



BioArray™
Terminal Labeling Kit
with Biotin-ddUTP
for DNA Probe Array Assays

25 labeling reactions

Cat. No. 42630

For Research Use Only

INTRODUCTION

The ENZO **BioArray™ Terminal Labeling Kit with Biotin-ddUTP** has been developed for DNA probe array assays. This method uses Biotin-ddUTP and terminal deoxynucleotide transferase to catalyze the addition of a single biotin-ddUMP to the 3'-OH terminus of an amplified and fragmented target DNA molecule. A standard reaction will label up to 100 picomoles (equivalent to 1 µg of a 30-nucleotide sequence). When DNA fragments are labeled at the 3'-OH terminus, the sequence bias that occurs with either nick translation or random priming is eliminated.

REAGENTS PROVIDED

The ENZO **BioArray™ Terminal Labeling Kit with Biotin-ddUTP** contains sufficient reagents for 25 labeling reactions of 100 picomoles of fragmented DNA, corresponding to 25 x 1µg of a 30-nucleotide sequence.

Biotin-ddUTP (100X), 25 µl

5X Reaction Buffer, 500 µl

Potassium cacodylate buffer, pH 7, containing β-mercaptoethanol and stabilizer

10X CoCl₂ Solution, 250 µl

Terminal Deoxynucleotide Transferase (50X), 50 µl
in storage buffer

WARNING: The **5X Reaction Buffer** and **Terminal Deoxynucleotide Transferase** contain potassium cacodylate which contains arsenic. It is toxic. Use gloves during the handling of these reagents. Dispose of these reagents and the individual labeling reaction waste materials according to local regulations.

REAGENTS AND MATERIALS REQUIRED

Preparation and Analysis of Labeled DNA Fragments

- 37°C Water Bath
- Agarose gel (e.g., 4%) - optional
- UV Transilluminator- optional

Termination of Labeling Reaction

- 0.2 M EDTA, pH 8

STORAGE

Upon receipt, store all reagents at -20°C, in a freezer that is not self-defrosting.

PROCEDURE FOR TARGET LABELING

1. To label amplified and fragmented target DNA, add the following reagents (in the indicated order) to a microcentrifuge tube, keeping the tube at **room temperature** while additions are made:

Reagent	Volume
5X Reaction Buffer	20 µl
10X CoCl ₂	10 µl
Amplified and fragmented target	variable (as required from target preparation protocol)
Biotin-ddUTP	1 µl
Terminal Deoxynucleotide Transferase	2 µl
Distilled or deionized water	variable (to give a final reaction volume of 100 µl)
Total Volume	100 µl

2. Carefully mix the reagents in the tube and collect the mixture in the bottom of the microcentrifuge tube by brief (5 seconds) microcentrifugation.
3. Incubate the tube for **15** minutes in a 37°C water bath. The reaction may be allowed to proceed for up to 1 hour.
4. If desired, transfer the tube to ice water (2-4°C), remove 2 µl and analyze on an agarose or acrylamide gel.
5. Stop the reaction by adding 5 µl of 0.2 M EDTA.
6. Store the labeled DNA fragments at -20°C. Do not freeze and thaw repeatedly.

For Research Use Only

This product is sold by ENZO LIFE SCIENCES, INC. for research purposes only by the end-user in the research market and is not intended for diagnostic or therapeutic use. Purchase does not include any right or license to use, develop or otherwise exploit this product commercially. Any commercial use, development or exploitation of this product or development using this product without the express prior written authorization of ENZO LIFE SCIENCES, INC. is strictly prohibited.

For Technical Assistance call ENZO:

Toll free from the U.S. and Canada: 1-800-221-7705

All others: 631-694-7070

Fax: 631-694-7501

e-mail: enzolifesciences.com

<http://www.enzolifesciences.com>

This product or the use of this product is covered by one or more Enzo patents, including, but not limited to, the following: U.S. Patent No. 4,994,373; Canadian Patent No. 1,309,672; and patents pending.

Rev. 1.0.4, October 2006