



EDA-ATPyS-5/6-TAMRA

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with 5/6-TAMRA, Triethylammonium salt

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

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: C₃₈H₄₂N₉O₁₇P₃S (free acid)

Molecular Weight: 1021.78 g/mol (free acid)

Purity: ≥ 90 % (HPLC)

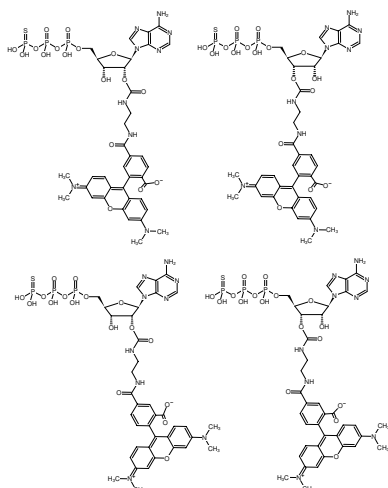
Form: solution in water

Color: pink to red

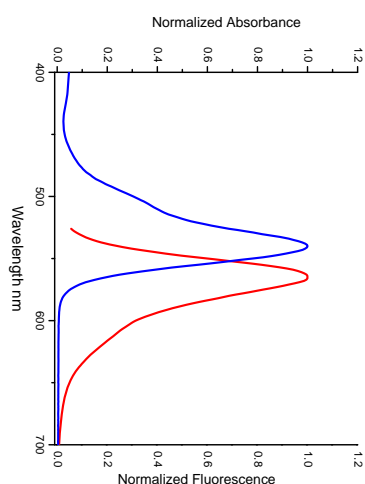
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 545 nm, λ_{em} 575 nm, ε 90.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-ATPyS-5/6-TAMRA



excitation and emission spectrum of 5/6-TAMRA