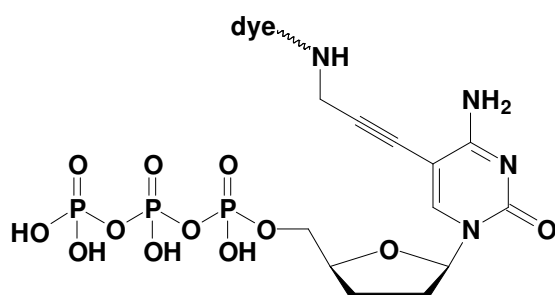


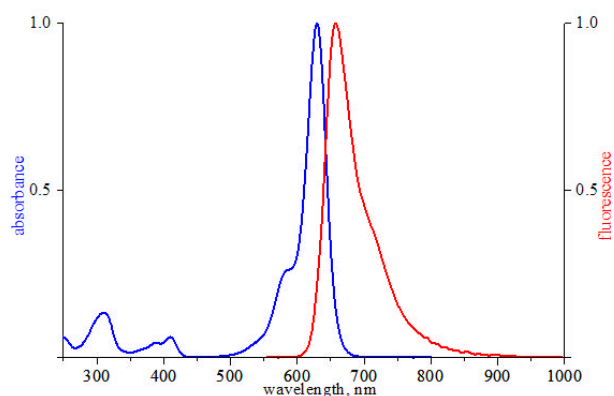
5-Propargylamino-ddCTP - ATTO 633

5-Propargylamino-2',3'-Dideoxy-cytidine-5'-triphosphate, labeled with ATTO 633,
 Triethylammonium salt

Cat. No.	Amount
NU-850-633	80 μ l (0,5 mM)



structural formula of 5-Propargylamino-ddCTP



excitation and emission spectrum of ATTO 633

Cat. No.: NU-850-633

Molecular Formula: C₁₂H₁₉N₄O₁₂P₃ - ATTO 633 (free acid)

Molecular Weight: 1037.22 (free acid)

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

Spectroscopic properties:

λ_{exc} 629 nm; λ_{em} 657 nm; ϵ 130,000 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!