



### 5-Propargylamino-ddUTP - DY-485XL

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with DY 485XL, Triethylammonium salt

Cat. No.	Amount
NU-1619-485XL	80 µl (0,5 mM)

**For research use only!**

**Shipping:** shipped on blue ice

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

Short term exposure (up to 1 week cumulative) to ambient temperature possible. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**Shelf Life:** 12 months

**Molecular Formula:** C<sub>37</sub>H<sub>46</sub>N<sub>5</sub>O<sub>19</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 989.77 g/mol (free acid)

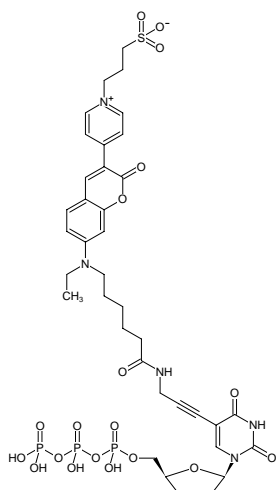
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution, pH 7.5 ±0.5

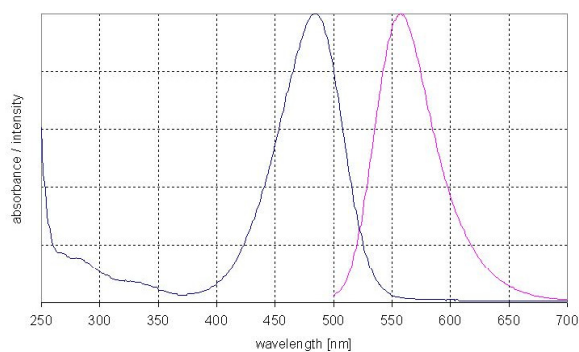
**Concentration:** 0.5 mM

**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 5-Propargylamino-ddUTP - DY-485XL



excitation and emission spectrum of DY 485XL