

CarrierScan™ Array Plates

Catalog Numbers 931930 and 951950

Doc. Part No. 703468 Pub. No. MAN0017674 Rev. A.0

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.

Product use

Applied Biosystems™ CarrierScan™ Array Plates are designed to enable the detection of both sequence and structural variants simultaneously. These variants including biallelic and multiallelic mutations such as single nucleotide variants (SNVs), insertion-deletion variants (indels), as well as structural genomic variants, such as small intragenic deletions and duplications (copy number variants) that are important in the field of carrier screening research. The array plates are available in 96-array format, and when used with the CarrierScan™ Reagent Kit 96 Reactions, they offer a unique opportunity to combine the number of assays and technologies that are required to perform expanded carrier screening research into one, to help save time and costs.

Array plates are run on the GeneTitan™ Multi-Channel Instrument for array washing, staining, and scanning.

The oligonucleotide probes on CarrierScan™ Array Plates are synthesized *in situ* using Thermo Fisher Scientific's photolithographic process, which avoids batch drop-outs that can cause relevant markers to be omitted.

Instructions for use

See the following documents for instructions on processing samples using the CarrierScan™ Assay.

- CarrierScan™ Assay 96-Array Format Manual Workflow User Guide (Pub. No. 703481, QR Pub. No. 703482)
- CarrierScan™ Assay 96-Array Format Automated Workflow on the Biomek™ FX^P User Guide (Pub. No. 703478, QR Pub. No. 703479)
- Axiom™ 2.0 gDNA Sample Preparation Quick Reference (Pub. No. 702987)
- GeneTitan™ MC Protocol for Axiom™ Array Plate Processing Quick Reference (Pub. No. 702988)

Library files

Library files contain information about the array design characteristics, probe use and content, and scanning and analysis parameters. These files are unique for each array. GeneTitan™ library files provide instructions for the fluidics and scanning activities on the GeneTitan™ MC Instrument. Analysis library files and annotation files are required for data analysis. Contact your field application scientist to get the appropriate library files for your array. The library files can also be found under the specific array product on our website.

Reagents, instruments, and software required

- CarrierScan™ Assay Kit (includes reagents, CarrierScan™ 96F Array Plate, and consumables) or CarrierScan™ 1S Assay Kit (includes reagents, CarrierScan™ 1S Array Plate, and consumables)
- GeneTitan™ Multi-Channel Instrument (GTMC)
- GeneChip™ Command Console™ software
- Axiom™ Analysis Suite software
- CarrierScan™ Reporter software

For a complete list of reagents and consumables that are required, see the appropriate site preparation guide.

Ordering information

Unless otherwise indicated, all materials are available through thermofisher.com.

Product name	Description	Cat. No.
CarrierScan™ Assay Kit	Includes reagents, an array plate, and GTMC consumables sufficient for 96 samples.	931931
CarrierScan™ 1S Assay Kit	Includes reagents, an array plate, and GTMC consumables sufficient for 96 samples.	951951

Storage, handling, and stability

The array plates should be stored at 2–8°C and must not be frozen.

The array plates must be protected at all times from damage or exposure to dust. Refer to the expiration date on the package label. Do not use array plates or reagents after the expiration date.

When handling the 96-array plate

Remove the array plate from the pouch with gloved hands. The array plate is packaged with a blue plastic base. Do not remove the protective blue plastic base from the array plate or touch the array plate directly. Keep the array plate in the protective base at all times, including when placed on the GeneTitan™ MC Instrument.

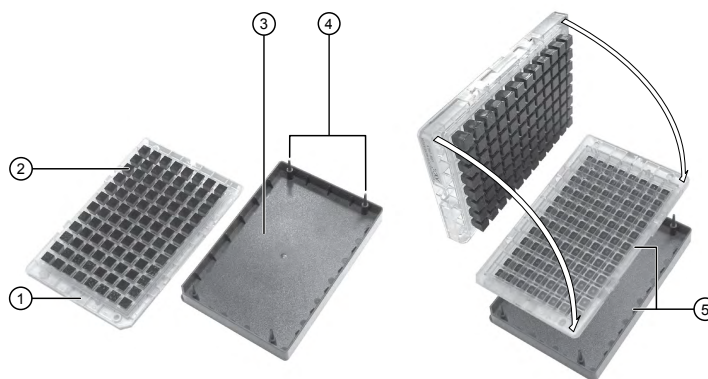


Fig. 1 Array plate assembly.

- Array plate
- Array on glass substrate that is mounted on a peg
- Blue plastic base
- Alignment pins
- Array plate and blue base assembly

Note: Displayed action is for demonstration purposes only.

When handling the 96-plate scan tray

Remove the scan tray from the pouch with gloved hands. The scan tray is packaged with a black plastic base. Do not remove the protective black plastic base from the scan tray or touch the scan tray directly. This protective base should stay with the scan tray at all times prior to loading into the GeneTitan™ MC Instrument.

CAUTION! The scan tray has protruding guiding posts that may be sharp and can stick out of the pouch if not handled carefully; therefore, take precaution to prevent unnecessary injury.

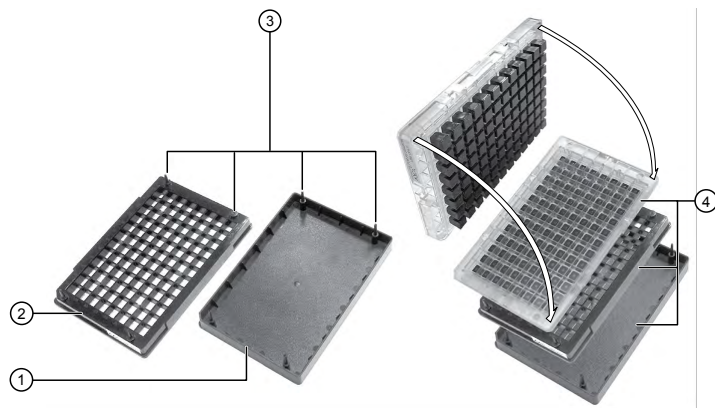


Fig. 2 Scan tray assembly.

- ① Black plastic base
- ② Scan tray
- ③ Alignment pins
- ④ Array plate, scan tray, and black plastic base assembly

Note: Displayed action is for demonstration purposes only. All movement of the array plate is performed during the fluidics protocol on the GeneTitan™ Instrument.

Manufacturer: 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, 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. MAN0017674

Revision	Date	Description
A.0	13 April 2018	Initial release in Thermo Fisher Scientific document control system. Supersedes legacy Affymetrix publication number 703468. Update document to add CarrierScan™ 1S product information.

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.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.