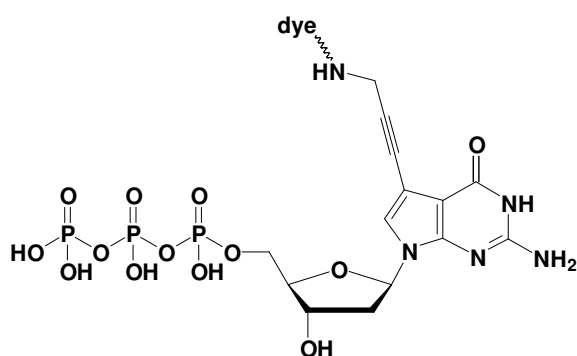


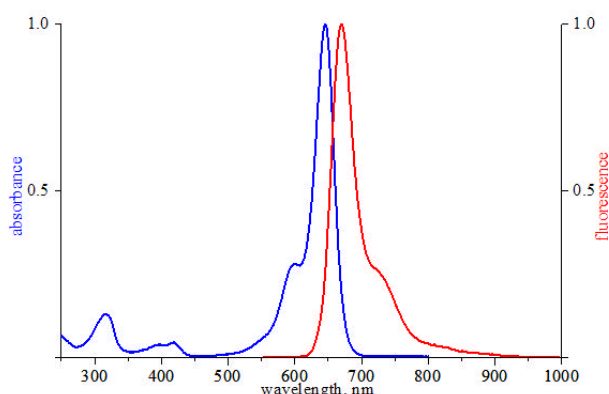
7-Propargylamino-7-deaza-dGTP - ATTO 647N

7-Deaza-7-propargylamino-2'-deoxy-guanosine-5'-triphosphate, labeled with ATTO 647N, Triethylammonium salt

Cat. No.	Amount
NU-1615-647N	40 μ l (1 mM)



structural formula of 7-Propargylamino-7-deaza-dGTP



excitation and emission spectrum of ATTO 647N

Cat. No.: NU-1615-647N

Molecular Formula: C₁₄H₂₀N₅O₁₃P₃ - ATTO 647N (free acid)

Molecular Weight: 1186.25 (free acid)

Purity: > 90%, contains approximately 3 – 7% labeled diphosphate, 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!