



7-Propargylamino-7-deaza-ATP-PEG₅-AZDye555

structural analog to Alexa Fluor® 555

7-Deaza-7-propargylamino-PEG₅-adenosine-5'-triphosphate, labeled with AZDye 555, Triethylammonium salt

| Cat. No. | Amount |
|-------------------|--------------|
| NU-957-PEG5-AZ555 | 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: C₆₂H₈₉N₈O₃₂P₃S₄ (free acid)

Molecular Weight: 1679.58 g/mol (free acid)

Exact Mass: 1678.37 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

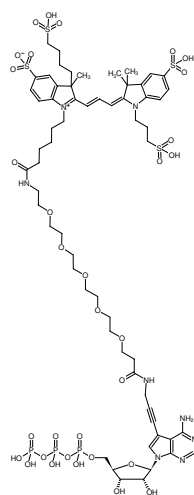
Form: solution in water

Color: red

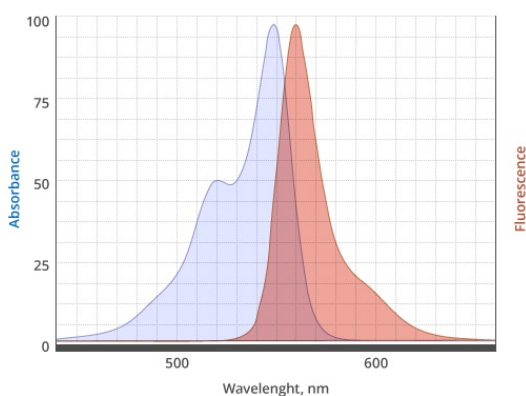
Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 555 nm, λ_{em} 572 nm, ε 155.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-ATP-PEG₅-AZDye555



excitation and emission spectrum of AZDye 555