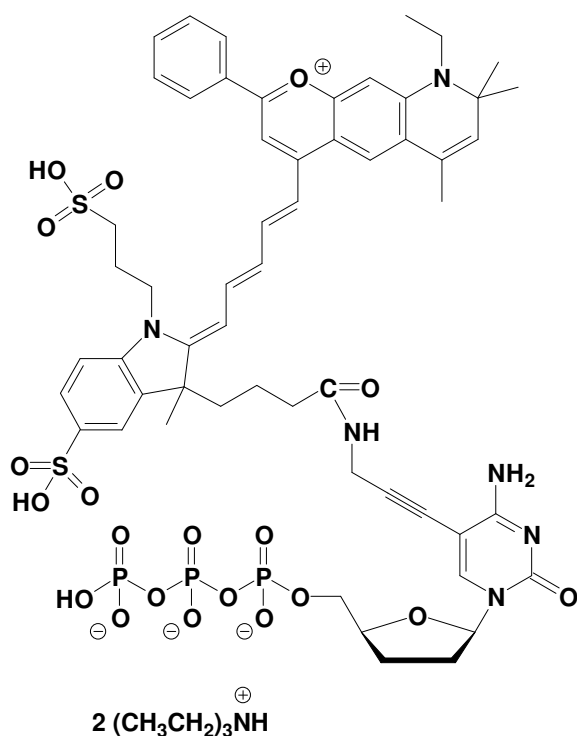


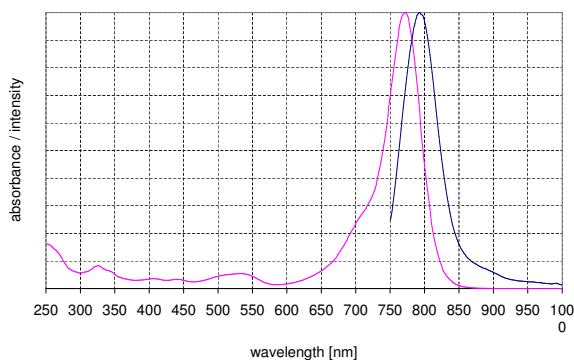
## 5-Propargylamino-ddCTP - DY 776

5-Propargylamino-2',3'-Dideoxy-cytidine-5'-triphosphate, labeled with DY 776,  
Triethylammonium salt

Cat. No.	Amount
NU-850-776	10 $\mu$ l / 1 mM



structural formula of 5-Propargylamino-ddCTP - DY 776



excitation and emission spectrum of DY 776

**Cat. No.:** NU-850-776

**Molecular Formula:**  $\text{C}_{56}\text{H}_{63}\text{N}_6\text{O}_{20}\text{P}_3\text{S}_2$  (Anion)

**Molecular Weight:** 1297.19 (Anion)

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

**Spectroscopic properties:**

$\lambda_{\text{exc}}$  771 nm;  $\lambda_{\text{em}}$  801 nm;  $\epsilon$  120,000  $\text{cm}^{-1} \text{M}^{-1}$

**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!**