



5-Propargylamino-ddUTP - DY-485XL

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with DY 485XL, Triethylammonium salt

Cat. No.	Amount
NU-1619-485XL	80 µl (0,5 mM)

For research use only!

Shipping: shipped on blue ice

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: C₃₇H₄₆N₅O₁₉P₃S (free acid)

Molecular Weight: 989.77 g/mol (free acid)

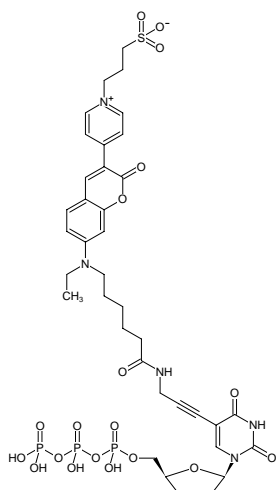
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

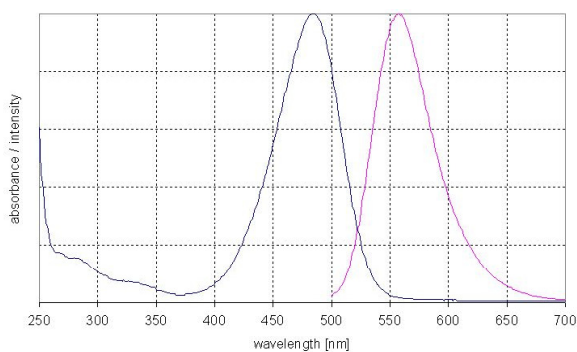
Concentration: 0.5 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 485 nm; λ_{em} 560 nm; ε 50.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 5-Propargylamino-ddUTP - DY-485XL



excitation and emission spectrum of DY 485XL