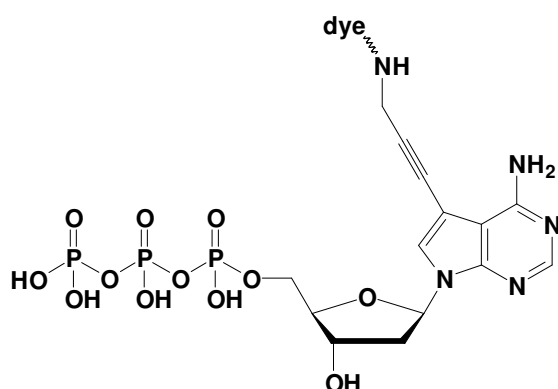


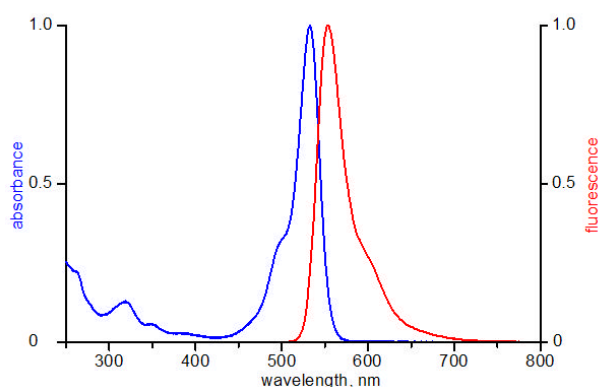
7-Propargylamino-7-deaza-dATP - ATTO 532

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with ATTO 532, Triethylammonium salt

Cat. No.	Amount
NU-1611-532	40 μ l (1 mM)



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



excitation and emission spectrum of ATTO 532

Cat. No.: NU-1611-532

Molecular Formula: $C_{14}H_{20}N_5O_{12}P_3$ - ATTO 532 (free acid)

Molecular Weight: 1170.26 (free acid)

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

Spectroscopic properties:

λ_{exc} 532 nm; λ_{em} 553 nm; ϵ 115,000 $cm^{-1} M^{-1}$

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at $-20^{\circ}C$. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!