



5-Propargylamino-ddUTP-DYQ-660

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with DYQ 660, Triethylammonium salt

Cat. No.	Amount
NU-1619-660Q	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: C₅₁H₆₀N₅O₁₈P₃S (free acid)

Molecular Weight: 1156.04 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

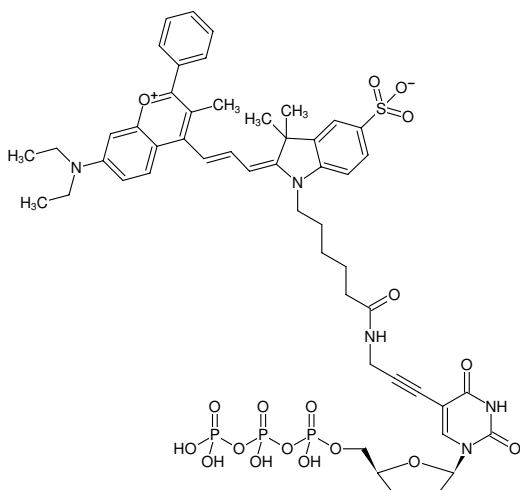
Color: blue

Concentration: 0.50 mM - 0.55 mM

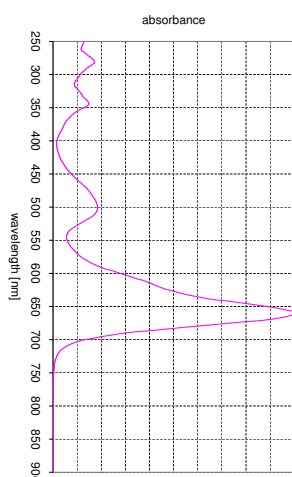
pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 660 nm, ε 109.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

λ_{exc} 660 nm, ε 140.0 L mmol⁻¹ cm⁻¹ (Ethanol)



Structural formula of 5-Propargylamino-ddUTP-DYQ-660



absorption spectrum of DYQ 660