

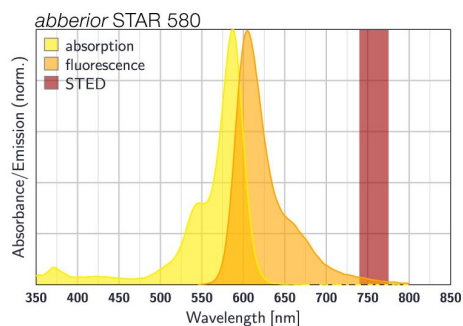
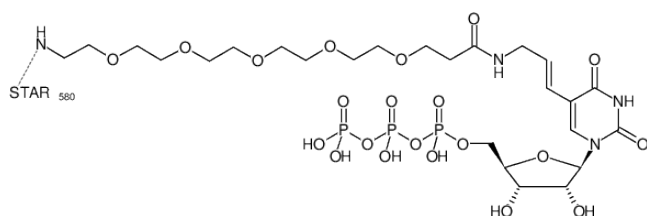


Aminoallyl-UTP-PEG₅-STAR 580

UTP-PEG₅-STAR 580

5-(3-Aminoallyl-PEG₅)-uridine-5'-triphosphate, labeled with abberior® STAR 580, Triethylammonium salt

Cat. No.	Amount
NU-821-PEG5-ST580	10 µl (5 mM)



excitation and emission spectrum of STAR 580

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: C₆₂H₈₂N₇O₂₄P₃

Molecular Weight: 1402.27 g/mol (free acid)

Exact Mass: 1401.47 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: purple

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 587 nm, λ_{em} 607 nm, ε 85.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

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