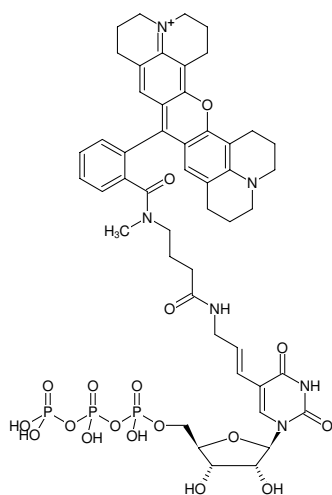




Aminoallyl-UTP-ATTO-Rho101

5-(3-Aminoallyl)-uridine-5'-triphosphate, labeled with ATTO Rho101, Triethylammonium salt

Cat. No.	Amount
NU-821-RHO101	40 μ l (1 mM)



Structural formula of Aminoallyl-UTP-ATTO-Rho101

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at $-20\text{ }^{\circ}\text{C}$

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: $\text{C}_{49}\text{H}_{57}\text{N}_6\text{O}_{18}\text{P}_3$ (free acid)

Molecular Weight: 1110.94 g/mol (free acid)

Exact Mass: 1110.29 g/mol (free acid)

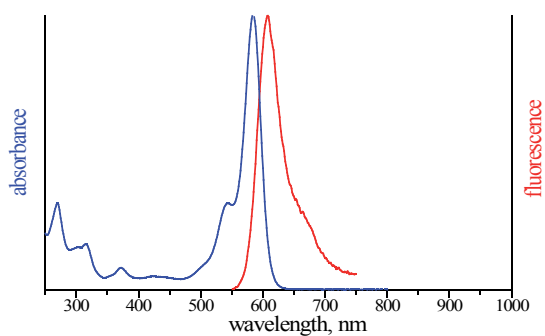
Purity: $\geq 95\%$ (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 587 nm, λ_{em} 609 nm, ϵ 120.0 L $\text{mmol}^{-1}\text{ cm}^{-1}$ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO Rho101