



EDA-m⁷GTP-ATTO-594

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate, labeled with ATTO 594, Triethylammonium salt

Cat. No.	Amount
NU-824-594	80 µl (0,5 mM)

For research use only!

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: C₅₅H₆₉N₁₀O₂₄P₃S₂ (free acid)

Molecular Weight: 1411.24 g/mol (free acid)

Exact Mass: 1410.31 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

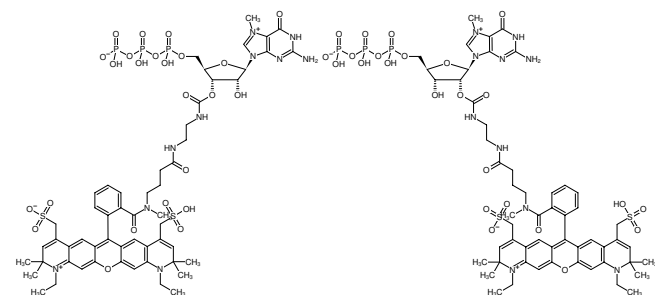
Form: solution in water

Color: red-violet

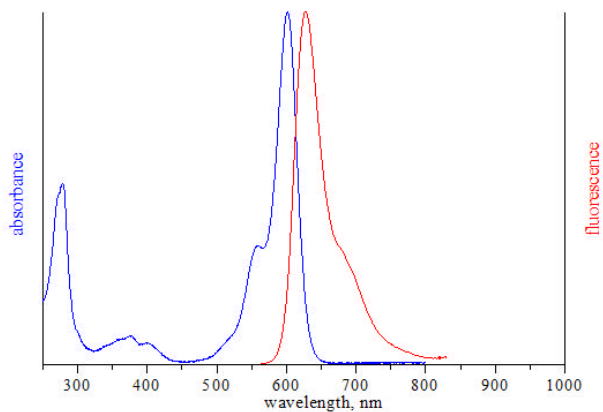
Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 603 nm, λ_{em} 626 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-m⁷GTP-ATTO-594



excitation and emission spectrum of ATTO 594