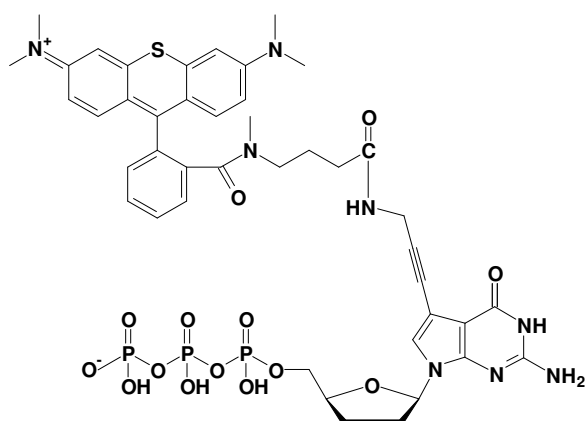


7-Propargylamino-7-deaza-ddGTP - ATTO Thio12

7-Deaza-7-propargylamino-2',3'-dideoxy-guanosine-5'-triphosphate, labeled with ATTO Thio12, Triethylammonium salt

Cat. No.	Amount
NU-1618-THIO12	80 μ l (0,5 mM)



structural formula of 7-Propargylamino-7-deaza-ddGTP - ATTO Thio12

Cat. No.: NU-1618-THIO12

Molecular Formula: C₄₃H₄₉N₈O₁₄P₃S (free acid)

Molecular Weight: 1026,88 (free acid)

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

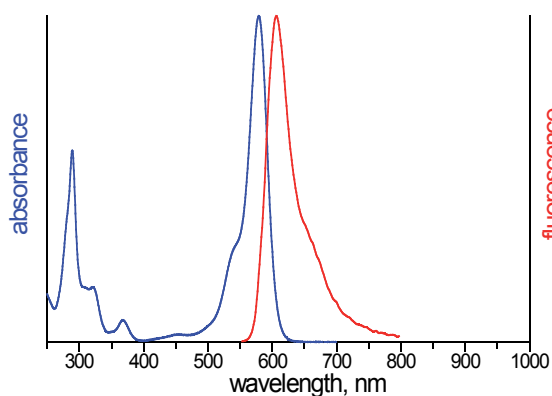
Spectroscopic properties:

λ_{exc} 579 nm; λ_{em} 609 nm; ϵ 1,1 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!



excitation and emission spectrum of ATTO Thio12