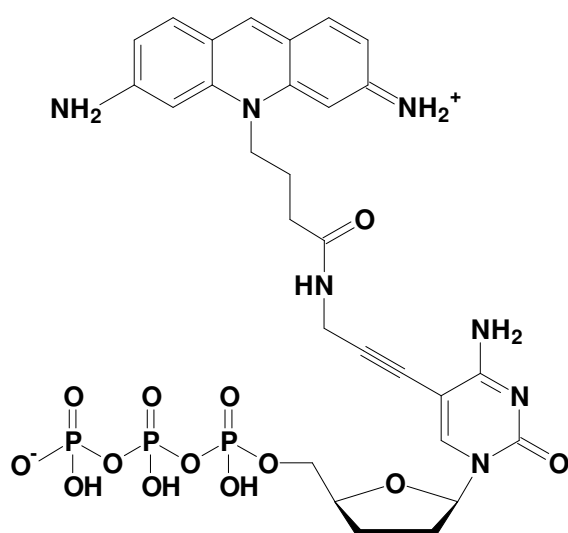


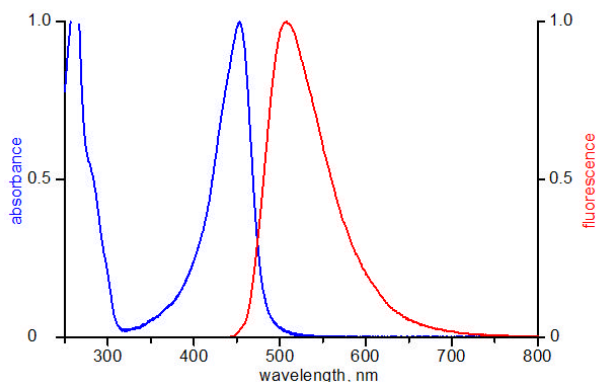
5-Propargylamino-ddCTP - ATTO 465

5-Propargylamino-2',3'-Dideoxy-cytidine-5'-triphosphate, labeled with ATTO 465,
 Triethylammonium salt

Cat. No.	Amount
NU-850-465	160 μ l (0,5 mM)



structural formula of 5-Propargylamino-ddCTP - ATTO 465



excitation and emission spectrum of ATTO 465

Cat. No.: NU-850-465

Molecular Formula: $C_{29}H_{34}N_7O_{13}P_3$ (free acid)

Molecular Weight: 781.54 (free acid)

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

Spectroscopic properties:

λ_{exc} 453 nm; λ_{em} 508 nm; ϵ 75,000 $cm^{-1} M^{-1}$

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!