

Fluorescein-12-ddUTP dideoxynucleotide pack for 3'- oligonucleotide

labeling

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 (1 µg 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 Dideoxynucleotide Pack contains a vial of modified dideoxynucleotide and a vial of control oligonucleotide labeled with the modified dideoxynucleotide. Terminal labeling reactions using dideoxynucleotides result in the addition of a single nucleotide at the 3'-OH terminus. For optimum results, the nucleotide contained in this Pack should be used with the 3'-Oligonucleotide Labeling System Reagent Pack.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-42733

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.
Handling	Avoid freeze/thaw cycles. Protect from light.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Contents

Fluorescein-12-ddUTP, 25µl 1mM Fluorescein-12-ddUTP, Prod. No. [ENZ-42833](#)
Labeled Control Oligonucleotide, 20µl 0.05µg/µl in TE Buffer (Fluorescein-12-ddUTP labeled Lac Z'- specific sequence identical to the unlabeled control oligonucleotide in the Reagent Pack, Prod. No. [ENZ-42730](#))

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.

For complete instructions for use, see Manual for Prod. No. ENZ-42730, ENZO BIO-PROBE[®] 3'-Oligonucleotide Labeling System Reagent Pack.

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)