

# BIO-PROBE® 3'- Oligonucleotide labeling system reagent pack

The procedure for labeling of DNA probes with a polynucleotide "tail" containing hapten-labeled nucleotides was developed by Enzo. In such terminal labeling reactions, terminal transferase catalyzes the addition of nucleotides to any 3'-OH terminus in a template independent manner. This rapid and convenient nonradioactive labeling procedure is free of any sequence bias that is normally observed in random priming or nick translation reactions. Terminal labeling is an ideal procedure for the labeling of oligonucleotides.

An oligonucleotide can be synthesized using standard, commercially available reagents and labeled after synthesis in a simple and reproducible enzyme reaction. The BIO-PROBE® 3'-Oligonucleotide Labeling System provides maximum flexibility for labeling oligonucleotides approximately 15 to 100 nucleotides in length. A standard reaction will label up to 100 picomoles (1µg of a 30-mer oligonucleotide sequence). A complete system combines a standard Reagent Pack (buffers and enzyme) with a Nucleotide Pack. All components contain sufficient reagents for 25 labeling reactions.

The system is available in two modular formats:

3'-Oligonucleotide Tailing (with a "tail" of nucleotides)

3'-Oligonucleotide Labeling (with a single nucleotide)

The BIO-PROBE® 3'-Oligonucleotide Labeling System Reagent Pack contains all reagents required for labeling except nucleotides. For optimum results, the reagents contained in this Pack should be used with the appropriate Nucleotide Pack.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-42730

25Reactions

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year after receipt when stored as recommended.
Long Term Storage	-20°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

### Application

Dot Blot, ISH (in situ hybridization), Northern Blot, Southern Blot

### Contents

**5X Reaction Buffer**, 100 µl Potassium cacodylate buffer, pH 7, containing β-mercaptoethanol  
**10X CoCl<sub>2</sub> Solution**, 50 µl 10mM solution  
**Terminal Deoxynucleotide Transferase**, 50 µl 30 units/µl in storage buffer  
**Control Oligonucleotide (unlabeled)**, 25 µl 0.2 µg/µl in TE Buffer (30-mer, 5'-TTG GGT AAC GCC AGG GTT TTC CCA GTC ACG-3', homologous to the lac Z' region of pUC-type vector DNAs)  
**Control Target DNA**, 25 µl 0.2 µg/µl in TE Buffer (Restriction enzyme digest of a pUC-type vector containing an Epstein-Barr Virus DNA insert)

### Technical Info / Product Notes

BIO-PROBE® is a trademark of Enzo Life Sciences, Inc. Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.

Last modified: May 29, 2024



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)