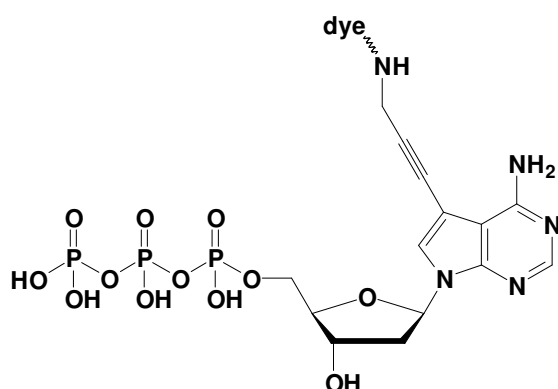


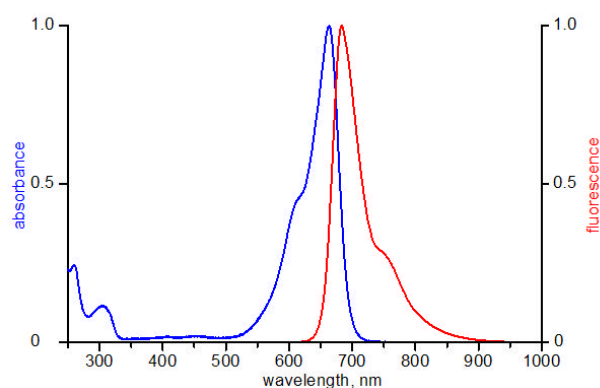
## 7-Propargylamino-7-deaza-dATP - ATTO 655

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with ATTO 655, Triethylammonium salt

Cat. No.	Amount
NU-1611-655	10 $\mu$ l / 1 mM



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



excitation and emission spectrum of ATTO 655

**Cat. No.:** NU-1611-655

**Molecular Formula:**  $C_{14}H_{20}N_5O_{12}P_3$  - ATTO 655 (free acid)

**Molecular Weight:** 1052.26 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  663 nm;  $\lambda_{em}$  684 nm;  $\epsilon$  125,000  $cm^{-1} M^{-1}$

**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!**