



GenetiSure Dx Wash Buffer Set

K1201-64300



Intended Use The GenetiSure Dx Wash Buffer 1, 4L is intended for use in the washing of oligonucleotide microarrays in laboratory defined, microarray-based, diagnostic assay workflows. The wash buffer is intended for use as part of molecular diagnostic test systems.

The GenetiSure Dx Wash Buffer 2, 4L is intended for use in the washing of oligonucleotide microarrays in laboratory defined, microarray-based, diagnostic assay workflows. The wash buffer is intended for use as part of molecular diagnostic test systems.

Rx Only

Summary The kit contains wash buffers for washing unhybridized labeled DNA from microarrays.

Materials Provided & Storage

Component	Part Number	Quantity	Storage Temperature
GenetiSure Dx Wash Buffer 1	K1201-64305	8 liters (2 boxes)	15° to 30°C
GenetiSure Dx Wash Buffer 2	K1201-64310	4 liters (1 box)	15° to 30°C

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

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 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.
- 9 Don't use products past their expiration dates.
- 10 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

Inspect packages upon arrival. If the tamper-evident label is opened at the perforations, 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.

Procedure

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



Technical Support

Visit 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


Europe: genomics_dx_tech_europe@agilent.com

Japan: 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

 REF	Catalog / code number	 IVD	In Vitro Diagnostic Medical Device
	Temperature limit		Manufacturer
	Consult Instructions for Use		Use-by date
	Contains sufficient for <n> tests		Caution
 UDI	Unique Device Identifier	 LOT	Batch code
	Do not reuse		Humidity limitation
	European Conformity	 RCNS	Reconstitute with
 EC REP	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

 EC REP Agilent Technologies Denmark ApS
 Produktionsvej 42
 2600 Glostrup, Denmark

© Agilent Technologies, Inc. 2016–2017, 2022



Revision C0, January 2022



K1201-90005