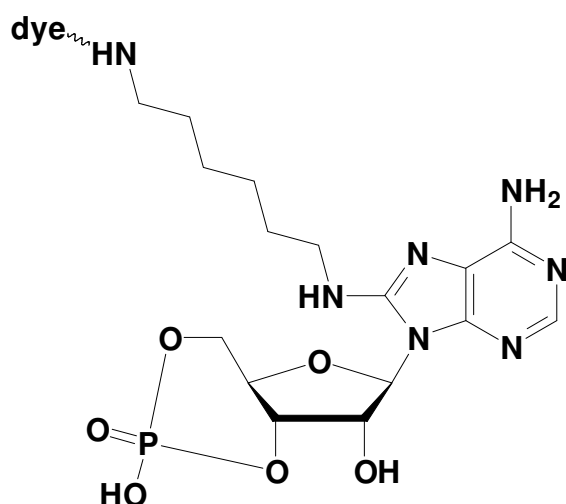


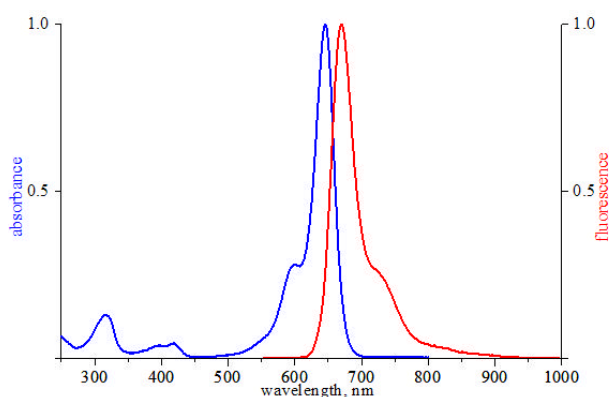
## 8-[(6-Amino)hexyl]-amino-cAMP - ATTO 647N

8-[(6-Amino)hexyl]-amino-adenosine-3',5'-cyclic monophosphate, labeled with ATTO 647N, Triethylammonium salt

Cat. No.	Amount
NU-851-647N	80 $\mu$ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-cAMP



excitation and emission spectrum of ATTO 647N

**Cat. No.:** NU-851-647N

**Molecular Formula:** C<sub>16</sub>H<sub>26</sub>N<sub>7</sub>O<sub>6</sub>P - ATTO 647N (free acid)

**Molecular Weight:** 1070.39 (free acid)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**

$\lambda_{exc}$  644 nm;  $\lambda_{em}$  669 nm;  $\epsilon$  150,000 cm<sup>-1</sup> M<sup>-1</sup>

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**