



## pCp-Desthiobiotin

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

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

### Applications:

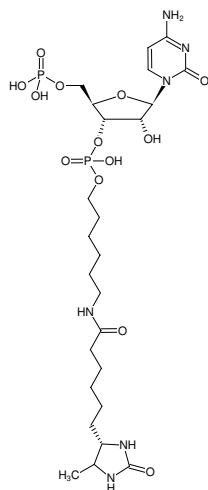
- Incorporation into RNA/miRNA by
- 3'-End Labeling with T4 RNA Ligase

### Description:

pCp-Desthiobiotin is recommended for 3'-End Labeling of RNA/miRNA with T4 RNA Ligase that conjugates a single Desthiobiotin molecule to the 3'-OH group of the RNA/miRNA template. The resulting Desthiobiotin-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 Desthiobiotin moiety can be performed using streptavidin conjugated with horseradish peroxidase (HRP), alkaline phosphatase (AP) or a fluorescent dye. Alternatively, Desthiobiotin-labeled RNA/miRNA probes can also be purified with Streptavidin-conjugated agarose/magnetic beads.

### Related Products:

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



Structural formula of pCp-Desthiobiotin

### For research use only!

**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:** C<sub>25</sub>H<sub>44</sub>N<sub>6</sub>O<sub>13</sub>P<sub>2</sub> (free acid)

**Molecular Weight:** 698.60 g/mol (free acid)

**Exact Mass:** 698.24 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in 10 mM Tris-HCl

**Color:** colorless to slightly yellow

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 271 nm, ε 8.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)