



7-Propargylamino-7-deaza-ATP-PEG₅-AZDye594

structural analog to Alexa Fluor® 594

7-Deaza-7-propargylamino-PEG₅-adenosine-5'-triphosphate, labeled with AZDye 594, Triethylammonium salt

Cat. No.	Amount
NU-957-PEG5-AZ594	10 µl (5 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

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: C₆₂H₇₇N₈O₂₉P₃S₂ (free acid)

Molecular Weight: 1555.37 g/mol (free acid)

Exact Mass: 1554.35 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

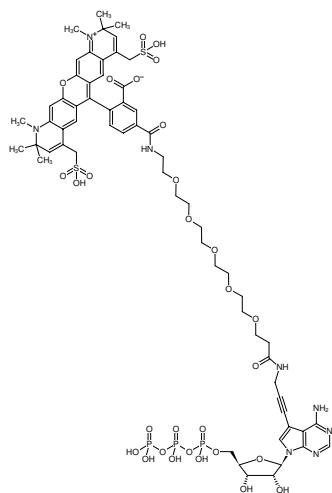
Form: solution in water

Color: red-violet

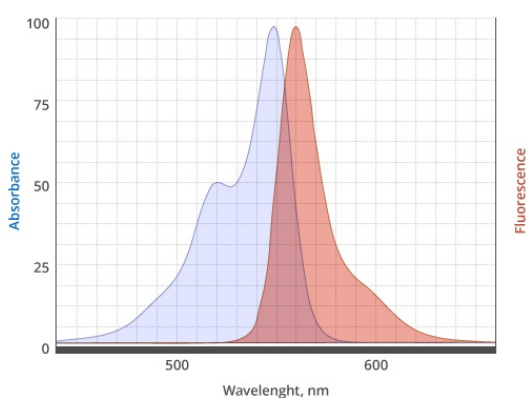
Concentration: 5.0 mM - 5.5 mM

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)



Structural formula of 7-Propargylamino-7-deaza-ATP-PEG₅-AZDye594



excitation and emission spectrum of AZDye 594