)
CLK-T07-XL	50 μl (100 mM)

Structural formula of 5-Ethynyl-dUTP (5-EdUTP)

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: C₁₁H₁₅N₂O₁₄P₃ (free acid)
Molecular Weight: 492.16 g/mol (free acid)
Exact Mass: 491.97 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow **Concentration:** 100 mM - 110 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 288 nm, ϵ 12.0 L mmol $^{\text{-1}}$ cm $^{\text{-1}}$ (Tris-HCl

pH 7.5)

Applications:

Incorporation into DNA by **PCR** with family B polymerases (Pwo, Deep Vent exo^- or KOD XL)^[1,2]

Incorporation into DNA by **Primer Extension** with family A polymerase Taq and family B polymerases (Pwo, Deep Vent exo^- or KOD $XL)^{[1]}$

The resulting ethynyl-functionalized DNA can subsequently be processed via Cu(I)-catalyzed click chemistry that offers the choice

- to introduce a Biotin group (via Azides of Biotin) for subsequent purification tasks
- to introduce fluorescent group (via Azides of fluorescent dyes) for subsequent microscopic imaging
- to crosslink the DNA to Azide-functionalized biomolecules e.g.proteins

Presolski et $al.^{[3]}$ and Hong et $al.^{[4]}$ provide a general protocol for Cu(I)-catalyzed click chemistry reactions that may be used as a starting point for the set up and optimization of individual assays.

Related Products:

Copper (II)-Sulphate (CuSO₄), #CLK-MI004 Tris(3-hydroxypropyltriazolylmethyl)amine (THPTA), #CLK-1010 Sodium Ascorbate (Na-Ascorbate), #CLK-MI005

Selected References:

[1] Gierlich *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.

[2] Burley *et al.* (2006) Directed DNA Metallization. *J. Am. Chem. Soc.* **128 (5)**:1398.

[3] Presolski et al. (2011) Copper-Catalyzed Azide-Alkyne Click Chemistry for Bioconjugation. Current Protocols in Chemical Biology **3**:153.

[4] Hong et al. (2011) Analysis and Optimization of Copper-Catalyzed Azide-Alkyne Cycloaddition for Bioconjugation. Angew. Chem. Int. Ed. **48**:9879.