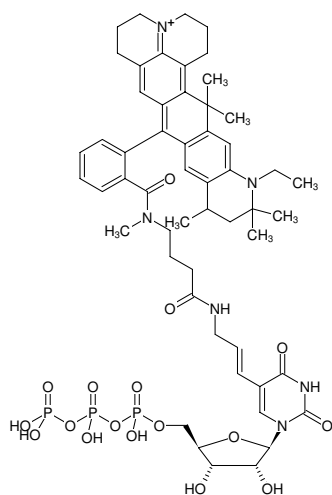




Aminoallyl-UTP-ATTO-647N

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with ATTO 647N, Triethylammonium salt

Cat. No.	Amount
NU-821-647N	10 µl (1 mM)



Structural formula of Aminoallyl-UTP-ATTO-647N

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: C₅₄H₆₉N₆O₁₇P₃ (free acid)

Molecular Weight: 1167.09 g/mol (free acid)

Exact Mass: 1166.39 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

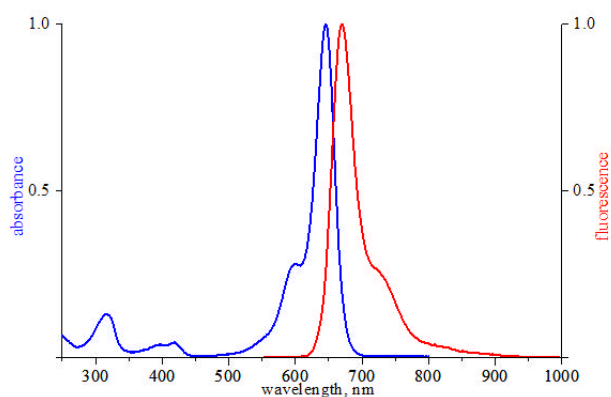
Form: solution in water

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 646 nm, λ_{em} 664 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 647N