

CytoScan™ 750K Accel Array

Catalog Numbers 952534

Pub. No. MAN1000007 Rev. A00

Product use

Chromosomal Microarray Analysis (CMA) has been adopted as a leading choice for the study of chromosomal aberrations and their impact on human disease. The Applied Biosystems™ CytoScan™ 750K Accel Assay and the CytoScan™ 750K Accel Array enables the detection of high-resolution copy number changes across the genome and provides allelic imbalance information from single nucleotide polymorphisms (SNPs).

The CytoScan™ 750K Accel Array is part of the CytoScan™ 750K Accel Cytogenetics Solution which includes the CytoScan™ Accel Reagent Kit, GeneChip™ Scanner 3000 instrument, and the Chromosome Analysis Suite software (ChAS).

The CytoScan™ 750K Accel Array is used in conjunction with the CytoScan™ Accel Reagent Kit (Cat. No. 952459) and the CytoScan™ Amplification Kit (Cat. No. 902975).

Instructions for use

See the *CytoScan™ 750K Accel Assay Manual Workflow User Guide* (Pub. No. [MAN1000001](#)) for procedures on DNA target preparation, target hybridization, fluidics setup, array scanning, and data analysis instructions.

Fluidics scripts

The fluidics script used depends on the array type, labeling protocol, and reagents used for cartridge processing. Refer to the GeneChip Fluidics Station Scripts support page to determine which fluidics script is appropriate for your application. The fluidics scripts can be downloaded from our website.

Library files

Library files contain information about the probe array design characteristics, probe use and content, and scanning and analysis parameters. These files are unique for each probe array. Additional information can be located under the specific array product on our website.

Instrumentation and software required

- GeneChip™ Scanner 3000 7G
- GeneChip™ Fluidics Station 450
- GeneChip™ Hybridization Oven 645
- GeneChip™ Data Collection Software (GCDC)
- Chromosome Analysis Suite (ChAS) software, version 4.3 or later

Critical specifications

Item	Specification
Array format	100
Feature size	5 µm
Hybridization volume	130 µL
Fluidics station protocol	CytoScan750K Accel_450
Library files	CytoScan 750K Accel Array
ChAS software	ChAS 4.3 and above
ChAS QC metric: MAPD	≤0.25
ChAS QC metric: SNPQC	≥15.0
ChAS QC metric: wSD	≤0.12

Ordering information

Unless otherwise indicated, all materials are available through [thermofisher.com](#).

Product name	Description	Cat. No.
CytoScan™ 750K Accel Array	6-pack	952534

Storage, handling, and stability

This cartridge array consists of a square glass substrate mounted in a plastic cartridge. The glass contains an array of oligonucleotides that, when mounted, is on the inner glass surface. A chamber in the plastic housing directly under the glass acts as a reservoir where hybridization and washing occur.

Although the inner glass surface of the probe array is protected, any contamination or scratches on the outer surface of the glass can compromise the accuracy of the scan. Avoid touching the surface of the glass with your fingers. Skin oils and other substances, such as lotions or ink, can fluoresce. If the surface of the glass is noticeably dirty, it can be carefully cleaned with a nonabrasive laboratory tissue.

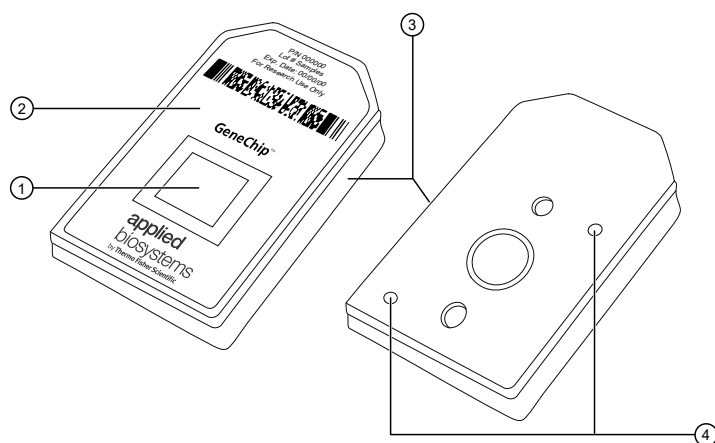


Figure 1 Cartridge array.

- | | |
|-------------------------------|---------|
| ① Probes on a glass substrate | ③ Notch |
| ② Plastic cartridge | ④ Septa |

The cartridge array should be stored at 2–8°C. Refer to the expiration date on the package label. Do not use arrays or reagents after the expiration date.

Customer and technical support

Visit [thermofisher.com/support](https://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.



Affymetrix Pte Ltd | 7 Gul Circle #2M-01 | Keppel Logistics Building | Singapore 629563

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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.

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.

©2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.