

Axiom™ PangenomiX Array Plate

Catalog Numbers 952518

Pub. No. MAN0029735 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™ PangenomiX Array Plate is designed for genotyping of single nucleotide polymorphisms (SNPs), simple insertion/deletions (indels), and copy number variations (CNV). This array plate, when used with the Axiom™ 2.0 Plus Reagent Kit 96F (or 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 solution that includes options for sample preparation, and instrumentation for array washing, staining, and scanning.

Markers are amplified using a whole-genome amplification step as part of the Axiom™ 2.0 Plus Reagent Kit 96F or the Axiom™ 2.0 Reagent Kit. A mPCR step is enabled with Axiom™ 2.0 Plus Reagent Kit 96F to offer accurate genotyping of markers in regions of high homology. Amplified DNA is fragmented into short segments of DNA (25–125 base pair) before array hybridization. The reagents for processing Axiom™ PangenomiX Array Plate are also included in these reagent kits.

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

Instructions for use

See the following documents for instructions on processing samples using the Axiom™ 2.0 Plus Assay or the Axiom™ 2.0 Assay.

Axiom™ 2.0 Plus Assay

- *Axiom™ 2.0 Plus Assay 96-Array Format Manual Workflow User Guide* (Pub. No. [MAN0026503](#), Site Preparation Guide Pub. No. [MAN0026504](#), Quick Reference Pub. No. [MAN0026505](#))
- *Axiom™ Propel Fast Wash Plus Workflow, 96-Array Format User Guide* (Pub. No. MAN0029733, Site Preparation Guide Pub. No. MAN0029734)

Axiom™ 2.0 Assay

- *Axiom™ 2.0 Assay 96-Array Format Manual Workflow User Guide* (Pub. No. [MAN0018000](#), Site Preparation Guide Pub. No. [MAN0018132](#), Quick Reference Pub. No. [MAN0018119](#))
- *Axiom™ Propel Fast Wash Workflow, 96-Array Format User Guide* (Pub. No. MAN0019450, Site Preparation Guide Pub. No. MAN0019451)

Related documentation

- *GeneTitan™ MC Protocol for Axiom™ Array Plate Processing Quick Reference* (Pub. No. [MAN0017718](#))
- *GeneTitan™ Multi-Channel Instrument User Guide* (Pub. No. [MAN0027694](#), Site Preparation Guide 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](#))
- *Axiom™ Copy Number Data Analysis User Guide* (Pub. No. [MAN0026736](#))

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 Plus Reagent Kit 96F (with reagents for mPCR step), or
 - Axiom™ 2.0 Reagent Kit
2. GeneTitan™ MC Fast Scan Instrument or the GeneTitan™ Multi-Channel Instrument
3. GeneChip™ Command Console™ (GCC) software
4. Axiom™ Analysis Suite software

For a complete list of reagents and consumables required, see the appropriate user guide for your assay.

Ordering information

Product name	Description	Source
Axiom™ PangenomiX Array Plate	One 96-array format plate	952518
Supporting products		
Axiom™ 2.0 Plus Reagent Kit 96F	Reagents sufficient for one 96-array plate (including a mPCR step)	951960
Axiom™ 2.0 Reagent Kit	Sufficient for one 96-array plate	901758
Axiom™ GeneTitan™ Consumables Kit	Sufficient consumables to process one 96-array plate on the GeneTitan™ MC Fast Scan Instrument or the GeneTitan™ MC Instrument	901606
Assay Kits		
Axiom™ PangenomiX Assay Kit	Sufficient for processing one 96-array plate. Consists of: <ul style="list-style-type: none"> • Axiom™ 2.0 Reagent Kit • Axiom™ GeneTitan™ Consumables Kit • Axiom™ PangenomiX Array Plate 	952519
Axiom™ Plus PangenomiX Assay Kit	Sufficient for processing one 96-array plate (with reagents for mPCR step). Consists of: <ul style="list-style-type: none"> • Axiom™ 2.0 Plus Reagent Kit 96F • Module A mPCR Primers • Axiom™ GeneTitan™ Consumables Kit • Axiom™ PangenomiX Array Plate 	952521
Axiom™ Propel PangenomiX Assay Kit, 4X	Sufficient for processing four 96-array plates. Consists of: <ul style="list-style-type: none"> • Axiom™ Propel Fast Reagent Kit, 4x96F • Axiom™ GeneTitan™ Barcoded Consumables Kit • Axiom™ PangenomiX Array Plate 	952528
Axiom™ Propel PangenomiX Assay Kit, 8X	Sufficient for processing eight 96-array plates. Consists of: <ul style="list-style-type: none"> • Axiom™ Propel Fast Reagent Kit, 8x96F • Axiom™ GeneTitan™ Barcoded Consumables Kit • Axiom™ PangenomiX Array Plate 	952529
Axiom™ Propel Plus PangenomiX Assay Kit, 4X	Sufficient for processing four 96-array plates (with reagents for mPCR step). Consists of: <ul style="list-style-type: none"> • Axiom™ Propel Fast Reagent Kit, 4x96F • Axiom™ GeneTitan™ Barcoded Consumables Kit • Module A mPCR Primers • Axiom™ PangenomiX Array Plate 	952530
Axiom™ Propel Plus PangenomiX Assay Kit, 8X	Sufficient for processing eight 96-array plates (with reagents for mPCR step). Consists of: <ul style="list-style-type: none"> • Axiom™ Propel Fast Reagent Kit, 8x96F • Axiom™ GeneTitan™ Barcoded Consumables Kit • Module A mPCR Primers • Axiom™ PangenomiX Array Plate 	952531

Product name	Description	Source
Training kits		
Axiom™ PangenomiX Training Kit	Sufficient for processing one 96-array plate. Consists of: <ul style="list-style-type: none"> • Axiom™ 2.0 Reagent Kit • Axiom™ GeneTitan™ Consumables Kit • Axiom™ DNA Training Plate, 96F (25 µL) • Axiom™ PangenomiX Array Plate 	952522
Axiom™ Plus PangenomiX Training Kit	Sufficient for processing one 96-array plate (with reagents for mPCR step). Consists of: <ul style="list-style-type: none"> • Axiom™ 2.0 Plus Reagent Kit 96F • Module A mPCR Primers • Axiom™ GeneTitan™ Consumables Kit • Axiom™ DNA Training Plate, 96F (25 µL) • Axiom™ PangenomiX Array Plate 	952523
Axiom™ PangenomiX Plus Array Training Kit with Propel Plus 4X Assay and Fast Wash	Sufficient for processing four 96-array plates (with reagents for mPCR step). Consists of: <ul style="list-style-type: none"> • Axiom™ Propel Fast Reagent Kit, 4x96F • Module A mPCR Primers • Axiom™ GeneTitan™ Barcoded Consumables Kit • Axiom™ DNA Training Plate, 96F (25 µL) • Axiom™ PangenomiX Array Plate 	952524
Axiom™ PangenomiX Plus Array Training Kit with Propel Plus 8X Assay and Fast Wash	Sufficient for processing eight 96-array plates (with reagents for mPCR step). Consists of: <ul style="list-style-type: none"> • Axiom™ Propel Fast Reagent Kit, 8x96F • Module A mPCR Primers • Axiom™ GeneTitan™ Barcoded Consumables Kit • Axiom™ DNA Training Plate, 96F (25 µL) • Axiom™ PangenomiX Array Plate 	952525

Storage, handling, and stability

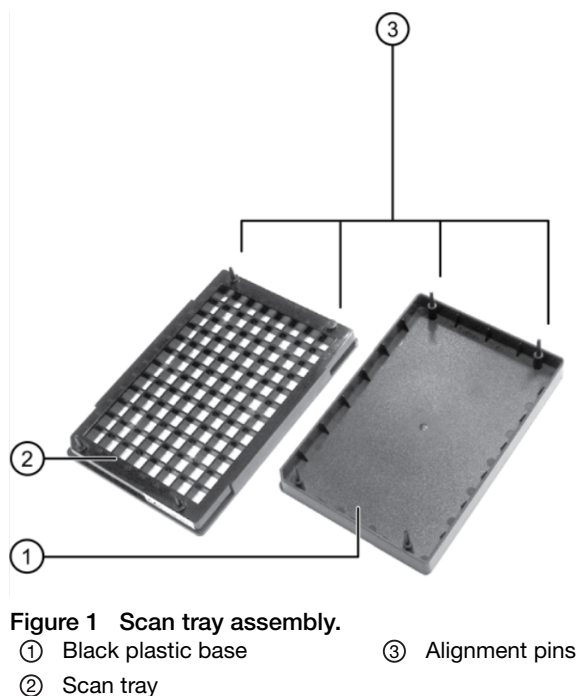
Store array plates at 2–8°C and do not allow to freeze. Do not open the packaging until ready for use—this is essential to protect the array plates from damage, exposure to dust, and excess light. Refer to the expiration date on the package label. Do not use array plates or reagents after the expiration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



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.

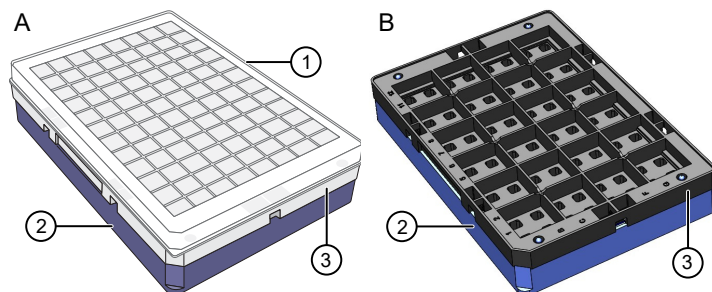


Figure 2 Array plate assembly.

- ① Clear plastic cover
- ② Blue base
- ③ Array plate

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

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

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.

©2023-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.