# GeneChip™ Human Genome U133A 2.0 Array

Catalog Numbers 900471, 900468, and 900469

Pub. No. 701501 Rev. 6



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from **thermofisher.com/support**.

#### Product use

The Applied Biosystems<sup>™</sup> GeneChip<sup>™</sup> Human Genome U133A 2.0 Array (HG-U133A 2.0) is one microarray comprised of over 500,000 unique oligonucleotide features covering over 18,000 transcripts and variants, which, in turn, represent approximately 14,500 of the best characterized human genes. All probe sets and normalization genes represented on the GeneChip™ Human Genome U133A Array are identically represented on the HG-U133A 2.0 Array. Sequences used in the design of the array were selected from GenBank $^{\scriptscriptstyle{\text{TM}}}$ , dbEST, and RefSeq. Sequence clusters were created from Build 133 of UniGene (April 20, 2001) and refined by analysis and comparison with a number of other publicly available databases, including the Washington University EST trace repository and the University of California, Santa Cruz, Golden-Path human genome database (April 2001 release). Sequences were further analyzed for correct orientation, false priming, false clustering, alternative splicing, and alternative polyadenylation.

Oligonucleotide probes are synthesized *in situ* complementary to each corresponding sequence. Eleven pairs of oligonucleotide probes are used to measure the level of transcription of each sequence represented.

Visit our website 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

- GeneChip<sup>™</sup> Scanner 3000 7G
- GeneChip Fluidics Station 450
- GeneChip<sup>™</sup> Hybridzation Oven 645
- GeneChip<sup>™</sup> Command Console<sup>™</sup> (GCC) software

### Critical specifications

Item	Specification
Feature size	11 µm
Probe pairs/sequence	11
Array format	100
Hybridization controls	bioB, bioC, bioD, and cre
Poly-A controls	dap, lys, phe, and thr
Normalization controls	100 probe sets
Housekeeping controls	GAPDH, β-Actin, ISGF-3 (STAT1)
Hybridization volume	130 μL.
	The total fill volume of the cartridge is 160 $\mu L$
Library files	HG_U133A 2.0

## **Accessory files**

### Fluidics scripts

The fluidics script used depends on the array type, labeling protocol, and reagents used for cartridge processing. Refer to the GeneChip Fluidics Station Scripts support page to determine which fluidics script is appropriate for your application. The fluidics scripts can be downloaded from our website.

### Library files

Library files contain information about the probe array design characteristics, probe use and content, and scanning and analysis parameters. These files are unique for each probe array. Additional information can be located under the specific array product on our website.

#### Mask files

The GeneChip™ Human Genome U133A 2.0 Array includes a set of human maintenance genes to facilitate the normalization and scaling of array experiments. This set of genes serves as a tool to normalize or scale your data prior to performing data comparison. This set of normalization genes shows consistent levels of expression over a diverse set of tissues. Mask files enabling the use of these probe sets for normalization and scaling are available on our website.

### Comparison spreadsheets

Comparison spreadsheets are designed to assist in understanding the relationship between the data generated using different, but related, GeneChip expression probe arrays. Because the content of this array is identical to the GeneChip Human Genome U133A Array, the same comparison spreadsheet applies to both arrays. Comparison spreadsheets are available on our website.

#### Ordering information

Unless otherwise indicated, all materials are available through **thermofisher.com**. MLS: Fisher Scientific (**fisherscientific.com**) or other major laboratory supplier.

Product	Description	Cat. No.	
GeneChip™ Human Genome U133A 2.0 Array	2 arrays	900471	
	6 arrays	900468	
	30 arrays	900469	
Supporting products			
GeneChip™ 3' IVT PLUS Reagent Kit	10 reactions	902415	
	30 reactions	902416	
GeneChip™ Hybridization, Wash, and Stain Kit <sup>[1]</sup>	30 reactions	900720	

<sup>[1]</sup> Each kit contains 1 Hybridization Module, 1 Stain Module, 3 bottles of Wash Buffer A, and 1 bottle of Wash Buffer B, sufficient for 30 reactions. Individual kit components may be ordered separately.



### Storage, handling, and stability

This cartridge array consists of a square glass substrate mounted in a plastic cartridge. 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 nonabrasive laboratory tissue.

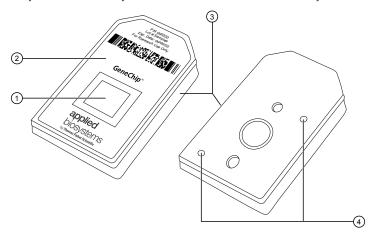


Fig. 1 Cartridge array.

- Probes on a glass substrate
- Plastic cartridge
- 3 Notch
- 4 Septa

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

# Customer and technical support

Visit **thermofisher.com/support** for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation, including:
  - User guides, manuals, and protocols
  - Certificates of Analysis
  - Safety Data Sheets (SDSs; also known as MSDSs)
    Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

### Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Affymetrix Pte Ltd | 7 Gul Circle #2M-01 | Keppel Logistics Building | Singapore 629563

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. 701501

Revision	Date	Description
6	17 November 2017	Update document to current template.
5	04 February 2009	Baseline for revision history.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are the property of their respective owners.

