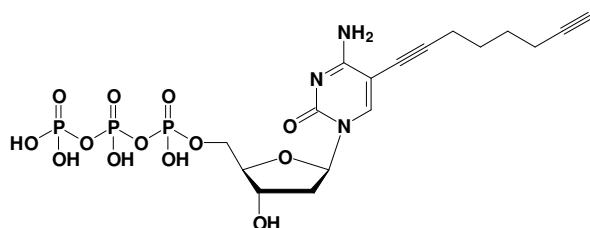


## C8-Alkyne-dCTP, 100 mM

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

Cat. No.	Amount
CLK-T06-S	5 $\mu$ l (0,5 $\mu$ mol)
CLK-T06-M	25 $\mu$ l (2,5 $\mu$ mol)
CLK-T06-L	50 $\mu$ l (5 $\mu$ mol)



**Cat. No.:** CLK-T06

**Molecular formula:** C<sub>17</sub>H<sub>24</sub>N<sub>3</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

**Molecular weight:** 571.31 (free acid)

**Spectroscopic properties:**

$\epsilon_{260\text{nm}}$  5000 cm<sup>-1</sup> M<sup>-1</sup> (in water)

$\epsilon_{292\text{nm}}$  9500 cm<sup>-1</sup> M<sup>-1</sup> (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.