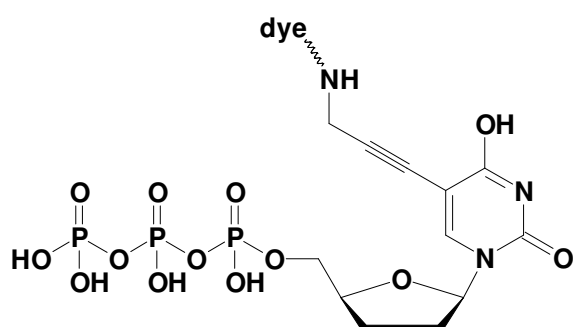


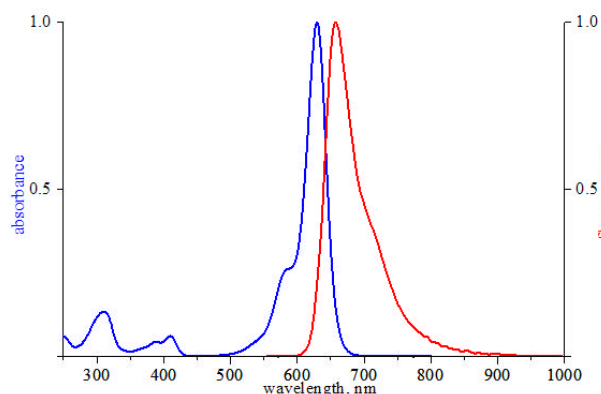
## 5-Propargylamino-ddUTP - ATTO 633

5-Propargylamino- 2', 3'-Dideoxy-uridine-5'-triphosphate, labeled with ATTO 633,  
 Triethylammonium salt

Cat. No.	Amount
NU-1619-633	80 $\mu$ l (0,5 mM)



structural formula of 5-Propargylamino-ddUTP



excitation and emission spectrum of ATTO 633

**Cat. No.:** NU-1619-633

**Molecular Formula:** C<sub>12</sub>H<sub>18</sub>N<sub>3</sub>O<sub>13</sub>P<sub>3</sub> - ATTO 633 (free acid)

**Molecular Weight:** 1037.20 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  629 nm;  $\lambda_{em}$  657 nm;  $\epsilon$  130,000 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!**