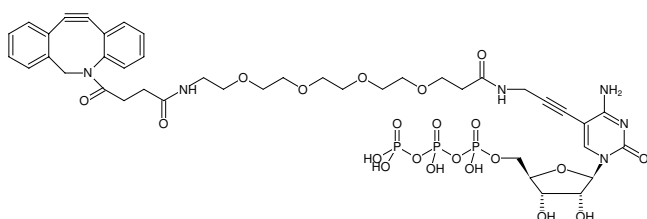


**5-DBCO-PEG₄-CTP**5-Dibenzylcyclooctyl-PEG₄-cytidine-5'-triphosphate, Triethylammonium salt

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

Structural formula of 5-DBCO-PEG₄-CTP**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:** C₄₂H₅₃N₆O₂₁P₃ (free acid)**Molecular Weight:** 1070.82 g/mol (free acid)**Exact Mass:** 1070.25 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** sterile solution in 100 mM Tris-HCl**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 308 nm, ε 13.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)