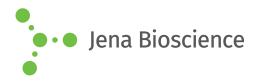
DATA SHEET

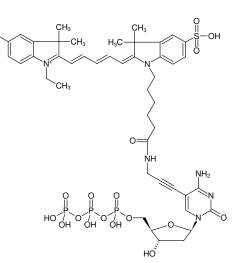




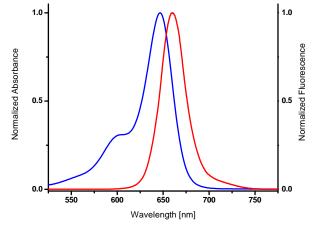
5-Propargylamino-dCTP-Cy5

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-809-CY5-S	10 μl (1 mM)
NU-809-CY5-L	5 x 10 μl (1 mM)



Structural formula of 5-Propargylamino-dCTP-Cy5



excitation and emission spectrum of Cy5

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: C₄₅H₅₇N₆O₂₀P₃S₂ (free acid)

Molecular Weight: 1159.01 g/mol (free acid)

Exact Mass: 1158.23 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 649 nm, λ_{em} 670 nm, ϵ 250.0 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:

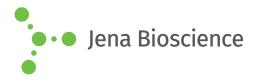
5-Propargylamino-dCTP-Cy5 is recommended for direct enzymatic labeling of DNA/cDNA e.g. by PCR and Nick Translation. It is incorporated as substitute for its natural counterpart dCTP. The resulting Dye-labeled DNA/cDNA probes are ideally suited for fluorescence hybridization applications such as FISH 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-Cy5/dCTP ratio for PCR and Nick Translation: 30-50% Propargylamino-dCTP-Cy5/ 70-50% dCTP

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 very 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.









5-Propargylamino-dCTP-Cy5

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Selected References:

Ramsay *et al.* (2010) CyDNA: Synthesis and Replication of Highly Cy-Dye Substituted DNA by an Evolved Polymerase. *J. Am. Chem. Soc.* **132 (14)**:5096.

