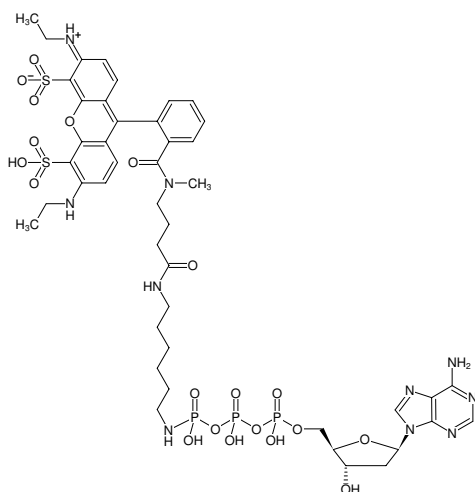




γ -[(6-Aminohexyl)-imido]-dATP-ATTO-532

γ -[(6-Aminohexyl)-imido]-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO-532, Triethylammonium salt

Cat. No.	Amount
NU-861-532	60 μ l (1 mM)



Structural formula of γ -[(6-Aminohexyl)-imido]-dATP-ATTO-532

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: C₄₅H₅₉N₁₀O₂₀P₃S₂ (free acid)

Molecular Weight: 1217.06 g/mol (free acid)

Exact Mass: 1216.26 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

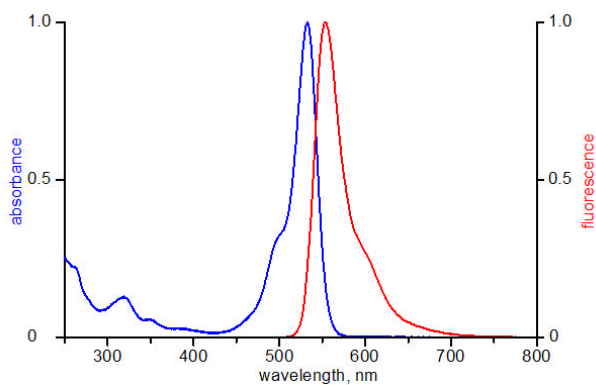
Color: orange

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 532 nm, λ_{em} 552 nm, ϵ 115.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Please note: This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.0.



excitation and emission spectrum of ATTO 532