



8-(6-Aminoethyl)-amino-GTP - ATTO-Rho13

8-(6-Aminoethyl)-amino-guanosine-5'-triphosphate, labeled with ATTO Rho13, Triethylammonium salt

| Cat. No. | Amount |
|--------------|--------------|
| NU-830-RHO13 | 80 µl (1 mM) |

For research use only!

Shipping: shipped on blue ice

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: C₅₇H₇₅N₁₀O₁₇P₃ (free acid)

Molecular Weight: 1264.45 g/mol (free acid)

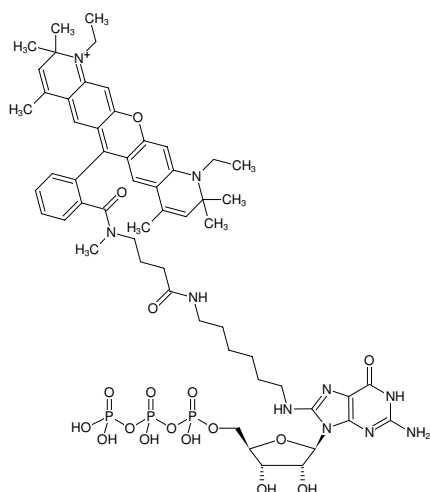
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

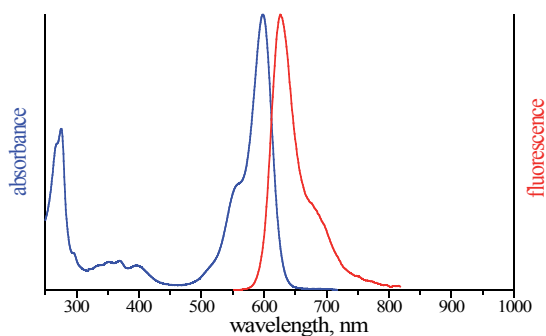
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

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



Structural formula of 8-(6-Aminoethyl)-amino-GTP - ATTO-Rho13



excitation and emission spectrum of ATTO Rho13