



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: C₃₇H₅₂N₁₀O₁₁P₂(free acid)

Molecular Weight: 874.82 g/mol (free acid)

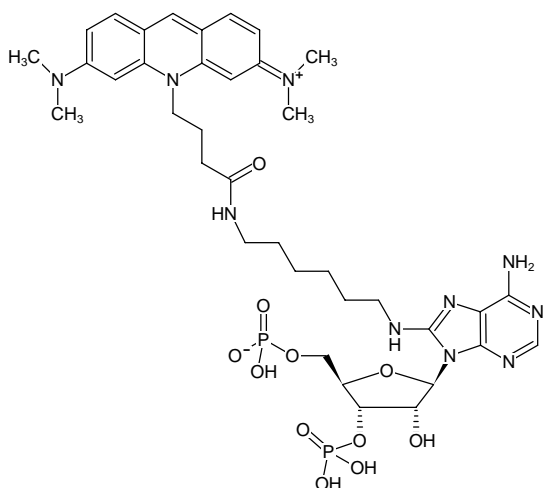
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution, pH 7.5 ± 0.5

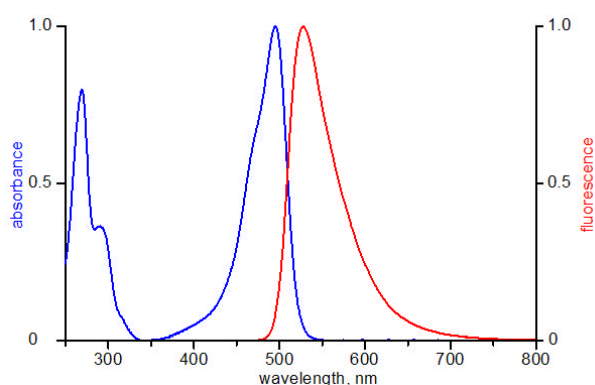
Concentration: 1 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 495 nm; λ_{em} 527 nm; ε 80.0 L mmol⁻¹ cm⁻¹ (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