



γ -(6-Aminoethyl)-N⁶-Benzyl-ATP-ATTO-647N

γ -(6-Aminoethyl)-N⁶-benzyl-adenosine-5'-triphosphate, labeled with ATTO 647N, Triethylammonium salt

Cat. No.	Amount
NU-911-647N	80 μ l (1 mM)

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: C₆₅H₈₄N₉O₁₅P₃ (free acid)

Molecular Weight: 1324.34 g/mol (free acid)

Exact Mass: 1323.53 g/mol (free acid)

Purity: \geq 95 % (HPLC)

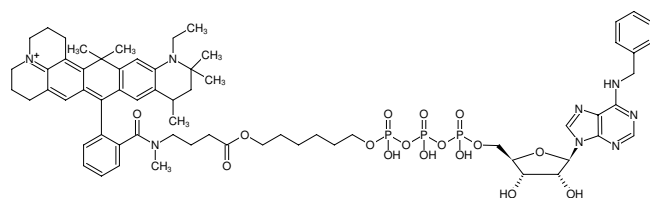
Form: solution in water

Color: blue

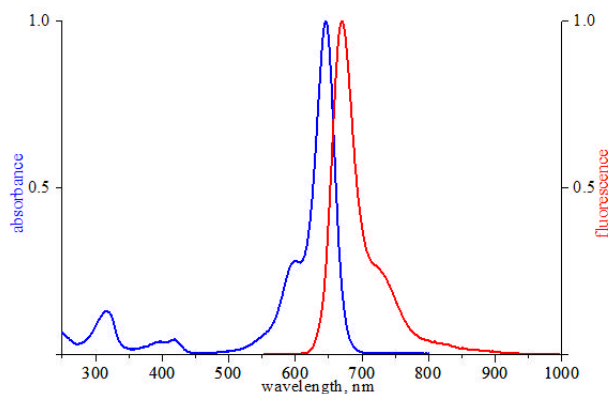
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 646 nm, λ_{em} 664 nm, ϵ 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of γ -(6-Aminoethyl)-N⁶-Benzyl-ATP-ATTO-647N



excitation and emission spectrum of ATTO 647N