



EDA-GTPyS-DY-776

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with DY 776, Triethylammonium salt

Cat. No.	Amount
NU-1610-776	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: C₅₇H₆₈N₉O₂₂P₃S₃ (free acid)

Molecular Weight: 1420.31 g/mol (free acid)

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

Form: solution in water

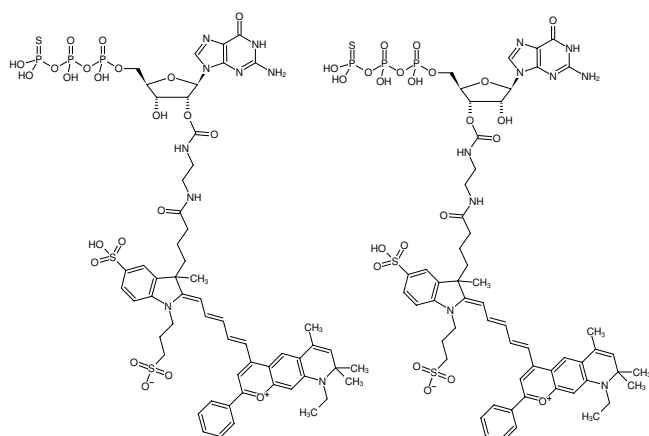
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

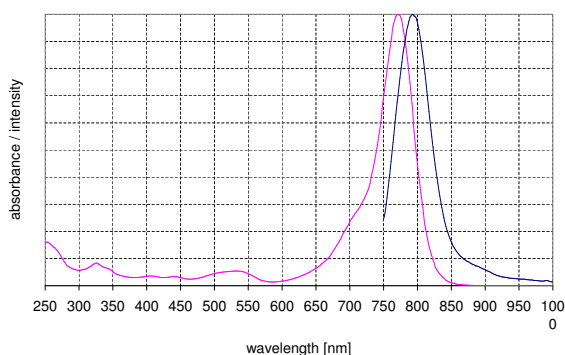
Spectroscopic Properties: λ_{exc} 771 nm, λ_{em} 793 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

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.



Structural formula of EDA-GTPyS-DY-776



excitation and emission spectrum of DY 776