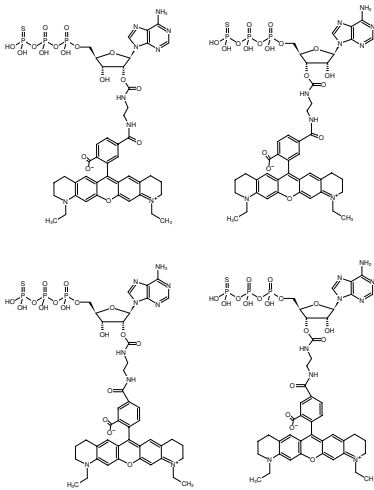




EDA-ATPyS-ATTO-565

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with ATTO 565, Triethylammonium salt

Cat. No.	Amount
NU-1609-565	50 µl (1 mM)



Structural formula of EDA-ATPyS-ATTO-565

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₄₄H₅₀N₉O₁₇P₃S (free acid)

Molecular Weight: 1101.91 g/mol (free acid)

Exact Mass: 1101.23 g/mol (free acid)

Purity: ≥ 90 % (HPLC)

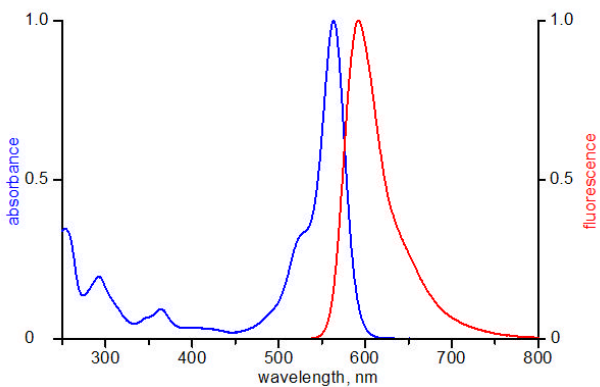
Form: solution in water

Color: red

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 564 nm, λ_{em} 590 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 565