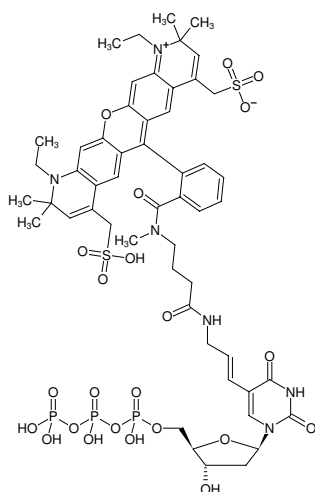




Aminoallyl-dUTP-ATTO-594

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO 594, Triethylammonium salt

Cat. No.	Amount
NU-803-594	40 µl (1 mM)



Structural formula of Aminoallyl-dUTP-ATTO-594

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: C₅₃H₆₅N₆O₂₃P₃S₂ (free acid)

Molecular Weight: 1311.16 g/mol (free acid)

Exact Mass: 1310.28 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

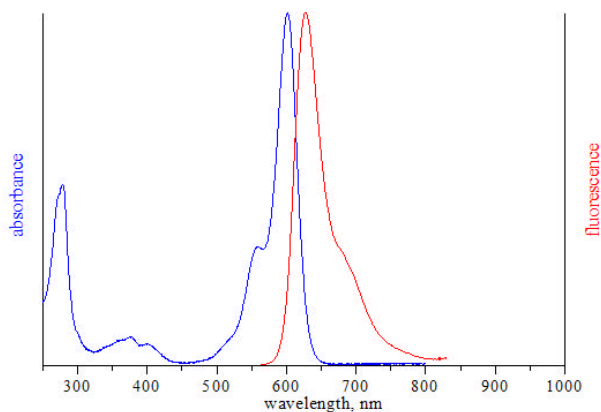
Form: solution in water

Color: red-violet

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 603 nm, λ_{em} 626 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 594