

BIOARRAY™ Terminal labeling kit with biotin-ddUTP for DNA probe array assays

The BIOARRAY™ 3'-OH Terminal Labeling System is the gold standard end-labeling system for use with Affymetrix® DNA SNP, Resequencing and Prokaryotic microarrays.

The BIOARRAY™ Terminal Labeling Kit with Biotin-ddUTP has been developed for DNA probe array assays. This method uses Bio-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 1µg 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.

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-42630	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

Contents

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

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)