



## 8-[(6-Amino)hexyl]-amino-adenosine-3',5'-bisphosphate - ATTO-495

8-[(6-Amino)hexyl]-amino-adenosine-3',5'-bisphosphate, labeled with ATTO 495, Triethylammonium salt

Cat. No.	Amount
NU-811-495	80 µl (1 mM)

**For research use only!**

**Shipping:** shipped on blue ice

**Storage Conditions:** store at -20 °C

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

**Shelf Life:** 12 months

**Molecular Formula:** C<sub>37</sub>H<sub>52</sub>N<sub>10</sub>O<sub>11</sub>P<sub>2</sub>(free acid)

**Molecular Weight:** 874.82 g/mol (free acid)

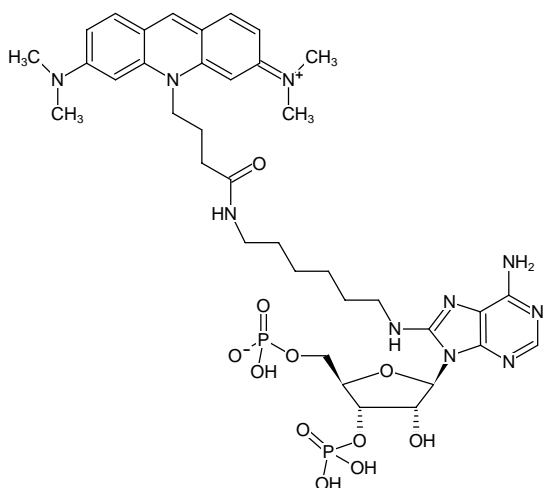
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution, pH 7.5 ± 0.5

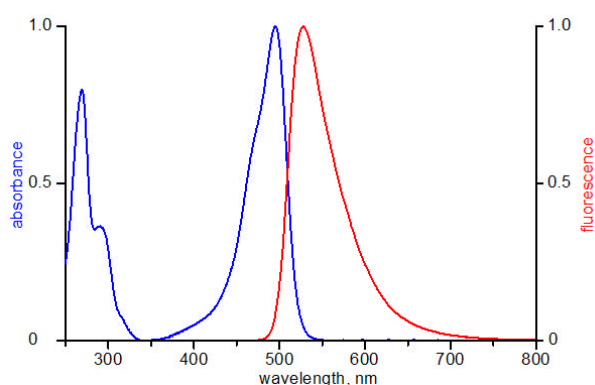
**Concentration:** 1 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 495 nm; λ<sub>em</sub> 527 nm; ε 80.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 8-[(6-Amino)hexyl]-amino-adenosine-3',5'-bisphosphate - ATTO-495



excitation and emission spectrum of ATTO 495