



EDA-GTPyS-DY-751

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with DY 751, Triethylammonium salt

Cat. No.	Amount
NU-1610-751	50 µl (1 mM)

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 6 months after date of delivery

Molecular Formula: C₅₅H₇₂N₉O₂₂P₃S₃ (free acid)

Molecular Weight: 1399.32 g/mol (free acid)

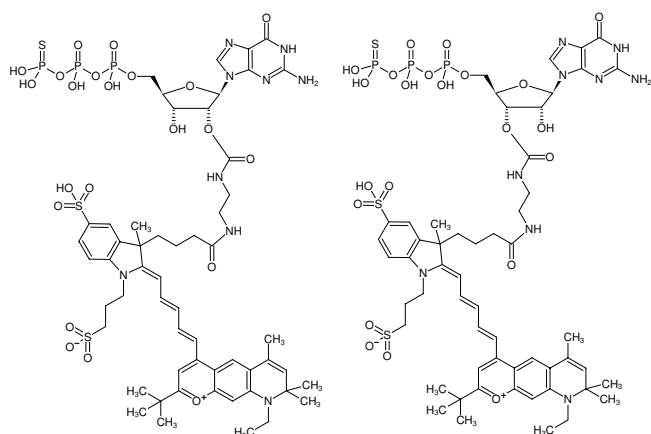
Purity: ≥ 90 % (HPLC), contains approx. 5 % EDA-GDP-dye

Form: solution in water

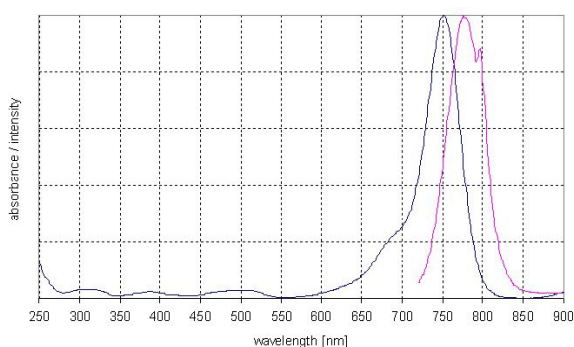
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 751 nm, λ_{em} 779 nm, ε 270.0 L mol⁻¹ cm⁻¹ (Ethanol)



Structural formula of EDA-GTPyS-DY-751



excitation and emission spectrum of DY 751

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.