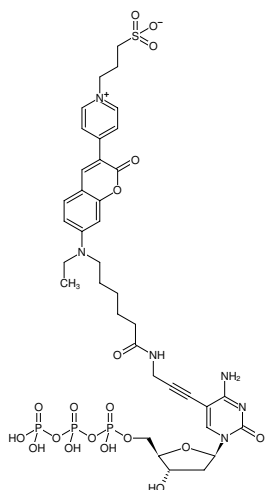




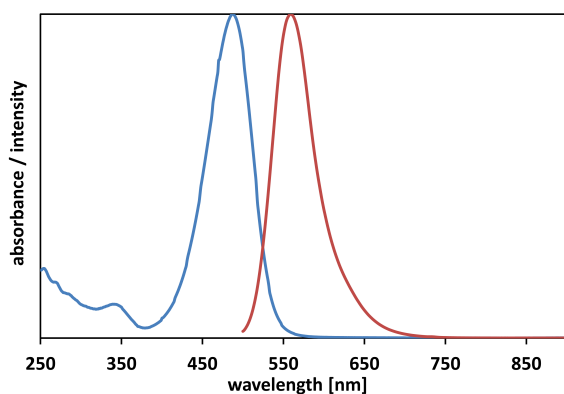
## 5-Propargylamino-dCTP-DY-485XL

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

| Cat. No.     | Amount       |
|--------------|--------------|
| NU-809-485XL | 40 µl (1 mM) |



Structural formula of 5-Propargylamino-dCTP-DY-485XL



excitation and emission spectrum of DY 485XL

**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>37</sub>H<sub>47</sub>N<sub>6</sub>O<sub>19</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1004.79 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)