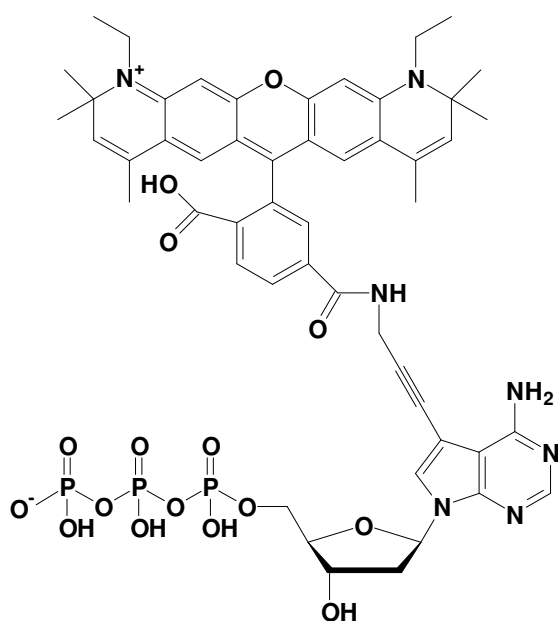


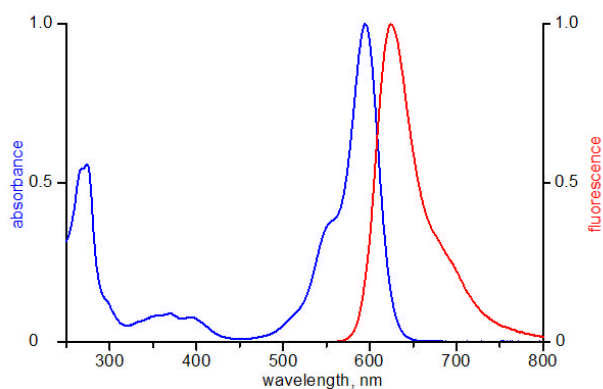
7-Propargylamino-7-deaza-dATP - ATTO 590

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with ATTO 590, Triethylammonium salt

Cat. No.	Amount
NU-1611-590	20 μ l / 1 mM



structural formula of 7-Propargylamino-7-deaza-dATP - ATTO 590



excitation and emission spectrum of ATTO 590

Cat. No.: NU-1611-590

Molecular Formula: C₅₁H₅₆N₇O₁₆P₃ (free acid)

Molecular Weight: 1115.96 (free acid)

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

Spectroscopic properties:

λ_{exc} 594 nm; λ_{em} 624 nm; ϵ 120,000 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!