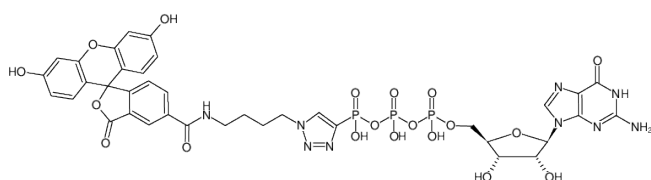




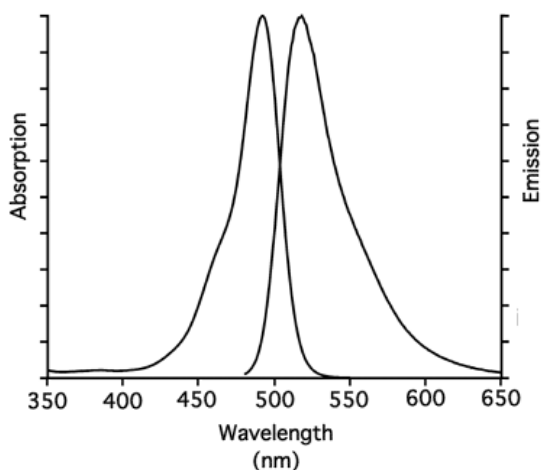
γ -[(4-Aminobutyl)triazolo]-GTP-5-FAM

γ -[(4-Aminobutyl)triazolo]- guanosine-5'-triphosphate , labeled with 5-FAM, Triethylammonium salt
Guanosine-5'-O-diphosphate[γ -((4-phosphono-1H-1,2,3-triazol-1-yl)butyl-4-amine-5-FAM)],
GTP γ -Triazolo-4-AB-5-FAM

Cat. No.	Amount
NU-1219-5FM	10 μ l (1 mM)



Structural formula of γ -[(4-Aminobutyl)triazolo]-GTP-5-FAM



excitation and emission spectrum of 5-FAM

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: C₃₇H₃₆N₉O₁₉P₃ (free acid)

Molecular Weight: 1003.66 g/mol (free acid)

Exact Mass: 1003.13 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: yellow to orange

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 492 nm, λ_{em} 517 nm, ϵ 83.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Pepanian *et al.* (2022) A fluorescence anisotropy assay with guanine nucleotides provides access to functional analysis of Gai1 proteins. *Anal Chem.* **94(41)**:14410.