



EDA-GTPyS-ATTO-740

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with ATTO 740, Triethylammonium salt

| Cat. No. | Amount |
|-------------|--------------|
| NU-1610-740 | 50 µl (1 mM) |

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Additional Storage Conditions: long time storage at pH 7.5

Shelf Life: 6 months after date of delivery

Molecular Formula: C₁₃H₂₂N₇O₁₄P₃S - ATTO 740 (free acid)

Molecular Weight: 1074.33 g/mol (free acid)

Purity: ≥ 90 % (HPLC)

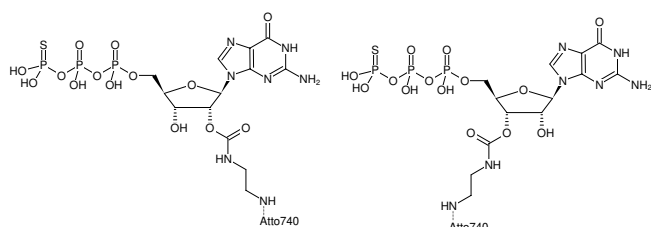
Form: solution in water

Color: green

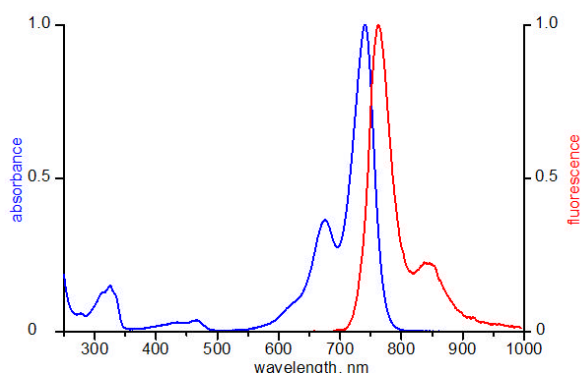
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 743 nm, λ_{em} 763 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-GTPyS-ATTO-740



excitation and emission spectrum of ATTO 740