



## $\gamma$ -(6-Aminoethyl)-AppCp-Cy3

gamma-(6-Aminoethyl)-AMPPCP-Cy3

$\gamma$ -(6-Aminoethyl)-adenosine-5'-[(beta,gamma)-methylene]triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-279-CY3	60 $\mu$ l (1 mM)

**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:** C<sub>48</sub>H<sub>67</sub>N<sub>8</sub>O<sub>19</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1217.14 g/mol (free acid)

**Exact Mass:** 1216.32 g/mol (free acid)

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

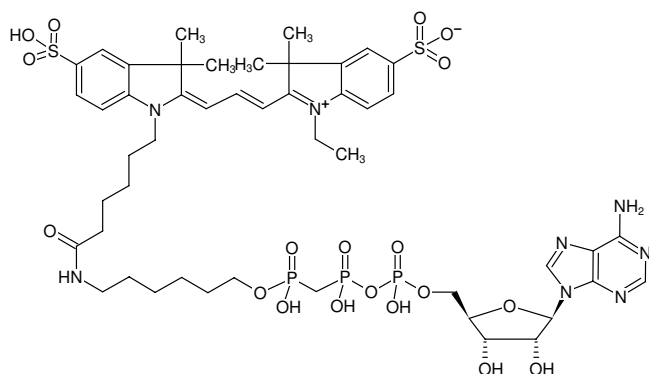
**Form:** solution in water

**Color:** pink

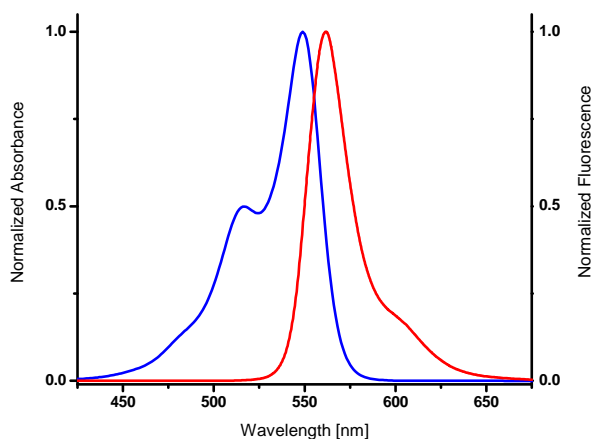
**Concentration:** 1.0 mM - 1.1 mM

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

**Spectroscopic Properties:**  $\lambda_{exc}$  550 nm,  $\lambda_{em}$  570 nm,  $\epsilon$  150.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of  $\gamma$ -(6-Aminoethyl)-AppCp-Cy3



excitation and emission spectrum of Cy3