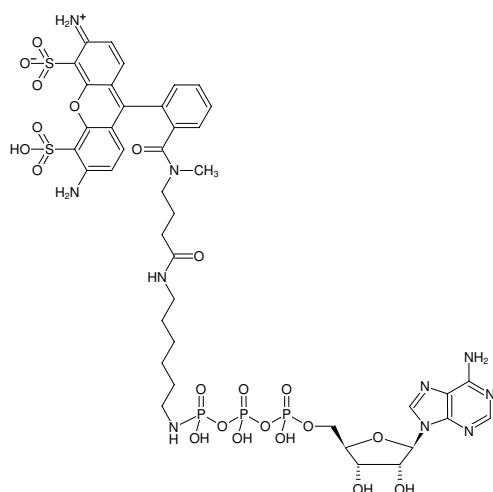




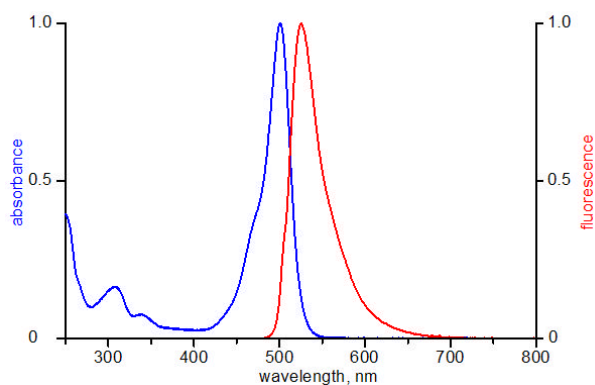
## $\gamma$ -[(6-Aminohexyl)-imido]-ATP-ATTO-488

$\gamma$ -[(6-Aminohexyl)-imido]-adenosine-5'-triphosphate, labeled with ATTO-488, Triethylammonium salt

Cat. No.	Amount
NU-814-488	120 $\mu$ l (1 mM)



Structural formula of  $\gamma$ -[(6-Aminohexyl)-imido]-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>51</sub>N<sub>10</sub>O<sub>21</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1176.19 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:**  $\lambda_{exc}$  500 nm,  $\lambda_{em}$  520 nm,  $\epsilon$  90.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Please note:** This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.0.