

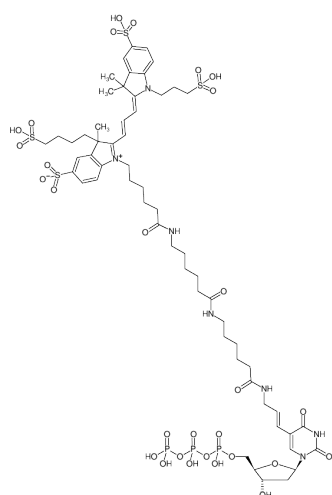


Aminoallyl-dUTP-XX-AZDye555

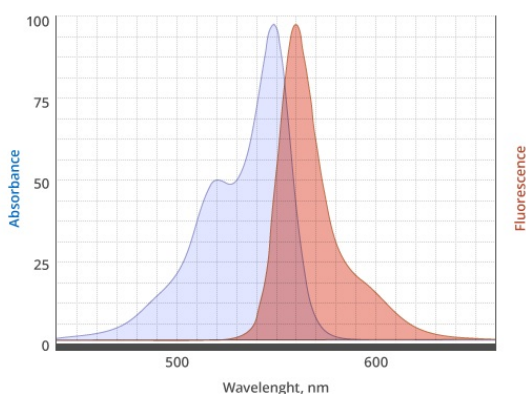
structural analog to Alexa Fluor® 555

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with AZDye 555, Triethylammonium salt

Cat. No.	Amount
NU-803-XX-AZ555-S	10 µl (1 mM)
NU-803-XX-AZ555-L	5 x 10 µl (1 mM)



Structural formula of Aminoallyl-dUTP-XX-AZDye555



excitation and emission spectrum of AZDye 555

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: C₅₉H₈₆N₇O₂₉P₃S₄ (free acid)

Molecular Weight: 1578.52 g/mol (free acid)

Exact Mass: 1577.36 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: red

Concentration: 1.0 mM - 1.1 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)

Applications:

Incorporation into DNA/cDNA by

- PCR with *Taq* polymerase in-house data

- Nick Translation with DNase I/ DNA Polymerase I in-house data

Description:

Aminoallyl-dUTP-XX-AZDye555 is recommended for direct enzymatic labeling of DNA/cDNA e.g. by PCR and Nick Translation. It is incorporated as substitute for its natural counterpart dTTP. The resulting Dye-labeled DNA/cDNA probes are ideally suited for fluorescence hybridization applications such as FISH or microarray-based gene expression profiling. Optimal substrate properties and thus labeling efficiency is ensured by an optimized linker attached to the C5 position of uridine. AZDye 555 (structural analog to Alexa Fluor 555®) is a hydrophilic dye with excellent photostability compared to fluorescein.

Recommended Aminoallyl-dUTP-XX-AZDye555/dTTP ratio for PCR and Nick Translation: 30-50% Aminoallyl-dUTP-XX-AZDye555/ 70-50% dTTP

Please note: Protect the Dye-labeled dUTP from exposure to light and carry out experimental procedures in low light conditions. The optimal final concentration of the Dye-labeled dUTP may vary depending on the application and assay conditions. For optimal



Aminoallyl-dUTP-XX-AZDye555

structural analog to Alexa Fluor® 555

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with AZDye 555, Triethylammonium salt

product yields and high incorporation rates an individual optimization of the Dye-labeled-dUTP/dTTP ratio is recommended.

Related Products:

HighFidelity AZDye555 PCR Labeling Kit, #APP-101-AZ555

HighFidelity YELLOW PCR Labeling Kit, #APP-101-YELLOW

AZDye555 NT Labeling Kit, #PP-305-AZ555