

BioArray[®] Eukaryotic Hybridization Controls

Catalog #:

ENZ-42661-30 30 reactions

ENZ-42661-150 150 reactions

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Please contact Enzo Life Sciences Technical Support if necessary.

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.

MATERIALS SUPPLIED

	ENZ-42661-30 (30 reactions)	ENZ-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. ENZ-42662-30 and ENZ-42662-150.*

STORAGE

Upon receipt, store all reagents at -20°C in a freezer that is not self-defrosting. For long-term storage, store the 20X Eukaryotic Hybridization Controls at -80°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.



Reagents require separate storage conditions.

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 (3nM)
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.

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com