

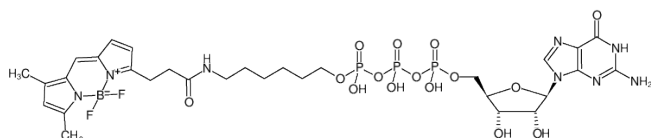


γ -(6-Aminoethyl)-GTP-BDP-FL

structural analog to BODIPY® FL

γ -(6-Aminoethyl)-guanosine-5'-triphosphate, labeled with BDP-FL, Triethylammonium salt

Cat. No.	Amount
NU-834-BDP	160 μ l (1 mM)



Structural formula of γ -(6-Aminoethyl)-GTP-BDP-FL

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: C₃₀H₄₂BF₂N₈O₁₅P₃ (free acid)

Molecular Weight: 896.44 g/mol (free acid)

Exact Mass: 896.20 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: orange

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 503 nm, λ_{em} 512 nm, ϵ 80.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)