

# BIO 3'-End Oligonucleotide Labeling Kit with Biotin-11-UTP

BIO 3'-End Oligonucleotide Labeling Kit with Biotin-11-UTP contains all reagents (except oligonucleotide template to be labeled and materials for Biotin detection) required for efficient 3'-End biotinylation 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 1-3 Biotin-11-UTPs to the 3'-OH group of ssDNA (e.g. an oligonucleotide) in the presence of CoCl<sub>2</sub>. It is similar to the principle of Biotin 3'-End DNA Labeling Kit (Pierce/Thermo Scientific). The resulting 3'-End biotinylated oligonucleotides are ideally suited for applications involving sequence-specific protein binding or hybridization such as EMSA, Northern or Southern blots. Compared to internal, random biotinylated probes, Biotin 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-001	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)