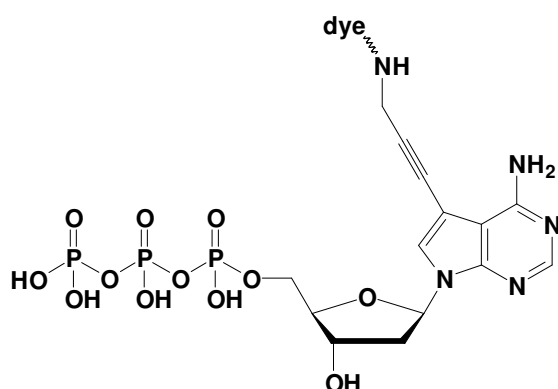


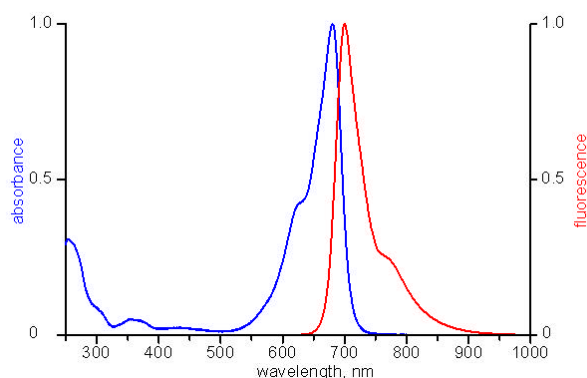
7-Propargylamino-7-deaza-dATP - ATTO 680

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with ATTO 680, Triethylammonium salt

Cat. No.	Amount
NU-1611-680	10 μ l / 1 mM



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



excitation and emission spectrum of ATTO 680

Cat. No.: NU-1611-680

Molecular Formula: C₁₄H₂₀N₅O₁₂P₃ - ATTO 680 (free acid)

Molecular Weight: 1050.26 (free acid)

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

Spectroscopic properties:

λ_{exc} 680 nm; λ_{em} 700 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!