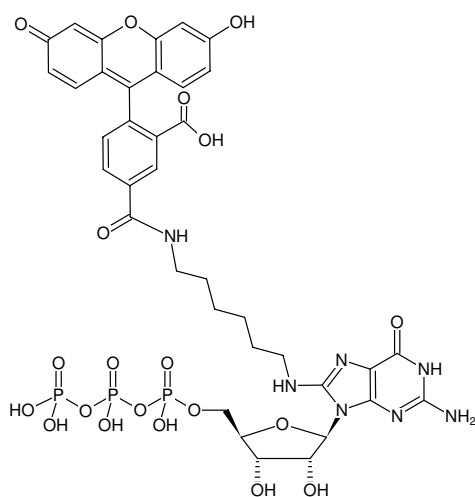




8-(6-Aminoethyl)-amino-GTP-5-FAM

8-(6-Aminoethyl)-amino-guanosine-5'-triphosphate, labeled with 5 FAM, Triethylammonium salt

Cat. No.	Amount
NU-830-5FM	160 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-GTP-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: C₃₇H₄₀N₇O₂₀P₃ (free acid)

Molecular Weight: 995.68 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

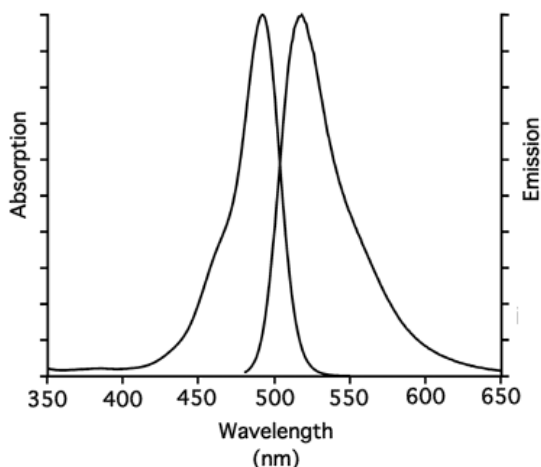
Form: solution in water

Color: yellow

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 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)



excitation and emission spectrum of 5-FAM