



EDA-ATPyS-ATTO-Rho101

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with ATTO Rho101, Triethylammonium salt

Cat. No.	Amount
NU-1609-RHO101	50 µl (1 mM)

For research use only!

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₅₀H₅₉N₁₀O₁₆P₃S (free acid)

Molecular Weight: 1180.30 g/mol (free acid)

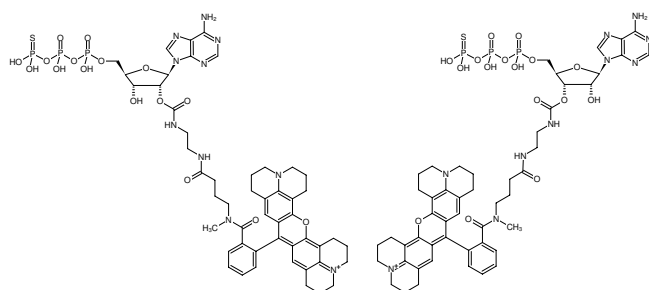
Purity: ≥ 90 % (HPLC), contains approx. 5 % EDA-ADP-dye

Form: solution in water

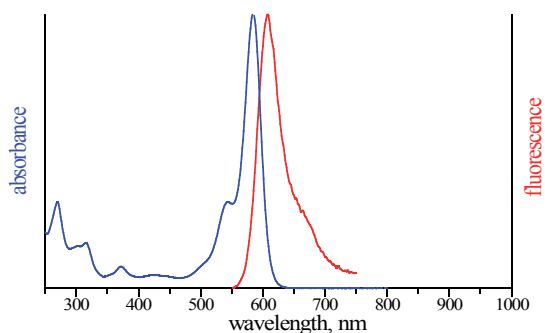
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 587 nm, λ_{em} 609 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-ATPyS-ATTO-Rho101



excitation and emission spectrum of ATTO Rho101