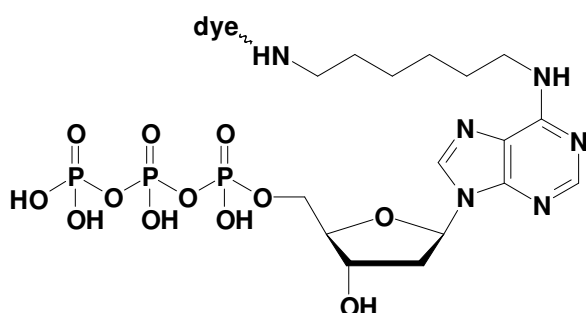


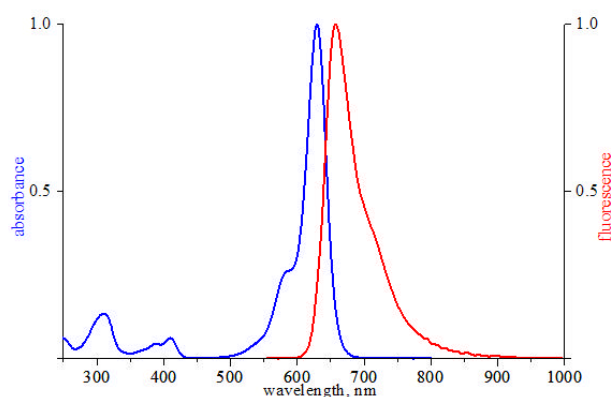
## N<sup>6</sup>-(6-Amino)hexyl-dATP - ATTO 633

N<sup>6</sup>-(6-Amino)hexyl-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO 633,  
 Triethylammonium salt

| Cat. No.   | Amount            |
|------------|-------------------|
| NU-835-633 | 40 $\mu$ l (1 mM) |



structural formula of N<sup>6</sup>-(6-Amino)hexyl-dATP



excitation and emission spectrum of ATTO 633

**Cat. No.:** NU-835-633

**Molecular Formula:** C<sub>16</sub>H<sub>29</sub>N<sub>6</sub>O<sub>12</sub>P<sub>3</sub> - ATTO 633 (free acid)

**Molecular Weight:** 1123.36 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  629 nm;  $\lambda_{em}$  657 nm;  $\epsilon$  130,000 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!**