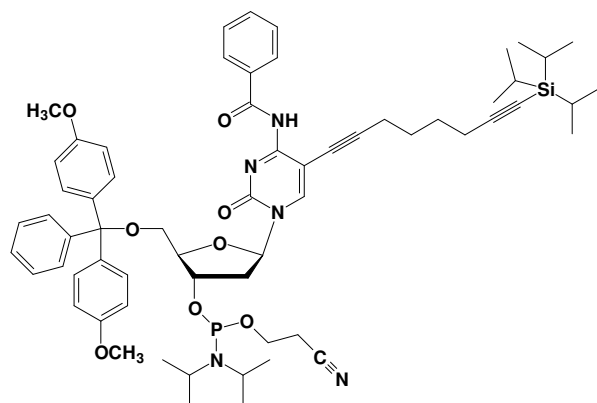


## C8-TIPS-dC-CEP

4-N-Benzoyl-5-(8-triisopropylsilyl-octa-1,7-diynyl)-3'-O-[(2-cyanoethoxy)(diisopropylamino)-phosphono]-5'-O-(4,4'-dimethoxytrityl)-2'-deoxycytidine

Cat. No.	Amount
CLK-A08-250MG	250 mg
CLK-A08-1G	1 g
CLK-A08-5G	5g



**Molecular formula:** C<sub>63</sub>H<sub>80</sub>N<sub>5</sub>O<sub>8</sub>P

**Molecular weight:** 1094.40 g/mol

**Storage conditions:** store at -20 °C, store dry and under inert gas

**Purity:** > 99 %

**Appearance:** off-white powder

**Shelf life:** 12 months

**Please note:**

**Commercial use of this product requires a license. Please contact us at [info@jenabioscience.com](mailto:info@jenabioscience.com) for more detailed information.**

***For research use only!***