



BIOARRAY™

Eukaryotic hybridization controls

Biotin-labeled transcripts for microarray internal controls

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.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-42661-150	150Reactions
---------------	--------------

Manuals, SDS & CofA

[View Online »](#)

- Provides known concentration of transcripts for microarray internal controls
- Use together with Deoxyoligo B2 (Prod No. [ENZ-42662](#)) to assay quality and performance of microarray

Handling & Storage

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Contents **20X Eukaryotic Hybridization Controls** (30 pM bioB, 100 pM bioC, 500 pM bioD and 2000 pM cre)
Control Deoxyoligo B2 (3 nM)



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)