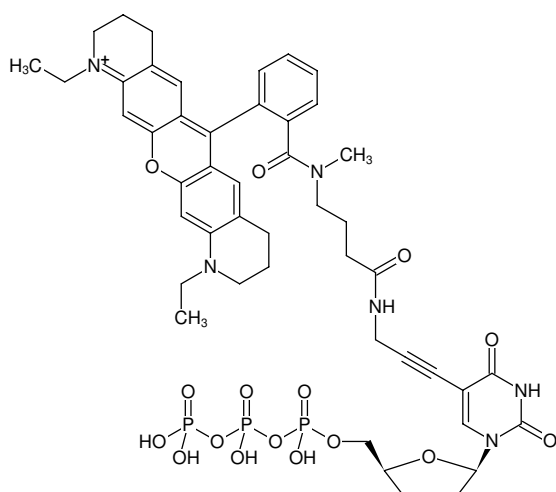




5-Propargylamino-ddUTP-ATTO-Rho11

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-1619-RHO11	80 μ l (0,5 mM)



Structural formula of 5-Propargylamino-ddUTP-ATTO-Rho11

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: C₄₇H₅₅N₆O₁₆P₃ (free acid)

Molecular Weight: 1052.90 g/mol (free acid)

Exact Mass: 1052.29 g/mol (free acid)

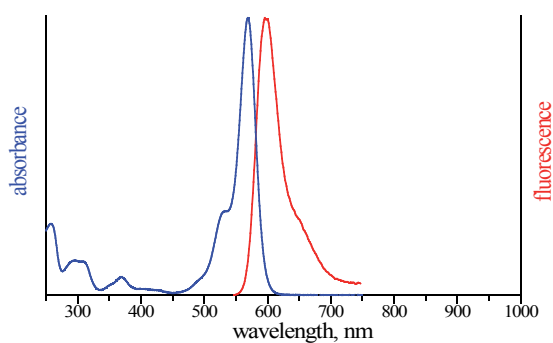
Purity: \geq 95 % (HPLC)

Form: solution in water

Concentration: 0.50 mM - 0.55 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 572 nm, λ_{em} 595 nm, ϵ 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO Rho11