



GeneChip® Expression 3'-Amplification Reagents for IVT Labeling

P/N 900449 (30 eukaryotic reactions)
In vitro transcription with MEGAscript™.

INTRODUCTION

The GeneChip® IVT Labeling Kit is optimized specifically for producing amplified and biotinylated targets to hybridize to eukaryotic GeneChip brand arrays for expression analysis. Each kit contains sufficient reagents for completing 30 *in vitro* transcription reactions.

The template DNA for this IVT reaction is typically double-stranded cDNA containing the T7 promoter sequence. In GeneChip target labeling experiments, the T7 promoter sequence is incorporated into the cDNA template by using a T7-Oligo(dT) Promoter Primer in the initial reverse transcription reaction (described in the *GeneChip® Expression Analysis Technical Manual*, available at www.affymetrix.com). Sufficient cDNA template can be obtained using one of the two protocols from high-quality starting materials:

- 1 to 15 µg of sample total RNA, or 0.2 to 2 µg of mRNA, using the One-Cycle cDNA Synthesis Procedure
- 10 to 100 ng of sample total RNA using the Two-Cycle cDNA Synthesis Procedure

Using this IVT Labeling Kit, over 30 µg of biotinylated cRNA targets are usually generated with the MEGAscript™ T7 Polymerase (**IVT Labeling Enzyme Mix**) in the presence of a mixture of the four natural ribonucleotides and one biotin-conjugated nucleotide analog at optimized ratio (**IVT Labeling NTP Mix**). The chemical structure of the analog is illustrated in Figure 1.

The synthetic analog is efficiently incorporated into the cRNA target during the *in vitro* transcription labeling reaction mediated by T7 RNA polymerase as a pseudouridine reagent. The biotinylated cRNA targets are then purified, fragmented, and hybridized to GeneChip expression arrays.

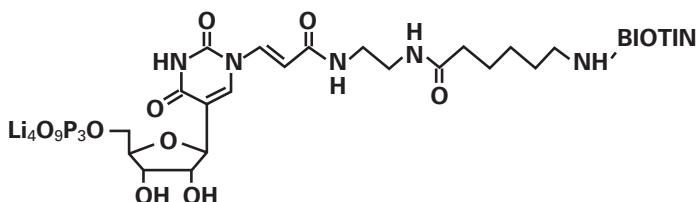


Figure 1. Structure of the biotin-conjugated nucleotide analog

A positive control, the **3'-Labeling Control**, is included in the kit and allows users to test the transcriptional activity of the kit components. Using 0.5 µg of the **3'-Labeling Control** as a template, routinely, over 30 µg of biotinylated IVT product is generated after 4 hours of incubation. The amplified, labeled

product can be visually inspected on an agarose gel or other electrophoretic system, and a distinct band at approximately 2 Kb should be observed.

NOTE Hybridization cocktail preparation and fluidics scripts have been optimized specifically for using this IVT Labeling Kit. Follow the instructions described below for best results.

CONTENTS

Volume	Component
120 µL	10X IVT Labeling Buffer
120 µL	IVT Labeling Enzyme Mix
360 µL	IVT Labeling NTP Mix
10 µL	3'-Labeling Control (0.5 µg/µL)
910 µL	RNase-free Water

STORAGE

Kit should be kept at -20°C in a non-frost-free freezer. Performance of the kit has been shown to be unaffected for up to eight freeze-thaw cycles.

GENECHIP EUKARYOTIC TARGET LABELING ASSAYS FOR EXPRESSION ANALYSIS

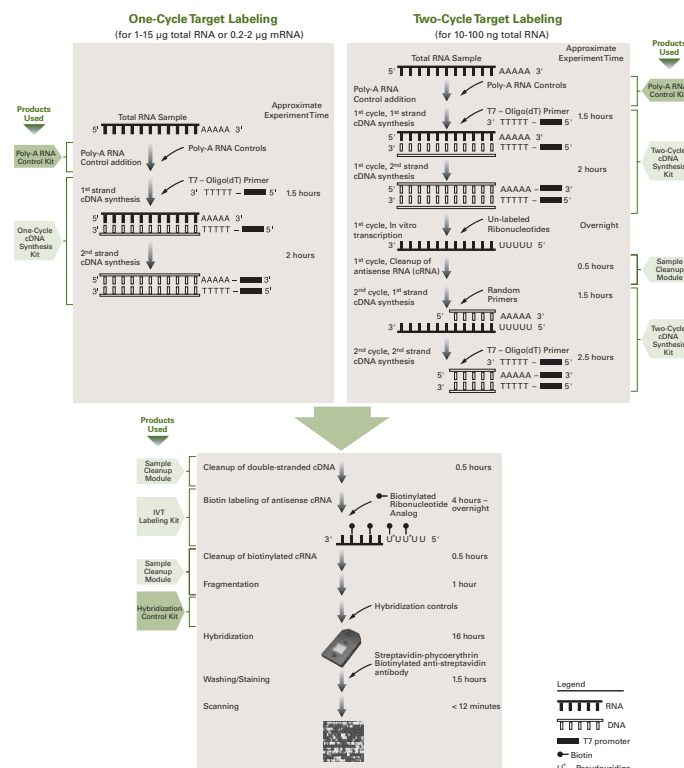


Figure 2. Schematic illustration of GeneChip® target labeling assays for expression analysis

OTHER NECESSARY REAGENTS NOT PROVIDED WITH THE KIT _____

Kit name	P/N	Size
One-Cycle cDNA Synthesis Kit	900431	30 rxns
Two-Cycle cDNA Synthesis Kit	900432	30 rxns
Sample Cleanup Module	900371	30 rxns
Eukaryotic Poly-A RNA Control Kit	900433	~ 100 rxns
Hybridization Controls	900454 and 900457	30 rxns and 150 rxns

IVT AMPLIFICATION AND LABELING PROCEDURE

NOTE This kit is only used for the IVT labeling step for generating biotin-labeled cRNA. For the IVT amplification step using unlabeled ribonucleotides in the First Cycle of the Two-Cycle cDNA Synthesis Procedure, a separate kit is recommended (MEGAscript® T7 Kit, Ambion, Inc.). Use only nuclease-free water, buffers, and pipette tips.

IMPORTANT Store all reagents in a -20°C freezer that is not self-defrosting. Prior to use, centrifuge all reagents briefly to ensure that the solution is collected at the bottom of the tube.

IMPORTANT The Target Hybridizations and Array Washing protocols have been optimized specifically for this IVT Labeling Protocol. Closely follow the recommendations described below for maximum array performance.

1. Use the following table to determine the amount of cDNA used for each IVT reaction following the cDNA cleanup step.

Starting Material	Volume of cDNA to use in IVT
Total RNA	
10 to 100 ng	all (~12 µL)
1.0 to 8.0 µg	all (~12 µL)
8.1 to 15 µg	6 µL
mRNA	
0.2 to 2 µg	all (~12 µL)

2. Transfer the needed amount of template cDNA to RNase-free microfuge tubes and add the following reaction components in the order indicated in the table below. If more than one IVT reaction is to be performed, a master mix can be prepared by multiplying the reagent volumes by the number of reactions. Do not assemble the reaction on ice, since spermidine in the **10X IVT Labeling Buffer** can lead to precipitation of the template cDNA.

Reagent	Volume
Template cDNA*	variable (see table above)
RNase-free Water	variable (to give a final reaction volume of 40 µL)
10X IVT Labeling Buffer	4 µL
IVT Labeling NTP Mix	12 µL
IVT Labeling Enzyme Mix	4 µL
Total Volume	40 µL

*1 µg of the 3'-Labeling Control can be used in place of the template cDNA sample in this reaction as a positive control for the IVT components in the kit.

3. Carefully mix the reagents and collect the mixture at the bottom of the tube by brief (5 second) microcentrifugation.
4. Incubate at 37°C for 16 hours. To prevent condensation that may result from water bath-style incubators, incubations are best performed in oven incubators for even temperature distribution, or in a thermal cycler.

NOTE Overnight IVT reaction time has been shown to maximize the labeled cRNA yield with high-quality array results. Alternatively, if a shorter incubation time (4 hours) is desired, 1 µL (200 units) of cloned T7 RNA polymerase (can be purchased directly from Ambion, P/N 2085) can be added to each reaction and has been shown to produce adequate labeled cRNA yield within 4 hours. The two different incubation protocols generate comparable array results, and users are encouraged to choose the procedure that best fits their experimental schedule and process flow.

5. Store labeled cRNA at -20°C, or -70°C if not purifying immediately. Alternatively, proceed to Cleanup and Quantification of Biotin-Labeled cRNA as described in the *GeneChip Expression Analysis Technical Manual*.

EUKARYOTIC TARGET HYBRIDIZATION _____

1. Mix the following for each target, scaling up volumes for hybridization to multiple probe arrays.

NOTE A final concentration of 10% DMSO is added in the Hybridization Cocktail for optimal results.

IMPORTANT It is imperative that the frozen 20X Hybridization Control Stock is heated to 65°C for 5 minutes to completely resuspend the cRNA before adding to the hybridization cocktail.

Component	49 Format/ 64 Format	100 Format	400 Format/ 169 Format	Final Concentration
Fragmented cRNA**	15 µg	10 µg	5 µg	0.05 µg/µL
Control Oligonucleotide B2 (3 nM)	5 µL	3.3 µL	1.7 µL	50 pM
20X Eukaryotic Hybridization Controls (<i>bioB</i> , <i>bioC</i> , <i>bioD</i> , <i>cre</i>)	15 µL	10 µL	5 µL	1.5, 5, 25 and 100 pM respectively
Herring Sperm DNA (10 mg/mL)	3 µL	2 µL	1 µL	0.1 mg/mL
Acetylated BSA (50 mg/mL)	3 µL	2 µL	1 µL	0.5 mg/mL
2X Hybridization Buffer	150 µL	100 µL	50 µL	1X
DMSO	30 µL	20 µL	10 µL	10%
H ₂ O	to final volume of 300 µL	to final volume of 200 µL	to final volume of 100 µL	
Final volume	300 µL	200 µL	100 µL	

*Please refer to specific probe array package insert for information on array format.

**Please see Section 2, Chapter 1 in the *GeneChip Expression Analysis Technical Manual* for amount of adjusted fragmented cRNA to use when starting from total RNA.

- Hybridize to expression probe arrays as described in the *Expression Analysis Technical Manual*.

WASHING AND SCANNING _____

- Prepare the SAPE solution and Antibody solution as described in Section 2, Chapter 4 of the *Expression Analysis Technical Manual* using the Antibody Amplification Stain Protocol for Eukaryotic Targets.
- For 49 Format and 64 Format arrays with 18 µm features**, use the existing EukGE-WS2v4 or EukGE-WS2v4_450 Fluidics Scripts.

For 49 Format and 64 Format arrays with 11 µm features, download the new EukGE-WS2v5 Fluidics Scripts from www.affymetrix.com. Alternatively, modify the FlexEukGE-WS2v4 or FlexEukGE-WS2v4_450 Fluidics Scripts as follows, and rename the scripts as EukGE-WS2v5 or EukGE-WS2v5_450.

For 100 Format arrays, use the existing Midi-euk2 or Midi-euk2_450 Fluidics Scripts.

	49 Format/ 64 Format EukGE-WS2v5	100 Format (Midi) Midi_euk2 ^a
Post Hyb Wash #1	10 cycles of 2 mixes/cycle with Wash Buffer A at 30°C	10 cycles of 2 mixes/cycle with Wash Buffer A at 30°C
Post Hyb Wash #2	6 cycles of 15 mixes/cycle with Wash Buffer B at 50°C	6 cycles of 15 mixes/cycle with Wash Buffer B at 50°C
Stain	Stain the probe array for 5 minutes in SAPE solution at 35°C	Stain the probe array for 5 minutes in SAPE solution at 35°C
Post Stain Wash	10 cycles of 4 mixes/cycle with Wash Buffer A at 30°C	10 cycles of 4 mixes/cycle with Wash Buffer A at 30°C
2nd Stain	Stain the probe array for 5 minutes in antibody solution at 35°C	Stain the probe array for 5 minutes in antibody solution at 35°C
3rd Stain	Stain the probe array for 5 minutes in SAPE solution at 35°C	Stain the probe array for 5 minutes in SAPE solution at 35°C
Final Wash	15 cycles of 4 mixes/cycle with Wash Buffer A at 35°C. The holding temperature is 25°C	15 cycles of 4 mixes/cycle with Wash Buffer A at 35°C. The holding temperature is 25°C

- Wash Buffer A = non-stringent wash buffer
- Wash Buffer B = stringent wash buffer

a. Midi_euk2_450 for the FS-450/250

FUNCTIONAL TESTING AND QUALITY CONTROL _____

GeneChip Array Functional Testing

Each lot is functionally tested in the GeneChip expression assay following the recommended protocol.

SAFETY INFORMATION _____

An MSDS is available from Affymetrix. The kit components are not believed to present any health hazard. However, the chemical, physical, and toxicological properties of these components have not been thoroughly investigated.

CONTACT INFORMATION _____

Affymetrix, Inc.

3380 Central Expressway
Santa Clara, CA 95051
USA

E-mail: support@affymetrix.com
Tel: 1-888-362-2447 (1-888-DNA-CHIP)
Fax: 1-408-731-5441

Affymetrix UK Ltd

Voyager, Mercury Park,
Wycombe Lane, Wooburn Green,
High Wycombe HP10 0HH
United Kingdom

E-mail: supporteurope@affymetrix.com
Tel: +44 (0) 1628 552550
Fax: +44 (0) 1628 552585



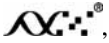
Affymetrix Japan, K.K.

Mita NN Bldg
16 Floor, 4-1-23 Shiba,
Minato-ku, Tokyo 108-0014
Japan

E-mail: supportjapan@affymetrix.com
Tel: +81-(0)3-5730-8200
Fax: +81-(0)3-5730-8201

www.affymetrix.com

TRADEMARKS _____

Affymetrix®, GeneChip®, , , ,
GenFlex®, HuSNP®, EASI™, MicroDB™, Jaguar™, Flying
Objective™, NetAffx™, CustomExpress™, CustomSeq™, 'Tools
to Take You As Far As Your Vision™', and 'The Way Ahead™'
are trademarks owned or used by Affymetrix, Inc.
MEGAscript™ is a trademark owned by Ambion, Inc.

LIMITED LICENSE _____

EXCEPT AS EXPRESSLY SET FORTH HEREIN, NO RIGHT TO COPY, MODIFY, DISTRIBUTE, MAKE DERIVATIVE WORKS OF, PUBLICLY DISPLAY, MAKE, HAVE MADE, OFFER TO SELL, SELL, USE, OR IMPORT PROBE ARRAYS OR ANY OTHER PRODUCT IS CONVEYED OR IMPLIED WITH THE PROBE ARRAYS, INSTRUMENTS, SOFTWARE, REAGENTS, OR ANY OTHER ITEMS PROVIDED HEREUNDER. EXCEPT FOR CERTAIN ARRAYS AND REAGENTS DESIGNATED AS "ANALYTE SPECIFIC REAGENTS" (SEE APPLICABLE PACKAGE INSERT) WHICH ARE LICENSED FOR USE AS ANALYTE SPECIFIC REAGENTS OR RESEARCH USE, ALL PRODUCTS (INCLUDING THE PROBE ARRAYS, INSTRUMENTS, SOFTWARE, AND REAGENTS) DELIVERED HEREUNDER ARE LICENSED TO BUYER FOR RESEARCH USE ONLY. THIS LIMITED LICENSE PERMITS ONLY THE USE BY BUYER OF THE PARTICULAR PRODUCT(S), IN ACCORDANCE WITH THE WRITTEN INSTRUCTIONS PROVIDED THEREWITH, THAT BUYER PURCHASES FROM AFFYMETRIX (AFX) OR ITS AUTHORIZED REPRESENTATIVE. THE PURCHASE OF ANY PRODUCT(S) DOES NOT BY ITSELF CONVEY OR IMPLY THE RIGHT TO USE SUCH PRODUCT(S) IN COMBINATION WITH ANY OTHER PRODUCT(S). IN PARTICULAR, (I) NO RIGHT TO MAKE, HAVE MADE, OR DISTRIBUTE OTHER PROBE ARRAYS IS CONVEYED OR IMPLIED BY THE PROBE ARRAYS, (II) NO RIGHT TO MAKE, HAVE MADE, IMPORT, DISTRIBUTE, OR USE PROBE ARRAYS IS CONVEYED OR IMPLIED BY THE INSTRUMENTS OR SOFTWARE, AND (III) NO RIGHT TO USE PROBE ARRAYS IN COMBINATION WITH INSTRUMENTS OR SOFTWARE IS CONVEYED UNLESS ALL COMPONENT PARTS HAVE BEEN PURCHASED FROM AFX OR ITS AUTHORIZED REPRESENTATIVE. FURTHERMORE, PROBE ARRAYS DELIVERED HEREUNDER ARE LICENSED FOR ONE (1) TIME USE ONLY AND MAY NOT BE REUSED. THE PRODUCTS DO NOT HAVE FDA APPROVAL. NO PATENT LICENSE IS CONVEYED TO BUYER TO USE, AND BUYER AGREES NOT TO USE, THE PRODUCTS IN ANY SETTING REQUIRING FDA OR SIMILAR REGULATORY APPROVAL OR EXPLOIT THE PRODUCTS IN ANY MANNER NOT EXPRESSLY AUTHORIZED IN WRITING BY AFX IN ADVANCE.

CERTAIN COMPONENTS OF THIS KIT OR COMBINATIONS THEREOF MAY BE COVERED BY U.S. PATENT NOS. 5,256,555, 6,586,218 AND 6,586,219, ASSIGNED TO AMBION, INC.

COPYRIGHT _____

©2003, 2004 Affymetrix, Inc. All rights reserved.