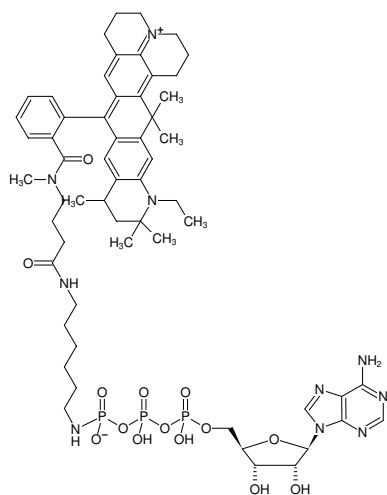




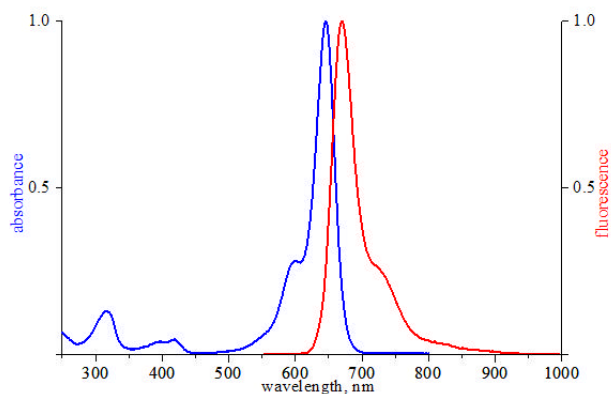
## $\gamma$ -[(6-Aminoethyl)-imido]-ATP-ATTO-647N

$\gamma$ -[(6-Aminoethyl)-imido]-adenosine-5'-triphosphate, labeled with ATTO-647N, Triethylammonium salt

Cat. No.	Amount
NU-814-647N	120 $\mu$ l (0,5 mM)



Structural formula of  $\gamma$ -[(6-Aminoethyl)-imido]-ATP-ATTO-647N



excitation and emission spectrum of ATTO 647N

**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>58</sub>H<sub>79</sub>N<sub>10</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1232.23 g/mol (free acid)

**Exact Mass:** 1232.50 g/mol (free acid)

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

**Form:** solution in 10 mM Tris-HCl

**Color:** blue

**Concentration:** 0.50 mM - 0.55 mM

**pH:** 8.0  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  646 nm,  $\lambda_{em}$  664 nm,  $\epsilon$  150.0 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.

**Selected References:**

Götz *et al.* (2016) A Multicolor Single-Molecule FRET Approach to Study Protein Dynamics and Interactions Simultaneously. *Methods Enzymol.* **581**:487.