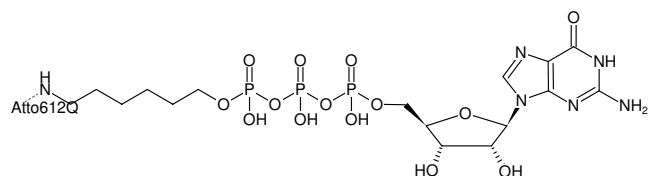




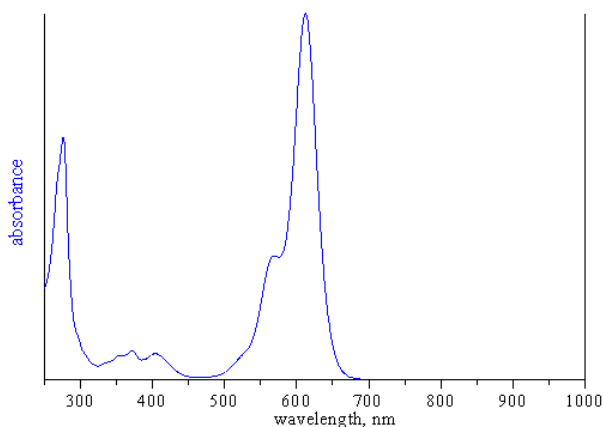
γ -(6-Aminoethyl)-GTP-ATTO-612Q

γ -(6-Aminoethyl)-guanosine-5'-triphosphate, labeled with ATTO 612Q, Triethylammonium salt

Cat. No.	Amount
NU-834-612Q	80 μ l (1 mM)



Structural formula of γ -(6-Aminoethyl)-GTP-ATTO-612Q



absorption spectrum of ATTO 612Q

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: C₁₆H₂₉N₆O₁₄P₃ - ATTO 612Q (free acid)

Molecular Weight: 1294.35 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 615 nm, ϵ 115.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)