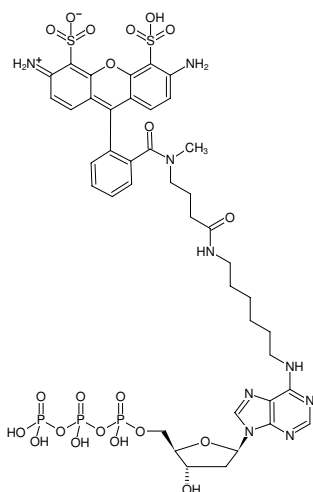




N⁶-(6-Aminoethyl)-dATP-ATTO-488

N⁶-(6-Aminoethyl)-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO 488, Triethylammonium salt

Cat. No.	Amount
NU-835-488	80 µl (1 mM)



Structural formula of N⁶-(6-Aminoethyl)-dATP-ATTO-488

For research use only!

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: C₄₁H₅₀N₉O₂₁P₃S₂ (free acid)

Molecular Weight: 1161.18 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

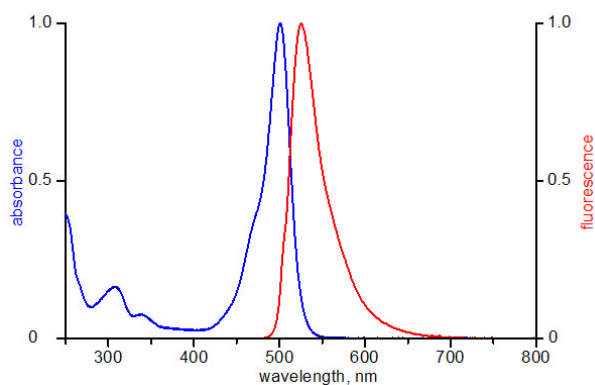
Form: solution in water

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