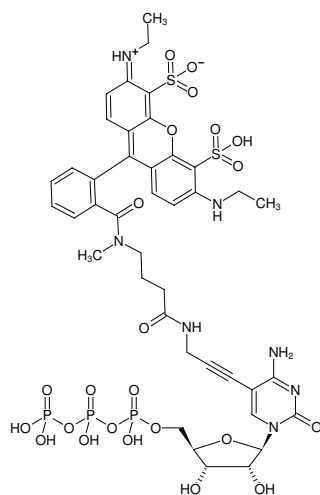




5-Propargylamino-CTP-ATTO-532

5-Propargylamino-cytidine-5'-triphosphate, labeled with ATTO 532, Triethylammonium salt

Cat. No.	Amount
NU-831-532	80 µl (1 mM)



Structural formula of 5-Propargylamino-CTP-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: C₄₁H₄₈N₇O₂₃P₃S₂ (free acid)

Molecular Weight: 1163.90 g/mol (free acid)

Exact Mass: 1163.15 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

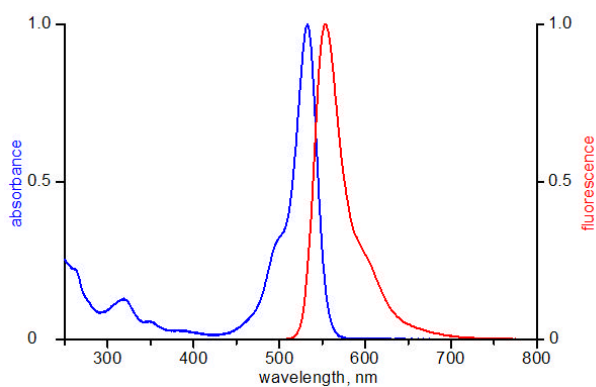
Form: solution in water

Color: orange

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 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)



excitation and emission spectrum of ATTO 532