

Whole Transcript Array Plates

Pub. No. MAN0019707 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](https://www.thermofisher.com/support).

Product description

The Applied Biosystems™ Whole Transcript Array Plates offer the most advanced and cost-effective gene expression profiling option for whole-transcript coverage. Probes are distributed across the full length of the gene, providing a more complete and accurate picture of overall gene expression.

Discovery content, such as transcript regions supported by more speculative sources including Expressed Sequence Tags (ESTs) and gene predictions, are not interrogated by the whole transcript arrays.

Oligonucleotide probes are synthesized *in situ* complementary to each corresponding sequence. For expression analysis, the samples need to be amplified, and the targets labeled in the sense orientation for hybridization to the arrays.

Visit our website for a complete list of available whole transcript arrays, supporting documentation, available reagent kits, and procedures regarding target preparation, target hybridization, washing, staining, and scanning.

Assay documentation

The following documentation is required for using Whole Transcript Array Plates:

- *GeneChip™ WT PLUS Reagent Kit User Guide* (Pub. No. MAN0018137)
- *GeneChip™ WT Pico Reagent Kit User Guide* (Pub. No. MAN0017893)
- *GeneChip™ 3' IVT Pico Kit Manual Workflow User Guide* (Pub. No. 703308)
- *GeneTitan™ Instrument User Guide for Expression Array Plates* (Pub. No. MAN0017794)

Instrumentation and software required

1. GeneTitan™ Multi-Channel (MC) Instrument
2. GeneChip™ Command Console™ (GCC) software
3. Expression Console™ software
4. Transcriptome Analysis Console software

For a complete list of reagents and consumables required, as well as a list of supporting documentation for array plates, refer to our website.

Reagents required

The following table lists the kits required to process whole transcript array plates.

Product name	Description	Cat. No.
GeneTitan™ Hybridization, Wash, and Stain Kit for WT Array Plates	96 reactions	901622

Accessory files

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.

Storage, handling, and stability

Store array plates at 2–8°C and do not allow to freeze. Do not open the packaging until ready for use—this is essential to protect the array plates from damage, exposure to dust, and excess light. 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.

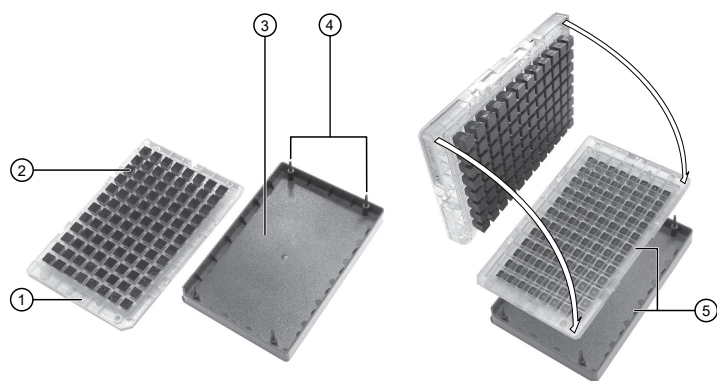


Figure 1 Array plate assembly.

- ① Array plate
- ② Array on glass substrate that is mounted on a peg
- ③ Blue plastic base
- ④ Alignment pins
- ⑤ 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.

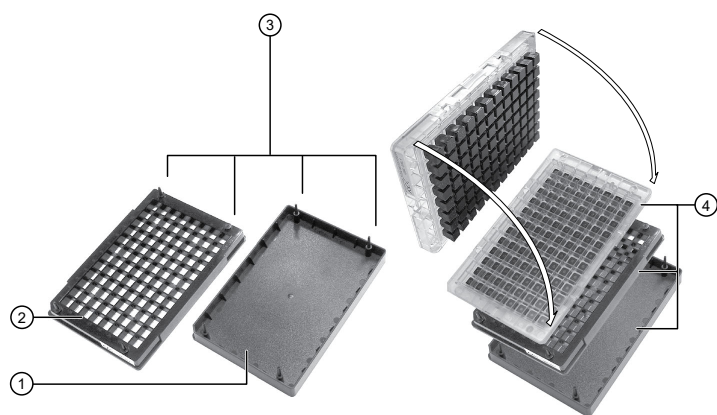


Figure 2 Scan tray assembly.

- 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™ MC Instrument.

Customer and technical support

Visit 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/or its affiliate(s) 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 any questions, please contact Life Technologies at www.thermofisher.com/support.



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.

Revision history: Pub. No. MAN0019707

Revision	Date	Description
A.0	21 December 2020	Initial release in Thermo Fisher Scientific document control system. Supersedes legacy Affymetrix publication number 703078. Updated to add a publication to the list of assay documents, and to correct a typo.

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.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are the property of their respective owners.