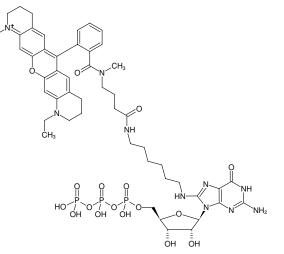




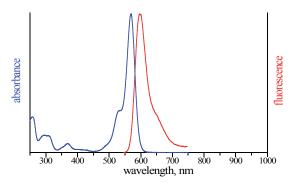
8-(6-Aminohexyl)-amino-GTP-ATTO-Rho11

8-(6-Aminohexyl)-amino-guanosine-5'-triphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-830-RHO11	80 μl (1 mM)



Structural formula of 8-(6-Aminohexyl)-amino-GTP-ATTO-Rho11



excitation and emission spectrum of ATTO Rho11

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: C₅₁H₆₇N₁₀O₁₇P₃ (free acid)

Molecular Weight: 1185.07 g/mol (free acid)

Exact Mass: 1184.39 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} 572 nm, λ_{em} 595 nm, ϵ 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

