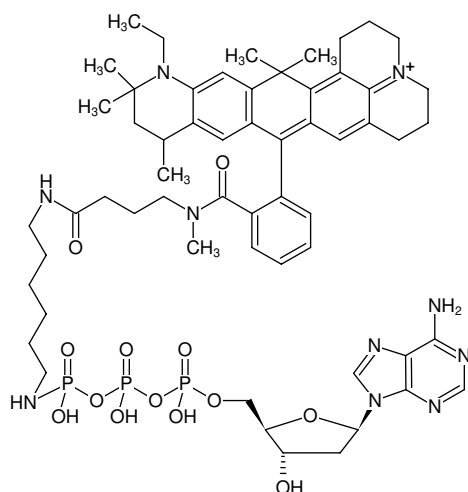




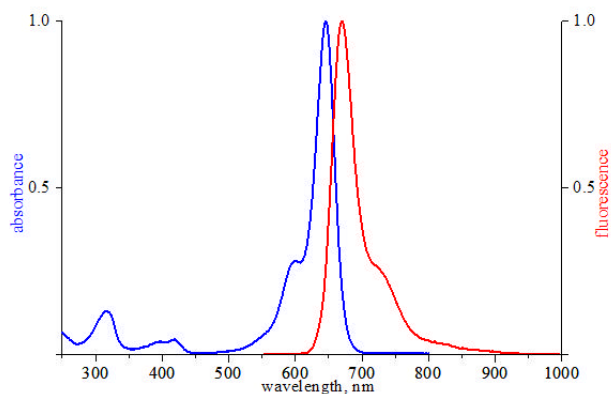
**$\gamma$ -[(6-Aminohexyl)-imido]-dATP-ATTO-647N**

$\gamma$ -[(6-Aminohexyl)-imido]-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO-647N, Triethylammonium salt

Cat. No.	Amount
NU-861-647N	120 $\mu$ l (0,5 mM)



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



excitation and emission spectrum of ATTO 647N

**For research use only!**

**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>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1217.23 g/mol (free acid)

**Exact Mass:** 1216.50 g/mol (free acid)

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

**Form:** solution in water

**Color:** blue

**Concentration:** 0.50 mM - 0.55 mM

**pH:** 7.5  $\pm$  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.