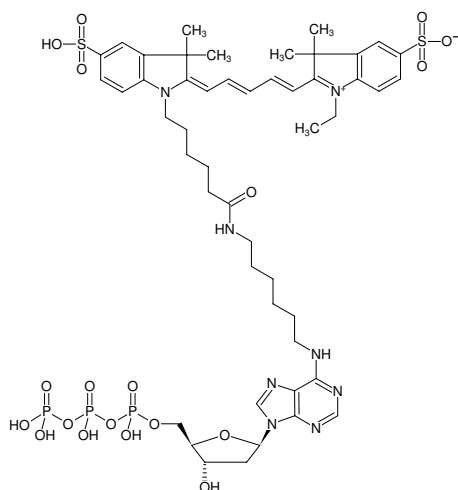




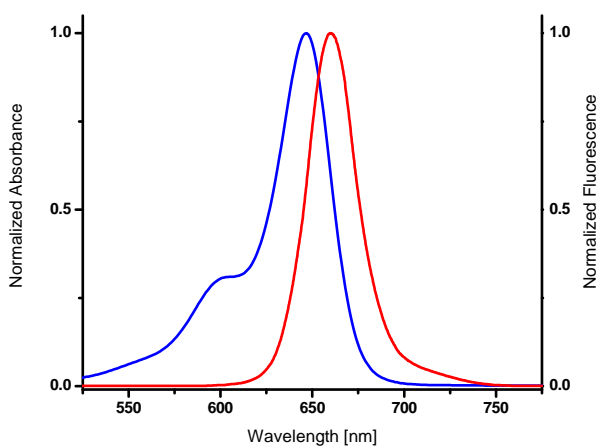
**N<sup>6</sup>-(6-Aminoethyl)-dATP-Cy5**

N<sup>6</sup>-(6-Aminoethyl)-2'-deoxyadenosine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-835-CY5	160 µl (1 mM)



Structural formula of N<sup>6</sup>-(6-Aminoethyl)-dATP-Cy5



excitation and emission spectrum of Cy5

**For research use only!**

**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>49</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:** 1229.15 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** blue

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 649 nm, λ<sub>em</sub> 670 nm, ε 250.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Selected References:**

Ando *et al.* (2019) ELTA: Enzymatic Labeling of Terminal ADP-Ribose. *Mol. Cell.* **73 (4):**845.

Evangelista *et al.* (1996) Characterization of fluorescent nucleoside triphosphates by capillary electrophoresis with laser-induced fluorescence detection: action of alkaline phosphatase and DNA polymerase. *Anal. Biochem.* **235 (1):**89.