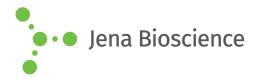
DATA SHEET

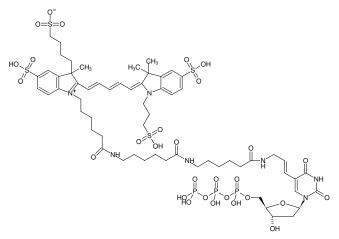




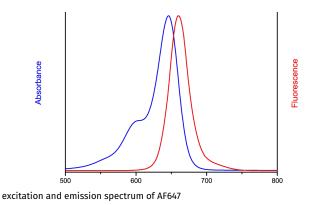
Aminoallyl-dUTP-XX-AF647

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with AF647, Triethylammonium salt

| Cat. No. | Amount |
|-------------------|------------------|
| NU-803-XX-AF647-S | 10 μl (1 mM) |
| NU-803-XX-AF647-L | 5 x 10 μl (1 mM) |



Structural formula of Aminoallyl-dUTP-XX-AF647



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: C₆₁H₈₈N₇O₂₉P₃S₂ (free acid)

Molecular Weight: 1604.56 g/mol (free acid)

Exact Mass: 1603.37 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

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

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 648 nm, λ_{em} 671 nm, ϵ 270.0 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-AF647 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. AF647 (structural analog to Alexa Fluor 647®) is a hydrophilic dye with excellent photostability compared to fluorescein.

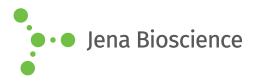
Recommended Aminoallyl-dUTP-XX-AF647/dTTP ratio for PCR and Nick Translation: 30-50% Aminoallyl-dUTP-XX-AF647/ 70-50% dTTP

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 very 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.



Jena Bioscience GmbH Löbstedter Str. 71 | 07749 Jena, Germany | Tel.:+49-3641-6285 000 https://www.jenabioscience.com







Aminoallyl-dUTP-XX-AF647

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with AF647, Triethylammonium salt

Related Products:

HighFidelity AF647 PCR Labeling Kit, #APP-101-AF647 HighFidelity RED PCR Labeling Testkit, #APP-101-RED AF647 NT Labeling Kit, #PP-305-AF647

