



EDA-ATPyS-ATTO-465

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with ATTO 465, Triethylammonium salt

Cat. No.	Amount
NU-1609-465	50 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤ 2 ml).

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₃₀H₃₇N₁₀O₁₄P₃S (free acid)

Molecular Weight: 886.66 g/mol (free acid)

Exact Mass: 886.14 g/mol (free acid)

Purity: ≥ 90 % (HPLC)

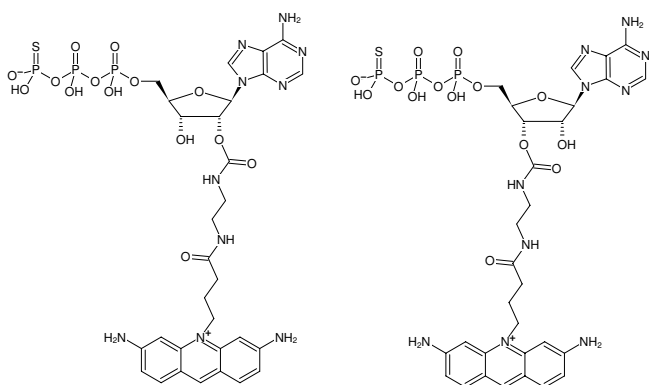
Form: solution in water

Color: yellow

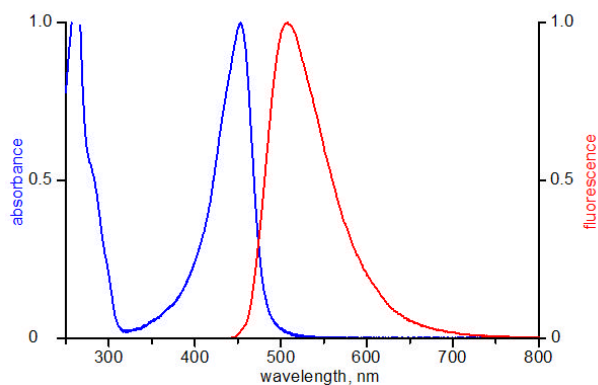
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 453 nm, λ_{em} 506 nm, ε 75.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-ATPyS-ATTO-465



excitation and emission spectrum of ATTO 465