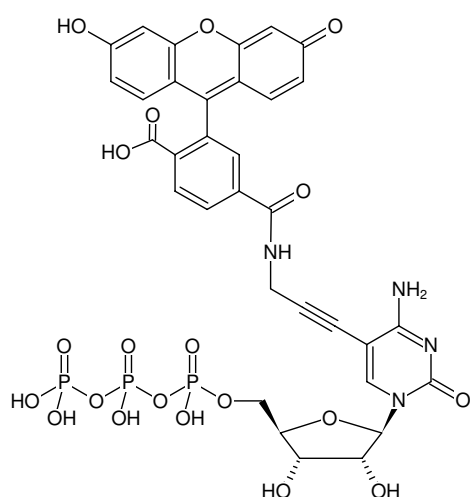




5-Propargylamino-CTP-6-FAM

5-Propargylamino-cytidine-5'-triphosphate, labeled with 6 FAM, Triethylammonium salt

Cat. No.	Amount
NU-831-6FM	30 µl (1 mM)



Structural formula of 5-Propargylamino-CTP-6-FAM

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: C₃₃H₂₉N₄O₂₀P₃ (free acid)

Molecular Weight: 894.52 g/mol (free acid)

Exact Mass: 894.06 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

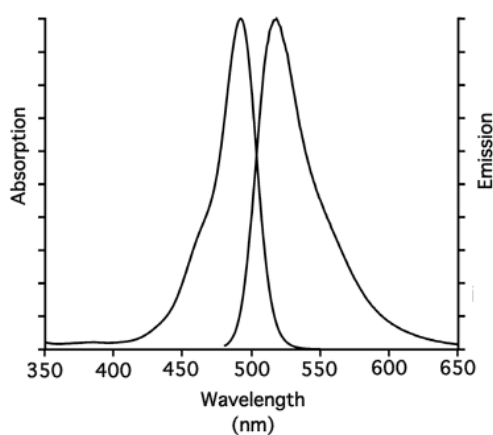
Form: solution in water

Color: yellow

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 492 nm, λ_{em} 517 nm, ε 83.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of 6-FAM