



EDA-m⁷GDP-ATTO-Rho13

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO Rho13, Triethylammonium salt

Cat. No.	Amount
NU-827-RHO13	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: C₅₅H₆₈N₁₀O₁₅P₂ (free acid)

Molecular Weight: 1171.15 g/mol (free acid)

Exact Mass: 1170.43 g/mol (free acid)

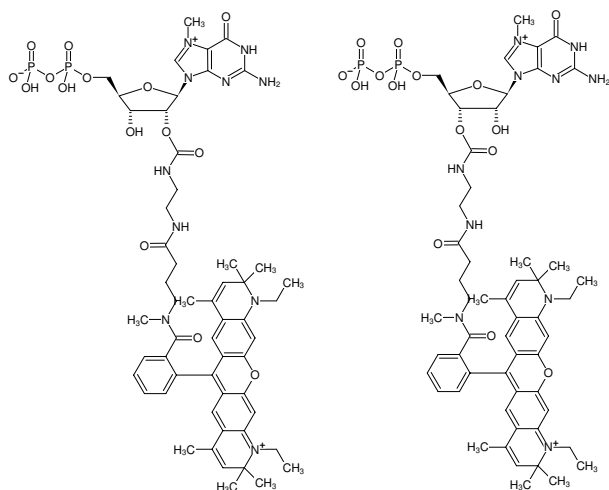
Purity: ≥ 95 % (HPLC)

Form: solution in water

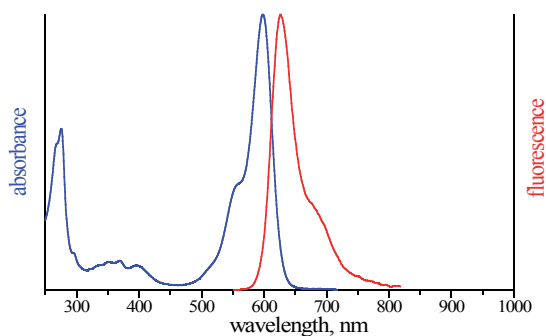
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 603 nm, λ_{em} 627 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-m⁷GDP-ATTO-Rho13



excitation and emission spectrum of ATTO Rho13