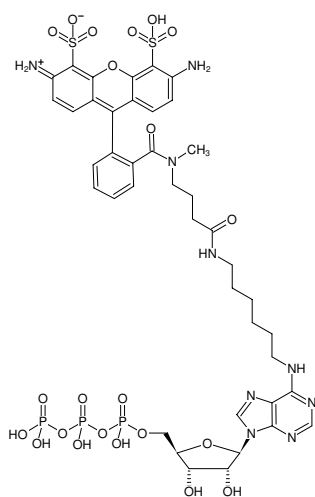




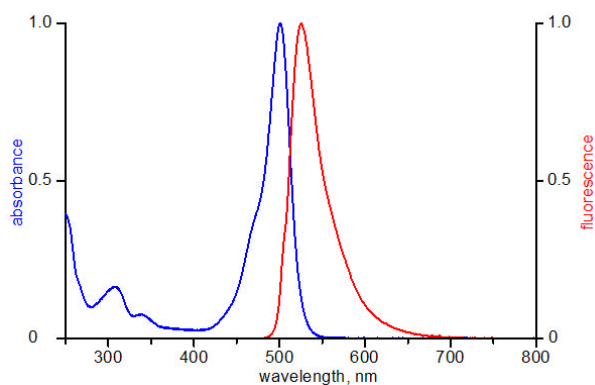
## N<sup>6</sup>-(6-Aminoheptyl)-ATP-ATTO-488

N<sup>6</sup>-(6-Aminoheptyl)-adenosine-5'-triphosphate, labeled with ATTO 488, Triethylammonium salt

Cat. No.	Amount
NU-805-488	30 µl (1 mM)



Structural formula of N<sup>6</sup>-(6-Aminoheptyl)-ATP-ATTO-488



excitation and emission spectrum of ATTO 488

**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>41</sub>H<sub>50</sub>N<sub>9</sub>O<sub>22</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1177.93 g/mol (free acid)

**Exact Mass:** 1177.17 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** yellow

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 500 nm, λ<sub>em</sub> 520 nm, ε 90.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Selected References:**

Pettinger *et al.* (2017) An Irreversible Inhibitor of HSP72 that Unexpectedly Targets Lysine-56. *Angew. Chem. Int. Ed. Engl.* **56** (13):3536.