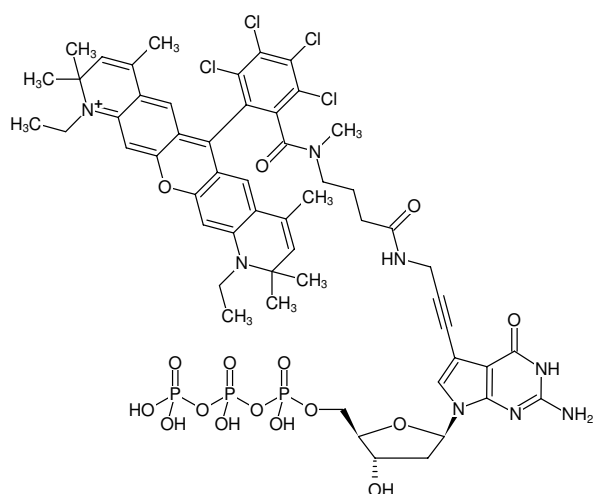




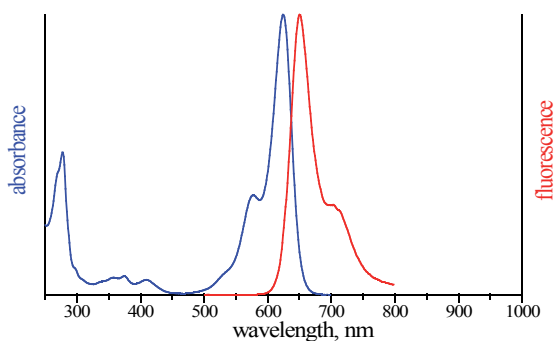
7-Propargylamino-7-deaza-dGTP-ATTO-Rho14

7-Deaza-7-propargylamino-2'-deoxyguanosine-5'-triphosphate, labeled with ATTO Rho14, Triethylammonium salt

Cat. No.	Amount
NU-1615-RHO14	40 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-dGTP-ATTO-Rho14



excitation and emission spectrum of ATTO Rho14

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: C₅₅H₆₁Cl₄N₈O₁₆P₃ (free acid)

Molecular Weight: 1324.85 g/mol (free acid)

Exact Mass: 1322.22 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} 626 nm, λ_{em} 646 nm, ε 140.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)