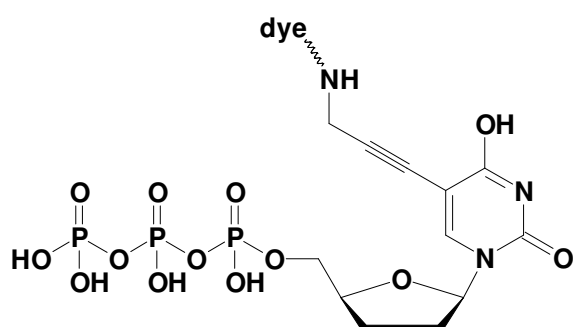


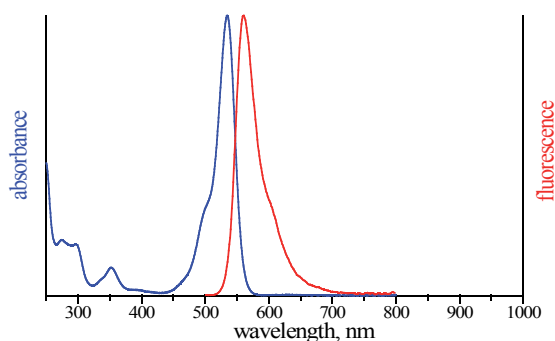
5-Propargylamino-ddUTP - ATTO Rho6G

5-Propargylamino- 2', 3'-Dideoxy-uridine-5'-triphosphate, labeled with ATTO Rho6G,
 Triethylammonium salt

Cat. No.	Amount
NU-1619-RHO6G	80 μ l (0,5 mM)



structural formula of 5-Propargylamino-ddUTP



excitation and emission spectrum of ATTO Rho6G

Cat. No.: NU-1619-RHO6G

Molecular Formula: C₁₂H₁₈N₃O₁₃P₃ - ATTO Rho6G
 (free acid)

Molecular Weight: 1000,20 (free acid)

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

Spectroscopic properties:

λ_{exc} 535 nm; λ_{em} 560 nm; ϵ 1,15 x 10⁵ 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!