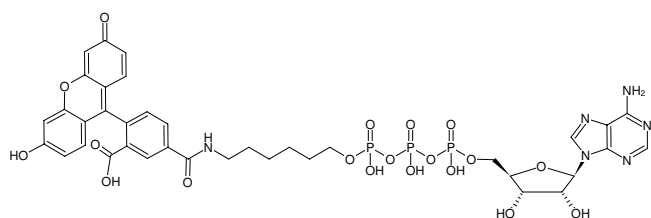




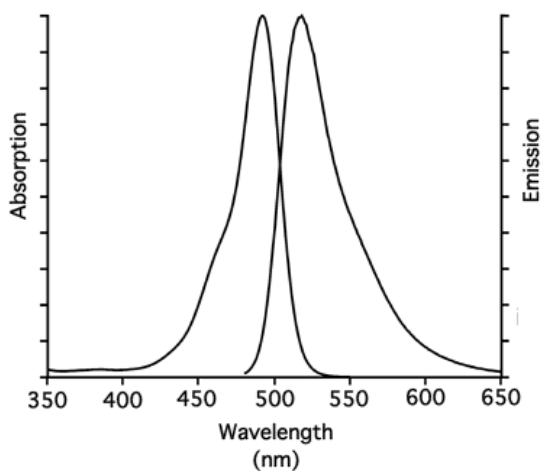
### $\gamma$ -(6-Aminoethyl)-ATP-5-FAM

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

Cat. No.	Amount
NU-833-5FM	160 $\mu$ l (1 mM)



Structural formula of  $\gamma$ -(6-Aminoethyl)-ATP-5-FAM



excitation and emission spectrum of 5-FAM

**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>37</sub>H<sub>39</sub>N<sub>6</sub>O<sub>19</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 964.66 g/mol (free acid)

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

**Form:** solution in water

**Color:** yellow

**Concentration:** 1.0 mM - 1.1 mM

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

**Spectroscopic Properties:**  $\lambda_{exc}$  492 nm,  $\lambda_{em}$  517 nm,  $\epsilon$  83.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Selected References:**

He *et al.* (2010) Aptamer based reversible DNA induced hydrogel system for molecular recognition and separation. *Chem. Commun.* **46** (34):6308.