

Digoxigenin NT Labeling Kit

Digoxigenin Nick Translation Labeling Kit contains all reagents (except template and materials for purification of the probe) required for nick translation-based Digoxigenin labeling of DNA providing a highly efficient, easy-to-perform and rapid labeling technology. The labeling principle is similar to the DIG-Nick Translation Mix (Roche). Nick translation labeling is based on the reverse activities of Polymerase I and DNase I. DNase I is able to introduce randomly distributed nicks to double stranded DNA at low enzyme concentrations. The 5'→3' exonuclease activity of Polymerase I removes nucleotides from the 3' side of the nick while synthesizing a partial new complementary strand using the 3'-OH termini as primer. In the presence of labeled dUTP Polymerase I incorporates labeled dUTP instead of dTTP. The DIGX NT labeling mix contains an optimized Digoxigenin-labeled dUTP for incorporation into DNA by nick translation using DNA Polymerase I. Digoxigenin-11-dUTP is enzymatically incorporated into DNA as substitute for its natural counterpart dTTP. Optimal substrate properties and thus labeling efficiency as well as an efficient detection of the Digoxigenin moiety is ensured by an 11-atom linker attached to the C5 position of uridine. The well balanced polymerase / nuclease activities of the enzyme mix ensure the generation of highly labeled double stranded DNA fragments. The resultant Biotin-labeled DNA is suitable for application in FISH and other nucleic acid hybridization assays.

Ordering Information

[Order Online »](#)

JBS-PP-310L-DIGX	75Reactions
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Quantity 20 µl each

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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-
be@enzolifesciences.com](mailto:info-be@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)