

# Applied Biosystems™ miRNA Array Plates

Pub. No. 703173 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™ miRNA Array Plates contain probe sets designed to interrogate small noncoding RNA transcripts. Information on the content including types and data sources used to design the array can be obtained from our website. All probe sets are PM (perfect match only).

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

## Instructions for use

Refer to the following guides for instructions on using the Applied Biosystems™ miRNA Array Plates:

- FlashTag™ Biotin HSR RNA Labeling Kit User Guide (Pub. No. 703095)
- GeneTitan™ Instrument User Guide for Expression Array Plates (Pub. No. 702933)

## 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

- GeneTitan™ Multi-Channel Instrument
- GeneChip™ Command Console™ (GCC) software
- Expression Console™ (EC) software
- Transcriptome Analysis Console (TAC) software

For a complete list of reagents and consumables required, refer to our website for a list of supporting manuals for miRNA array plates.

## Reagents required

Unless otherwise indicated, all materials are available through [thermofisher.com](http://thermofisher.com). MLS: Fisher Scientific ([fisherscientific.com](http://fisherscientific.com)) or other major laboratory supplier.

Product name	Description	Cat. No.
FlashTag™ Biotin HSR RNA Labeling Kit	10 reactions	901910
	30 reactions	901911
GeneTitan™ Hybridization, Wash, and Stain Kit for miRNA Array Plates <sup>[1]</sup>	96 reactions	902276
GeneTitan™ Hybridization Module for miRNA Plates	96 reactions	902275
Array Holding Buffer (110 mL)	96 reactions	901582
GeneTitan™ Wash Buffers A and B Module	96 reactions	901583

<sup>[1]</sup> Each kit contains 1 Hybridization Module, Array Holding Buffer, and GeneTitan™ Wash Buffers A and B Module, sufficient for 96-reactions. Individual kit components may be ordered separately.

## 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 all times, including when placed on the GeneTitan™ MC Instrument.

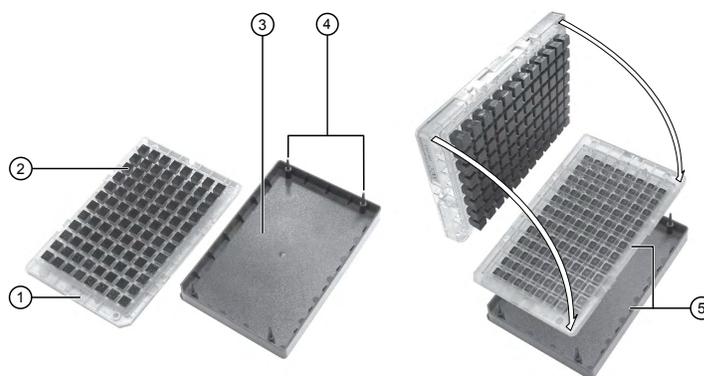


Fig. 1 Array plate assembly.

- Array plate
- Probe array on glass substrate mounted on 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.

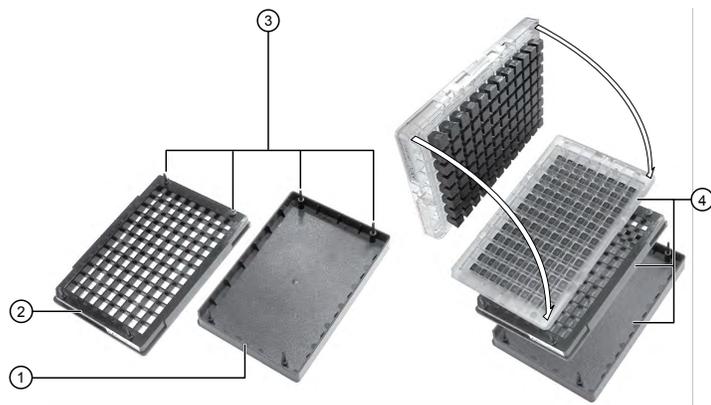


Fig. 2 Scan tray assembly.

- 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.

## Customer and technical support

Visit [thermofisher.com/support](http://thermofisher.com/support) for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation, including:
  - 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 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).



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. 703173

Revision	Date	Description
3	10 November 2017	Update document to current template.
2	December 2013	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. All other trademarks are the property of their respective owners.