



## GeneTitan® Wash Buffer A and B Module

### Intended Use

This package insert provides information for the GeneTitan® Wash Buffer A and B Module.

### Introduction

During the laboratory procedure for various target labeling assays, biotin-labeled fragments (referred to as the "target") are generated. After target preparation, arrays are processed in preparation for scanning. A hybridization cocktail is prepared including the labeled, fragmented target and hybridization controls. The target is then hybridized to the GeneChip array in array plate format during a 16-hour incubation. Immediately following hybridization, the array is washed and stained with streptavidin phycoerythrin conjugate using an automated protocol, followed by scanning on a GeneTitan Instrument.

Wash Buffer A and Wash Buffer B are required for the processing of array plates on the GeneTitan Instrument and the GeneTitan Wash Buffer A and B Module is sufficient for processing one array plate.

The GeneTitan Hybridization, Wash, and Stain Kit reagents are formulated into the fewest number of individual components possible, reducing preparation and handling steps. Stain Cocktail 1 & 3, Stain Cocktail 2, Array Holding Buffer and Wash Buffers A and B are ready-to-use solutions. The hybridization cocktail is prepared by combining the Hybridization Mix, and Nuclease-free Water with the target prepared by the user during the specific target labeling assay.

**NOTE:** Hybridization controls are required and must be purchased separately. The GeneChip® Hybridization Control Kit includes the Control Oligonucleotide B2. The Control Oligonucleotide B2 is also available for individual purchase.

### Kit Components

Component	Volume	Storage
Wash Buffer A	1,045 mL	2°C to 8°C
Wash Buffer B	450 mL	2°C to 8°C

### Shipping and Storage

The GeneTitan Wash Buffers A and B Module will be sent at ambient temperature and should be stored at 2° to 8°C.

### Functional Testing

Each lot of reagents is rigorously tested for performance.

### Safety Information

A Material Safety Data Sheet(s) (MSDS) is available at [www.affymetrix.com](http://www.affymetrix.com). If the product is a kit or is supplied with more than one material, please refer to the MSDS for each component for hazard information.

### Ordering Information

P/N	Product Name	Description/Size
901583	GeneTitan® Wash Buffers A and B Module	Wash Buffer A = 1,045 mL Wash Buffer B = 450 mL
<b>Related Products</b>		
901530	GeneTitan® Hybridization, Wash, and Stain Kit for 3' IVT Arrays	96 Reactions
901622	GeneTitan® Hybridization, Wash, and Stain Kit for WT Arrays	96 Reactions
900454	GeneChip® Hybridization Control Kit	30 Reactions
900457		150 Reactions
900301	GeneChip® Control Oligo B2, 3 nM	30 Reactions

Additional Affymetrix® reagents are available for purchase separately. Please refer to [www.affymetrix.com](http://www.affymetrix.com) for related product ordering information.

Affymetrix products can be purchased directly from Affymetrix in the United States, many European countries, and many 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).

### 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  
E-mail: [supportjapan@affymetrix.com](mailto:supportjapan@affymetrix.com)  
Tel: +81-3-6430-4020  
Fax: +81-3-6430-4021

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

### Copyright

© 2010 Affymetrix, Inc. All rights reserved.