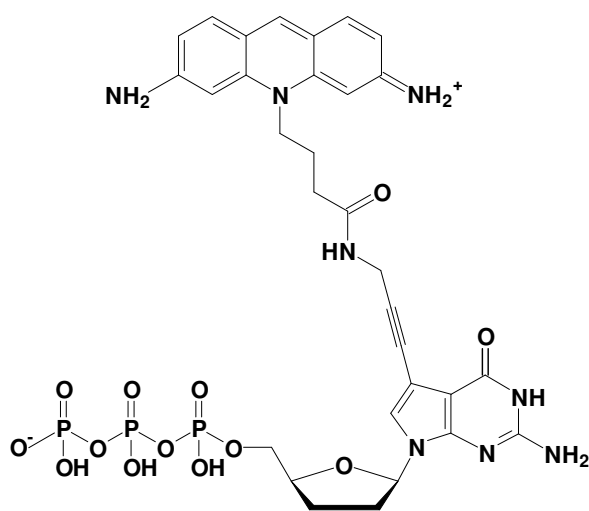


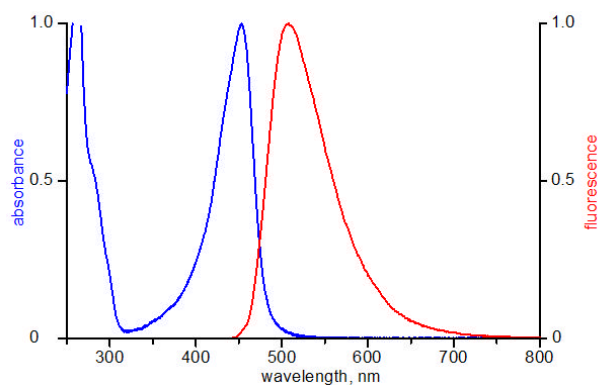
## 7-Propargylamino-7-deaza-ddGTP - ATTO 465

7-Deaza-7-propargylamino-2',3'-dideoxy-guanosine-5'-triphosphate, labeled with ATTO 465, Triethylammonium salt

Cat. No.	Amount
NU-1618-465	160 $\mu$ l (0,5 mM)



structural formula of 7-Propargylamino-7-deaza-ddGTP - ATTO 465



excitation and emission spectrum of ATTO 465

**Cat. No.:** NU-1618-465

**Molecular Formula:** C<sub>31</sub>H<sub>33</sub>N<sub>8</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 820.58 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  453 nm;  $\lambda_{em}$  508 nm;  $\epsilon$  75,000 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!**