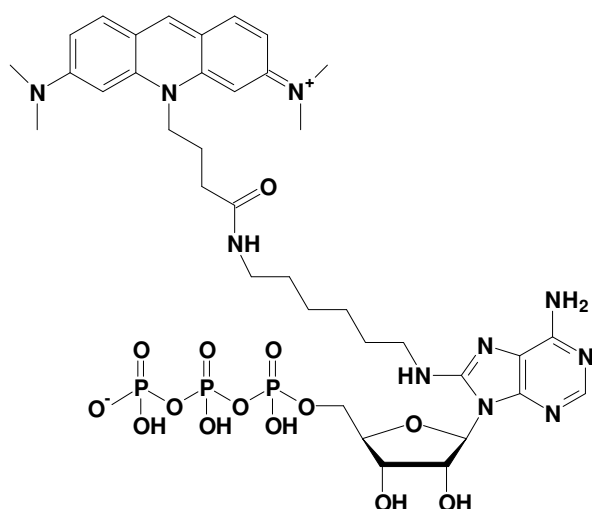


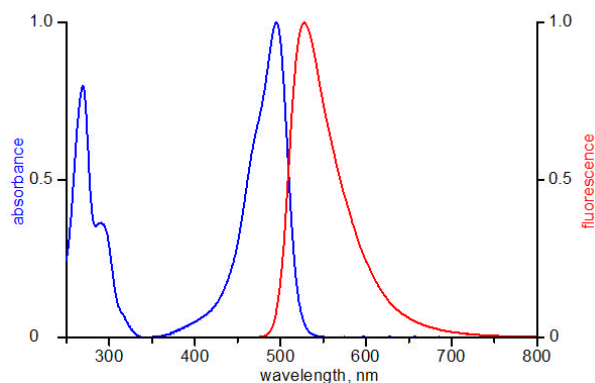
## 8-[(6-Amino)hexyl]-amino-ATP - ATTO 495

8-[(6-Amino)hexyl]-amino-adenosine-5'-triphosphate, labeled with ATTO 495,  
 Triethylammonium salt

Cat. No.	Amount
NU-807-495	120 $\mu$ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-ATP - ATTO 495



excitation and emission spectrum of ATTO 495

**Cat. No.:** NU-807-495

**Molecular Formula:** C<sub>37</sub>H<sub>53</sub>N<sub>10</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 954.80 (free acid)

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

**Spectroscopic properties:**

$\lambda_{\text{exc}}$  495 nm;  $\lambda_{\text{em}}$  527 nm;  $\epsilon$  80,000 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!**