



5-Propargylamino-CTP-ATTO-Rho12

5-Propargylamino-cytidine-5'-triphosphate, labeled with ATTO Rho12, Triethylammonium salt

Cat. No.	Amount
NU-831-RHO12	40 μ l (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: C₅₃H₆₈N₇O₁₇P₃ (free acid)

Molecular Weight: 1168.08 g/mol (free acid)

Exact Mass: 1167.39 g/mol (free acid)

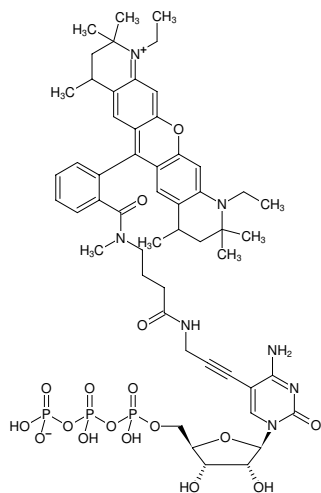
Purity: \geq 95 % (HPLC)

Form: solution in water

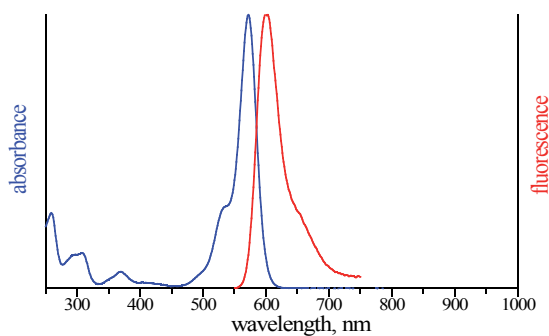
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 576 nm, λ_{em} 601 nm, ϵ 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 5-Propargylamino-CTP-ATTO-Rho12



excitation and emission spectrum of ATTO Rho12