

## GeneChip® Globin-Reduction RNA Controls

P/N 900586  
MANUFACTURED BY QIAGEN®  
ON BEHALF OF PREANALYTIX™  
FOR AFFYMETRIX®



### INTRODUCTION

The GeneChip® Globin-Reduction RNA Controls are designed to provide both positive and negative control RNA samples that can be used to validate the globin reduction procedure and are intended for use in conjunction with the GeneChip® Blood RNA Concentration Kit (P/N 900585).

These kits seamlessly integrate the globin reduction procedure with the GeneChip® Target Labeling Assay described in the *GeneChip® Expression Analysis Technical Manual* and require no additional steps or time. The kits are intended for use with methods that isolate RNA from unfractionated human whole blood only, such as samples isolated using the PAXgene™ Blood RNA System from PreAnalytiX.

The Globin-Reduction RNA Controls box contains 2 vials of control RNA:

- **Jurkat RNA:** This control RNA is purified from the Jurkat cell line that originated from human peripheral blood and contains negligible amounts of globin mRNA. The Jurkat RNA control, therefore, represents RNA purified from the leukocyte fraction of whole blood (i.e., RNA purified from whole blood that had been subjected to erythrocyte lysis before RNA purification).
- **Jurkat + Globin RNA:** This control is the Jurkat RNA described above but spiked with *in vitro* synthesized polyadenylated globin transcripts. The Jurkat + Globin RNA control mimics RNA purified from unfractionated whole blood. The amount of globin *in vitro* transcripts in the Jurkat + Globin RNA control has been adjusted so that the derived cRNAs will provide GeneChip array analysis that are similar to those obtained from human whole blood.

The kit provides sufficient Jurkat RNA and Jurkat + Globin RNA for 6 reactions each and, therefore, a total of 12 reactions. The protocol described in the handbook of the GeneChip® Blood RNA Concentration Kit recommends preparing both Jurkat RNA and Jurkat + Globin RNA with and without globin reduction treatment each in three technical replicates.

### ORDERING INFORMATION

#### From Affymetrix

P/N	Product Name	Description
<b>KIT</b>		
900586	GeneChip® Globin Reduction RNA Controls	
	• Jurkat RNA	6 rxns
	• Jurkat + Globin RNA	6 rxns
<b>SUPPORTING PRODUCTS</b>		
900585	GeneChip® Blood RNA Concentration Kit	30 rxns
900493	One-Cycle Target Labeling and Control Reagents	30 rxns
<i>These reagents are necessary for performing the globin reduction procedure and GeneChip Eukaryotic Target Labeling Assay.</i>		

#### From QIAGEN

The following products are recommended for blood collection, storage, and RNA isolation.

P/N	Product Name	Description
<b>PAXGENE BLOOD RNA SYSTEM - FOR BLOOD COLLECTION, RNA STABILIZATION, AND PURIFICATION</b>		
762115 <sup>a</sup> 762105 <sup>b</sup> 762125 <sup>c</sup>	PAXgene™ Blood RNA Tubes	100 tubes
<i>a: Canada and USA, b: Japan, c: All other countries.</i>		
762134	PAXgene™ Blood RNA Kit	50 preps
762132	PAXgene™ Blood RNA Validation Kit	10 preps and 10 tubes
762331	PAXgene™ 96 Blood RNA Kit	4 x 96 preps
<b>RNASE-FREE DNASE SET - FOR DNASE DIGESTION DURING RNA PURIFICATION</b>		
79254	RNase-Free DNase Set	50 preps

### STORAGE CONDITIONS

The GeneChip® Globin-Reduction RNA Controls are shipped frozen on dry ice. All components should be stored immediately upon receipt at -20°C. When stored at -20°C and handled correctly, the RNA samples can be kept for at least nine months and subjected to up to ten freeze-thaw cycles without any compromise in performance.

### 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.

### 24-Hour Emergency Information

Please call CHEMTREC at 1-800-424-9300 within the United States, Canada, Puerto Rico, and the Virgin Islands. All other callers can use CHEMTREC's international number (emergency collect calls are accepted) at 1-703-527-3887.

### PRODUCT USE LIMITATIONS

For research use only. Not for use in diagnostic procedures.

---

## 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

Please visit our web site for international distributor contact information.

[www.affymetrix.com](http://www.affymetrix.com)

---

## TRADEMARKS

Affymetrix<sup>®</sup>, GeneChip<sup>®</sup>, <sup>®</sup>, <sup>®</sup>, <sup>®</sup>, GenFlex<sup>®</sup>, HuSNP<sup>®</sup>, EASI<sup>™</sup>, MicroDB<sup>™</sup>, Jaguar<sup>™</sup>, Flying Objective<sup>™</sup>, NetAffx<sup>™</sup>, NimbleExpress<sup>™</sup>, CustomExpress<sup>®</sup>, CustomSeq<sup>™</sup>, "Tools to Take You As Far As Your Vision<sup>®</sup>", and The Way Ahead<sup>™</sup> are trademarks owned or used by Affymetrix, Inc.

PAXgene<sup>™</sup> and PreAnalytiX<sup>™</sup> are trademarks of PreAnalytiX GmbH. QIAGEN<sup>™</sup> is a registered trademark of QIAGEN Group. Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are not to be considered unprotected by law.

---

## 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.

---

## COPYRIGHT

©2004 PreAnalytiX and Affymetrix, Inc. All rights reserved.