

Digoxigenin-UTP, alkali-stable

**Digoxigenin-labeled nucleotide suitable for
non-radioactive labeling of RNA probes**

Digoxigenin-UTP can replace UTP in reactions catalyzed by T3, T7 or SP6 RNA polymerases. The digoxigenin-labeled RNA transcripts produced by these reactions are suitable for a wide range of applications such as nucleic acid hybridization, sequencing, and genome analysis. The transcription reaction produces multiple RNA copies of the DNA template(s) during a short incubation period. RNA probes offer higher target specificity and greater sensitivity than the corresponding DNA-DNA hybrids. The single-stranded RNA probes offer selectivity unavailable with double-stranded DNA probes, because they are strand-specific. Furthermore, RNA probes hybridize much more efficiently to target molecules than DNA probes because there is no self-hybridization. The hybridized digoxigenin-labeled DNA probes can be detected by their interaction with antibodies coupled to fluorescent dyes or color-producing enzymes.

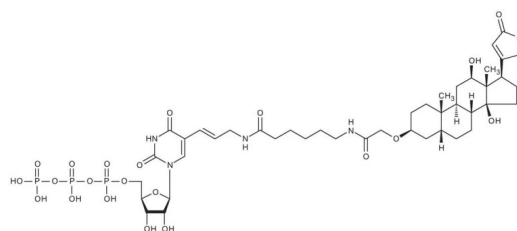
Ordering Information

[Order Online »](#)

| | |
|-----------------|---------|
| ENZ-NUC114-0250 | 250nmol |
|-----------------|---------|

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-UTP

Appearance Clear, colorless liquid.

Concentration 10mM

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

Formula $C_{43}H_{65}N_4O_{22}P_3$

Formulation Liquid. Solution in water.

MW 1082.92 (free acid)

Purity ≥93% (HPLC)

Quantity 25µl



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)