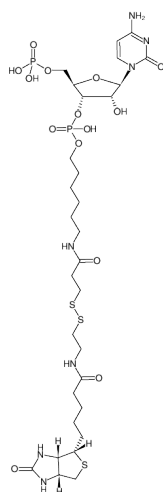


**pCp-SS-Biotin**

Cytidine-5'-phosphate-3'-(6-aminohexyl)phosphate, labeled with a cleavable disulfide linker containing Biotin, Triethylammonium salt

Cat. No.	Amount
NU-1706-SS-BIO	20 µl (1 mM)



Structural formula of pCp-SS-Biotin

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: C₃₀H₅₁N₇O₁₄P₂S₃ (free acid)

Molecular Weight: 891.90 g/mol (free acid)

Exact Mass: 891.21 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

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)