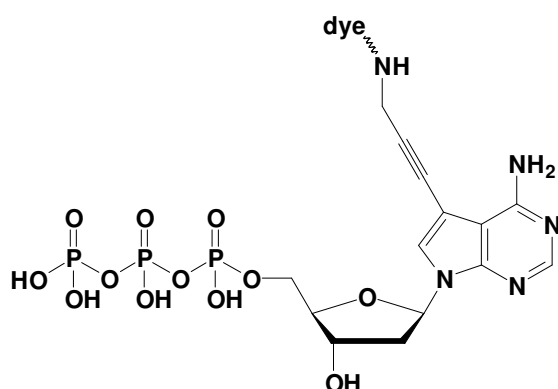


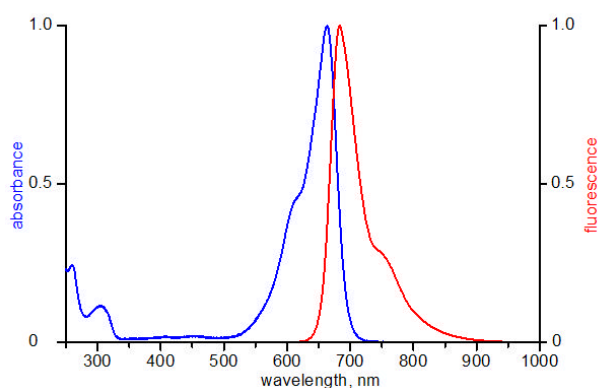
7-Propargylamino-7-deaza-dATP - ATTO 655

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with ATTO 655, Triethylammonium salt

Cat. No.	Amount
NU-1611-655	40 μ l (1 mM)



structural formula of 7-Propargylamino-7-deaza-dATP



excitation and emission spectrum of ATTO 655

Cat. No.: NU-1611-655

Molecular Formula: C₁₄H₂₀N₅O₁₂P₃ - ATTO 655 (free acid)

Molecular Weight: 1052.26 (free acid)

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

Spectroscopic properties:

λ_{exc} 663 nm; λ_{em} 684 nm; ϵ 125,000 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!