



## BioArray™ Eukaryotic Hybridization Controls

Biotinylated control RNA sequences for assaying quality and performance of GeneChip® Eukaryotic Expression Assays

Cat. No. 42661-30 30 reactions

Cat. No. 42661-150 150 reactions

For Research Use Only

### INTRODUCTION

The ENZO **BioArray™ Eukaryotic Hybridization Controls** contain biotin-labeled transcripts prepared from three *E.coli* genes, *bioB*, *bioC* and *bioD*, and one from bacteriophage P1, *cre*. These together with the synthetic biotin-labeled deoxyoligonucleotide control (Deoxyoligo B2) are added to biotin-labeled target cRNA before hybridization to probe sequences on the array. These sequences of known concentration provide an internal control for assaying quality and performance. Deoxyoligo B2 aids software programs such as the Affymetrix Microarray Suite by aligning the grid with the array image.

### REAGENTS PROVIDED

	Cat. 42661-30 (30 reactions)	Cat. 42661-150 (150 reactions)
20X Eukaryotic Hybridization Controls (30 pM <i>bioB</i> , 100 pM <i>bioC</i> , 500 pM <i>bioD</i> and 2000 pM <i>cre</i> )	6 x 75 µL	2 x 1125 µL
Control Deoxyoligo B2 (3 nM)*	1 x 150 µL	1 x 750 µL

\*This reagent is also available as a separate product, Cat. Nos. 42662-30 and 43662-150.

### STORAGE

Upon receipt, store all reagents at -20°C or -70°C in a freezer that is not self-defrosting. For long-term storage, store the 20X Eukaryotic Hybridization Controls at -70°C.

In order to prevent any deterioration due to repeated freezing and thawing, aliquot the reagents in volumes suitable for future use.

### PROCEDURE

1. Thaw one vial of the 20X Eukaryotic Hybridization Controls and the Control Deoxyoligo B2 vial at 65°C for 5 minutes.
2. When the vials are completely thawed, gently vortex. Centrifuge the tubes for about 5 seconds in a microcentrifuge.

**NOTE:** When the reagents are thawed the first time, they should be distributed into smaller volumes and refrozen for future use. This will avoid deterioration of the probe that results from freezing and thawing more than three times.

3. Add the Eukaryotic Hybridization Controls and Control Deoxyoligo B2 according to the following table.

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

### 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™ Control Deoxyoligo B2**, Cat. No. 42662-30 (30 reactions) and Cat. No. 42662-150 (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™ High Yield™ RNA Transcript Labeling Kits (T7)**, Cat. Nos. 42655-10 and 42655-20
- **BioArray™ High Yield™ 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*, *High Yield* and *GeneBeam* are trademarks of Enzo Life Sciences, Inc. GeneChip is a registered trademark of Affymetrix, 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: [enzolifesciences.com](mailto:enzolifesciences.com)

<http://www.enzolifesciences.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.

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 No. 4,994,373; Canadian Patent No. 1,309,672; and patents pending.

Version 4.0.2, May 2006