

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.

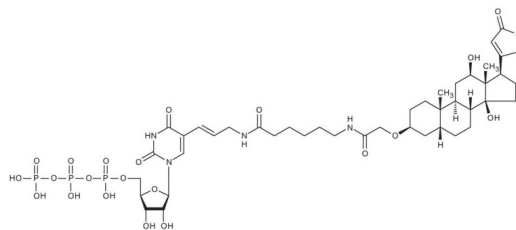
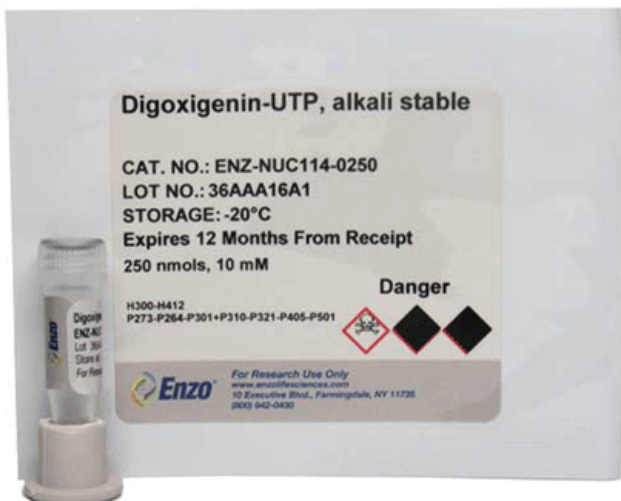
Ordering Information

[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 ⁻¹ cm ⁻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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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