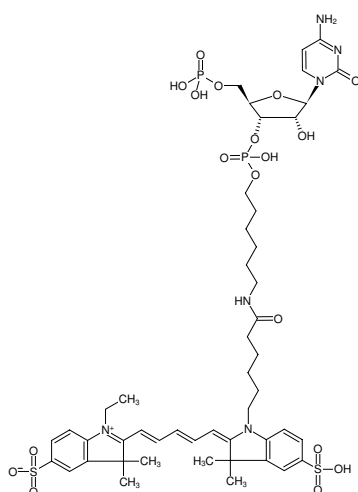




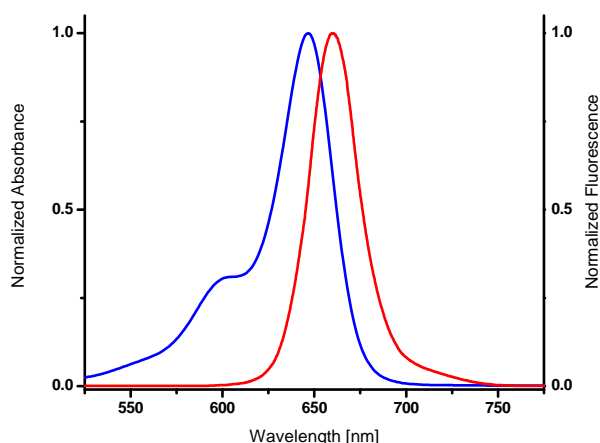
pCp-Cy5

Cytidine-5'-phosphate-3'-(6-aminohexyl)phosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-1706-CY5	20 µl (1 mM)



Structural formula of pCp-Cy5



Excitation and Emission spectrum of Cy5

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: C₄₈H₆₆N₆O₁₈P₂S₂ (free acid)

Molecular Weight: 1141.14 g/mol (free acid)

Exact Mass: 1140.34 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 649 nm, λ_{em} 670 nm, ε 250.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Incorporation into RNA/miRNA by

- 3'-End Labeling with T4 RNA Ligase^[1]

Description:

pCp-Cy5 is recommended for 3'-End Labeling of RNA/miRNA with T4 RNA Ligase that conjugates a single dye molecule to the 3'-OH group of the RNA/miRNA template. The resulting Dye-labeled RNA/miRNA probes are ideally suited for hybridization experiments such as microarray-based expression profiling or electro mobility shift assays (EMSA).

Please note: Protect the Dye-labeled pCp from exposure to light and carry out experimental procedures under low light conditions.

Related Products:

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

Selected References:

[1] Wang *et al.* (2007) Direct and sensitive miRNA profiling from low-input total RNA. *RNA* **13** (1):151.