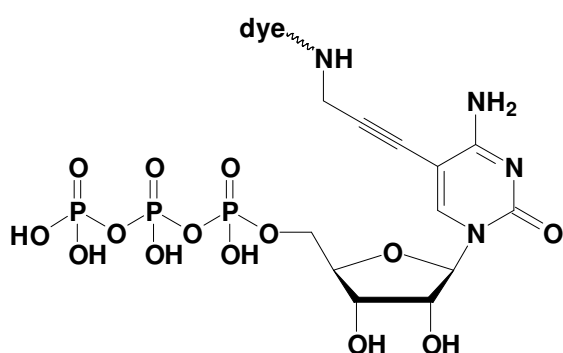


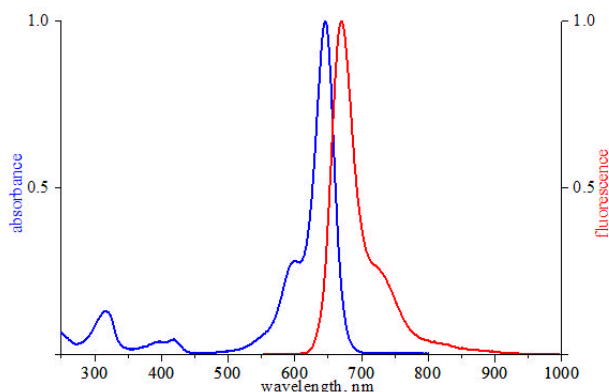
5-Propargylamino-CTP - ATTO 647N

5-Propargylamino-cytidine-5'-triphosphate, labeled with ATTO 647N, Triethylammonium salt

Cat. No.	Amount
NU-831-647N	40 μ l (1 mM)



structural formula of 5-Propargylamino-CTP



excitation and emission spectrum of ATTO 647N

Cat. No.: NU-831-647N

Molecular Formula: $C_{12}H_{19}N_4O_{14}P_3$ - ATTO 647N (free acid)

Molecular Weight: 1163.22 (free acid)

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

Spectroscopic properties:

λ_{exc} 644 nm; λ_{em} 669 nm; ϵ 150,000 $cm^{-1} M^{-1}$

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!