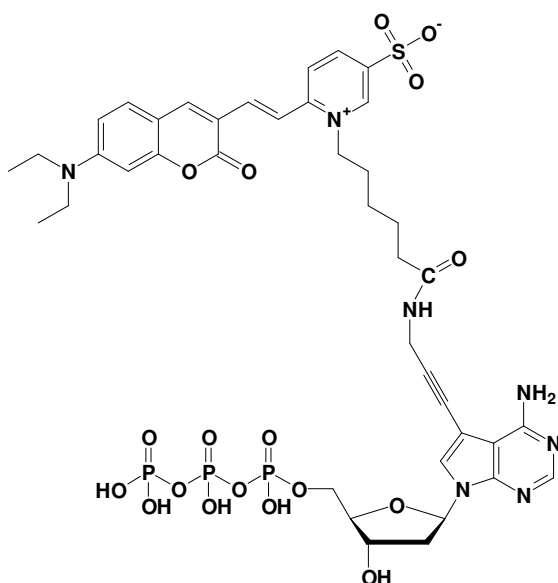


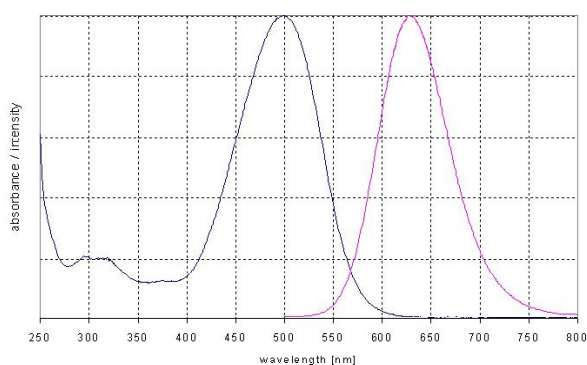
## 7-Propargylamino-7-deaza-dATP - DY 480XL

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with DY 480XL, Triethylammonium salt

Cat. No.	Amount
NU-1611-480XL	40 $\mu$ l (1 mM)



structural formula of 7-Propargylamino-7-deaza-dATP - DY 480XL



excitation and emission spectrum of DY 480XL

**Cat. No.:** NU-1611-480XL

**Molecular Formula:** C<sub>40</sub>H<sub>48</sub>N<sub>7</sub>O<sub>18</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1039,83 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  500 nm;  $\lambda_{em}$  630 nm;  $\epsilon$  50,000 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!**