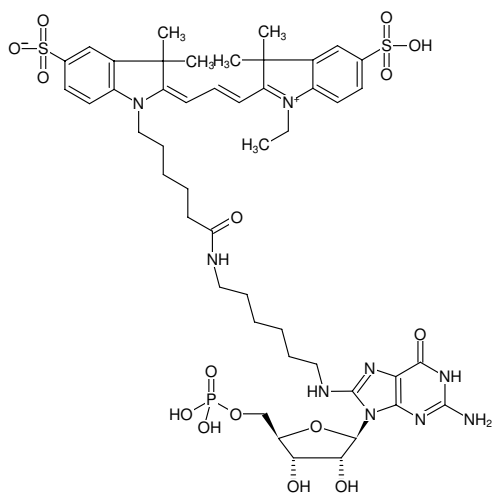




8-(6-Aminoethyl)-amino-GMP-Cy3

8-(6-Aminoethyl)-amino-guanosine-5'-monophosphate, labeled with Cy3, Triethylammonium salt

| Cat. No. | Amount |
|------------|--------------------|
| NU-829-CY3 | 160 μ l (1 mM) |



Structural formula of 8-(6-Aminoethyl)-amino-GMP-Cy3

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: C₄₇H₆₄N₉O₁₅PS₂ (free acid)

Molecular Weight: 1090.17 g/mol (free acid)

Purity: \geq 95 % (HPLC)

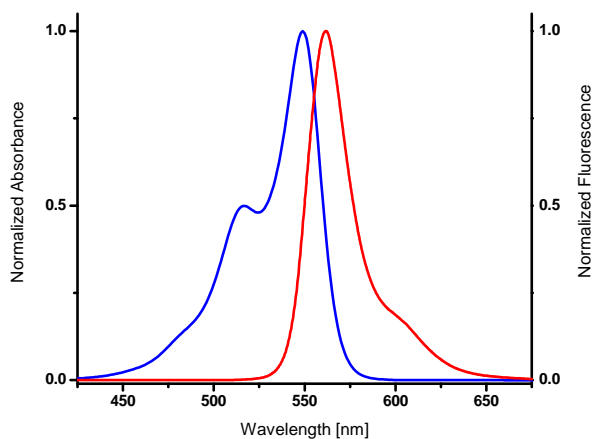
Form: solution in water

Color: pink

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 550 nm, λ_{em} 570 nm, ϵ 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of Cy3