



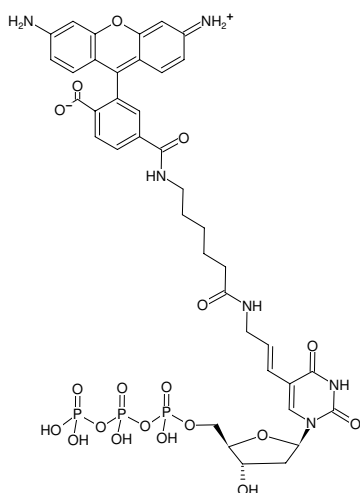
Rhodamine-12-dUTP

Rhodamine-X-(5-aminoallyl)-dUTP

Rho-green-X-dUTP, RhG-dUTP, R110-dUTP

5/6-Rhodamine-X-(5-aminoallyl)-2'-deoxyuridine-5'-triphosphate, Triethylammonium salt

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



Structural formula of Rhodamine-12-dUTP

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: C₃₉H₄₃N₆O₁₉P₃ (free acid)

Molecular Weight: 992.71 g/mol (free acid)

Exact Mass: 992.18 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: orange

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 505 nm, λ_{em} 530 nm, ε 85.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), CF_{280 nm} 0.27, CF_{260 nm} 0.39

Applications:

Incorporation into DNA/cDNA by

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

Description:

Rhodamine-12-dUTP is recommended for direct enzymatic labeling of DNA/cDNA by 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.

Recommended Rhodamine-12-dUTP/dTTP ratio for Nick Translation: 30-50% Rhodamine-12-dUTP/ 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.