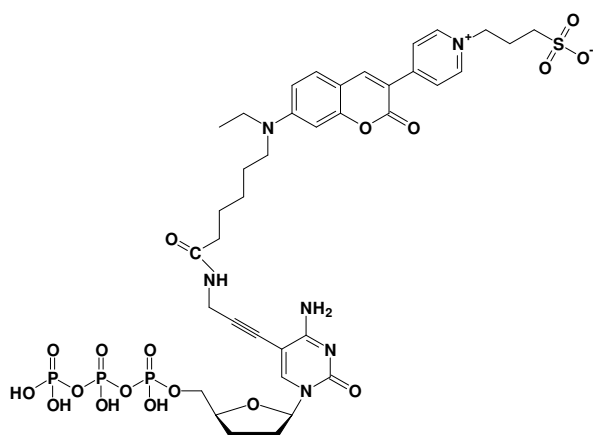


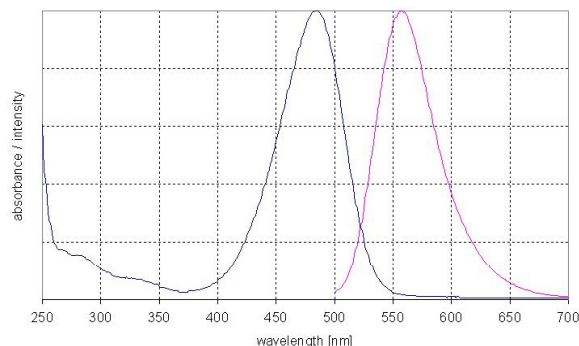
5-Propargylamino-ddCTP - DY 485XL

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

Cat. No.	Amount
NU-850-485XL	160 μ l (0,5 mM)



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



excitation and emission spectrum of DY 485XL

Cat. No.: NU-850-485XL

Molecular Formula: C₃₇H₄₇N₆O₁₈P₃S (free acid)

Molecular Weight: 988,79 (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!