



EDA-GTPyS-ATTO-594

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with ATTO 594, Triethylammonium salt

Cat. No.	Amount
NU-1610-594	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: C₅₄H₆₇N₁₀O₂₃P₃S₃ (free acid)

Molecular Weight: 1413.28 g/mol (free acid)

Exact Mass: 1412.28 g/mol (free acid)

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

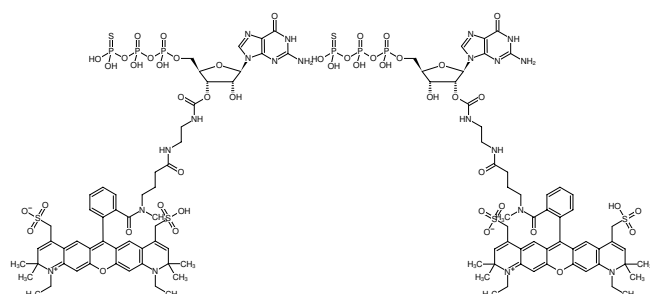
Form: solution in water

Color: red-violet

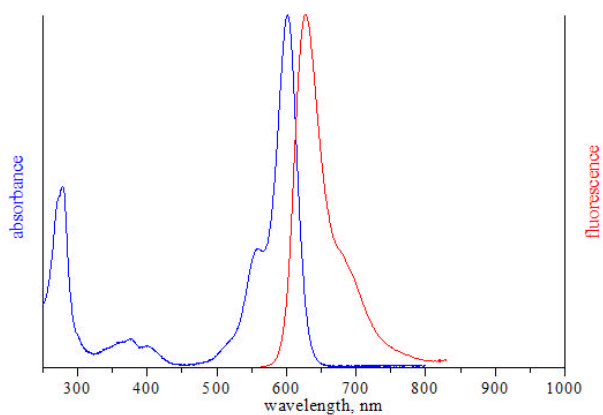
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 603 nm, λ_{em} 626 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-GTPyS-ATTO-594



excitation and emission spectrum of ATTO 594