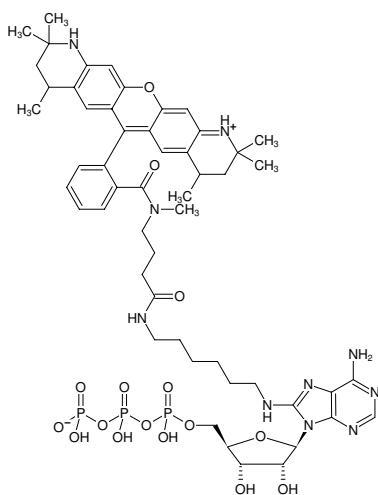




8-(6-Aminoethyl)-amino-ATP-ATTO-550

8-(6-Aminoethyl)-amino-adenosine-5'-triphosphate, labeled with ATTO 550, Triethylammonium salt

Cat. No.	Amount
NU-807-550	120 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-ATP-ATTO-550

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: C₅₃H₇₁N₁₀O₁₆P₃ (free acid)

Molecular Weight: 1197.13 g/mol (free acid)

Exact Mass: 1196.43 g/mol (free acid)

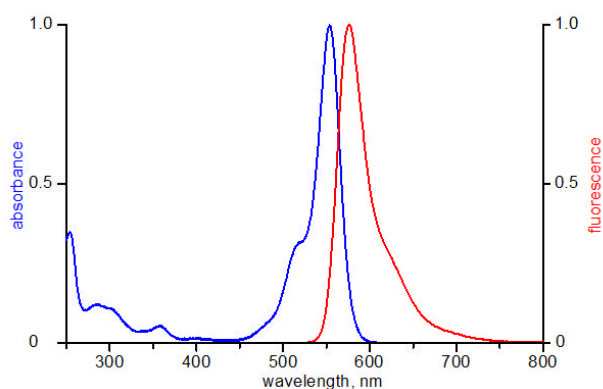
Purity: ≥ 95 % (HPLC)

Form: red-violet solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 554 nm, λ_{em} 576 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 550