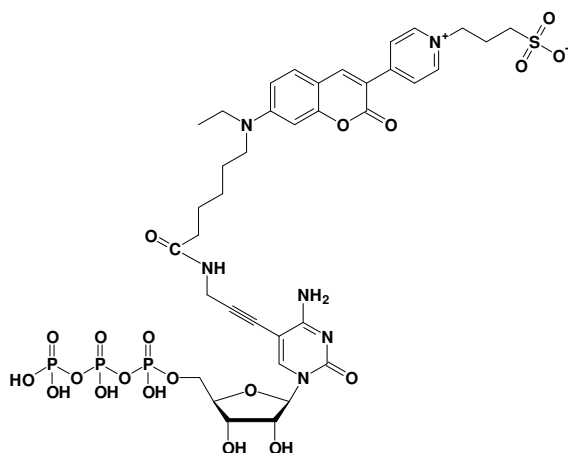


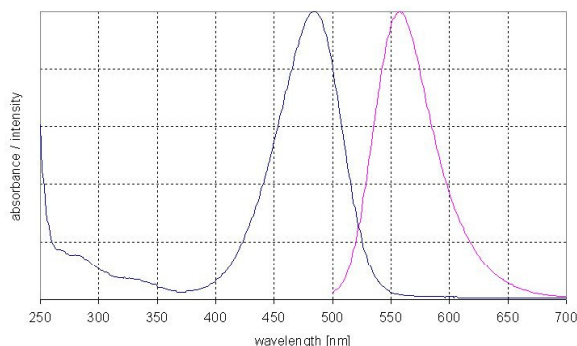
5-Propargylamino-CTP - DY 485XL

5-Propargylamino-cytidine-5'-triphosphate, labeled with DY 485XL, Triethylammonium salt

Cat. No.	Amount
NU-831-485XL	40 μ l (1 mM)



structural formula of 5-Propargylamino-CTP-DY 485XL



excitation and emission spectrum of DY 485XL

Cat. No.: NU-831-485XL

Molecular Formula: C₃₇H₄₇N₆O₂₀P₃S (free acid)

Molecular Weight: 1020,78 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{exc} 485 nm; λ_{em} 560 nm; ϵ 50,000 cm⁻¹ M⁻¹

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!