



EDA-m⁷GTP-5/6-TAMRA

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-triphosphate, labeled with 5/6-TAMRA, Triethylammonium salt

| Cat. No. | Amount |
|------------|---------------|
| NU-824-TAM | 160 µl (1 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: C₃₉H₄₄N₉O₁₉P₃ (free acid)

Molecular Weight: 1035.74 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

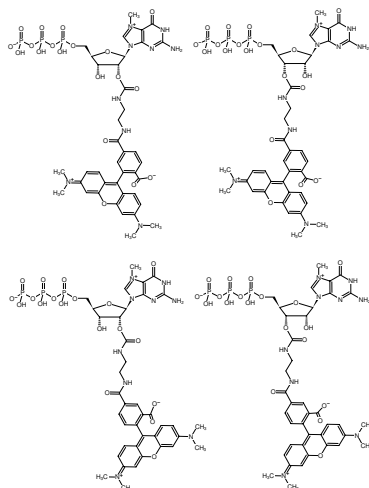
Form: solution in water

Color: pink to red

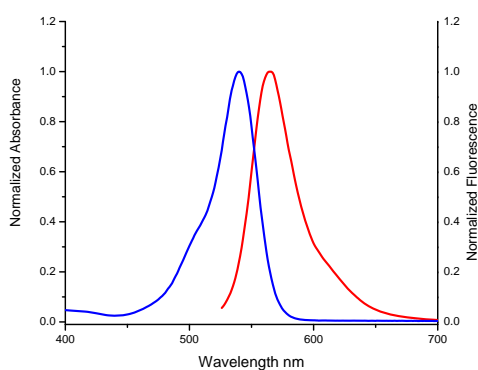
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 545 nm, λ_{em} 575 nm, ε 90.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-m⁷GTP-5/6-TAMRA



excitation and emission spectrum of 5/6-TAMRA