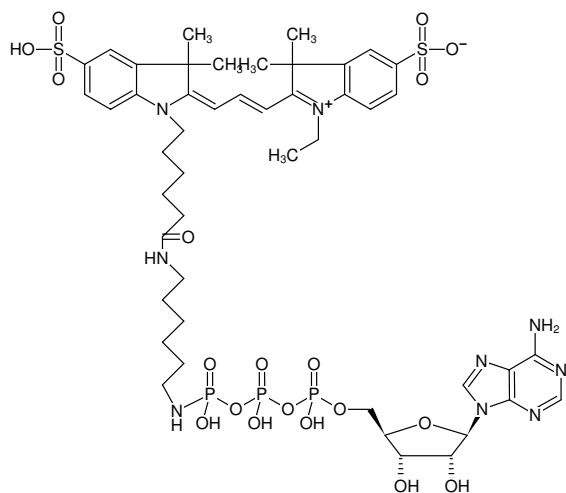




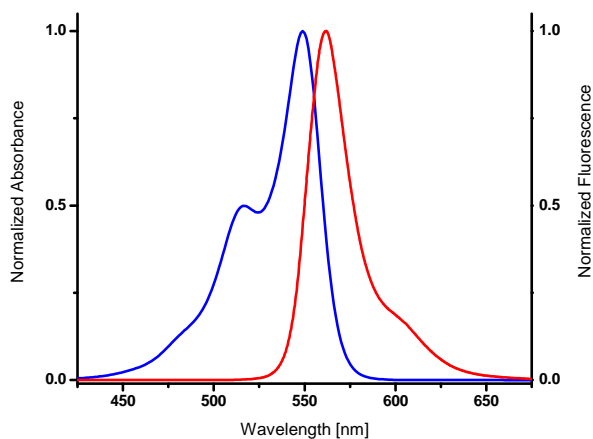
### $\gamma$ -[[6-Aminohexyl)-imido]-ATP-Cy3

$\gamma$ -[[6-Aminohexyl)-imido]-adenosine-5'-triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-814-CY3	120 $\mu$ l (1 mM)



Structural formula of  $\gamma$ -[[6-Aminohexyl)-imido]-ATP-Cy3



excitation and emission spectrum of Cy3

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at  $-20\text{ }^{\circ}\text{C}$

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:**  $\text{C}_{47}\text{H}_{66}\text{N}_9\text{O}_{19}\text{P}_3\text{S}_2$  (free acid)

**Molecular Weight:** 1218.13 g/mol (free acid)

**Exact Mass:** 1217.31 g/mol (free acid)

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

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

**Color:** pink

**Concentration:** 1.0 mM - 1.1 mM

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

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

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