


Axiom™ Reagent Kits

Catalog Number Multiple

Pub. No. MAN0017595 Rev. D.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

The Applied Biosystems™ Axiom™ 2.0 Reagent Kit enables genotyping of single nucleotide polymorphisms (SNPs) and insertions/deletions (indels) (on Axiom™ genome-wide array plates and Axiom™ myDesign™ Array Plates), the detection of microbial organisms in a sample (Axiom™ Microbiome Array Plates), and targeted structural genomic variants, such as small intragenic deletions and duplications (copy number variants, CarrierScan™ Assay, PharmacoFocus™ and PharmacoScan™ Solutions). The kit includes the reagents required to:

- Produce hybridization-ready targets for Axiom™ genotyping array plates, Axiom™ Microbiome Array Plates, Axiom™ myDesign™ Array Plates, Axiom™ PMD Array Plates, PharmacoScan™ Array Plates, and Axiom™ PharmacoFocus™ Array Plates.
- Process array plates on the Applied Biosystems™ GeneTitan™ Multi-Channel (MC) Instrument.

Ordering information

Product name	Description	Source
Axiom™ 2.0 Reagent Kit	Sufficient for 1 Axiom™ Array Plate (96-array format).	901758
Axiom™ 2.0 384HT Reagent Kit	Sufficient for 1 Axiom™ 384HT Array Plate.	902245
Axiom™ 2.0 Reagent Kit 4x24 Reactions ^[1]	Sufficient for 4 Axiom™ Array Plates(24-array format).	902798
Axiom™ 2.0 Assay Mini 96 Reagent Kit	Sufficient for 2 Axiom™ Mini 96-Array Plates.	903013
Axiom™ Microbiome Reagent Kit 4x24 Reactions	Sufficient for 4 Axiom™ Array Plates(24-array format).	902910
Axiom™ 2.0 Plus Reagent Kit 96F	Sufficient for 1 Axiom™ Array Plate with mPCR step.	951960
PharmacoScan™ Reagent Kit 4x24 Reactions ^[1]	Sufficient for 4 PharmacoScan™ 24F Array Plates.	902908TS
PharmacoScan™ Reagent Kit 96 Reactions	Sufficient for 1 PharmacoScan™ 96F Array Plate.	913025
Axiom™ PharmacoFocus™ Assay Mini 96 Reagent Kit	Sufficient for 1 Axiom™ PharmacoFocus™ Mini 96-Array Plate.	952389
Axiom™ PharmacoFocus™ 384HT Reagent Kit	Sufficient for 1 Axiom™ PharmacoFocus™ 384HT Array Plate.	952390

^[1] The PharmacoScan™ Reagent Kit 4x24 Reactions and the Axiom™ 2.0 Reagent Kit 4x24 Reactions contain extra ligase and hold buffer. These extra buffers are not required when running a single 96-array format plate.

Instructions for use

For instructions on using this reagent kit, Axiom™ genotyping array plates, Axiom™ Microbiome Array Plates, Axiom™ myDesign™ Array Plates, Axiom™ PMD Array Plates, PharmacoScan™ Array Plates, and Axiom™ PharmacoFocus™ Array Plates, see the following documentation:

Axiom™ 2.0 Assay

- *Axiom™ 2.0 Assay 96-Array Format Automated Workflow User Guide—Biomek™ FX^P*, Pub. No. 702963 (QR Pub. No. MAN0018118).
- *Axiom™ 2.0 Assay 96-Array Format Automated Workflow User Guide—Applied Biosystems™ NIMBUS™ Target Preparation Instrument*, Pub. No. MAN0017740 (QR Pub. No. MAN0017742).
- *Axiom™ 2.0 Assay 96-Array Format Manual Workflow User Guide*, Pub. No. MAN0018000 (QR Pub. No. MAN0018119).
- *Axiom™ 2.0 Assay 24-Array Format Manual Workflow User Guide*, Pub. No. 703335 (QR Pub. No. MAN0018147).
- *Axiom™ 2.0 Assay 384HT Array Format Automated Workflow User Guide—Biomek™ FX^P*, Pub. No. 703268 (QR Pub. No. 703270).
- *Axiom™ 2.0 Assay 384HT Array Format Automated Workflow User Guide—Biomek™ i7*, Pub. No. MAN0017555 (QR Pub. No. MAN0017557 and MAN0019790).
- *Axiom™ 2.0 Assay Mini 96-Array Format Manual Workflow User Guide*, Pub. No. 703434 (QR Pub. No. 703436).
- *Axiom™ Microbiome Solution User Guide*, Pub. No. 703408.

Axiom™ 2.0 Plus Assay

- *Axiom™ 2.0 Plus Assay 96-Array Format Automated Workflow User Guide—Applied Biosystems™ NIMBUS™ Target Preparation Instrument*, Pub. No. MAN0018331 (QR Pub. No. MAN0018333).

PharmacoScan™ Assay

- *PharmacoScan™ Assay 96-Array Format Automated Workflow User Guide—Biomek™ FX^P*, Pub. No. 703472 (QR Pub. No. 703474).
- *PharmacoScan™ Assay 96-Array Format Automated Workflow User Guide—Applied Biosystems™ NIMBUS™ Target Preparation Instrument*, Pub. No. MAN0018101 (QR, Pub. No. MAN0018103).
- *PharmacoScan™ Assay 96-Array Format Manual Workflow User Guide*, Pub. No. 703459 (QR Pub. No. 703461).
- *PharmacoScan™ Assay 24-Array Format Manual Workflow User Guide*, Pub. No. 703286 (QR Pub. No. 703288).

Axiom™ PharmacoFocus™ Assay

- *Axiom™ PharmacoFocus™ Assay Mini 96-Array Format Manual Workflow User Guide*, Pub. No. MAN0019172 (QR Pub. No. MAN0019174).
- *Axiom™ PharmacoFocus™ Assay Mini 96-Array Format Automated Workflow User Guide—Applied Biosystems™ NIMBUS™ Target Preparation Instrument*, Pub. No. MAN0019222 (QR Pub. No. MAN0019224).
- *Axiom™ PharmacoFocus™ Assay 384HT Array Format Automated Workflow User Guide—Applied Biosystems™ NIMBUS™ Target Preparation Instrument*, Pub. No. MAN0019212 (QR Pub. No. MAN0019214).

Contents and storage

Module	Storage	Axiom™ 96 901758	Axiom™ 384HT 902245	Axiom™ 4x24 902798	Axiom™ Mini 96 903013	Axiom™ Microbiome 902910	Axiom™ 2.0 Plus 951960	PS 4x24 902908TS	PS 96 913025	PF Mini 96 952389	PF 384HT 952390
Module 1 (Part No. 901711)	-25°C to -15°C	1	2	1	1	1	1	1	1	1	4
Axiom™ 2.0 Denat Soln 10X		1	2	1	1	1	1	1	1	1	4
Axiom™ 2.0 Neutral Soln		1	2	1	1	1	1	1	1	1	4
Axiom™ Water		1	2	1	1	1	1	1	1	1	4
Axiom™ 2.0 Amp Soln		1	2	1	1	1	1	1	1	1	4
Axiom™ 2.0 Amp Enzyme		1	2	1	1	1	1	1	1	1	4
Module 2—Pouch 1 of 2 (Part No. 901528)	-25°C to -15°C	1	2	1	1	1	1	1	1	1	4
Axiom™ Frag Enzyme		1	2	1	1	1	1	1	1	1	4
Axiom™ 10X Frag Buffer		1	2	1	1	1	1	1	1	1	4
Axiom™ Precip Soln 2		1	2	1	1	1	1	1	1	1	4
Axiom™ Hyb Buffer		1	2	1	1	1	1	1	1	1	4
Axiom™ Hyb Soln 1		1	2	1	1	1	1	1	1	1	4
Module 2—Pouch 2 of 2 (Part No. 901529)	2°C to 8°C	1	2	1	1	1	1	1	1	1	4
Axiom™ Frag Diluent		1	2	1	1	1	1	1	1	1	4
Axiom™ Frag Rxn Stop		1	2	1	1	1	1	1	1	1	4
Axiom™ Precip Soln 1		1	2	1	1	1	1	1	1	1	4
Axiom™ Resusp Buffer		1	2	1	1	1	1	1	1	1	4
Axiom™ Hyb Soln 2		1	2	1	1	1	1	1	1	1	4
Module 3	Room temperature	1	1	4	2	4	1	4	1	1	1
Axiom™ Wash Buffer A (Part No. 901446)		2	2	8	4	8	2	8	2	2	2
Axiom™ Wash Buffer B (Part No. 901447)		1	1	4	2	4	1	4	1	1	1
Axiom™ Water (Part No. 901578)		1	1	4	2	4	1	4	1	1	1

Module	Storage	Axiom™ 96 901758	Axiom™ 384HT 902245	Axiom™ 4x24 902798	Axiom™ Mini 96 903013	Axiom™ Microbiome 902910	Axiom™ 2.0 Plus 951960	PS 4x24 902908TS	PS 96 913025	PF Mini 96 952389	PF 384HT 952390
Module 4—Pouch 1 of 2 (Part No. 901278)	-25°C to -15°C	1	2	1	1	1	1	1	1	1	2
Axiom™ Ligate Buffer		1	2	1	1	1	1	1	1	1	2
Axiom™ Ligate Enzyme		1	2	1	1	1	1	1	1	1	2
Axiom™ Ligate Soln 1		1	2	1	1	1	1	1	1	1	2
Axiom™ Probe Mix 1		1	2	1	1	1	1	1	1	1	2
Axiom™ Stain Buffer		1	2	1	1	1	1	1	1	1	2
Axiom™ Stabilize Soln		1	2	1	1	1	1	1	1	1	2
Module 4—Pouch 2 of 2 (Part No. 901276)	2°C to 8°C	1	2	1	1	1	1	1	1	1	2
Axiom™ Ligate Soln 2		1	2	1	1	1	1	1	1	1	2
Axiom™ Probe Mix 2		1	2	1	1	1	1	1	1	1	2
Axiom™ Wash A		1	2	1	1	1	1	1	1	1	2
Axiom™ Stain 1-A		1	2	1	1	1	1	1	1	1	2
Axiom™ Stain 1-B		1	2	1	1	1	1	1	1	1	2
Axiom™ Stain 2-A		1	2	1	1	1	1	1	1	1	2
Axiom™ Stain 2-B		1	2	1	1	1	1	1	1	1	2
Axiom™ Stabilize Diluent		1	2	1	1	1	1	1	1	1	2
Axiom™ Water		1	2	1	1	1	1	1	1	1	2
Axiom™ Hold Buffer	1	2	1	1	1	1	1	1	1	2	
Module 5—Pouch 1 of 2 (Part No. 902796)	-25°C to -15°C	—	—	1	—	1	1	1	—	—	—
Axiom™ Ligate Enzyme				1		1	1	1			
Module 5—Pouch 2 of 2 (Part No. 902797)	2°C to 8°C	—	—	1	—	1	1	1	—	—	—
Axiom™ Hold Buffer				3		3	3	3			
Genomic DNA Standard (Part No. 901012)	-25°C to -15°C	—	—	—	—	3	—	—	—	—	—
Axiom™ Reference gDNA 103											

Module	Storage	Axiom™ 96 901758	Axiom™ 384HT 902245	Axiom™ 4x24 902798	Axiom™ Mini 96 903013	Axiom™ Microbiome 902910	Axiom™ 2.0 Plus 951960	PS 4x24 902908TS	PS 96 913025	PF Mini 96 952389	PF 384HT 952390
Module A—mPCR Primers (Part No. 912896) 10X Primer Mix	-25°C to -15°C	—	—	—	—	—	1 1	1 4	1 ^[1] 1	1	4 4
Module B—Control DNA (Part No. 912897) Control DNA 1, 5 ng/μL Control DNA 2, 5 ng/μL	-25°C to -15°C	—	—	—	—	—	—	1 4 4	1 ^[2] 1 1	1 1 1	4 4 4
Axiom™ Hold Buffer (Part No. 903012)	2°C to 8°C	—	—	—	1	—	—	—	—	—	—

^[1] The Part No. for Module A of the PharmacoScan™ Reagent Kit 96 Reactions is 912896. This differs from the PharmacoScan™ Reagent Kit 4x24 Reactions Module A Part No. listed.

^[2] The Part No. for Module B of the PharmacoScan™ Reagent Kit 96 Reactions is 912897. This differs from the PharmacoScan™ Reagent Kit 4x24 Reactions Module B Part No. listed.

Related documentation

- *GeneTitan™ MC Protocol for Axiom™ Array Plate Processing Quick Reference* (Pub. No. MAN0017718)
- *GeneTitan™ MC Protocol for Axiom™ 384HT Array Plate Processing Quick Reference* (Pub. No. MAN0017596)
- *Axiom™ gDNA Sample Preparation Quick Reference* (Pub. No. MAN0017720)
- *Axiom™ 384HT and Mini 96 gDNA Sample Preparation Quick Reference* (Pub. No. MAN0017719)

Customer and technical support

Visit [thermofisher.com/support](https://www.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.



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 | Vilnius, Lithuania

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.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.

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