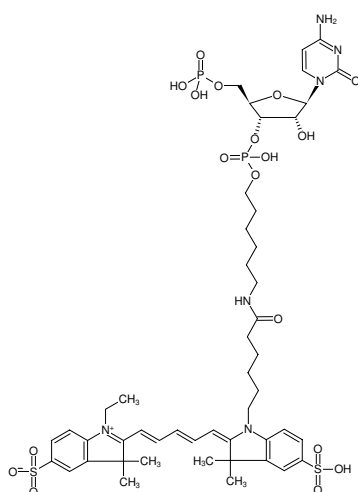




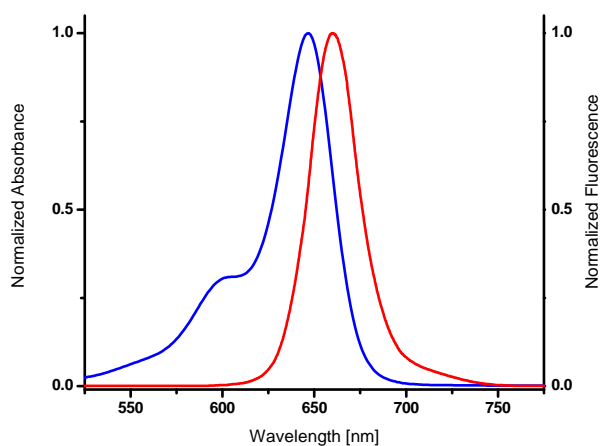
## pCp-Cy5

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 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:** C<sub>48</sub>H<sub>66</sub>N<sub>6</sub>O<sub>18</sub>P<sub>2</sub>S<sub>2</sub> (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<sub>exc</sub> 649 nm, λ<sub>em</sub> 670 nm, ε 250.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

### Applications:

- Incorporation into RNA/miRNA by
- 3'-End Labeling with T4 RNA Ligase<sup>[1]</sup>

### 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.