



BioArray™ Control Deoxyoligo B2 (3 nM)

Biotin-labeled deoxyoligonucleotide sequence
for alignment of Eukaryotic Expression Arrays

Cat. No. 42662-30	30 reactions
Cat. No. 42662-150	150 reactions

For Research Use Only

DESCRIPTION

The ENZO **BioArray™ Control Deoxyoligo B2** is a synthetic biotin-labeled sequence that can be used in eukaryotic expression assays and in aiding software programs such as the Affymetrix Microarray Suite by aligning the grid with the array image. The **Control Deoxyoligo B2**, together with the **Eukaryotic Hybridization Controls** (included in Part Nos. 42661-30 and 42661-150), provides an internal control for assaying quality and performance. This is done by adding the controls to biotin-labeled target cRNA before hybridization to probe sequences on the array. The **Control Deoxyoligo B2** is supplied as a 3 nM solution.

REAGENTS PROVIDED

	Cat. 42662-30 (30 reactions)	Cat. 42662-150 (150 reactions)
Control Deoxyoligo B2 (3 nM)	1 x 150 µL	1 x 750 µL

STORAGE

Upon receipt, store the reagent at -20°C or -70°C in a freezer that is not self-defrosting. In order to prevent any deterioration due to repeated freezing and thawing, aliquot the reagent in volumes suitable for future use.

PREPARATION AND USE

1. Thaw the Control Deoxyoligo B2 at 65 °C for 5 minutes.
2. When completely thawed, gently vortex. Centrifuge the tube for about 5 seconds in a microcentrifuge.

NOTE: When thawed the first time, the Control Deoxyoligo B2 should be distributed into smaller volumes and refrozen for future use. This will prevent deterioration that results from freezing and thawing more than three times.

3. Add the Control Deoxyoligo B2 according to the following table. The required amounts for the 20X Eukaryotic Hybridization Controls are also given below.

Array	Hybridization Mix Volume	Control Deoxyoligo B2 (3 nM)	20X Eukaryotic Hybridization Controls
Mini or Micro	100 µL	1.67 µL	5 µL
Midi	200 µL	3.33 µL	10 µL
Standard	300 µL	5.00 µL	15 µL

This product or the use of this product is covered by one or more claims of Enzo patents, including, but not limited to, the following: U.S. Patent Nos. 4,711,955; 5,328,824; 5,449,767; 5,476,928; 4,994,373 and 5,328,824; Canadian Patent Nos. 1,309,672; and patents pending.

IMPORTANT NOTES

1. Take precautions to prevent microbial contamination.
2. Use RNase-free tubes and pipet tips.
3. Discoloration or precipitation indicates reagent deterioration.
4. Do not use reagents after three freeze-thaw cycles.

RELATED PRODUCTS

- **BioArray™** Eukaryotic Hybridization Control, Cat. No. 42661-30 (30 reactions) and Cat. No. 42661-150 (150 reactions)
- **BioArray™** 3'-OH Terminal Labeling Kit with Biotin-ddUTP, Cat. No. 42630
- **BioArray™** 3'-OH Terminal Labeling Kit with Fluorescein-ddCTP, Cat. No. 42631
- **BioArray™** RNA Transcript Labeling Kit (T3/T7) with Fluorescein-UTP, Cat. No. 42650
- **BioArray™ HighYield™** RNA Transcript Labeling Kits (T7), Cat. Nos. 42655-10 and 42655-20
- **BioArray™ HighYield™** RNA Transcript Labeling Kit (T3), Cat. No. 42656
- **BioArray™** RNA Amplification Kits, Cat. Nos. 42410S, 42410X, 42410S-B, 42410S-BB and 42410X-BB
- **BioArray™ GeneBeam™** First-Strand cDNA Labeling Kit, Cat. No. 42530

BioArray™, HighYield™ and GeneBeam™ are trademarks of Enzo Life Sciences, Inc.

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: custserv@enzobio.com

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.

Version 1, June 2004