



## 7-Propargylamino-7-deaza-dGTP-DY-485XL

7-Deaza-7-propargylamino-2'-deoxyguanosine-5'-triphosphate, labeled with DY 485XL, Triethylammonium salt

Cat. No.	Amount
NU-1615-485XL	40 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>39</sub>H<sub>48</sub>N<sub>7</sub>O<sub>19</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1043.82 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

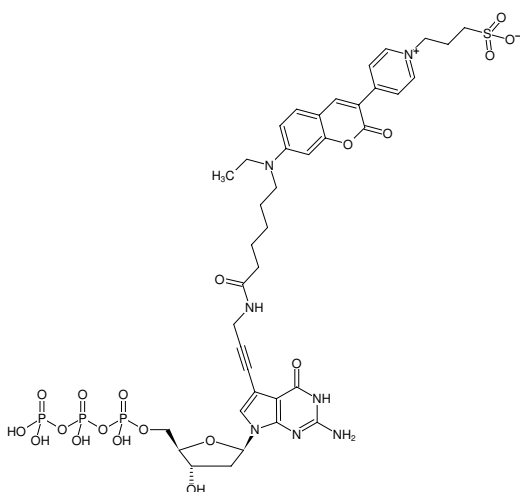
**Form:** solution in water

**Color:** orange

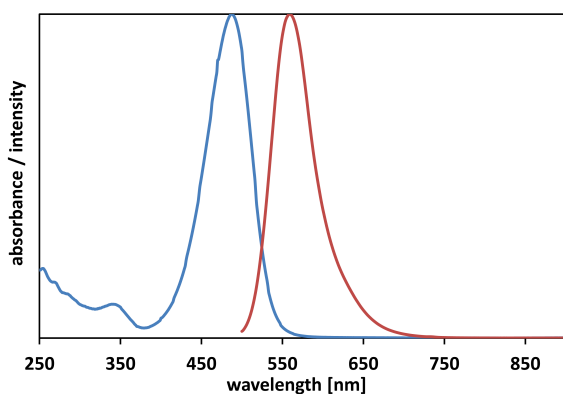
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 485 nm, λ<sub>em</sub> 560 nm, ε 50.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-dGTP-DY-485XL



excitation and emission spectrum of DY 485XL