



EDA-m⁷GTP-DY-751

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate, labeled with DY 751, Triethylammonium salt

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate labeled with DY 751, Triethylammonium salt

Cat. No.	Amount
NU-824-751	40 µl (1 mM)

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

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

Shelf Life: 12 months after date of delivery

Molecular Formula: C₅₆H₇₄N₉O₂₃P₃S₂ (free acid)

Molecular Weight: 1397.36 g/mol (free acid)

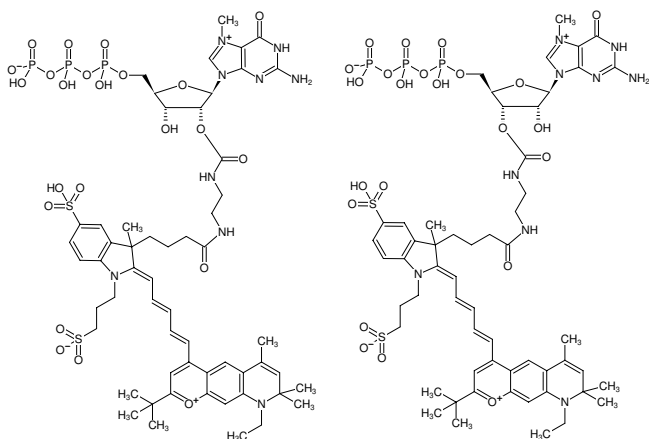
Purity: ≥ 95 % (HPLC)

Form: solution in water

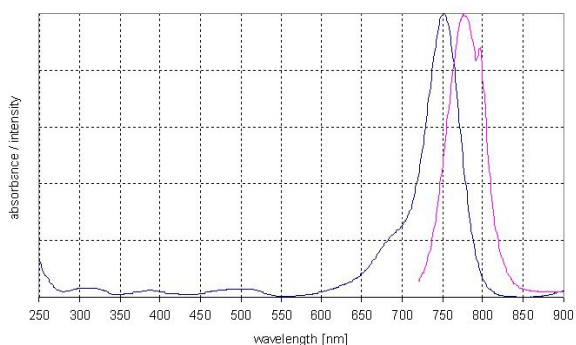
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

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



Structural formula of EDA-m⁷GTP-DY-751



excitation and emission spectrum of DY 751