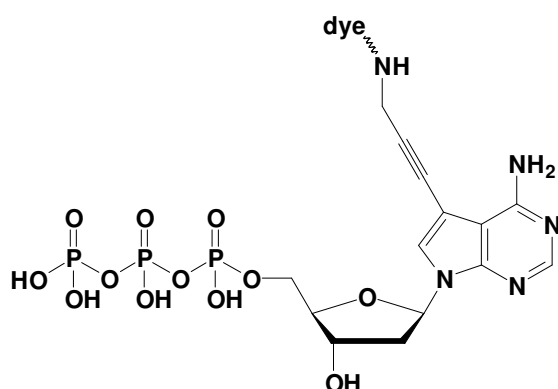


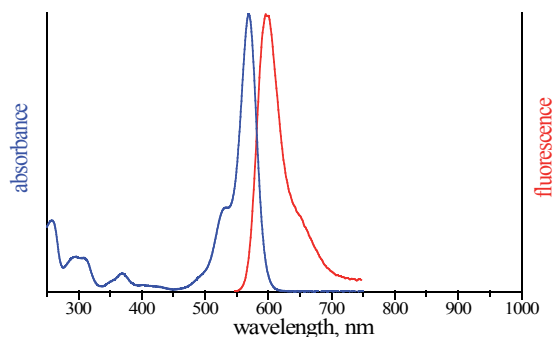
7-Propargylamino-7-deaza-dATP - ATTO Rho 11

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with ATTO Rho 11, Triethylammonium salt

Cat. No.	Amount
NU-1611-RHO11	40 μ l (1 mM)



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



excitation and emission spectrum of ATTO Rho 11

Cat. No.: NU-1611-RHO11

Molecular Formula: C₁₄H₂₀N₅O₁₂P₃ - ATTO Rho11 (free acid)

Molecular Weight: 1091,26 (free acid)

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

Spectroscopic properties:

λ_{exc} 571 nm; λ_{em} 595 nm; ϵ 1,2 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!