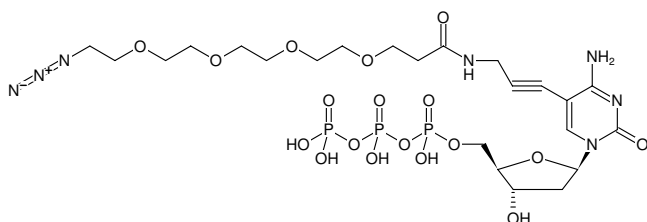


**5-Azido-PEG<sub>4</sub>-dCTP**5-Azido-PEG<sub>4</sub>-2'-deoxycytidine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
CLK-070S	50 µl (10 mM)
CLK-070L	5 x 50 µl (10 mM)

Structural formula of 5-Azido-PEG<sub>4</sub>-dCTP**For research use only!****Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** avoid freeze/thaw cycles

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>23</sub>H<sub>38</sub>N<sub>7</sub>O<sub>18</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 793.51 g/mol (free acid)**Exact Mass:** 793.15 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** sterile solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 294 nm, ε 9.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)