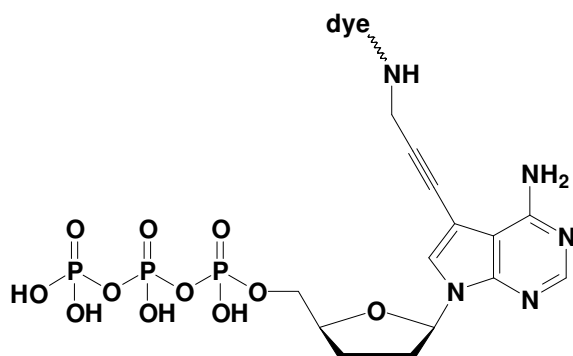


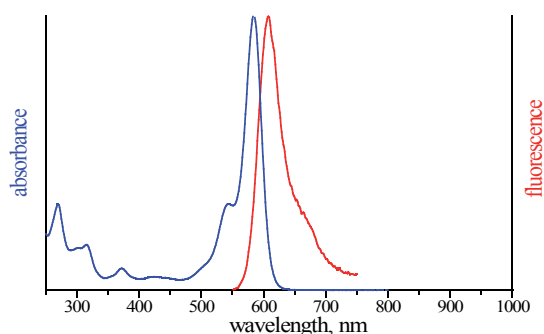
## 7-Propargylamino-7-deaza-ddATP - ATTO Rho101

7-Deaza-7-propargylamino-2'/3'-dideoxy-adenosine-5'-triphosphate, labeled with ATTO Rho101, Triethylammonium salt

Cat. No.	Amount
NU-1612-RHO101	80 $\mu$ l (0,5 mM)



structural formula of 7-Propargylamino-7-deaza-ddATP



excitation and emission spectrum of ATTO Rho101

**Cat. No.:** NU-1612-RHO101

**Molecular Formula:** C<sub>14</sub>H<sub>20</sub>N<sub>5</sub>O<sub>11</sub>P<sub>3</sub> - ATTO Rho101 (free acid)

**Molecular Weight:** 1098,26 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  586 nm;  $\lambda_{em}$  610 nm;  $\epsilon$  1,2 x 10<sup>5</sup> cm<sup>-1</sup> M<sup>-1</sup>

**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!**