



EDA-GTPyS-ATTO-Rho6G

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with ATTO Rho6G, Triethylammonium salt

Cat. No.	Amount
NU-1610-RHO6G	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: C₄₄H₅₅N₁₀O₁₇P₃S (free acid)

Molecular Weight: 1120.27 g/mol (free acid)

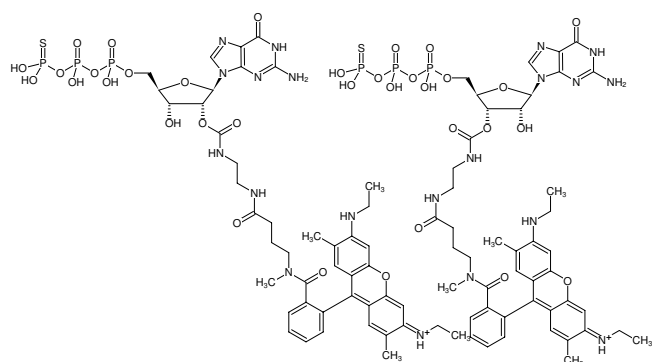
Purity: ≥ 90 % (HPLC), contains approx. 5 % EDA-GDP-dye

Form: solution in water

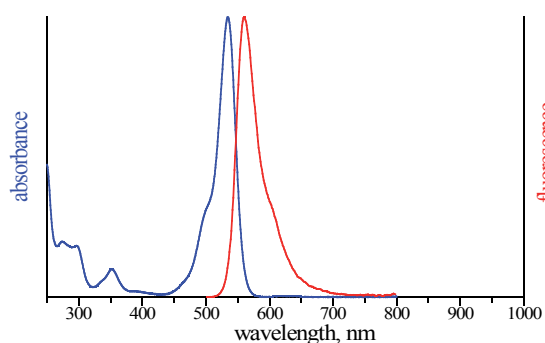
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 533 nm, λ_{em} 557 nm, ε 115.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-GTPyS-ATTO-Rho6G



excitation and emission spectrum of ATTO Rho6G

Please note: For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0.

The limited stability of labeled ATPyS and GTPyS nucleotides makes it impossible to keep stocks of them and to offer them as catalog products that are available immediately. Therefore, any labeled ATPyS or GTPyS is freshly prepared per order. For this reason the price is higher than for stable nucleotide-dye conjugates which can be synthesized and kept in stock.