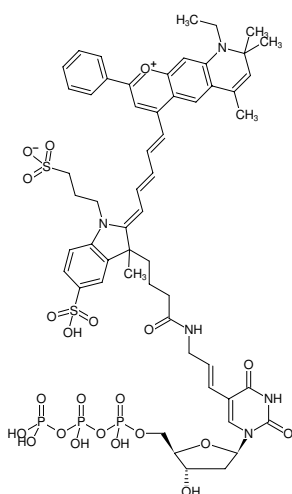




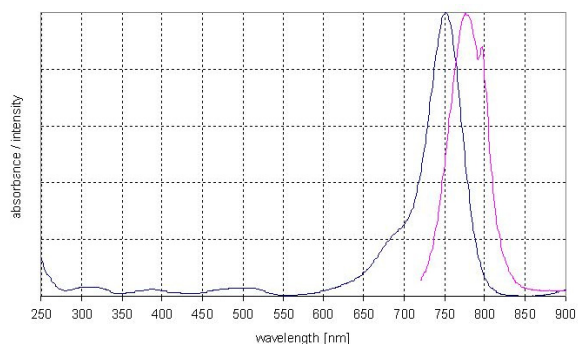
Aminoallyl-dUTP-DY-776

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with DY 776, Triethylammonium salt

Cat. No.	Amount
NU-803-776	80 µl (0.5 mM)



Structural formula of Aminoallyl-dUTP-DY-776



excitation and emission spectrum of DY 776

For general laboratory use.

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: C₅₆H₆₇N₅O₂₂P₃S₂ (free acid)

Molecular Weight: 1318.20 g/mol (free acid)

Exact Mass: 1317.29 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: dark blue

Concentration: 0.5 mM - 0.55 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 771 nm, λ_{em} 793 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)