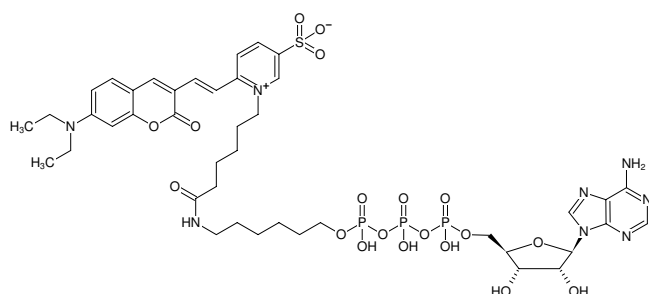




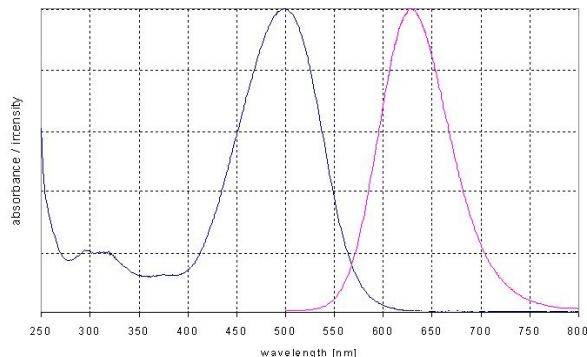
### $\gamma$ -(6-Aminoethyl)-ATP-DY-480XL

$\gamma$ -(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with DY 480XL, Triethylammonium salt

Cat. No.	Amount
NU-833-480XL	80 $\mu$ l (1 mM)



Structural formula of  $\gamma$ -(6-Aminoethyl)-ATP-DY-480XL



excitation and emission spectrum of DY 480XL

**For general laboratory use.**

**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>42</sub>H<sub>57</sub>N<sub>8</sub>O<sub>19</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1102.93 g/mol (free acid)

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

**Form:** solution in water

**Concentration:** 1.0 mM - 1.1 mM

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

**Spectroscopic Properties:**  $\lambda_{exc}$  500 nm,  $\lambda_{em}$  630 nm,  $\epsilon$  50.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)