

DIG 3'-End Oligonucleotide Labeling Kit with Digoxigenin-11-ddUTP

DIG 3'-End Oligonucleotide Labeling Kit with Digoxigenin-11-ddUTP contains all reagents (except oligonucleotide template to be labeled and materials for Digoxigenin detection) required for efficient 3'-End digoxigenylation of DNA oligonucleotides (length: 20 -100 bp, 5 pmol per reaction). The labeling principle is based on Terminal deoxynucleotidyl Transferase (TdT) that template-independently transfers 1 Digoxigenin-11-ddUTP to the 3'-OH group of ssDNA (e.g. an oligonucleotide) in the presence of CoCl₂. It is similar to the principle of DIG Oligonucleotide 3'-End DNA Labeling Kit, 2nd Generation (Roche). The resulting 3'-End digoxigenylated oligonucleotides are ideally suited for applications involving sequence-specific protein binding or hybridization such as EMSA, Northern or Southern blots. Compared to internal, random digoxigenylated probes, Digoxigenin is located at the 3'-End only and less likely interferes with probe binding. TdT possesses a preference for single-stranded DNA (ssDNA) over dsDNA with 3'-overhangs or blunt ends. For the preparation of labeled dsDNA complexes, label each complementary oligonucleotide separately and anneal them before use.

Ordering Information

[Order Online »](#)

JBS-APP-002	25Reactions
-------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Notes



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)