



GeneChip® Human Genome U95 Set

Intended Use

The GeneChip® Human Genome U95 Set consists of 5 probe arrays (HG-U95Av2, HG-U95B, HG-U95C, HG-U95D, and HG-U95E) which allows one to monitor the relative abundance of greater than 60,000 mRNA transcripts from genes and ESTs. We selected these sequences from Build 95 of the UniGene Database (derived from GenBank®113, dbEST/10-02-99).

GeneChip probe arrays are made by synthesizing oligonucleotide probes directly onto a glass surface. Each 25-mer oligonucleotide probe is uniquely complementary to a gene, with approximately 16 pairs of oligonucleotide probes used to measure the transcript level of each of the sequences represented on the Human Genome U95 Set.

All sequences represented on the GeneChip® HuGeneFL Array are also represented on the HG-U95Av2 Array. Sequences represented on the GeneChip® Hu35K Set are represented on the HG-U95B and HG-U95C Arrays. Due to the evolving nature of public databases, probe sets for these sequences will not be identical and in some cases will be represented by a completely new probe set. As a result, data generated with different versions of the human arrays may not always produce concordant results.

GeneChip brand probe arrays are for research use only and not intended for use in diagnosis of disease. Please visit www.affymetrix.com for a complete list of supporting documentation including procedures regarding target preparation, target hybridization, fluidics station setup, probe array scan, and data analysis.

Instrumentation and Software Required

1. GeneChip® Scanner 3000
2. Affymetrix® GeneChip® Command Console® Software (AGCC)
3. GeneChip® Fluidics Station

Precautions

1. GENECHIP PROBE ARRAYS ARE FOR RESEARCH USE ONLY; NOT FOR DIAGNOSTIC PROCEDURES.
2. Avoid microbial contamination, which may cause erroneous results.
3. **WARNING: All biological specimens and materials with which they come into contact should be handled as if capable of transmitting infection and disposed of with proper precautions in accordance with federal, state, and local regulations. This includes adherence to the OSHA Bloodborne Pathogens Standard (29 CFR 1910.1030) for blood-derived and other samples governed by this act. Never pipet by mouth. Avoid specimen contact with skin and mucous membranes.**
4. **CAUTION:** Exercise standard precautions when obtaining, handling, and disposing of potentially carcinogenic reagents.
5. Exercise care to avoid cross-contamination of samples during all steps of this procedure, as this may lead to erroneous results.
6. Use powder-free gloves whenever possible to minimize introduction of powder particles into sample or probe array cartridges.

Critical Specifications

| | |
|----------------------------|--|
| Number of Arrays/Set | 5 |
| Feature Size | 20 µm |
| Probe Pairs/Sequence | ~16 |
| Array Format | 49 |
| Hybridization Controls | <i>bioB</i> , <i>bioC</i> , <i>bioD</i> , and <i>cre</i> |
| Poly-A Controls | <i>dap</i> , <i>lys</i> , <i>phe</i> , and <i>thr</i> |
| Normalization Controls | 100 probe sets replicated on both A and B arrays |
| Housekeeping/Control Genes | GAPDH, beta-Actin, ISGF-3 (STAT1) |
| Fluidics Protocol | EukGE-WS2v4 <i>Fluidics Station 450/250 scripts are designated by the suffix "_450"</i> |
| Hybridization Volume | 200 µL. The total fill volume of the cartridge is 250 µL. |
| Library Files | HG_U95A, HG_U95B, HG_U95C, HG_U95D, and HG_U95E |

Ordering Information

| P/N | Product Name | Description |
|---|---------------------|---------------|
| Arrays | | |
| 510559 | GeneChip® HG-U95Av2 | 40 ± 5 Arrays |
| 510462 | GeneChip® HG-U95B | 40 ± 5 Arrays |
| 510463 | GeneChip® HG-U95C | 40 ± 5 Arrays |
| 510464 | GeneChip® HG-U95D | 40 ± 5 Arrays |
| 510465 | GeneChip® HG-U95E | 40 ± 5 Arrays |
| The arrays in the GeneChip Human Genome U95 Set are part of the Made-to-Order Program. Made-to-Order Arrays are ordered in whole lots. The lot size for each array in the GeneChip Human Genome U95 Set is 40 ± 5 arrays. | | |

Supporting Products

| | | |
|------------------|---|--|
| 901228 901229 | GeneChip® 3' IVT Express Kit | Sufficient for 10 Reactions Sufficient for 30 Reactions |
| 900720 | GeneChip® Hybridization, Wash, and Stain Kit ¹ | Sufficient for 30 Reactions |

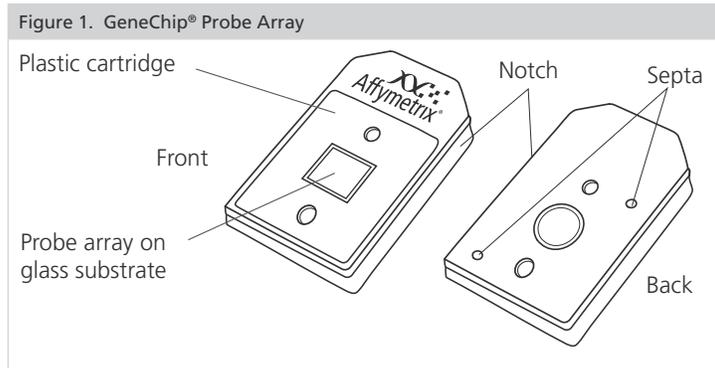
1. Each kit contains one (1) Hybridization Module, one (1) Stain Module, three (3) bottles of Wash Buffer A and one (1) bottle of Wash Buffer B, sufficient for 30 reactions. Individual kit components may be ordered separately.

Affymetrix® products can be purchased directly from Affymetrix in the United States, and many European and Asian countries. For all other territories, please view a list of our distribution partners, which can be located at: www.affymetrix.com/site/contact/index.affx.

Storage, Handling and Stability

The probe array consists of a square glass substrate mounted in a plastic cartridge (Figure 1). The glass contains an array of oligonucleotides that, when mounted, is on the inner glass surface. A chamber in the plastic housing directly under the glass acts as a reservoir where hybridization and washing occur.

Although the inner glass surface of the probe array is protected, any contamination or scratches on the outer surface of the glass can compromise the accuracy of the scan. Avoid touching the surface of the glass with your fingers. Skin oils and other substances, such as lotions or ink, can fluoresce. If the surface of the glass is noticeably dirty, it can be carefully cleaned with a non-abrasive laboratory tissue.



The GeneChip® probe arrays should be stored at 2° to 8°C. Refer to the expiration date on the package label. Do not use probe arrays or reagents after the expiration date.

Limited License

Subject to the Affymetrix terms and conditions that govern your use of Affymetrix products, Affymetrix grants you a non-exclusive, non-transferable, non-sub-licensable license to use this Affymetrix product only in accordance with the manual and written instructions provided by Affymetrix. You understand and agree that except as expressly set forth in the Affymetrix terms and conditions, that no right or license to any patent or other intellectual property owned or licensable by Affymetrix is conveyed or implied by this Affymetrix product. In particular, no right or license is conveyed or implied to use this Affymetrix product in combination with a product not provided, licensed or specifically recommended by Affymetrix for such use.

Contact Information

Affymetrix, Inc.

3420 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
UK and Others Tel: +44 (0) 1628 552550
France Tel: 0800919505
Germany Tel: 01803001334
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

Patents

Products may be covered by one or more of the following patents: U.S. Patent Nos. 5,445,934; 5,744,305; 5,945,334; 6,140,044; 6,261,776; 6,291,183; 6,346,413; 6,399,365; 6,420,169; 6,551,817; 6,610,482; 6,733,977; 6,955,915 and D430,024 and other U.S. or foreign patents. Products are manufactured and sold under license from OGT under 5,700,637 and 6,054,270.

Trademarks

Affymetrix®, GeneChip®, NetAffx®, Command Console®, Genotyping Console™, DMET™, Powered by Affymetrix™, GeneChip-compatible™, and GeneTitan™ are trademarks or registered trademarks of Affymetrix, Inc. All other trademarks are the property of their respective owners.

Copyright

© 2000-2009 Affymetrix, Inc. All rights reserved.