

## Affymetrix GeneChip® SNP Kit



### INTENDED USE

The GeneChip® SNP Kit is intended for research use only for SNP genotyping using the Affymetrix GeneChip® Scanner 3000 Targeted Genotyping System. The kit and MIP Assay Protocol are based on patented proprietary Molecular Inversion Probe (MIP) technology. This technology enables users to genotype tens of thousands of SNPs in a single assay. It eliminates the need to amplify the DNA segment around each SNP prior to analysis by using circularizing intra-molecular interactions. The use of a single probe per SNP increases the accuracy of the assay by ensuring that the test only reports information about a known target sequence among the billions of possible sequences found in the human genome.

The contents of this kit are sufficient to process 24 reactions. This insert describes the kit contents and storage conditions.

### REAGENT AND OTHER KIT COMPONENTS

Important! Do not thaw and refreeze reagents. Use immediately once thawed.

#### Affymetrix GeneChip® SNP Kit

<b>Assay Panel</b> (Store at -20°C)	Assay Panel 160 µL Kit Control DNA 70 µL
<b>Pre-Amp Kit</b> (Store at -20 °C)	Enzyme A 60 µL Buffer A 200 µL Water 2 mL Gap Fill Mix 130 µL Exo Mix 530 µL Cleavage Tube 3000 µL Cleavage Enzyme 30 µL Amp Mix 3000 µL dNTP Mix 1 plate; 20 µL/well
<b>Stain Cocktail</b> (Store at 2-8 °C even if frozen upon arrival).	Stain Cocktail 5 mL
<b>Post-Amp Kit</b> (Store at -20 °C)	HY A Allele Tube 950 µL HY C Allele Tube 950 µL HY G Allele Tube 950 µL HYT Allele Tube 950 µL HY Digest Mix 260 µL Hyb Cocktail 900 µL Buffer H 400 µL
<b>Wash Solution Kit</b> (Store at room temperature.)	Wash Soln A 500 mL Wash Soln B 500 mL

### INSTRUCTIONS FOR USE

Please refer to the Affymetrix GeneChip® Scanner 3000 Targeted Genotyping System User Guide for procedures on processing genomic DNA samples and controls.

---

**STORAGE, HANDLING, AND STABILITY**

Do not thaw and refreeze reagents. Use immediately once thawed. Store as directed on the box. Reagents are stable until the expiration date indicated.

---

**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 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

**Affymetrix UK Ltd**

Voyager, Mercury Park,  
Wycombe Lane, Wooburn Green,  
High Wycombe HP10 0HH  
United Kingdom  
supporteurope@affymetrix.com  
UK and Others Tel:  
+44 (0) 1628 552550  
France Tel: 0800919505  
Germany Tel: 01803001334  
Fax: +44 (0) 1628 552585

Please visit our web site for international distributor contact information.

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


---

**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.

---

**TRADEMARKS**

Affymetrix®, GeneChip®, ®, HuSNP®, GenFlex®, NetAffx™, Flying Objective™, CustomExpress®, CustomSeq®, Tools to Take You As Far As Your Vision®, The Way Ahead™, Powered by Affymetrix™, and GeneChip-compatible™, are trademarks of Affymetrix, Inc.

Qdot is a trademark of Quantum Dot Corporations.

---

**PATENTS**

Reagent Products and their use may be covered by U.S. Patent Nos. 5,871,921, 5,866,337, and 6,858,412 and other U.S. or foreign patents that are owned or licensed by Affymetrix.

---

**COPYRIGHT**

©2005 Affymetrix, Inc. All rights reserved.

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