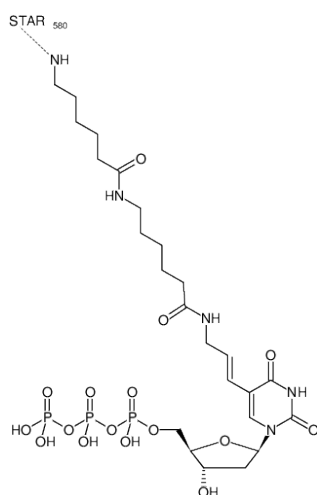




Aminoallyl-dUTP-XX-STAR 580

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with abberior® STAR 580, Triethylammonium salt

Cat. No.	Amount
NU-803-XX-ST580-S	10 µl (1 mM)
NU-803-XX-ST580-L	5 x 10 µl (1 mM)



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: C₆₁H₇₉N₈O₁₉P₃ (free acid)

Molecular Weight: 1321.25 g/mol (free acid)

Exact Mass: 1320.47 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

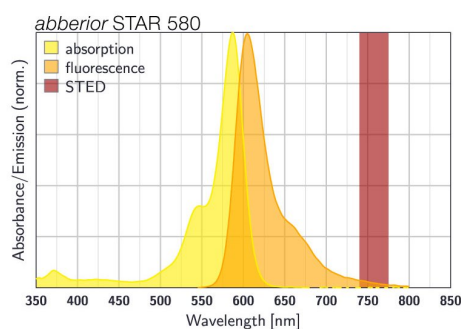
Form: solution in 10 mM Tris-HCl

Color: pink

Concentration: 1.0 mM - 1.1 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)



excitation and emission spectrum of STAR 580