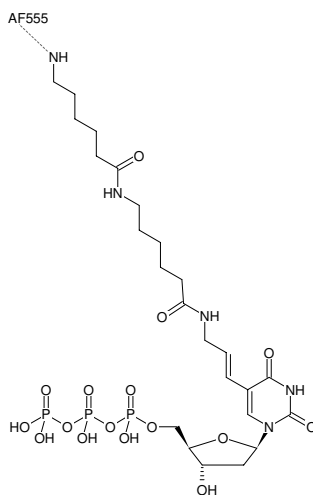




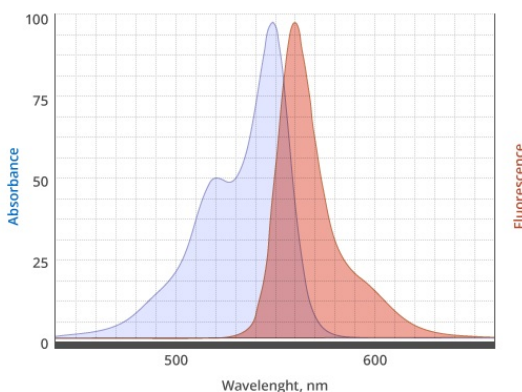
Aminoallyl-dUTP-XX-AF555

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with AF555, Triethylammonium salt

Cat. No.	Amount
NU-803-XX-AF555-S	10 µl (1 mM)
NU-803-XX-AF555-L	5 x 10 µl (1 mM)



Structural formula of Aminoallyl-dUTP-XX-AF555



excitation and emission spectrum of AF555

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: C₂₄H₄₁N₅O₁₆P₃ - AF555(free acid)

Molecular Weight: 1577.50 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: filtered solution (30 kDa) in 10 mM Tris-HCl

Color: pink

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-AF555 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. AF555 (structural analog to Alexa Fluor 555®) is a hydrophilic dye with excellent photostability compared to fluorescein.

Recommended Aminoallyl-dUTP-XX-AF555/dTTP ratio for PCR and Nick Translation: 30-50% Aminoallyl-dUTP-XX-AF555/ 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 product yields and high incorporation rates an individual optimization of the Dye-labeled-dUTP/dTTP ratio is recommended.



Aminoallyl-dUTP-XX-AF555

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with AF555, Triethylammonium salt

Related Products:

HighFidelity AF555 PCR Labeling Kit, #APP-101-AF555

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

AF555 NT Labeling Kit, #PP-305-AF555