

# BIOARRAY HIGHYIELD<sup>®</sup> RNA Transcript Labeling Kit (T7)

The BIOARRAY HIGHYIELD<sup>®</sup> RNA Transcript Labeling System is the gold standard transcript labeling system and is intended for use with Affymetrix<sup>®</sup> GeneChip<sup>®</sup> brand eukaryotic expression microarrays.

**The gold standard in transcript labeling, for use with Affymetrix GeneChip microarrays.**

The BIOARRAY HIGHYIELD<sup>®</sup> RNA Transcript Labeling Kit (T7) has been developed for the production of large amounts of hybridizable biotin-labeled RNA targets by *in vitro* transcription from bacteriophage T7 RNA polymerase promoters. Using T7 RNA polymerase and biotin-labeled nucleotides, large amounts of single stranded nonradioactive RNA molecules can be produced *in vitro*. Because of the nature of transcription reactions, many RNA copies of the template DNA are produced. Because RNA-DNA hybrids have a higher melting temperature than corresponding DNA-DNA hybrids, single-stranded RNA targets offer higher target avidity and greater sensitivity than DNA probes. RNA targets offer selectivity unavailable with DNA targets—being single stranded, they are strand specific and hybridize more effectively to probes because the target population does not self-hybridize.

Citations: 67

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-42655-40

40Reactions

## Manuals, SDS & CofA

[View Online »](#)

- Produces large amounts of RNA from template DNA
- Incorporates biotin during transcription

## 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

10X HY Reaction Buffer  
10X Biotin-Labeled Ribonucleotides  
10X DTT  
10X RNase Inhibitor Mix  
20X T7 RNA Polymerase

### Specificity

**Yield:** Each lot is shown to generate 50 µg of cRNA when transcribed from a double stranded cDNA library generated from 2 µg total RNA (mixed tissue) in a single round of amplification.

**Incorporation:** Each lot is tested to ensure at least 10 % of incorporated nucleotides are biotin labeled.

**RNase activity:** Insignificant amounts of RNase (less than 100 fg RNase A activity in 1 µL of reagent) are detected.

**% P value and 3'/5' ratio:** The cRNA synthesized by the kit is tested on Affymetrix HU133A 2.0 chip and shown to give > 35% P-calls, with 3'/5' ratios for both β-actin and GAPDH of < 2.5.

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)