



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

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

Cat. No.	Amount
NU-829-5FM	160 µl (1 mM)

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: C₃₇H₃₈N₇O₁₄P (free acid)

Molecular Weight: 835.72 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

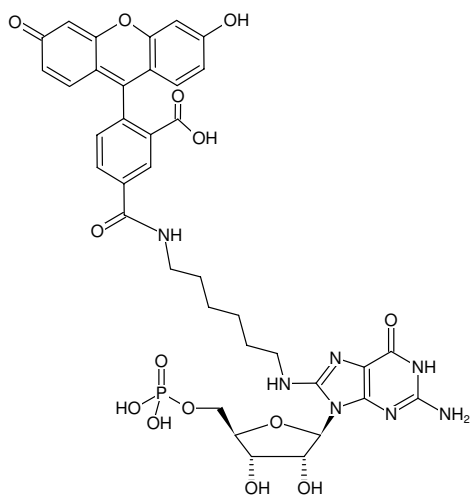
Form: solution in water

Color: yellow

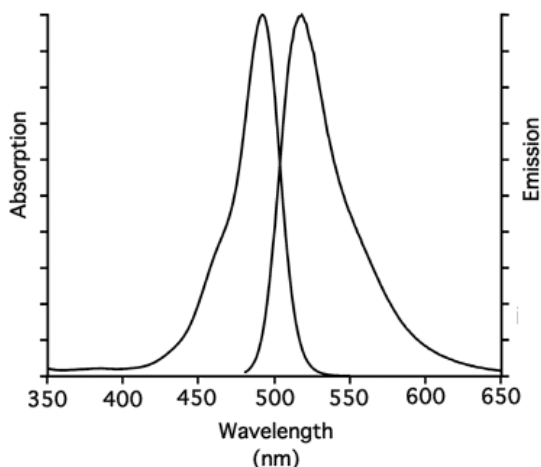
Concentration: 1.0 mM - 1.1 mM

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)



Structural formula of 8-(6-Aminoethyl)-amino-GMP-5-FAM



excitation and emission spectrum of 5-FAM