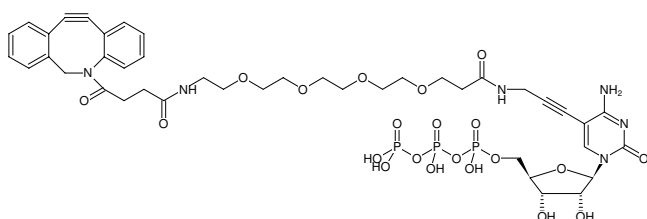


**5-DBCO-PEG<sub>4</sub>-CTP**5-Dibenzylcyclooctyl-PEG<sub>4</sub>-cytidine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
CLK-056S	20 µl (10 mM)
CLK-056L	5 x 20 µl (10 mM)

Structural formula of 5-DBCO-PEG<sub>4</sub>-CTP**For general laboratory use.****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>42</sub>H<sub>53</sub>N<sub>6</sub>O<sub>21</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 1070.82 g/mol (free acid)**Exact Mass:** 1070.25 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** filtered solution in 100 mM Tris-HCl**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 308 nm, ε 13.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)