



Package Insert

Axiom® Genome-Wide EUR 1 Array Plate

Intended Use

The Axiom® Genome-Wide EUR 1 Array Plate consists of array plates designed for the genotyping of single nucleotide polymorphisms (SNPs) and simple insertion/deletions (in/dels) in populations with European ancestry. This array plate, when used with the Axiom® 2.0 Reagent Kit (see **Ordering Information** for further details) enables researchers to perform large-scale genotyping studies with minimal hands-on processing per plate.

Each plate consists of 96 microarrays and is offered as a part of a complete automated solution from Affymetrix that includes instrumentation for sample preparation, and array washing, staining and scanning.

Each microarray on the Axiom Genome-Wide EUR 1 Array Plate enables the genotyping of up to 668,602 SNPs and 5,915 simple indels. SNPs on this array are chosen from the Axiom Genomic Database and were selected based on their MAF and their ability to tag European haplotypes. For detailed info on the recommended analysis workflow for this array, see the *Advanced Analysis Workflow for the Axiom Genome-Wide EUR 1, EAS 1, LAT 1 and AFR 1 Arrays Tech Note* (P/N DNA01006-1).

SNPs were chosen and represented on the array in part based upon genotyping reproducibility and call rate, with accuracy determined using genotypes from the International HapMap Project. These SNPs have been validated across multiple samples, including the original 270 HapMap Phase I samples as well as additional extended HapMap Phase 3 populations.

These markers are amplified using the Axiom 2.0 Reagent Kit and are present on short randomly generated fragments of genomic DNA (25 to 125 basepair) prior to array hybridization. The reagents for processing Axiom arrays are also included in this reagent kit.

Identical to the cartridge array manufacturing process, the oligonucleotide probes on Axiom Genome-Wide EUR 1 Array Plates are synthesized *in situ* using Affymetrix' photolithographic process.

Instructions for Use

Refer to the following documents for instructions on processing samples using the Axiom® 2.0 Assay.

- *Advanced Analysis Workflow for the Axiom® Genome-Wide EUR 1, EAS 1, LAT 1 and AFR 1 Arrays Tech Note*, P/N DNA01006-1
- *Axiom® 2.0 Assay Automated Workflow User Guide*, P/N 702963
- *Axiom® 2.0 Assay Automated Workflow Site Prep Guide*, P/N 702984
- *Axiom® 2.0 Assay Automated Target Prep Protocol QRC*, P/N 702962
- *Axiom® 2.0 Assay Manual Workflow User Guide*, P/N 702990
- *Axiom® 2.0 Assay Manual Workflow Site Prep Guide*, P/N 702991
- *Axiom® 2.0 Assay Manual Target Prep Protocol QRC*, P/N 702989
- *Axiom® 2.0 gDNA Sample Prep Protocol QRC*, P/N 702987
- *GeneTitan® MC Protocol for Axiom 2.0 Array Plate Processing QRC*, P/N 702988

Related Documents

The following documents contain information related to the instruments and software required to perform the Axiom 2.0 Genotyping Assay:

- *GeneTitan® Multichannel Instrument User's Manual*, P/N 08-0306
- *GeneTitan® Multichannel Instrument Site Preparation Guide*, P/N 08-0305
- *Affymetrix® GeneChip® Command Console® Software User Manual*, P/N 702569
- *Affymetrix® Genotyping Console™ 4.1 User Manual*, P/N 702982
- *Biomek® Liquid Handler User's Manual*, Beckman Coulter P/N 987834
- *Biomek® Software User's Manual*, Beckman Coulter P/N 987835

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 type. Additional information can be located under the specific array product on the Affymetrix web site at www.affymetrix.com/support/technical/libraryfilesmain.affx.

Reagents, Instrumentation and Software Required

1. Axiom® 2.0 Reagent Kit
2. Axiom® Genome-Wide EUR 1 Array Plate
3. Biomek FX[®] Target Prep Express by Beckman Coulter
4. GeneTitan® Multichannel Instrument
5. Affymetrix GeneChip® Command Console® Software
6. Affymetrix Genotyping Console

For a complete list of reagents and consumables required, please refer to the *Axiom® 2.0 Assay Automated Workflow Site Prep Guide*, P/N 702984, or the *Axiom® 2.0 Assay Manual Workflow Site Prep Guide*, P/N 702991.

Ordering Information

Product Name	Description	Part Number
Axiom® Genome-Wide EUR 1 96-Array Plate	One 96-Array Plate	901787
Axiom® GeneTitan® Consumables Kit ¹	Contains all GeneTitan consumables required to process an Axiom array plate	901606
Axiom® 2.0 Reagent Kit	One 96-Reaction Kit	901758

1. One Hyb Tray, one Scan Tray, and five Stain Trays with covers are included for use with each array plate. These consumables are required for processing Axiom array plates on the GeneTitan Multichannel Instrument.

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

Axiom Genome-Wide EUR 1 Array Plates should be stored at 2 to 8 °C and must not be frozen. Refer to the expiration date on the package label. Do not use arrays or reagents after the expiration date.

When Handling the Axiom Array Plate

Remove the Array Plate from the pouch with gloved hands. The plate is packaged with a blue plastic base. Do not remove the Array Plate from this protective base, or touch the Array Plate directly. Keep the Array Plate in the protective base at all times, including when placed on the GeneTitan Multichannel Instrument.

When Handling the 96 Plate Scan Tray

Remove the 96 Plate Scan Tray (Scan Tray) from the pouch with gloved hands. The Scan Tray is packaged with a black protective base and cover. Keep the Scan Tray in the protective base with the cover at all times prior to loading into the GeneTitan Multichannel Instrument. Do not touch the bottom of the Scan Tray directly.

The Scan Tray has protruding guide posts that may be sharp and can puncture the pouch if not handled carefully. Take the necessary precautions to avoid injury.

Precautions

1. AXIOM GENOME-WIDE EUR 1 ARRAY PLATES ARE FOR RESEARCH USE ONLY. NOT FOR USE IN 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 plates.
7. CAUTION: Use care when handling the Scan Tray as it has protruding guiding posts that may be sharp and can stick out of the pouch if not handled carefully.

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-sublicensable 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.

Patents

Products may be covered by one or more of the following patents: U.S. Patent Nos. 5,445,934; 5,744,305; 6,399,365; 6,733,977; 7,790,389 and other U.S. or foreign patents. Products are manufactured and sold under license from OGT under 5,700,637.

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.

ORIX Hamamatsucho Bldg, 7F
1-24-8 Hamamatsucho, Minato-ku
Tokyo 105-0013, Japan
Tel: +81-3-6430-4020
Fax: +81-3-6430-4021
E-mail: supportjapan@affymetrix.com

Please visit our web site for international distributor contact information
www.affymetrix.com

Trademarks

Affymetrix®, Axiom®, Command Console®, CytoScan™, DMET™, GeneAtlas®, GeneChip®, GeneChip-compatible™, GeneTitan®, Genotyping Console™, myDesign™, NetAffx®, OncoScan™, Powered by Affymetrix™, PrimeView™, Procarta®, and QuantiGene® are trademarks or registered trademarks of Affymetrix, Inc. All other trademarks are the property of their respective owners.

Copyright

© 2011 Affymetrix, Inc. All rights reserved.