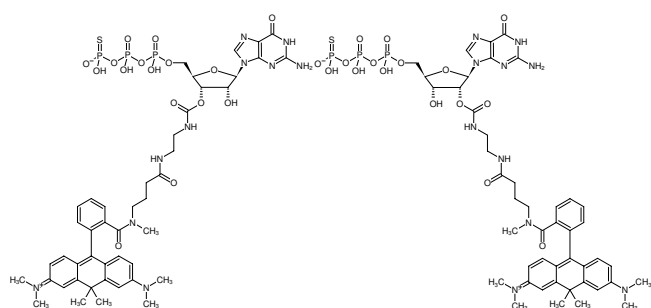




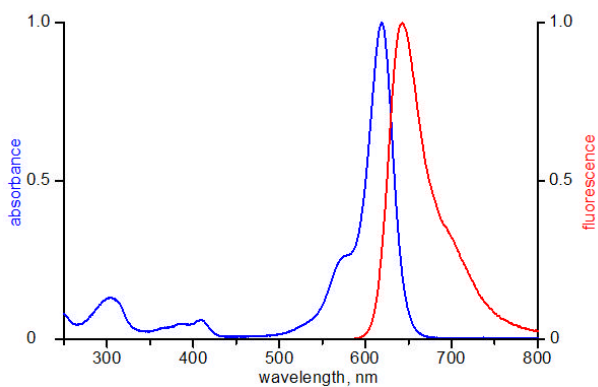
EDA-GTPyS-ATTO-620

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with ATTO 620, Triethylammonium salt

Cat. No.	Amount
NU-1610-620	50 µl (1 mM)



Structural formula of EDA-GTPyS-ATTO-620



excitation and emission spectrum of ATTO 620

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: C₄₅H₅₇N₁₀O₁₆P₃S (free acid)

Molecular Weight: 1118.99 g/mol (free acid)

Exact Mass: 1118.29 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} 620 nm, λ_{em} 642 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

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.