

5-Propargylamino-ddUTP - DY 776

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

Cat. No.	Amount
NU-1619-776	80 μ l (0,5 mM)

Cat. No.: NU-1619-776

Molecular Formula: C₅₆H₆₄N₅O₂₁P₃S₂ (free acid)

Molecular Weight: 1300.18 (free acid)

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

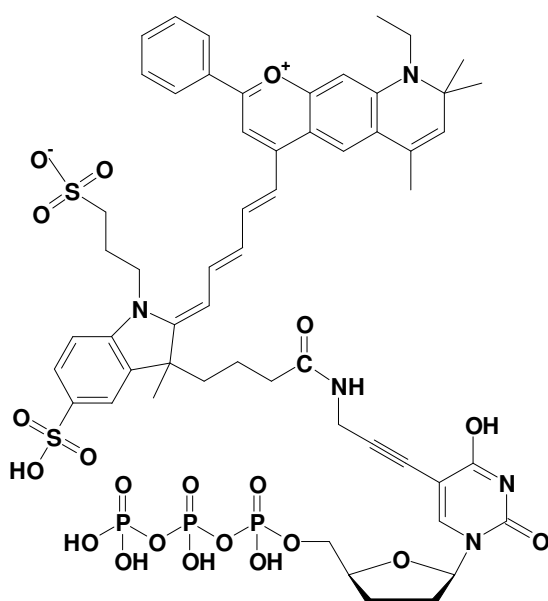
Spectroscopic properties:

λ_{exc} 771 nm; λ_{em} 801 nm; ϵ 120,000 cm⁻¹ M⁻¹

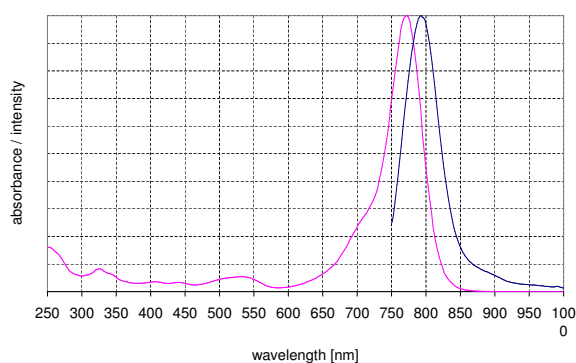
Storage conditions:

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!



structural formula of 5-Propargylamino-ddUTP - DY 776



excitation and emission spectrum of DY 776