

Digoxigenin-dUTP, alkali-stable

Digoxigenin-labeled nucleotide suitable for
non-radioactive labeling of DNA probes

Digoxigenin-dUTP replaces dTTP in reactions in which it serves as a substrate for *E. coli* DNA polymerase (holoenzyme and Klenow fragment), T4 and Taq DNA polymerases, reverse transcriptase (from AMV and M-MuLV) and terminal transferase. Digoxigenin-dUTP can be used to produce digoxigenin-labeled DNA probes in a variety of labeling reactions including nick translation, random prime labeling, cDNA labeling and 3'-end labeling. The resulting digoxigenin-labeled probe can be used in a variety of hybridization applications including Southern blots, Northern blots, or dot blots. The probes can also be used for *in situ* hybridization procedures on fixed cells and tissues. Digoxigenin-labeled probes have been shown to hybridize to homologous nucleic acid at the same rate and to the same extent as non-labeled probes. The hybridized digoxigenin-labeled DNA probes can be detected by their interaction with antibodies coupled to fluorescent dyes or color-producing enzymes.

This labeled dUTP can be used with the Nick Translation DNA Labeling System 2.0 (Prod. No. ENZ-GEN111).

Citations: 3

[View Online »](#)

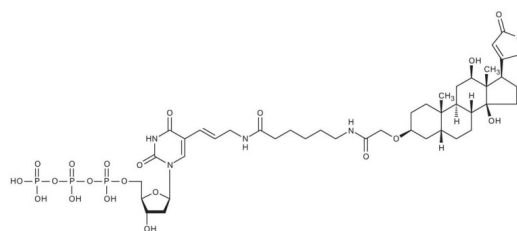
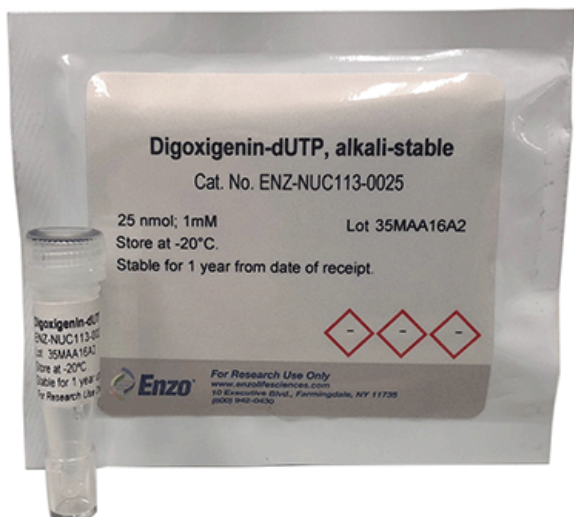
Ordering Information

[Order Online »](#)

ENZ-NUC113-0025	25nmol
-----------------	--------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Digoxigenin-11-dUTP

Appearance Clear, colorless liquid.

Concentration 1mM

Extinction Coefficient 22,600 M-1cm-1 (220nm, pH 7)

Formula $C_{43}H_{65}N_4O_{21}P_3$

Formulation Liquid. Solution in water.

MW 1066.92 (free acid)

Purity ≥93% (HPLC)

Quantity Sufficient for approximately 25 reactions, following the recommended protocol of Prod. No. ENZ-GEN111.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)