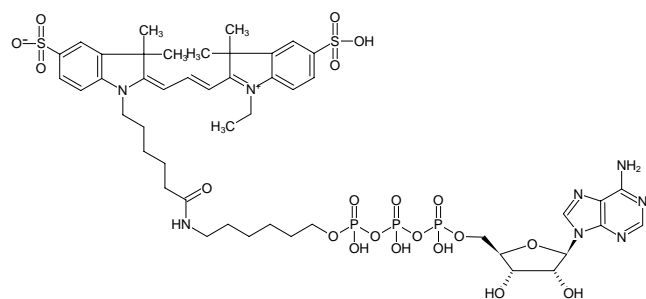




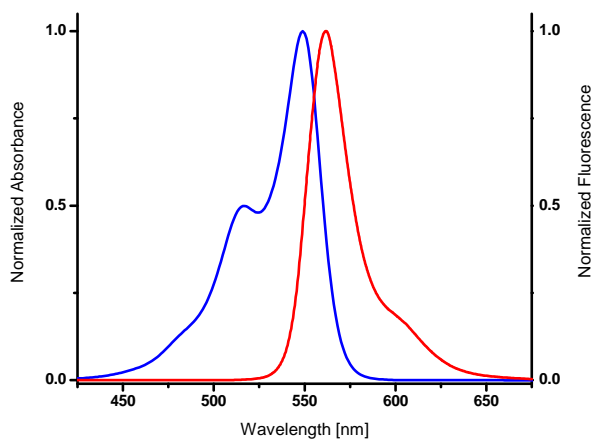
**$\gamma$ -(6-Aminoethyl)-ATP - Cy3**

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

Cat. No.	Amount
NU-833-CY3	120 $\mu$ l (1 mM)



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



excitation and emission spectrum of Cy3

**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>47</sub>H<sub>65</sub>N<sub>8</sub>O<sub>20</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1219.11 g/mol (free acid)

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

**Form:** pink solution in water, pH 7.5  $\pm$  0.5

**Concentration:** 1 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)

**Selected References:**

Schax *et al.* (2014) Microarray-based screening of heat shock protein inhibitors. *J. Biotechnol.* **180**:1.