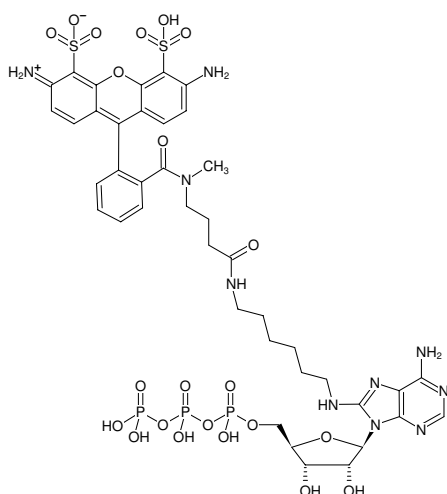




8-(6-Aminoethyl)-amino-ATP-ATTO-488

8-(6-Aminoethyl)-amino-adenosine-5'-triphosphate, labeled with ATTO 488, Triethylammonium salt

Cat. No.	Amount
NU-807-488	120 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-ATP-ATTO-488

For general laboratory use.

Shipping: shipped on gel packs

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: C₄₁H₅₁N₁₀O₂₂P₃S₃ (free acid)

Molecular Weight: 1192.95 g/mol (free acid)

Exact Mass: 1192.18 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

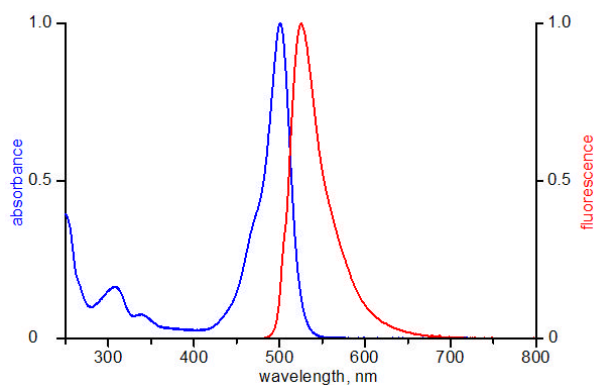
Form: solution in water

Color: yellow

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 500 nm, λ_{em} 520 nm, ε 90.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 488