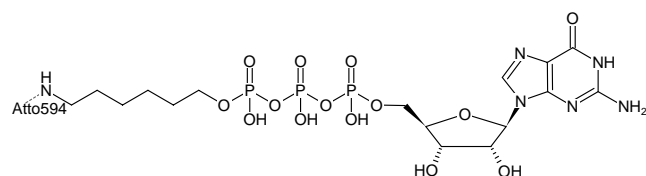




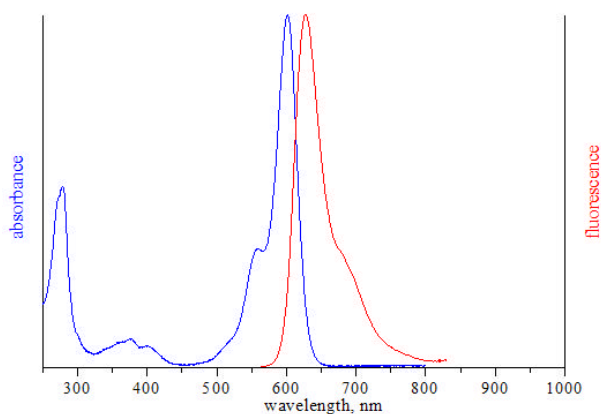
γ -(6-Aminoethyl)-GTP - ATTO-594

γ -(6-Aminoethyl)-guanosine-5'-triphosphate, labeled with ATTO 594, Triethylammonium salt

Cat. No.	Amount
NU-834-594	80 μ l (1 mM)



Structural formula of γ -(6-Aminoethyl)-GTP - ATTO-594



excitation and emission spectrum of ATTO 594

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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months

Molecular Formula: C₁₆H₂₉N₆O₁₄P₃ - ATTO 594 (free acid)

Molecular Weight: 1409.35 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: clear aqueous solution, pH 7.5 \pm 0.5

Concentration: 1 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 601 nm; λ_{em} 627 nm;
 ϵ 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)