

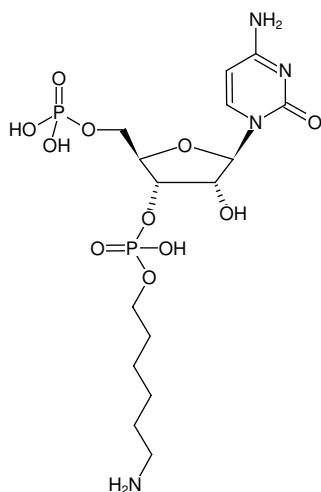
**pCp-Amine**

Cytidine-5'-phosphate-3'-(6-aminohexyl)phosphate, Sodium salt

Cat. No.	Amount
NU-1706	20 µl (1 mM)

Related Products:

pCp-BIO, #NU-1706-BIO
 pCp-Desthiobiotin, #NU-1706-Desthiobio
 pCp-Cy3, #NU-1706-CY3
 pCp-Cy5, #NU-1706-CY5
 pCp-Azide, #NU-1708
 pCp-Alkyne, #NU-1709



Structural formula of pCp-Amine

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:** C₁₅H₂₈N₄O₁₁P₂ (free acid)**Molecular Weight:** 502.35 g/mol (free acid)**Exact Mass:** 502.12 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 1.0 mM - 1.1 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)