



N⁶-(6-Aminohexyl)-ATP-ATTO-495

N⁶-(6-Aminohexyl)-adenosine-5'-triphosphate, labeled with ATTO 495, Triethylammonium salt

Cat. No.	Amount
NU-805-495	120 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.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₃₇H₅₂N₉O₁₄P₃ (free acid)

Molecular Weight: 939.78 g/mol (free acid)

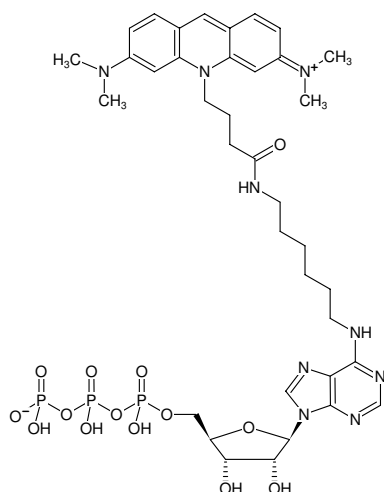
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

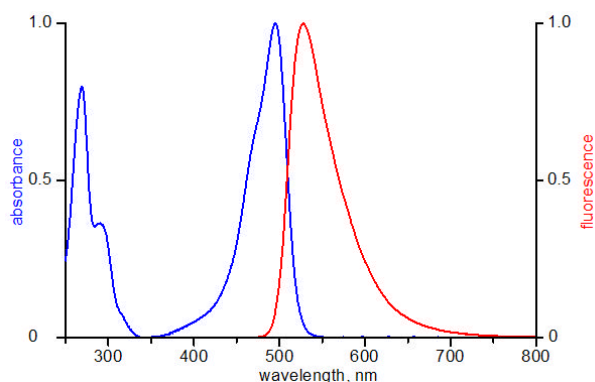
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 498 nm, λ_{em} 526 nm, ε 80.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of N⁶-(6-Aminohexyl)-ATP-ATTO-495



excitation and emission spectrum of ATTO 495