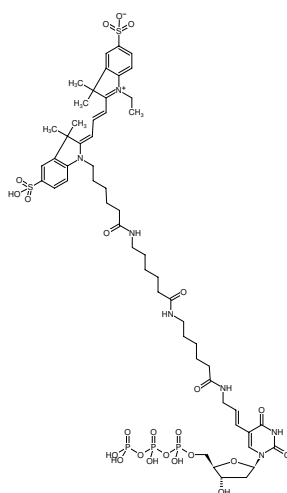




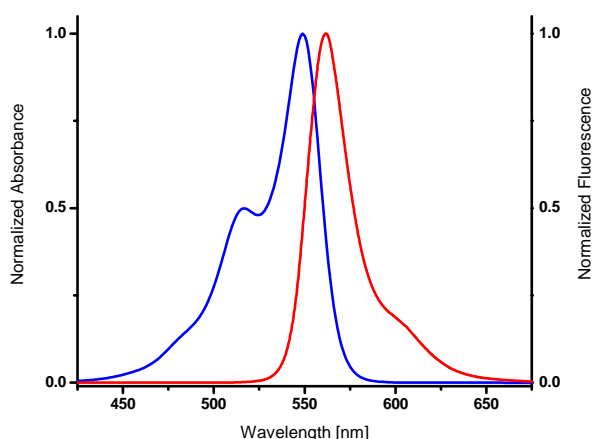
Aminoallyl-dUTP-XX-Cy3

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-803-XX-CY3-S	10 µl (1 mM)
NU-803-XX-CY3-L	5 x 10 µl (1 mM)



Structural formula of Aminoallyl-dUTP-XX-Cy3



Excitation and Emission spectrum of Cy3

For research use only!

Shipping: shipped on blue ice

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: C₅₅H₇₈N₇O₂₃P₃S₂ (free acid)

Molecular Weight: 1362.29 g/mol (free acid)

Exact Mass: 1361.38 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: sterile solution in 10 mM Tris-HCl

Color: pink

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 550 nm, λ_{em} 570 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), CF₂₆₀ 0.08

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-Cy3 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. **Higher incorporation and thus labeling rates can be achieved compared to Aminoallyl-dUTP-Cy3 (NU-803-Cy3).**

Recommended Aminoallyl-dUTP-XX-Cy3/dTTP ratio for PCR and Nick Translation: 30-50% Aminoallyl-dUTP-XX-Cy3/ 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.

Related Products:



Aminoallyl-dUTP-XX-Cy3

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with Cy3, Triethylammonium salt

HighFidelity Cy3 PCR Labeling Kit, #APP-101-Cy3
HighFidelity YELLOW PCR Labeling Testkit, #APP-101-YELLOW
Cy3 NT Labeling Kit, #PP-305-Cy3