

Axiom™ Array Plate

Pub. No. MAN0017712 Rev. B00



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 use

The Applied Biosystems™ Axiom™ Array Plate is designed for the genotyping of single nucleotide polymorphisms (SNPs) and simple insertion/deletions (indels). This array plate, when used with the Axiom™ 2.0 Reagent Kit, enables researchers to perform large-scale genotyping studies with minimal hands-on processing per plate.

Each plate consists of 96 microarrays and is offered as a part of a complete automated solution that includes instrumentation for sample preparation, and array washing, staining, and scanning.

These markers are amplified using the Axiom™ 2.0 Reagent Kit and are present on short randomly generated fragments of genomic DNA (25–125 base pair) before array hybridization. The reagents for processing Axiom™ arrays are also included in this reagent kit.

Identical to the cartridge array manufacturing process, the oligonucleotide probes on Axiom™ Array Plates are synthesized *in situ* using the photolithographic process.

Instructions for use

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

Axiom 2.0 Assay

- *Axiom™ 2.0 Assay 96-Array Format Automated Workflow User Guide—Biomek™ FX^P* (Pub. No. [MAN0027801](#), SPG Pub. No. [MAN0019897](#), QR Pub. No. [MAN0018118](#))
- *Axiom™ 2.0 Assay 96-Array Format Automated Workflow User Guide—Applied Biosystems™ NIMBUS® Target Preparation Instrument* (Pub. No. [MAN0017740](#), SPG Pub. No. [MAN0017741](#), QR Pub. No. [MAN0017742](#))
- *Axiom™ 2.0 Assay 96-Array Format Manual Workflow User Guide* (Pub. No. [MAN0018000](#), SPG Pub. No. [MAN0018132](#), QR Pub. No. [MAN0018119](#))
- *Axiom™ 2.0 Assay 24-Array Format Manual Workflow User Guide* (Pub. No. [MAN1000233](#), SPG Pub. No. [MAN0018141](#), QR Pub. No. [MAN0018147](#))

Related documentation

- *Axiom™ 2.0 gDNA Sample Preparation Quick Reference* (Pub. No. [MAN0017720](#))
- *GeneTitan™ MC Protocol for Axiom™ Array Plate Processing Quick Reference* (Pub. No. [MAN0017718](#))
- *GeneTitan™ Multi-Channel Instrument User Guide* (Pub. No. [MAN0027694](#), SPG Pub. No. [MAN0025571](#))
- *GeneChip™ Command Console™ User Guide* (Pub. No. [MAN0027771](#))
- *Axiom™ Analysis Suite User Guide* (Pub. No. [MAN0027928](#))
- *Axiom™ Genotyping Solution Data Analysis User Guide* (Pub. No. [MAN0018363](#))

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.

Reagents, instrumentation, and software required

1. Reagent kit:
 - Axiom™ 2.0 Reagent Kit
 - Axiom™ 2.0 Reagent Kit 4x24 Reactions
2. Automated workstation (optional):
 - Beckman Coulter™ Biomek™ FX^P Target Prep Express
 - Applied Biosystems™ NIMBUS® Target Preparation Instrument
3. GeneTitan™ Multi-Channel Instrument or GeneTitan™ MC Fast Scan Instrument
4. GeneChip™ Command Console™ (GCC) software
5. Axiom™ Analysis Suite software

For a complete list of reagents and consumables required, see to the appropriate Axiom™ 2.0 Assay site preparation guide.

Ordering information

Product name	Description	Cat. No.
Axiom™ Array Plate	One array plate, 96-array format or 24-array format	Varies
Axiom™ 2.0 Reagent Kit	Sufficient for one 96-array plate	901758
Axiom™ 2.0 Reagent Kit 4x24 Reactions	Sufficient for four 24-array plates	902798
Axiom™ GeneTitan™ Consumables Kit ^[1]	Contains all GeneTitan™ consumables required to process one Axiom™ array plate.	901606


^[1] One hybridization tray, one scan tray, and five stain trays with covers are included for use with each array plate. These consumables are required for processing Axiom™ Array Plates on the GeneTitan™ Multi-Channel Instrument.

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-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.

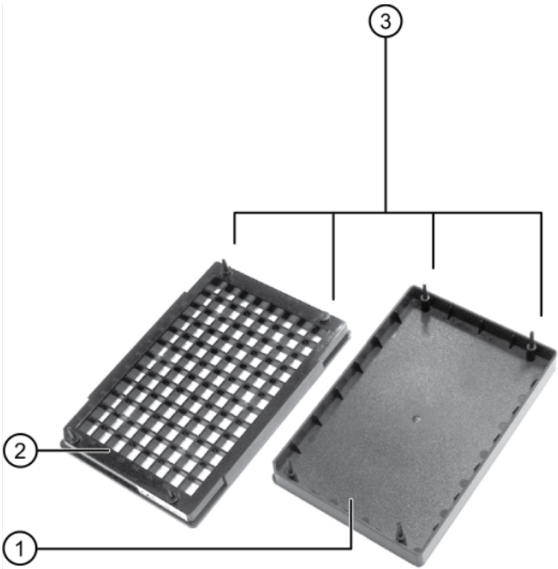


Figure 1 Scan tray assembly.
① Black plastic base ③ Alignment pins
② Scan tray

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. The 96-format array plate is shipped in 1 of 2 different configurations: A or B, below. Configuration A includes a clear shipping cover and a clear array plate. Configuration B does not include a cover and has a black array plate. Both configurations are acceptable. Keep the array plate in the protective base at all times.

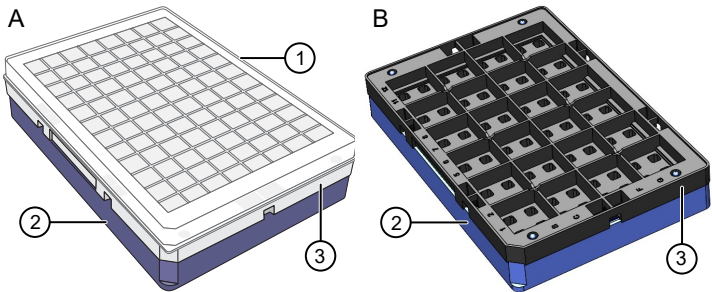


Figure 2 Array plate assembly.
① Clear plastic cover ③ Array plate
② Blue base

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.



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.

©2018-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.