



## $\gamma$ -(6-Aminoethyl)-AppCp-Cy5

gamma-(6-Aminoethyl)-AMPPCP-Cy5

$\gamma$ -(6-Aminoethyl)-adenosine-5'-[(beta,gamma)-methylene]triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-279-CY5	60 $\mu$ l (1 mM)

**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>50</sub>H<sub>68</sub>N<sub>7</sub>O<sub>20</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1244.16 g/mol (free acid)

**Exact Mass:** 1243.32 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

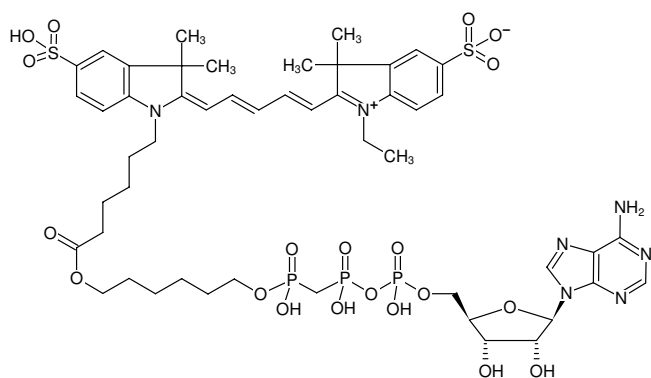
**Form:** solution in water

**Color:** blue

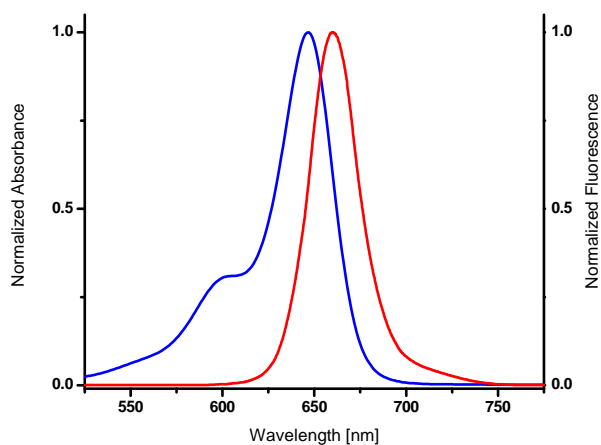
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  649 nm,  $\lambda_{em}$  670 nm,  $\epsilon$  250.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of  $\gamma$ -(6-Aminoethyl)-AppCp-Cy5



excitation and emission spectrum of Cy5