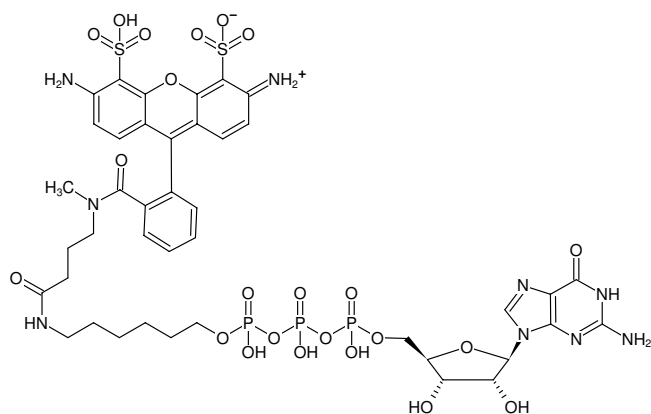




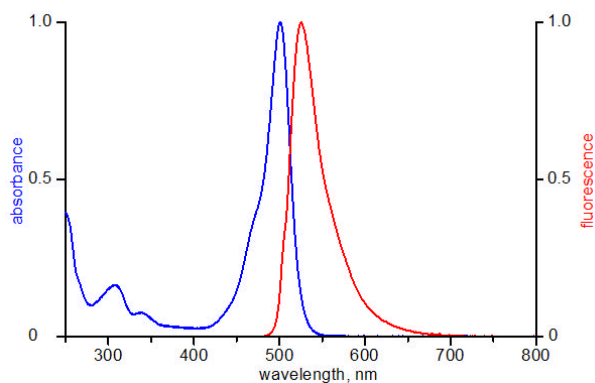
γ -(6-Aminoethyl)-GTP-ATTO-488

γ -(6-Aminoethyl)-guanosine-5'-triphosphate, labeled with ATTO 488, Triethylammonium salt

Cat. No.	Amount
NU-834-488	120 μ l (1 mM)



Structural formula of γ -(6-Aminoethyl)-GTP-ATTO-488



excitation and emission spectrum of ATTO 488

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: C₄₁H₅₀N₉O₂₃P₃S₂ (free acid)

Molecular Weight: 1193.93 g/mol (free acid)

Exact Mass: 1193.17 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: yellow solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 500 nm, λ_{em} 520 nm, ϵ 90.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Furness *et al.* (2016) Ligand-Dependent Modulation of G Protein Conformation Alters Drug Efficacy. *Cell*. **167** (3):739.