



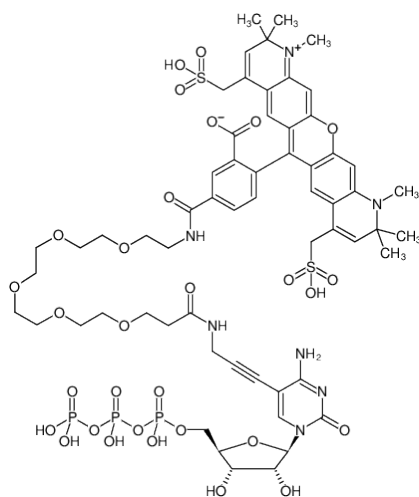
5-Propargylamino-CTP-PEG₅-AZDye594

structural analog to Alexa Fluor® 594

CTP-PEG₅-AZDye594

5-(PEG₅-Propargylamino)-cytidine-5'-triphosphate, labeled with AZDye 594, Triethylammonium salt

Cat. No.	Amount
NU-831-PEG5-AZ594	10 µl (5 mM)



Structural formula of 5-Propargylamino-CTP-PEG₅-AZDye594

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: C₆₀H₇₆N₇O₃₀P₃S₂ (free acid)

Molecular Weight: 1532.33 g/mol (free acid)

Exact Mass: 1531.33 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: red-violet

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ± 0.5

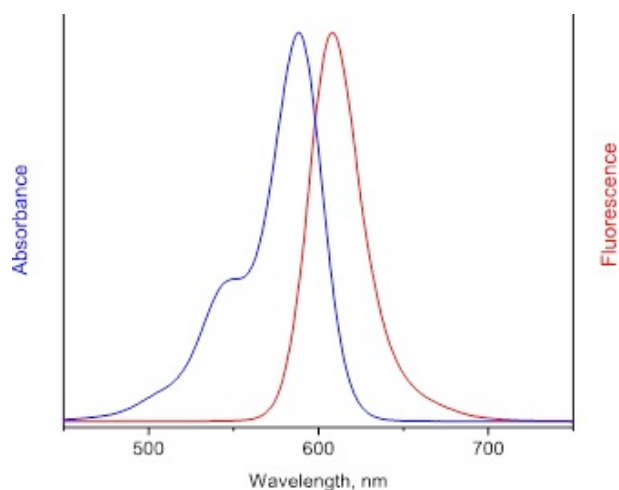
Spectroscopic Properties: λ_{exc} 590 nm, λ_{em} 617 nm, ε 92.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

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

Related Products:

HighYield T7 RNA Synthesis Kit, #RNT-101



excitation and emission spectrum of AZDye 594