



## $\gamma$ -[(6-Aminohexyl)-imido]-dGTP-Cy5

$\gamma$ -[(6-Aminohexyl)-imido]-2'-deoxyguanosine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-849-CY5	120 $\mu$ 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>49</sub>H<sub>68</sub>N<sub>9</sub>O<sub>19</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

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

**Exact Mass:** 1243.33 g/mol (free acid)

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

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

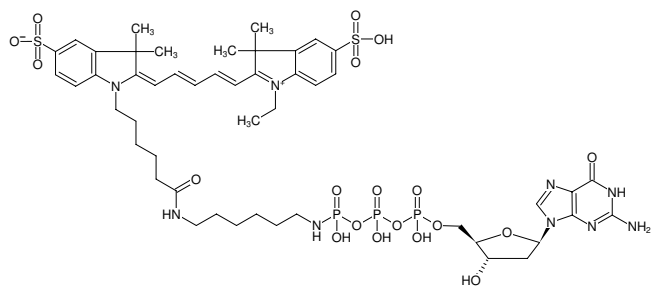
**Color:** blue

**Concentration:** 1.0 mM - 1.1 mM

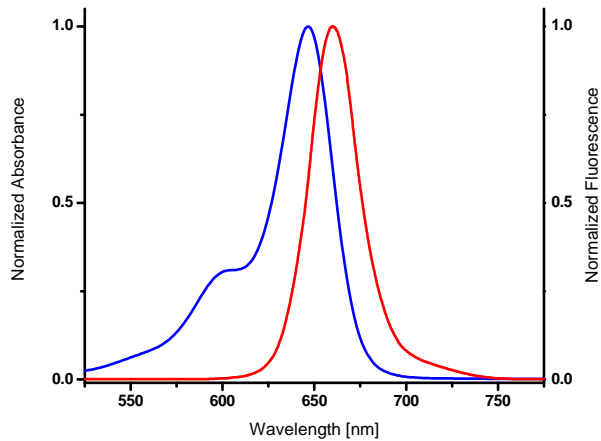
**pH:** 8.0  $\pm$  0.5

**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)

**Please note:** This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.0.



Structural formula of  $\gamma$ -[(6-Aminohexyl)-imido]-dGTP-Cy5



excitation and emission spectrum of Cy5