

# Genome-Wide Human SNP Nsp/Sty Assay Kit 5.0/6.0

Catalog Numbers 901013, 901152, and 901015

Pub. No. 702508 Rev. 3

**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](http://thermofisher.com/support).

## Product use

The Applied Biosystems™ Genome-Wide Human SNP Nsp/Sty Assay Kit 5.0/6.0 enables genotyping and DNA copy number analysis on the Applied Biosystems™ Genome-Wide Human SNP Array 5.0 and Genome-Wide Human SNP Nsp/Sty 6.0 Array.

## Kit components

Components and storage
<b>Box 1 of 3</b> (store in a room free of DNA templates) Adaptor, Nsp (50 µM) PCR Primer, 002 (100 µM)
<b>Box 2 of 3</b> (store in a room free of DNA templates) Adaptor, Sty (50 µM) PCR Primer, 002 (100 µM)
<b>Box 3 of 3</b> (can store in a room with PCR amplicons) GeneChip™ Fragmentation Reagent (see tube label for unit concentration) 10X Fragmentation Buffer GeneChip™ DNA Labeling Reagent (30 mM) Terminal Deoxynucleotidyl Transferase (30 U/µL) 5X Terminal Deoxynucleotidyl Transferase Buffer Oligonucleotide Control Reagent 0100
<b>Box 1 of 1</b> (store in a room free of PCR amplicons) Reference Genomic DNA, 103 (50 ng/µL)

## Instructions for use

See the Applied Biosystems™ *Genome-Wide Human SNP Nsp/Sty 6.0 Assay User Guide* (Pub. No. 702504), or the Applied Biosystems™ *Cytogenetics Copy Number Assay User Guide* (Pub. No. 702607) for procedures on DNA target and oligonucleotide control preparation, target hybridization, fluidics setup, array scanning, and data analysis instructions.

## Ordering information

Cat. No.	Product name	Description
901013	Genome-Wide Human SNP Nsp/Sty Assay Kit 5.0/6.0	30 reactions
901152	Genome-Wide Human SNP Nsp/Sty Assay Kit 5.0/6.0	50 reactions
901015	Genome-Wide Human SNP Nsp/Sty Assay Kit 5.0/6.0	100 reactions
Supporting products		
901182	Genome-Wide Human SNP Nsp/Sty 6.0 Array	5 arrays
901153	Genome-Wide Human SNP Nsp/Sty 6.0 Array	50 arrays
901150	Genome-Wide Human SNP Nsp/Sty 6.0 Array	100 arrays
901012	Reference Genomic DNA, 103	2 reactions

## Storage, handling, and stability

The reagents that are contained in the Genome-Wide Human SNP Nsp/Sty Assay Kit 5.0/6.0 are shipped on dry ice and must be stored at -20°C. See the expiration date on the package label. Do not use reagent components after the expiration date.

The GeneChip™ Fragmentation Reagent and the Terminal Deoxynucleotidyl Transferase contain active enzymes. The following precautions are recommended to ensure optimal activity:

- Store at -20°C (do not store at -80°C).
- Minimize repeated exposure to temperatures >0°C.
- Use a benchtop cooler when handling the reagent.

**IMPORTANT!** Store the unopened Genome-Wide Human SNP Nsp/Sty Assay Kit 5.0/6.0 in a room that is free of PCR amplicons. The reagent components are separated into four boxes to support the recommended workflow, and minimize the chance of contamination during use. Each of these boxes has different recommended storage locations: Box 1 of 3—store in a room free of DNA templates. Box 2 of 3—store in a room free of DNA templates. Box 3 of 3—can store in a room with PCR amplicons. Box 1 of 1 (Reference Genomic DNA)—store in a room free of PCR amplicons.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and

Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 | Vilnius, Lithuania

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Revision history:** Pub. No. 702508

Revision	Date	Description
3	10 August 2017	Updated document to current template.
2	07 Jan 2008	Baseline for revision history.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.