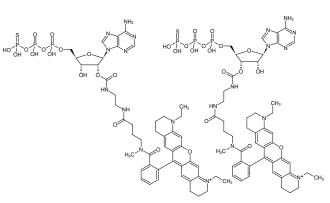




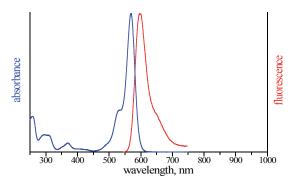
EDA-ATPyS-ATTO-Rho11

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-1609-RHO11	50 µl (1 mM)



Structural formula of EDA-ATPyS-ATTO-Rho11



excitation and emission spectrum of ATTO Rho11

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: $C_{48}H_{59}N_{10}O_{16}P_3S$ (free acid) Molecular Weight: 1156.30 g/mol (free acid) Exact Mass: 1156.30 g/mol (free acid) Purity: ≥ 90 % (HPLC) Form: solution in water Concentration: 1.0 mM - 1.1 mM pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 572 nm, λ_{em} 595 nm, ϵ 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

