

Axiom™ 2.0 Assay for 96 Samples Manual Workflow

Site Preparation Guide

Information in this document is subject to change without notice.

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC 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.

Corporate entity

Life Technologies | Carlsbad, CA 92008 USA | Toll free in USA 1.800.955.6288

Trademarks

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.

© 2017 Thermo Fisher Scientific Inc. All rights reserved.

P/N 702991

Contents

Chapter 1 Equipment and Supplies for the Axiom™ 2.0 Assay for 96 Samples Manual Workflow.....	3
Equipment, Software, Reagents and Arrays Required	3
Related Documentation.....	4
Equipment Required for Manual Target Prep	5
Pre-Amplification Area/Amplification Staging Area	5
Oven Requirements	5
Spectrophotometer	6
Thermal Cycler Recommendations.....	7
Thermal Cycler Consumables.....	7
Plate Centrifuge	7
Sample Plates.....	9
Pipettes and Tips.....	9
Other Labware Required.....	10
Other Reagents and Gels Required	10
Other Lab Equipment Required	11
Appendix A Supplier Contact List	12

Chapter 1

Equipment and Supplies for the Axiom™ 2.0 Assay for 96 Samples Manual Workflow

This site preparation guide includes the supplier and part number information for the equipment, software, reagents, arrays, labware and other consumables that have been validated for use with the Axiom™ 2.0 Assay for 96 Samples Manual Workflow.

Information on suppliers of the equipment and supplies is listed in Appendix A, *Supplier Contact List*.

Equipment, Software, Reagents and Arrays Required

Table 1.1 Equipment, Software, Reagents and Arrays Required

✓ Item	Part Number
Equipment	
<ul style="list-style-type: none"> GeneTitan™ Multi-Channel Instrument* 	Contact Thermo Fisher, Inc.
Software	
<ul style="list-style-type: none"> GeneChip™ Command Console™ (AGCC) 	version 3.1 or later
<ul style="list-style-type: none"> Axiom™ Analysis Suite 	version 1.1 or later
Reagents	
<ul style="list-style-type: none"> Axiom™ 2.0 Reagent Kit 	901758 (96 reactions)
Arrays	
<ul style="list-style-type: none"> Axiom™ Genotyping Solution Array Plates <p>Please visit our web site for a complete list of array plates available for use with the Axiom™ 2.0 Assay.</p>	Varies, depending on array design
<ul style="list-style-type: none"> Axiom™ Microbiome 96-Array Plates 	550666/902904
Consumables	
<ul style="list-style-type: none"> Each Axiom™ GeneTitan™ Consumables Kit contains the following labware for each array plate: <ul style="list-style-type: none"> 1 Hyb Tray 1 Scan Tray with top cover and protective base 5 Stain Trays with top covers <p>These trays are required for processing Axiom array plates on the GeneTitan™ Multichannel Instrument.</p> <p>NOTE: Consumables are included with the Axiom Genome-Wide BOS and Pan-African Genotyping Bundles.</p> 	901606

* For a complete list of all equipment and supplies required for GeneTitan Instrument installation and operation, please consult the GeneTitan Multi-Channel Instrument Site Preparation Guide (P/N 08-0305).

Table 1.1 Equipment, Software, Reagents and Arrays Required (Continued)

✓ Item	Part Number
Training Module	
<ul style="list-style-type: none"> • Axiom™ 2.0 Assay Training Module for CEU, EUR, LAT, and AFR Arrays • Axiom™ 2.0 Assay Training Module for ASI, CHB, and EAS Arrays • Axiom™ 2.0 Assay Training Module for BOS Arrays <p>Each Kit Includes:</p> <ul style="list-style-type: none"> - Two Genome-Wide 96-Array Plates for specific array plate types listed above, except for BOS which includes 1 CEU Array Plate and 1 BOS Array Plate. - Two Axiom™ 2.0 Reagent Kits (96 reactions) - Two Axiom™ GeneTitan™ Consumables Kits - Two full plates of 96 HapMap DNA samples (except for BOS training modules which only include one 96 HapMap DNA sample plate) - One Zerostat Anti-Static Gun 	<p>901767</p> <p>901768</p> <p>901750</p>
<ul style="list-style-type: none"> • Axiom™ 2.0 Pan-African Training Module <p>Each Kit Includes:</p> <ul style="list-style-type: none"> - One Axiom Genome-Wide Pan-African Genotyping Bundle (1 set of 96-array plates, 3 Axiom 2.0 Reagent Kits, and 3 Axiom GeneTitan Consumables Kits) - Three identical 96 HapMap DNA plates - One Zerostat Anti-Static Gun 	<p>901769</p>
<ul style="list-style-type: none"> • Axiom™ 2.0 Assay Training Module for Microbiome 96-Array Plates <p>Each Kit Includes:</p> <ul style="list-style-type: none"> - One Axiom™ CEU 96-Array Plate - One Axiom™ Microbiome 96-Array Plate - Two Axiom™ 2.0 Reagent Kits (96 reactions) - Two Axiom™ GeneTitan™ Consumables Kits - One full plate of 96 HapMap DNA samples - One Zerostat Anti-Static Gun 	<p>902988</p>

Related Documentation

- *Axiom™ 2.0 Assay for 96 Samples Manual Workflow User Guide*, P/N 702990
- *Axiom™ 2.0 Assay for 96 Samples Manual Workflow Site Preparation Guide*, P/N 702991
- *Axiom™ 2.0 Assay for 96 Samples Manual Target Prep Protocol QRC*, P/N 702989
- *Axiom™ Microbiome Solution User Guide*, P/N 703408
- *Axiom™ 2.0 Assay for 24 Samples User Guide*, P/N 703335
- *Axiom™ 2.0 Assay for 24 Samples Site Preparation Guide*, P/N 703336
- *Axiom™ 2.0 Assay for 24 Samples Manual Target Preparation Protocol QRC*, P/N 703337
- *Axiom™ 2.0 Assay Automated Workflow on NIMBUS™ User Guide*, P/N 703349
- *Axiom™ 2.0 Assay Automated Workflow on NIMBUS™ Site Preparation Guide*, P/N 703349
- *Axiom™ 2.0 Automated Target Prep Protocol on NIMBUS™ QRC*, P/N 703351
- *Axiom™ 2.0 gDNA Sample Preparation QRC*, P/N 702987

- *Axiom™ gDNA Sample Prep for Axiom Genome-Wide BOS 1 Array Plate QRC*, P/N 702975
- *Axiom™ Genotyping Solution Data Analysis Guide*, P/N 702961
- *GeneTitan™ MC Protocol for Axiom 2.0 Array Plate Processing QRC*, P/N 702988
- *GeneChip™ Command Console™ (AGCC) 4.0 User Guide*, P/N 702569
- *Axiom™ Analysis Suite User Guide*, P/N 703307
- *Axiom™ 2.0 Assay Automated Workflow User Guide*, P/N 702963
- *Axiom™ 2.0 Assay Automated Target Prep Protocol QRC*, P/N 702962
- *GeneTitan™ Multi-Channel Instrument User Guide*, P/N 08-0306
- *GeneTitan™ Multi-Channel Instrument Site Preparation Guide*, P/N 08-0305

Equipment Required for Manual Target Prep

Table 1.2 Additional Instruments Required for the Axiom™ 2.0 Assay Manual Workflow

✓ Equipment	Manufacturer/ Distributor	Part Number
<ul style="list-style-type: none"> • Two or three ovens (see <i>Oven Requirements</i>) 		
<ul style="list-style-type: none"> • Fume Hood (Strongly recommended: see the <i>Axiom 2.0 Assay Manual Workflow User Guide</i> for more information) 		
<ul style="list-style-type: none"> • One of the following Thermal Cyclers: <ul style="list-style-type: none"> - Bio-Rad DNA Engine PTC-200/MJ PTC-200 Thermal Cycler - Bio-Rad / MJ Tetrad™ 2 PTC-0240G - ABI 9700 - ABI 2720 • (see <i>Thermal Cycler Recommendations</i>, below) 	<ul style="list-style-type: none"> - Bio-Rad - Bio-Rad - ABI - ABI 	
<ul style="list-style-type: none"> • One of the following shakers: <ul style="list-style-type: none"> • Compact Digital Microplate Shaker • Jitterbug™ 	<ul style="list-style-type: none"> - Thermo Scientific - Boekel Scientific 	<ul style="list-style-type: none"> - 88880023 - Model 130 000

Pre-Amplification Area/Amplification Staging Area

Precautions are required when manipulating genomic DNA or setting up amplification reactions to avoid contamination with foreign DNA amplified in other reactions and procedures. It is recommended that genomic DNA manipulations and amplification reaction set up are performed in a dedicated rooms or areas separate from the main laboratory.

These areas should have dedicated sets of pipettes and plasticware. If no dedicated area is available, use of a dedicated bench or a dedicated biosafety hood and dedicated pipettes is suggested. If no dedicated bench or biosafety hood is available, a set of dedicated pipettes is recommended.

Ideally, the pre-amplification and amplification staging areas would be separate; however these areas may be combined due to space and equipment limitations.

Oven Requirements

Multiple ovens are required for manual target preparation. The exact number depends upon whether you are running only a single sample plate and array plate through the workflow, or if you are trying to run the three plate/week manual target preparation workflow.

- If you are running individual plates, you will need two ovens for the workflow.
- If you are running the three plate / week workflow, a third oven is highly recommended.

Refer to the chapter, *Processing Three Axiom Array Plates per Week with Manual Target Preparation in the Axiom™ 2.0 Assay Manual Workflow User Guide*(P/N 702990) for more information.

Table 1.3 Suggested Settings for Ovens When Performing Three Plate/week Manual Target Prep Workflow

Day of Workflow	Oven 1	Oven 2	Oven 3
Day 1	37°C	N/A	N/A
Day 2	37°C	65°C	37°C
Day 3	48°C	65°C	37°C
Day 4	48°C	65°C	37°C
Day 5	N/A	N/A	N/A

Table 1.4 Ovens Required for the Axiom™ 2.0 Assay Manual Workflow

✓ Equipment	Manufacturer/ Distributor	Part Number
<ul style="list-style-type: none"> • Two to three ovens, any combination of the following types: - ED 53 drying oven by Binder 	VWR	50/60 Hz, 115 V, 1200W VWR P/N 47746-744 Mfg. P/N 9010-0131
		50/60 Hz, 230 V, 1200 W VWR P/N 47746-690 Mfg. P/N 9010-0078
GeneChip™ Hyb Oven 645*	Thermo Fisher	00-0331

*The GeneChip™ Hybridization Oven 640 is currently not supported with the Axiom 2.0 Assay.

Spectrophotometer

We recommend using one of the spectrophotometers listed in Table 1.5.

Table 1.5 Spectrophotometers

✓ Item	Supplier	Part Number
<ul style="list-style-type: none"> • One of the following spectrophotometers: - DTX 880 Multimode Detector, with: <ul style="list-style-type: none"> • Genomic Filter Slide or	Beckman Coulter	987921 – detector
	Beckman Coulter	A30184 – filter slide
<ul style="list-style-type: none"> • SpectraMax High throughput Microplate Spectrophotometer 	Molecular Devices	Plus384

Thermal Cycler Recommendations

We have verified the performance of this assay using the following thermal cyclers:

- Bio-Rad DNA Engine PTC-200/MJ PTC-200, or
- Bio-Rad / MJ Tetrad™ 2 PTC-0240G, or
- ABI 9700 (with gold, sliver, or aluminum block), or
- ABI 2720

The performance of this assay has not been verified with other thermal cyclers. Use of other thermal cyclers may result in assay failure and may violate the Axiom Array and Reagent replacement policy. The thermal cycler needs to be programmed with the “Axiom 2.0 Denature” protocol:

1. 95°C 10 min
2. 48°C 3 min
3. 48°C hold

Use the heated lid option when setting up or running the protocol.

Thermal Cycler Consumables

Table 1.6 provides details into the consumables to be used with each thermal cycler.

Table 1.6 Thermal Cycler Consumables for the Axiom 2.0 Assay Manual Workflow

Thermal Cycler Model	PCR Plate Type	Seal*
Bio-Rad PTC-200	Bio-Rad Hard-Shell Thin-Wall 96-Well Skirted PCR Plates, P/N HSP-9631	MicroAmp Clear Adhesive Film from Applied Biosystems (P/N 4306311)
Bio-Rad Tetrad 2 PTC-0240	Bio-Rad Hard-Shell Thin-Wall 96-Well Skirted PCR Plates, P/N HSP-9631	MicroAmp Clear Adhesive Film from Applied Biosystems (P/N 4306311)
ABI 9700	Bio-Rad P/N HSS-9601 (half skirtedplate)	MicroAmp Clear Adhesive Film from Applied Biosystems (P/N 4306311)
ABI 2720	Bio-Rad P/N HSS-9601 (half skirted plate)	MicroAmp Clear Adhesive Film from Applied Biosystems (P/N 4306311)

**Microseal “B” film from Bio-Rad (P/N MSB-1001) may be used in place of MicroAmp Clear Adhesive Film for the Bio-Rad and ABI thermal cyclers. For Axiom™Microbiome applications, only use the MicroAmp Clear Adhesive Film (P/N 4306311). The use of the Bio-Rad Microseal ‘B’ Film has not been validated for Axiom Microbiome use.*

Plate Centrifuge

One plate centrifuge is required for the Axiom™ 2.0 Assay Manual Workflow. We recommend the plate centrifuges listed in Table 1.7. When centrifuging and drying pellets, the centrifuge must be able to spin down plates at:

- rcf: 3200 xg with an appropriate rotor - bucket combination (4000 RPM for the Eppendorf 5810R configuration described below)
- temperature: 4°C

In addition, the bottom of the rotor buckets should be soft rubber to ensure that the deep well plates do not crack. Do not use buckets where the plates sit directly on a metal or hard plastic bottom. For the Eppendorf 5810R, do not use the A-4-62 rotor with a WO-15 plate carrier (hard bottom).

Table 1.7 Plate Centrifuges Recommended for the Axiom™ 2.0 Assay Manual Workflow

✓ Item	Supplier	Part Number
<ul style="list-style-type: none"> One of the following centrifuges: 		
<ul style="list-style-type: none"> Allegra™ 25R Refrigerated Benchtop Centrifuge 	Beckman Coulter	369434 (60 Hz, 280 V) 369435 (50/60 Hz, 200 V) 369436 (50 Hz, 230 V) 368954 S5700 Microplate rotor
<ul style="list-style-type: none"> Centrifuge 5810R (refrigerated) A-4-81 MTP/Flex swinging bucket rotor with 4 microtest-plate buckets 	Eppendorf	022625501 (60 Hz, 120 V) 022625101 (50 Hz, 120 V, 20 AMP version) 022625551 (50 Hz, 230 V) 022638807
<ul style="list-style-type: none"> Sorvall™ Legend™ XTR (refrigerated), with: - TX-750 high visibility swinging rotor bucket - Carrier for microplates (including plate trays and neoprene pads) 	Thermo Scientific	75004521 (60 Hz, 120 V) 75004520 (50 Hz, 230 V) 75004523 (50/60 Hz, 230 V USA and Canada) 75003607 One of either: <ul style="list-style-type: none"> Set of two Carriers (75003795) Set of four carriers (75003617)

Relative centrifugal force (*rcf*) can be calculated as follows:

$$rcf = (1.118 \times 10^{-5}) R S^2$$

where R is the radius of the rotor in centimeters, and S is the speed of the centrifuge in revolutions per minute.

Sample Plates

Table 1.8 Plates Required for Axiom 2.0 Assay Manual Workflow

✓ Equipment	Manufacturer/ Distributor	Part Number
<ul style="list-style-type: none"> ABgene™ 96 Square Well Storage Plate Mark II, 2.2 mL (for use with Axiom Genotyping Assay) 	Thermo Scientific	AB-0932
<ul style="list-style-type: none"> Eppendorf 96 Deep-well Plate, 2000 µL (for use with Axiom Microbiome Assay) 	Eppendorf	951033481
<ul style="list-style-type: none"> Bio-Rad Hard Shell 96-well plate 	Bio-Rad	HSP-9631
<ul style="list-style-type: none"> 96-well UV Star Plates, 370 µL/well 	E&K Scientific	25801
<ul style="list-style-type: none"> Hard-Shell Full-Height 96-Well Semi- Skirted PCR Plate (required only if using ABI 2720 or ABI 9700 Thermal Cyclers) 	Bio-Rad	HSS-9601

Pipettes and Tips

Pipettes and tips recommended for manual target prep are listed in Table 1.9.

Table 1.9 Recommended Pipettes and Tips for Axiom 2.0 Assay Manual Workflow

✓ Equipment	Manufacturer/ Distributor	Part Number
<ul style="list-style-type: none"> Pipet-Lite™, Magnetic-Assist single channel P20 	Rainin	L-20
<ul style="list-style-type: none"> Pipet-Lite™, Magnetic-Assist single channel P200 	Rainin	L-200
<ul style="list-style-type: none"> Pipet-Lite™, Magnetic-Assist single channel P1000 	Rainin	L-1000
<ul style="list-style-type: none"> Pipette, 12-channel P20 	Rainin	L12-20
<ul style="list-style-type: none"> Pipette, 12-channel P50 (optional) 	Rainin	L12-50
<ul style="list-style-type: none"> Pipette, 12-channel P200 	Rainin	L12-200
<ul style="list-style-type: none"> Pipette, 12-channel P1200 	Rainin	L12-1200
<ul style="list-style-type: none"> Pipette tips GP = refill 	Rainin	GP-L10F
<ul style="list-style-type: none"> Pipette tips GP = refill 	Rainin	GP-L200F
<ul style="list-style-type: none"> Pipette tips GP = refill 	Rainin	GP-L1000F
<ul style="list-style-type: none"> Pipette tips RT = with rack 	Rainin	RT-L10F
<ul style="list-style-type: none"> Pipette tips RT = with rack 	Rainin	RT-L200F
<ul style="list-style-type: none"> Pipette tips RT = with rack 	Rainin	RT-L1000F

Other Labware Required

Table 1.10 Other Labware Required for the Axiom 2.0 Assay Manual Workflow

✓ Item	Manufacturer/ Distributor	Part Number
• Conical tube, 50 mL	VWR	21008-178
• Conical tube, 15 mL	VWR	21008-202
• Tube Holder, Falcon, 50 mL	various	various
• Solution Basin, 100 mL sterile, multichannel	VWR	89092-836
• 96 well plate metal chamber	Diversified Biotech	CHAM-1000
• Serological Pipettes (10 mL)	VWR	89130-898
• Serological Pipettes (5 mL)	VWR	89130-896
• Adhesive film for 96-well plates – use one of the following:		
• MicroAmp™ Clear Adhesive Film	Applied Biosystems	4306311
• Microseal 'B' Film*	Bio-Rad	MSB1001
• Kimwipes™	Your choice	—

* For Axiom™ Microbiome applications, only use the MicroAmp Clear Adhesive Film (P/N 4306311). The use of the Bio-Rad Microseal 'B' Film has not been validated for Axiom Microbiome use.

Other Reagents and Gels Required

Table 1.11 Other Reagents Required

✓ Item	Supplier	Part Number
• 2-Propanol, anhydrous, 99.5% (Isopropanol)	Sigma-Aldrich	278475
• Reduced EDTA TE Buffer (10 mM Tris-HCl pH 8.0, 0.1 mM EDTA)	Thermo Fisher	75793

Table 1.12 Reagents and Gels Required to Run QC Steps

✓ Item	Supplier	Part Number
• Mother E-Base™ Device	Thermo Fisher Scientific (formerly Life Technologies™ / Invitrogen)	EB-M03
• Daughter E-Base™ Device		EB-D03
• E-Gel™48 4% agarose gels (for Axiom QC)	Thermo Fisher Scientific (formerly Life Technologies™ / Invitrogen)	G8008-04
• TrackIt™ 25 bp DNA Ladder (for Axiom QC)		10488-022
• TrackIt™ Cyan/Orange Loading Buffer (for Axiom QC)		10482-028
• E-Gel™48 1% agarose gels (for gDNA QC)		G8008-01
• RediLoad™ (for gDNA QC)		750026
• E-Gel™96 High Range DNA Marker (for gDNA QC)		12352-019
• Water, nuclease-free, ultrapure MB grade (only required for performing rehybridization protocols and gel QC)		Thermo Fisher

Other Lab Equipment Required

The remaining equipment required is listed in Table 1.13.

Table 1.13 Other Equipment Required

✓ Item	Supplier	Part Number
• Freezer, -20°C	Any vendor	—
• Refrigerator, 2 to 8°C	Any vendor	—
• Vortex-Genie™ (for plates and microtubes)*	Scientific Industries	SI-0236 (120 V/60 Hz) SI-0246 (230 V/50 Hz)
• Mini Microcentrifuge, for 2 mL tubes*	VWR	93000-196 (120V) 93000-196 (230V)
• Bel-Art Cryo-Safe Mini Cooler, -15°C*	VWR	47751-730
• Ice bucket, 4 to 9 liters	Any vendor	—
• Markers, permanent, fine point	Your choice	—
• Pipet aid*	VWR	53106-220

*Equivalent items from other manufacturers are acceptable.

Appendix A Supplier Contact List

Table A.1 Supplier Contact List

Supplier	Website Address
Thermo Fisher	www.thermofisher.com
Applied Biosystems	www.appliedbiosystems.com
Beckman Coulter	www.beckmancoulter.com
Bio-Rad	bio-rad.com
E&K Scientific	eandkscientific.com
Eppendorf	eppendorf.com
Rainin	www.rainin.com
Scientific Industries	www.scientificindustries.com
Sigma-Aldrich	www.sigmaaldrich.com
Thermo Fisher Scientific (Thermo Scientific, Applied Biosystems, Invitrogen, Fisher Scientific)	www.thermofisher.com
USB Corporation (Thermo Fisher)	www.usb.thermofisher.com
VWR	vwr.com

Documentation and support

Obtaining support

Technical support For the latest services and support information for all locations, visit www.thermofisher.com.

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support (thermofisher.com/support)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

Safety Data Sheets (SDS) Safety Data Sheets (SDSs) are available at thermofisher.com/support.

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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

For support visit thermofisher.com/support or emailtechsupport@lifetech.com

thermofisher.com

23 January 2017

ThermoFisher
SCIENTIFIC