



EDA-GTP-DY-776

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

Cat. No.	Amount
NU-820-776	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: C₅₇H₆₈N₉O₂₃P₃S₂ (free acid)

Molecular Weight: 1404.25 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

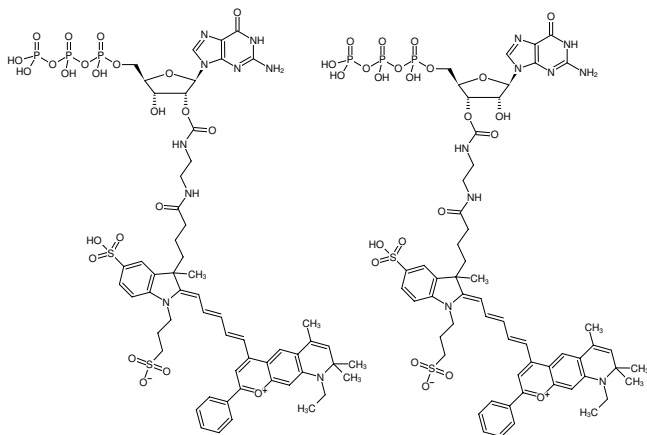
Form: solution in water

Color: dark blue

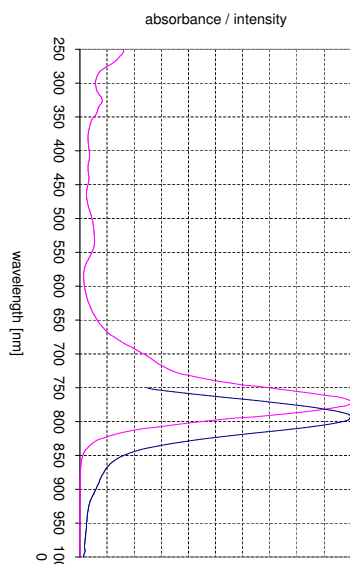
Concentration: 1.0 mM - 1.1 mM

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)



Structural formula of EDA-GTP-DY-776



excitation and emission spectrum of DY 776