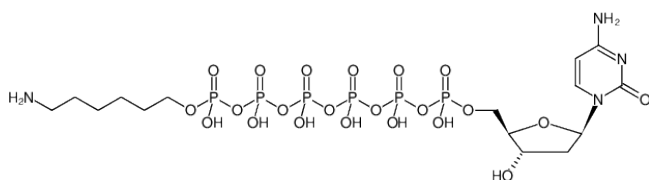




zeta-(6-Aminohexyl)-dC6P

zeta-(6-Aminohexyl)-2'-deoxycytidine-5'-hexaphosphate, Sodium salt

Cat. No.	Amount
NU-283	50 µl (10 mM)



Structural formula of zeta-(6-Aminohexyl)-dC6P

For general laboratory use.

Shipping: shipped on gel packs

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: C₁₅H₃₂N₄O₂₂P₆ (free acid)

Molecular Weight: 806.27 g/mol (free acid)

Exact Mass: 805.99 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)