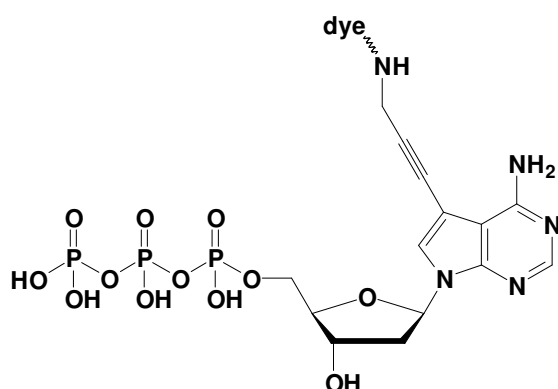


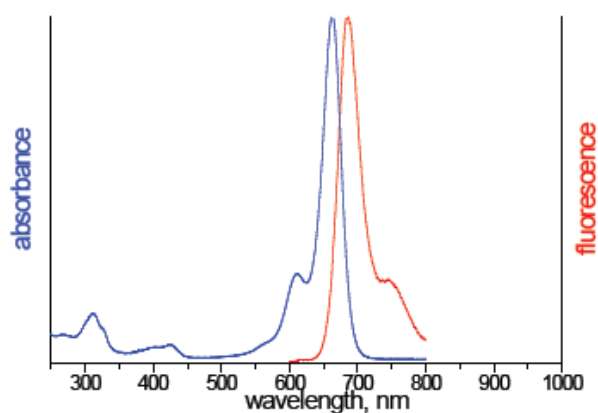
7-Propargylamino-7-deaza-dATP - ATTO 665

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with ATTO 665, Triethylammonium salt

Cat. No.	Amount
NU-1611-665	400 μ l (0,1 mM)



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



excitation and emission spectrum of ATTO 665

Cat. No.: NU-1611-665

Molecular Formula: $C_{14}H_{20}N_5O_{12}P_3$ - ATTO 665 (free acid)

Molecular Weight: 1147,26 (free acid)

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

Spectroscopic properties:

λ_{exc} 663 nm; λ_{em} 684 nm; ϵ 160,000 $cm^{-1} M^{-1}$

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!