

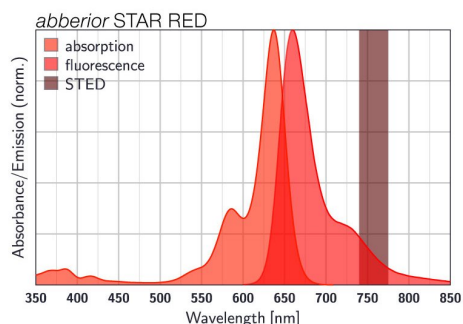
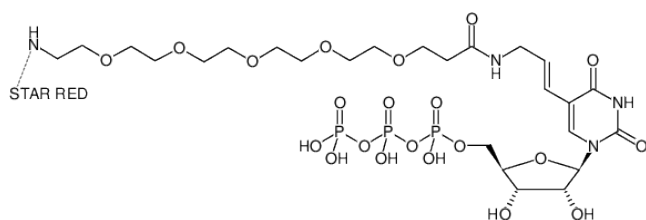


Aminoallyl-UTP-PEG₅-STAR RED

UTP-PEG₅-STAR RED

5-(3-Aminoallyl-PEG₅)-uridine-5'-triphosphate, labeled with abberior® STAR RED, Triethylammonium salt

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



excitation and emission spectrum of STAR RED

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: C₆₉H₈₈F₄N₇O₃₁P₃S₂

Molecular Weight: 1744.51 g/mol (free acid)

Exact Mass: 1743.41 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: dark blue

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 638 nm, λ_{em} 655 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

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