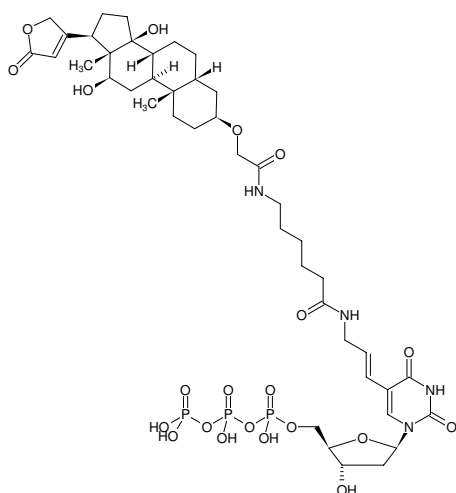


**DIG-11-dUTP**

alkali-stable

Digoxigenin-X-(5-aminoallyl)-2'-deoxyuridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-803-DIGXS	25 µl (1 mM)
NU-803-DIGXL	5 x 25 µl (1 mM)



Structural formula of DIG-11-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:** C₄₃H₆₅N₄O₂₁P₃ (free acid)**Molecular Weight:** 1066.91 g/mol (free acid)**Exact Mass:** 1066.34 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** sterile solution in 10 mM Tris-HCl**Color:** colorless to slightly yellow**Concentration:** 1.0 mM - 1.1 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 290 nm, ε 8.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Applications:**

Incorporation into DNA/cDNA by

- PCR with *Taq* polymerase [1,2] & in-house data
- Nick Translation with DNase I/ DNA Polymerase I [3] and in-house data
- Primer Extension with Klenow *exo*⁻ [4]
- 3'-End Labeling with Terminal deoxynucleotidyl Transferase (TdT) [5]
- Reverse Transcription with MMLV Reverse Transcriptase [6]

Incorporation into RNA by

- 3'-End Labeling with Terminal deoxynucleotidyl Transferase (TdT) [7]

Description:

Digoxigenin-11-dUTP is enzymatically incorporated into DNA/cDNA as substitute for its natural counterpart dTTP. The resulting Digoxigenin-labeled DNA/cDNA probes are subsequently detected using Digoxigenin-antibodies conjugated with horseradish peroxidase (HRP), alkaline phosphatase (AP) or a fluorescent dye. Optimal substrate properties and thus labeling efficiency is ensured by a 11-atom linker attached to the C5 position of uridine.

Recommended Digoxigenin-11-dUTP/dTTP ratio for PCR and Nick Translation: 35% Digoxigenin-11-dUTP/ 65% dTTP

Please note: The optimal final concentration of Digoxigenin-11-dUTP may very depending on the application and assay conditions. For optimal product yields and high incorporation rates an individual optimization of the Digoxigenin-11-dUTP/dTTP ratio is recommended.

Related Products:

HighFidelity Digoxigenin PCR Labeling Kit, #APP-101-DIGX
 Digoxigenin NT Labeling Kit, #PP-310-DIGX

Selected References:

- [1] Anderson *et al.* (2005) Incorporation of reporter-labeled nucleotides by DNA polymerases. *Biotechniques* **38**:257.
- [2] Jackson *et al.* (1991) Detection of Shiga Toxin-Producing *Shigella dysenteriae* Type 1 and *Escherichia coli* by Using Polymerase Chain Reaction with Incorporation of Digoxigenin-11-dUTP. *J Clin Microbiol.* **29** (9):1910.
- [3] Dauwerse *et al.* (1999) Two-colour FISH detection of the *inv* (16) in interphase nuclei of patients with acute myeloid leukemia. *Br J Haematol* **106**:111.
- [4] Wiegant *et al.* (2008) Probe Labeling and Fluorescence In Situ Hybridization. *Current Protocols in Cytometry*: Unit 8.3.
- [5] Schmitz *et al.* (2008) Nonradioactive labeling of oligonucleotides in vitro with the haptent digoxigenin by tailing with terminal transferase. *Anal Biochem* **199**:222.
- [6] Grimmond *et al.* (2001) Expression Profiling with cDNA Microarray's: A User's Perspective and Guide. In: *DNA Microarrays: Gene Expression*



DIG-11-dUTP

alkali-stable

Digoxigenin-X-(5-aminoallyl)-2'-deoxyuridine-5'-triphosphate, Triethylammonium salt

Applications (Jordan). Springer Verlag Berlin Heidelberg.

[7] Rosemeyer *et al.* (1995) Nonradioactive 3'-end-labeling of RNA molecules of different lengths by terminal deoxynucleotidyltransferase *Anal Biochem* **224**:446.