



γ -(6-Aminoethyl)-dGTP - Cy3

γ -(6-Aminoethyl)-2'-deoxyguanosine-5'-triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-916-CY3	120 μ l (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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months

Molecular Formula: C₄₇H₆₅N₈O₂₀P₃S₂ (free acid)

Molecular Weight: 1219.11 g/mol (free acid)

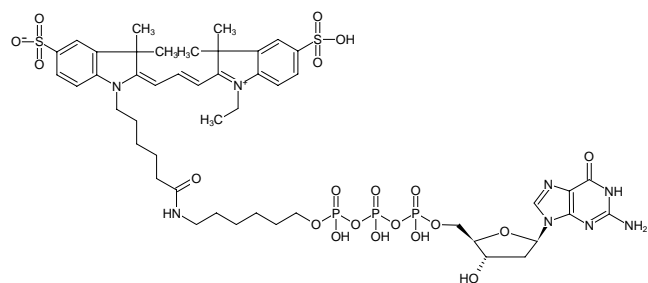
Purity: \geq 95 % (HPLC)

Form: pink solution in water, pH 7.5 \pm 0.5

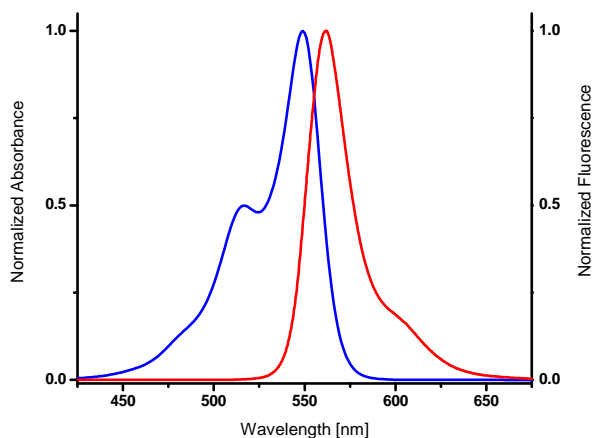
Concentration: 1 mM - 1.1 mM

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)



Structural formula of γ -(6-Aminoethyl)-dGTP - Cy3



excitation and emission spectrum of Cy3