



5-Propargylamino-ddCTP-ATTO-540Q

5-Propargylamino-2',3'-dideoxycytidine-5'-triphosphate, labeled with ATTO 540Q, Triethylammonium salt

Cat. No.	Amount
NU-850-540Q	80 µl (0,5 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

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: C₁₂H₁₉N₄O₁₂P₃ - ATTO 540Q (free acid)

Molecular Weight: 1044.22 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

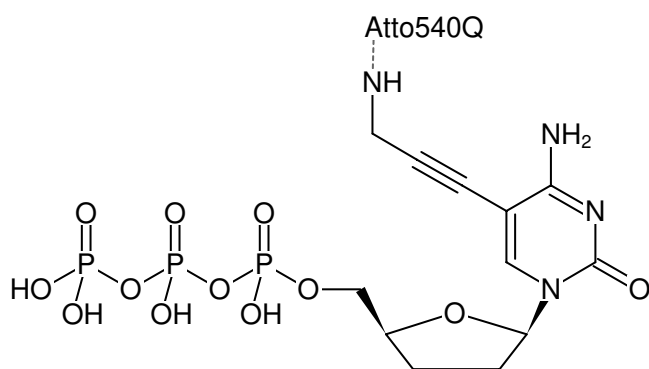
Form: solution in water

Color: red

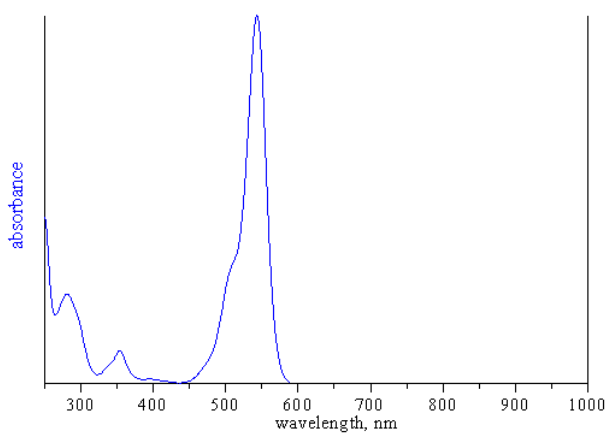
Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 543 nm, ε 105.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 5-Propargylamino-ddCTP-ATTO-540Q



absorption spectrum of ATTO 540Q