



N⁶-(6-Aminoethyl)-dATP-DYQ-660

N⁶-(6-Aminoethyl)-2'-deoxyadenosine-5'-triphosphate, labeled with DYQ 660, Triethylammonium salt

Cat. No.	Amount
NU-835-660Q	40 µl (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: C₅₅H₇₁N₈O₁₇P₃S (free acid)

Molecular Weight: 1241.19 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

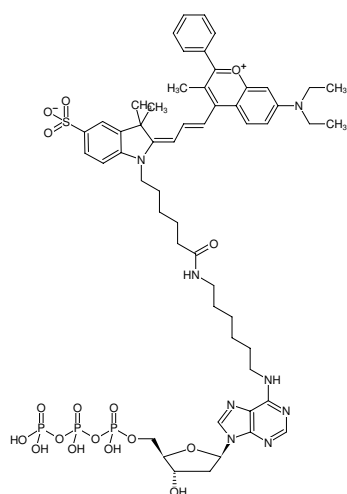
Form: solution in water

Concentration: 1.0 mM - 1.1 mM

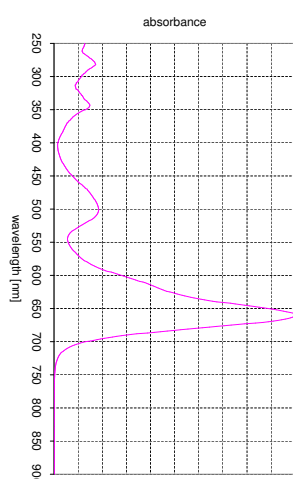
pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 660 nm, ε 109.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

λ_{exc} 660 nm, ε 140.0 L mmol⁻¹ cm⁻¹ (Ethanol)



Structural formula of N⁶-(6-Aminoethyl)-dATP-DYQ-660



absorption spectrum of DYQ 660