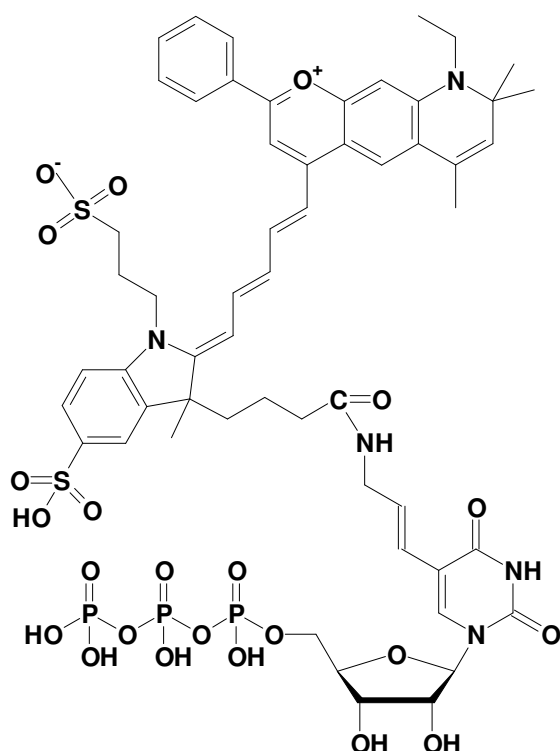


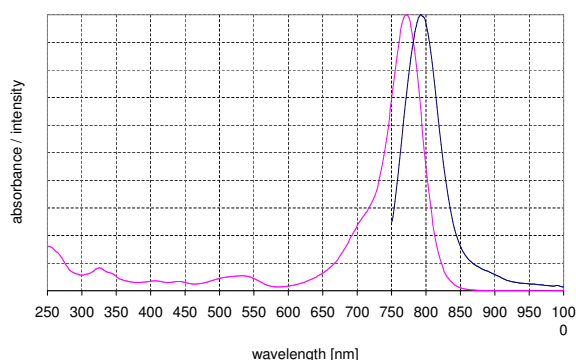
Aminoallyl-UTP - DY 776

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with DY 776, Triethylammonium salt

Cat. No.	Amount
NU-821-776	10 μ l / 1 mM



structural formula of Aminoallyl-UTP - DY 776



excitation and emission spectrum of DY 776

Cat. No.: NU-821-776

Molecular Formula: C₅₆H₆₆N₅O₂₃P₃S₂ (free acid)

Molecular Weight: 1334.20 (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!

Selected References:

Köhn *et al.* (2010) Near-infrared (NIR) dye-labeled RNAs identify binding of ZBP1 to the noncoding Y3-RNA. *RNA* **16**(7):1420.