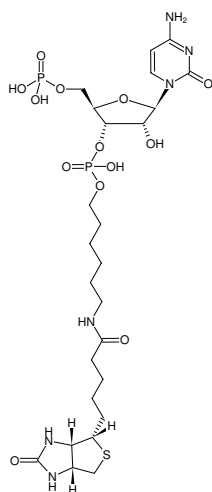




pCp-Biotin

Cytidine-5'-phosphate-3'-(6-aminohexyl)phosphate, labeled with Biotin, Triethylammonium salt

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



Structural formula of pCp-Biotin

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: C₂₅H₄₂N₆O₁₃P₂S₁ (free acid)

Molecular Weight: 728.65 g/mol (free acid)

Exact Mass: 728.20 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution in 10 mM Tris-HCl

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)

Applications:

Incorporation into RNA/miRNA by

- 3'-End Labeling with T4 RNA Ligase

Description:

pCp-Biotin is recommended for 3'-End Labeling of RNA/miRNA with T4 RNA Ligase that conjugates a single Biotin molecule to the 3'-OH group of the RNA/miRNA template. The resulting Biotin-labeled RNA/miRNA probes are ideally suited for hybridization experiments such as microarray-based expression profiling or electro mobility shift assays (EMSA). Detection of the Biotin moiety can be performed using streptavidin conjugated with horseradish peroxidase (HRP), alkaline phosphatase (AP) or a fluorescent dye. Alternatively, Biotin-labeled RNA/miRNA probes can also be purified with Streptavidin-conjugated agarose/magnetic beads.

Related Products:

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

Selected References:

Kashida *et al.* (2018) Nanoparticle-based local translation reveals mRNA as translation-coupled scaffold with anchoring function. [doi:10.1101/483727](https://doi.org/10.1101/483727).

Booy *et al.* (2016) RNA Helicase Associated with AU-rich Element (RHAU/DHX36) Interacts with the 3'-Tail of the Long Non-coding RNA BC200 (BCYRN1). *J. Biol. Chem.* **291**(10):5355.