



Aminoallyl-UTP-5/6-TAMRA

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with 5/6-TAMRA, Triethylammonium salt

Cat. No.	Amount
NU-821-TAM	30 µl (1 mM)

For research use only!

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: C₃₇H₄₀N₅O₁₉P₃ (free acid)

Molecular Weight: 951.66 g/mol (free acid)

Exact Mass: 951.15 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

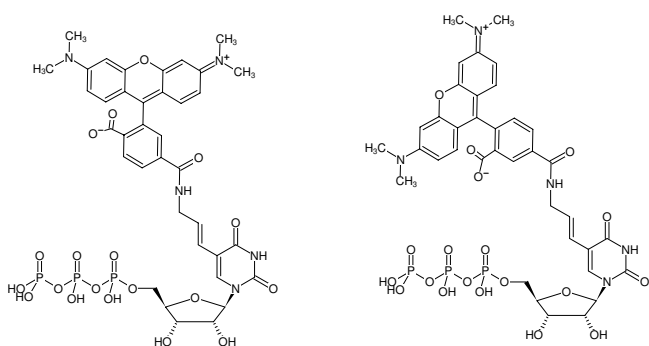
Form: solution in water

Color: pink to red

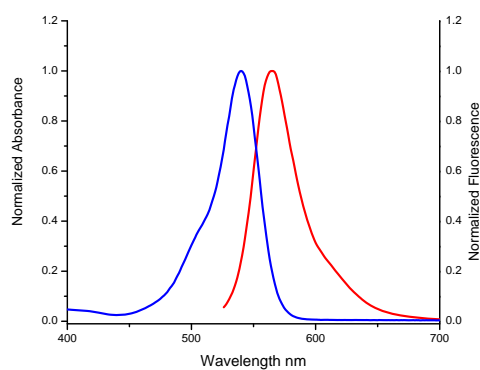
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 545 nm, λ_{em} 575 nm, ε 90.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of Aminoallyl-UTP-5/6-TAMRA



excitation and emission spectrum of 5/6-TAMRA