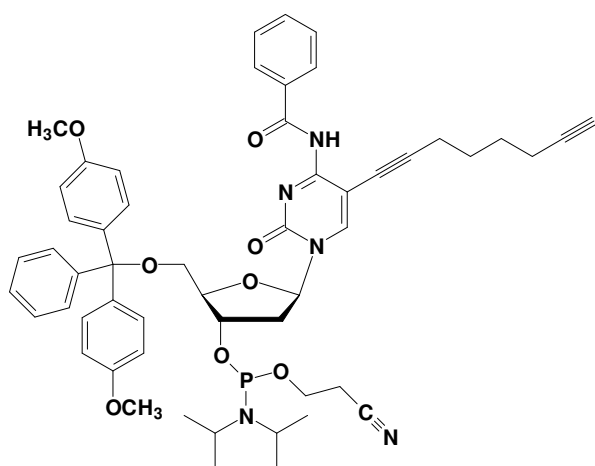


C8-Alkyne-dC-CEP

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

Cat. No.	Amount
CLK-A07-250MG	250 mg
CLK-A07-1G	1 g
CLK-A07-5G	5g



Molecular formula: C₅₄H₆₀N₅O₈P

Molecular weight: 938.06 g/mol

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

Purity: > 99.5 %

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 for more detailed information.

For research use only!