

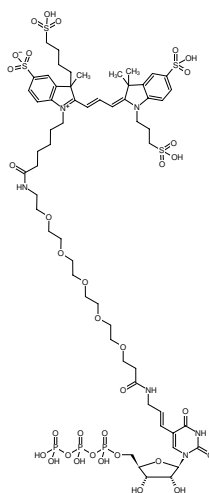


Aminoallyl-UTP-PEG₅-AF555

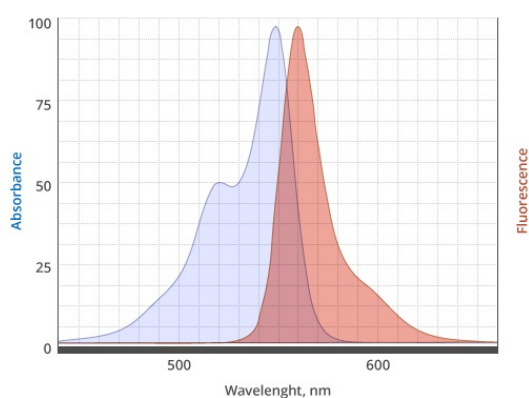
UTP-PEG₅-AF555

5-(3-Aminoallyl-PEG₅)-uridine-5'-triphosphate, labeled with AF555, Triethylammonium salt

Cat. No.	Amount
NU-821-PEG5-AF555	10 µl (5 mM)



Structural formula of Aminoallyl-UTP-PEG₅-AF555



excitation and emission spectrum of AF555

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: C₆₀H₈₉N₆O₃₄P₃S₄

Molecular Weight: 1659.54 g/mol (free acid)

Exact Mass: 1658.35 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: red

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 555 nm, λ_{em} 572 nm, ε 155.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

RNA Labeling by in vitro Transcription with T7 RNA Polymerase (in-house tested)

Related Products:

HighYield T7 RNA Synthesis Kit, #RNT-101

HighYield T7 AF555 RNA Labeling Kit, #RNT-101-AF555