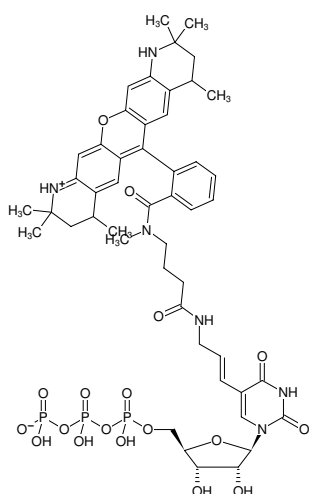




Aminoallyl-UTP-ATTO-550

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with ATTO 550, Triethylammonium salt

Cat. No.	Amount
NU-821-550	20 µl (5 mM)



Structural formula of Aminoallyl-UTP-ATTO-550

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: C₄₉H₆₁N₆O₁₈P₃ (free acid)

Molecular Weight: 1114.97 g/mol (free acid)

Exact Mass: 1114.33 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

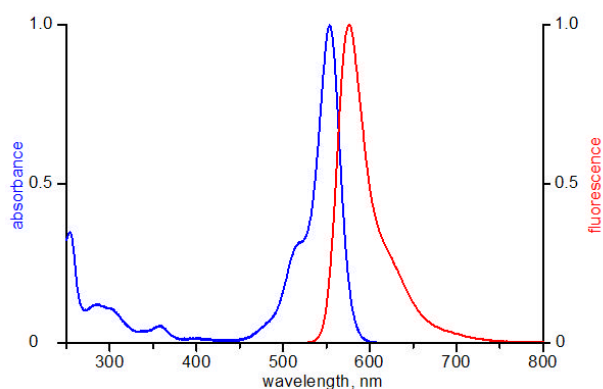
Form: filtered solution (30 kDa) in 10 mM Tris-HCl

Color: red-violet

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 554 nm, λ_{em} 576 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 550