

# BIO-PROBE<sup>®</sup> 3'- Oligonucleotide labeling kit with bio- 16-ddUTP

Terminal labeling is an ideal procedure for the labeling of oligonucleotides with haptens such as biotin or fluorescein. An oligonucleotide can be synthesized using standard, commercially available reagents and labeled after synthesis in a simple and reproducible enzyme reaction. The BIO-PROBE<sup>®</sup> 3'-Oligonucleotide Labeling Kit with Bio-16-ddUTP contains all reagents required for terminal labeling of oligonucleotides with biotin except the oligonucleotides to be labeled.

The BIO-PROBE<sup>®</sup> 3'-Oligonucleotide Labeling Kit provides maximum flexibility for labeling oligonucleotides approximately 15 to 100 nucleotides in length. A standard reaction will label up to 100 picomoles (1 µg of a 30-mer oligonucleotide sequence). A complete kit contains two component packs that are available separately: a Reagent Pack and a Bio-16-ddUTP Dideoxynucleotide Pack. Both packs contain sufficient reagents for 25 labeling reactions.

## Ordering Information

[Order Online »](#)

ENZ-42730-31	25Reactions
--------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

**Use/Stability** 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** BIO-PROBE® 3'-Oligonucleotide Labeling System Reagent Pack (Prod. No. ENZ-42730)  
Bio-16-ddUTP Deoxynucleotide Pack for 3'-Oligonucleotide Labeling (Prod. No. ENZ-42731)

**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-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)