



N⁶-(6-Aminohexyl)-dATP-ATTO-565

N⁶-(6-Aminohexyl)-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO 565, Triethylammonium salt

Cat. No.	Amount
NU-835-565	120 µl (1 mM)

For research use only!

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: C₄₇H₅₇N₈O₁₆P₃ (free acid)

Molecular Weight: 1082.93 g/mol (free acid)

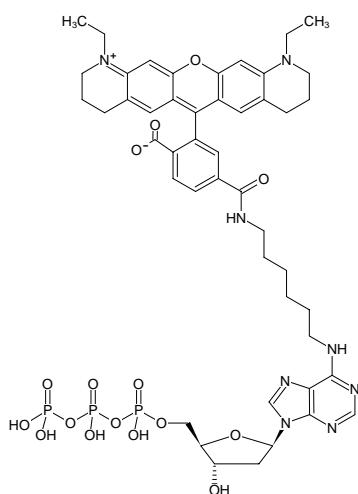
Purity: ≥ 95 % (HPLC)

Form: solution in water

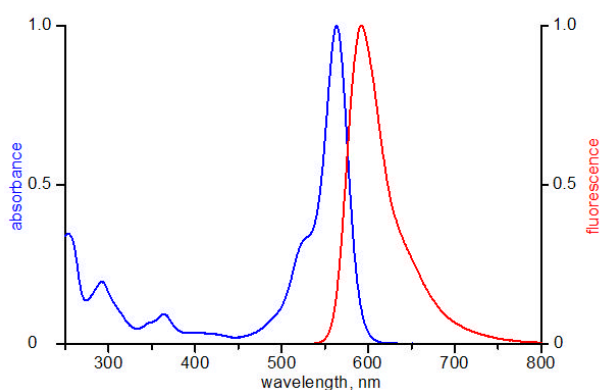
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 564 nm, λ_{em} 590 nm, ε 120.0 L mol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of N⁶-(6-Aminohexyl)-dATP-ATTO-565



excitation and emission spectrum of ATTO 565