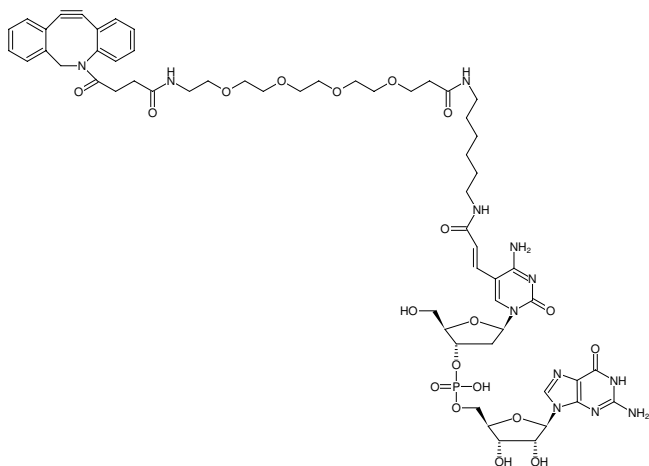


**5-DBCO-PEG₄-dCpG**

Triethylammonium salt

Cat. No.	Amount
CLK-062	1 μ mol

Structural formula of 5-DBCO-PEG₄-dCpG**For research use only!****Shipping:** shipped on blue ice**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:** C₅₈H₇₅N₁₂O₁₉P (free acid)**Molecular Weight:** 1275.26 g/mol (free acid)**Purity:** \geq 95 % (HPLC)**Form:** solid**Solubility:** DMSO, very slightly soluble in water and MeOH**Spectroscopic Properties:** λ_{max} 308 nm; ϵ 13.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)