



# GenetiSure Dx DNA Labeling Kit

## K1201-64100



**Intended Use** The GenetiSure Dx DNA Labeling Kit is for use in the labeling of samples to be hybridized to oligonucleotide microarrays in laboratory defined, microarray-based, diagnostic assay workflows. The kit is intended for use as part of molecular diagnostic test systems.

### Rx Only

**Summary** The kit contains Random Primers and the Exo (-) Klenow fragment to differentially label genomic DNA (gDNA) samples with fluorescently labeled nucleotides for the purpose of labeling samples prior to microarray hybridization.

### Materials Provided & Storage

Component	Part Number	Quantity	Storage Temperature
Exo (-) Klenow	5190-7306	55 µL	-25° to -15°C
5X gDNA Reaction Buffer	5190-7310	550 µL	-25° to -15°C
Nuclease Free Water	5190-7311	1.5 mL	-25° to -15°C
Alu I Restriction Enzyme	5190-7312	28 µL	-25° to -15°C
Rsa I Restriction Enzyme	5190-7313	28 µL	-25° to -15°C
10X Restriction Enzyme Buffer	5190-7314	142 µL	-25° to -15°C
BSA	5190-7315	15 µL	-25° to -15°C
10X dNTP Mix	5190-7316	265 µL	-25° to -15°C
Cyanine 3-dUTP	5190-7308	78 µL	-25° to -15°C
Cyanine 5-dUTP	5190-7309	78 µL	-25° to -15°C
Random Primers	5190-7307	265 µL	-25° to -15°C
Human Reference DNA Female	5190-7317	125 µL; 0.2 µg/µL	2° to 8°C
Human Reference DNA Male	5190-7318	125 µL; 0.2 µg/µL	2° to 8°C
Purification Columns	—	50	15° to 30°C

Store the kit components at the recommended storage temperature until the expiration date.

Cyanine 3-dUTP and Cyanine 5-dUTP are light sensitive and are subject to degradation by multiple freeze-thaw cycles. To avoid prolonged exposure to light, keep the tubes of Cyanine 3-dUTP and Cyanine 5-dUTP covered with foil or a darkened lid.

### Warnings and Precautions

- 1 For In Vitro Diagnostic Use
- 2 Avoid microbial contamination, which may cause erroneous results.
- 3 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. Refer to the appropriate blood borne pathogen precautions indicated by your facility or local regulations. Never pipet by mouth. Avoid specimen contact with skin and mucous membranes.
- 4 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 kit materials.
- 7 This product is stable until the expiration date when stored at the recommended storage temperature.
- 8 Performance of the kit has been shown to be unaffected for up to 8 freeze-thaw cycles.
- 9 The clinical interpretation of any test results should be evaluated within the context of the patient's medical history and other diagnostic laboratory test results by qualified personnel.

**10** Don't use products past their expiration dates.

**11** Proper storage and handling of reagents and samples are essential for the performance. All laboratory equipment used to prepare the target during this procedure should be calibrated and maintained to ensure accuracy. Incorrect measurement of reagents may affect the outcome of the procedure.

**Indications of  
Instability or  
Deterioration**

This product contains components that ship at different temperature conditions (room temperature and dry ice). Product may arrive as separate packages or in a single box that contains different shipping areas. Inspect packages upon arrival. If the tamper-evident label is opened at the perforations, or if the dry ice used for shipping has evaporated, do not use the contents of the package. For customer service or technical support, please contact Agilent Technical Support.

**Safety Information**

A Material Safety Data Sheet(s) (MSDS) is available at [www.agilent.com](http://www.agilent.com).

**Procedure**

Refer to the GenetiSure Dx Postnatal Assay for instructions for use of this product and limitations of the procedure.

**Limitations of the  
Procedure**

Proper storage and handling of reagents and samples are essential for the performance. All laboratory equipment used to prepare the target during this procedure should be calibrated and maintained to ensure accuracy. Incorrect measurement of reagents may affect the outcome of the procedure.



**Technical Support** Visit [www.agilent.com/en/contact-us/page](http://www.agilent.com/en/contact-us/page) to find country-specific phone numbers, or send an email to one of the email addresses below.
















**United States:** [clinical.support@agilent.com](mailto:clinical.support@agilent.com)

**Europe:** [genomics\\_dx\\_tech\\_europe@agilent.com](mailto:genomics_dx_tech_europe@agilent.com)

**Japan:** [email\\_japan@agilent.com](mailto:email_japan@agilent.com)

Any serious incident that has occurred in relation to the device shall be reported to the manufacturer and the competent authority of the country in which the user and/or the patient is established.

**Symbol Table**

 Catalog / code number	 In Vitro Diagnostic Medical Device
 Temperature limit	 Manufacturer
 Consult Instructions for Use	 Use-by date
 Contains sufficient for <n> tests	 Caution
 Unique Device Identifier	 Batch code
 Do not reuse	 Humidity limitation
 European Conformity	 Reconstitute with
 Authorized representative in the European Community	

 Agilent Technologies, Inc.  
5301 Stevens Creek Blvd, Santa Clara, CA 95051, USA  
Manufactured at:  
1834 State Hwy 71 W, Cedar Creek, TX 78612, USA  
[www.agilent.com](http://www.agilent.com)

 Agilent Technologies Denmark ApS  
Produktionsvej 42  
2600 Glostrup, Denmark



© Agilent Technologies, Inc. 2016–2017, 2022

Revision C0, January 2022



K1201-90003