



5-Propargylamino-dCTP-Texas Red

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with Texas Red, Triethylammonium salt

Cat. No.	Amount
NU-809-TXR-S	10 µl (1 mM)
NU-809-TXR-L	5 x 10 µl (1 mM)

Applications:

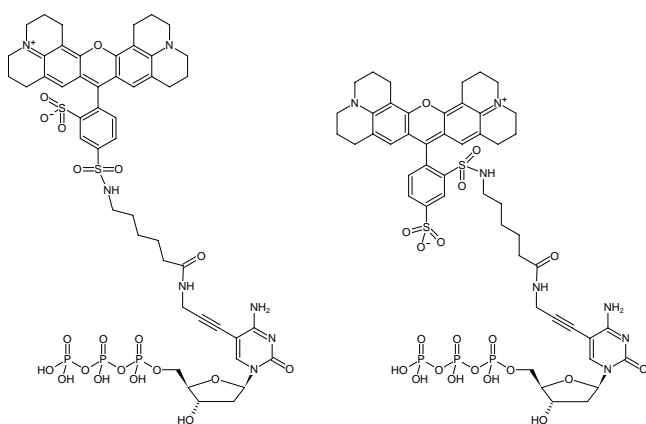
- Incorporation into DNA/cDNA by
- PCR with *Taq* polymerase in-house data
 - Nick Translation with DNase I/ DNA Polymerase I in-house data

Description:

5-Propargylamino-dCTP-Texas Red 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 dCTP. 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 cytidine.

Recommended Propargylamino-dCTP-TexasRed/dCTP ratio for PCR and Nick Translation: 20-30% Propargylamino-dCTP-TexasRed/70-70% dCTP (PCR), 30-50% Propargylamino-dCTP-TexasRed/70-50% dCTP (Nick Translation)

Please note: Protect the Dye-labeled dCTP from exposure to light and carry out experimental procedures in low light conditions. The optimal final concentration of the Dye-labeled dCTP may vary depending on the application and assay conditions. For optimal product yields and high incorporation rates an individual optimization of the Dye-labeled-dCTP/dCTP ratio is recommended.



Structural formula of 5-Propargylamino-dCTP-Texas Red

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: C₄₉H₅₈N₇O₂₀P₃S₂ (free acid)

Molecular Weight: 1222.07 g/mol (free acid)

Exact Mass: 1221.24 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: red-violet

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 588 nm, λ_{em} 609 nm, ε 80.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)