

# Package Insert

## OncoScan® CNV Array

**For Research Use Only**  
**Not for use in diagnostic procedures.**

### Introduction

The OncoScan® CNV FFPE Assay Kit includes the OncoScan® CNV Arrays and the OncoScan® CNV Reagent Kit. The OncoScan CNV FFPE Assay Kit detects genome wide copy number gains and losses, as well as loss of heterozygosity (LOH) including copy neutral LOH using limited DNA from Formalin-Fixed Paraffin-Embedded (FFPE) samples of various ages. The OncoScan CNV Array contains 328,000 tags, in triplicate, that hybridize to amplified DNA target generated during the assay.

OncoScan® CNV FFPE Assay Kit is for research use only and not intended for use in diagnosis of disease. Please refer to [www.affymetrix.com](http://www.affymetrix.com) for a list of supporting manuals for assay and data analysis procedures.

### Instrumentation and Software Required

1. GeneChip® Scanner 3000 7G
2. GeneChip® Fluidics Station 450
3. GeneChip® Hybridization Oven 645
4. GeneChip® Command Console® Software (AGCC)
5. OncoScan® Console Software (analysis) (AGCC)

### Ordering Information

P/N	Product Name	Description
902695	OncoScan® CNV FFPE Assay Kit Array and Reagent Kit Bundle	Sufficient for 24 samples and includes 48 OncoScan® CNV Arrays
902693	OncoScan® CNV Training Kit Array and Reagent Kit Bundle	Sufficient for 18 samples and includes 36 OncoScan® CNV Arrays

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](http://www.affymetrix.com/site/contact/index.affx).

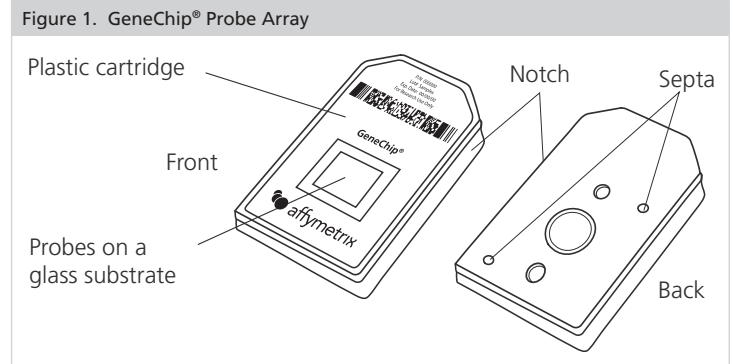
### Precautions

1. ONCOSCAN CNV 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.

### 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 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® 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

Affymetrix hereby grants to buyer a non-exclusive, non-transferable, non-sublicensable license to Affymetrix' Core Product IP to use the product(s), but only in accordance with the product labels, inserts, manuals and written instructions provided by Affymetrix. "Core Product IP" is the intellectual property owned or controlled by Affymetrix as of the shipment date of a product that covers one or more features of the product that are applicable in all applications of the product that are in accordance with the product labels, inserts, manuals and written instructions provided by Affymetrix. The license granted herein to buyer to the Core Product IP expressly excludes any use that: (i) is not in accordance with the product labels, inserts, manuals and written instructions provided by Affymetrix, (ii) requires a license to intellectual property that covers one or more features of a product that are only applicable within particular fields of use or specific applications, (iii) involves reverse engineering, disassembly, or unauthorized analysis of the product and/or its methods of use, or (iv) involves the re-use of a consumable product. Buyer understands and agrees that except as expressly set forth, no right or license to any patent or other intellectual property owned or controlled by Affymetrix is granted upon purchase of any product, whether by implication, estoppel or otherwise. In particular, no right or license is conveyed or implied to use any product provided hereunder in combination with a product or service not provided, licensed or specifically recommended by Affymetrix for such use. Furthermore, buyer understands and agrees that buyer is solely responsible for determining whether buyer possesses all intellectual property rights that may be necessary for buyer's specific use of the product, including any rights from third parties.

---

## Contact Information

### Affymetrix, Inc.

3420 Central Expressway  
Santa Clara, CA 95051 USA  
E-mail: [support@affymetrix.com](mailto: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](mailto: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](mailto:supportjapan@affymetrix.com)

Please visit our web site for international distributor contact information  
[www.affymetrix.com](http://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; 6,399,365; 6,551,817; 6,733,977; 7,629,164; 7,790,389 and D430,024 and other U.S. or foreign patents. Products are manufactured and sold under license from OGT under 5,700,637.

## Trademarks

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

## Copyright

©2015 Affymetrix, Inc. All rights reserved.