



Product Information

Automated Protocol for GenomePlex® Whole Genome Amplification (WGA) Kit Using the Biomek® FX Workstation (Beckman Coulter)

WGA Product Codes **WGA1** and **WGA2**

Automation Guide	2
I. Description	2
II. Product Components	3
III. Storage	3
IV. Materials to Be Supplied by the User	3
V. Instrument Requirements for the Biomek FX Workstation	3
VI. Deck Setup	4
VII. Sample Requirements	4
VIII. Reagent Preparation	5
1. <i>Genomic DNA Samples</i>	5
2. <i>5× Fragmentation Buffer</i>	5
3. <i>Library Preparation and Stabilization Solution Mixture</i>	5
4. <i>Library Preparation Buffer and Enzyme Mixture</i>	5
5. <i>Amplification Master Mix</i>	5
IX. Automated Method Description	5
A. <i>Getting Started</i>	5
B. <i>Method Overview</i>	6
X. Recommended Parameters for PCR Amplification:	6
XI. Method Customization	7
A. <i>Use of a different PCR plate for Amplification</i>	7
B. <i>Use of a different PCR plate or Reservoir for the Reagents</i>	7
C. <i>Use of a different Reservoir for the Amplification Master Mix</i>	7
XII. Performance Characteristics	8
XIII. Troubleshooting	10
XIV. Contact Information	11

Automation Guide

I. Description

The GenomePlex Whole Genome Amplification (WGA) kit has been developed for use as a high-throughput system for the rapid and highly representative, 500-fold amplification of genomic DNA from trace amounts of sample. The kit utilizes a proprietary amplification technology based upon random fragmentation of genomic DNA and conversion of the resulting small fragments to PCR-amplifiable OmniPlex[®] Library molecules flanked by universal priming sites. The OmniPlex library is then amplified using universal oligonucleotide primers and a limited number of cycles.

The GenomePlex kit is suitable for use with purified genomic DNA from a variety of sources including buccal swabs, whole blood, blood card, and formalin-fixed, paraffin-embedded (FFPE) tissue. GenomePlex WGA kits require 10 ng of starting DNA, and yield 5–10 µg of WGA product after amplification.

A semi-automated method is available for use on the Biomek FX Liquid Handling Workstation from Beckman Coulter. Amplification of genomic DNA from various sources is accomplished in a few simple steps:

1. Fragmentation Buffer is added to genomic DNA in a 96-well plate and incubated off-line at 95 °C for 4 minutes.
2. A mixture of Library Preparation and Stabilization Solution is added to the samples followed by an off-line incubation at 95 °C for 2 minutes.
3. A mixture of Library Preparation Enzyme and Library Preparation Buffer is added to the samples. Samples are then subjected to a series of stepped isothermal reactions in a thermal cycler to generate the OmniPlex library.
4. An amplification master mix with a thermal stable DNA Polymerase is added to the samples followed by 14 cycles of thermal cycling to amplify the OmniPlex library.

All 96 DNA samples can be amplified in less than 4 hours.

II. Product Components

Reagents Provided	Product Code	GenomePlex Whole Genome Amplification Kit WGA1*	GenomePlex Complete Whole Genome Amplification Kit WGA2*
		50 reactions	50 reactions
10× Fragmentation Buffer	F4304	55 µl	55 µl
Library Preparation Buffer	L7167	110 µl	110 µl
Library Stabilization Solution	L7292	55 µl	55 µl
Library Preparation Enzyme	E0531	55 µl	55 µl
10× Amplification Master Mix	A5604	410 µl	410 µl
Nuclease-Free Water	W4765	2 × 1.5 ml	2 × 1.5 ml
Control Human Genomic DNA (5 ng/µl)	D7192	10 µl	10 µl
WGA DNA Polymerase	W3891	Not Included	275 µl

* To perform WGA in a 96-well plate of samples, 3 kits must be ordered. Alternatively, custom packaging is available. Contact your local Sigma sales representative for more information.

III. Storage

All components should be stored at –20 °C. Do not store in a frost-free freezer.

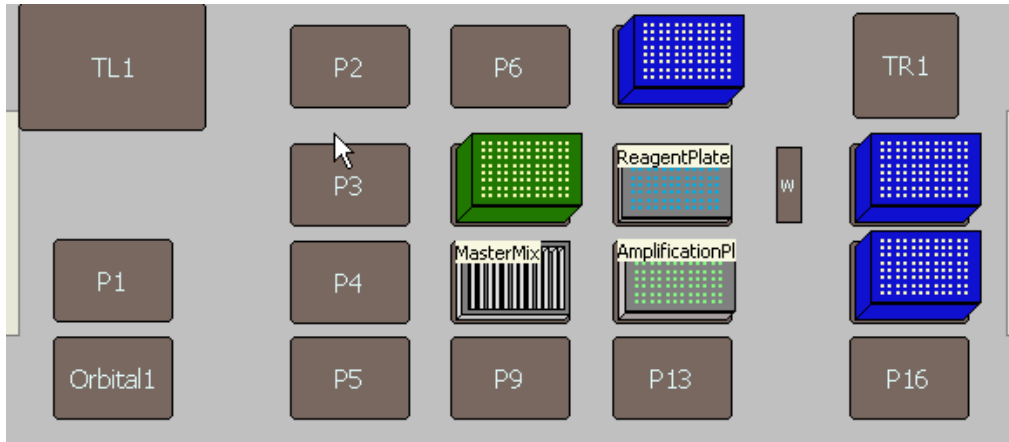
IV. Materials to Be Supplied by the User

1. Genomic DNA samples
2. Water, molecular biology reagent (Sigma, W4502)
3. (Optional) DNA Polymerase for use with WGA1 (Sigma, W3891 or WGA3)
4. 96-well PCR plates, with full skirt (Sigma, P4616)
5. 96-well PCR plates, with half skirt (ABgene, AB-1100)
6. Thermowell® sealing mat (Fisher, 07200614)
7. Corning plate holder (Corning, 6525)
8. Sealing film, SealPlate (Sigma, Z369659)
9. 12 column reagent reservoir with low profile (Innovative Microplates, S30028)
10. (Optional) 12 column reagent reservoir with high profile (Innovative Microplates, S30019)
11. Thermal Cycler

V. Instrument Requirements for the Biomek FX Workstation

Part Description	Qty	Ordering Information
Span-8 Pod	1	Contact Beckman Coulter
Span-8 Tip Trash	1	Contact Beckman Coulter
Span-8 Tip Wash	1	Contact Beckman Coulter
Standard Passive ALPs (One by Three)	3	Contact Beckman Coulter
Standard Passive ALPs (One by One)	1	Contact Beckman Coulter
Span-8 P250 Barrier Tips, Sterile	1	BK379503 (Beckman Coulter)
Span-8 P20 Barrier Tips, Sterile	3	BK379506 (Beckman Coulter)

VI. Deck Setup



Deck Position	Equipment
P7	Span-8 P250 Barrier Tips
P8	12 column reservoir for amplification master mix
P10	Span-8 P20 Barrier Tips
P11	96-well PCR plate with full skirt containing reagents
P12	96-well PCR plate with genomic DNA samples to be amplified (seated into a plate holder)
P14	Span-8 P20 Barrier Tips
P15	Span-8 P20 Barrier Tips

VII. Sample Requirements

1. Protocols for purifying genomic DNA from different biological sources may be downloaded from: www.sigmaaldrich.com/wgaadvisor
2. A minimum of 10 ng of genomic DNA is recommended for amplification with the GenomePlex WGA kit. DNA concentration is critical for successful whole genome amplification and needs to be determined by absorbance. Other methods tend to underestimate single stranded DNA present in a sample.
3. DNA isolated from formalin-fixed, paraffin-embedded (FFPE) tissue may contain impurities and should be quantified using PicoGreen[®] Quantitation Reagent. In this and other samples where DNA damage may have occurred, the starting concentration of genomic DNA may need to be increased. Successful WGA amplification has been performed with degraded samples using 10 ng of starting template, but may require increasing the starting template to 25–100 ng.

VIII. Reagent Preparation

1. Genomic DNA Samples

The GenomePlex kit requires 10 ng of starting genomic DNA. Prepare DNA solutions at a concentration of 1.25 ng/ μ l and aliquot 8 μ l of this solution into each well of a 96-well PCR plate located at position P12. Positive or negative controls can be added to the same plate.

2. 5 \times Fragmentation Buffer

To process a single plate of 96 samples, dilute the 10 \times Fragmentation Buffer 2-fold by adding 140 μ l of water to 140 μ l of Fragmentation Buffer. Aliquot 35 μ l of the solution into each well the first column of the 96-well PCR plate located at position P11.

3. Library Preparation and Stabilization Solution Mixture

To process a single plate of 96 samples, combine the Library Preparation and Stabilization Solution at a 1:1 ratio in a total volume of 280 μ l. Aliquot 35 μ l of the mixture into each well of the third column of 96-well PCR plate located at position P11.

4. Library Preparation Buffer and Enzyme Mixture

To process a single plate of 96 samples, combine the Library Preparation Buffer and Enzyme at a 1:1 ratio in a total volume of 280 μ l. Aliquot 35 μ l of the mixture into each well of the fifth column of 96-well PCR plate located at position P11.

5. Amplification Master Mix

To prepare the Master Mix, add water and DNA Polymerase to the 10 \times Amplification Master Mix as described in the table below.

	Water	Amplification Master Mix	DNA Polymerase
Stock		10 \times	50 \times
Working (6 ml)	5.175 ml	750 μ l	500 μ l

To process one 96-well plate, a total of 6 ml of Amplification Master Mix needs to be added to the first column of the 12-column low profile reservoir (S30028) located at position P8. If setting up more than one 96-well plate of the samples, it will be necessary to use a 12-column high profile reservoir (S30019).

IX. Automated Method Description

This overview describes the general liquid handling steps required to execute the automated WGA method and can be customized to a variety of applications. For custom applications see Section XI.

A. Getting Started

1. Set up the deck layout by placing the tip boxes, plates and reservoirs at the appropriate positions on the deck as described in section VI.
2. Add reagents to the appropriate reservoirs as described in section VIII.
3. Run the method using Biomek Software Version 3.1.
4. When prompted, place the sealing mat onto the amplification plate and move the plate to a thermal cycler for incubations or amplifications as indicated in the dialog box.
5. Store the amplified DNA samples at -20 °C until ready for analysis or purification.

B. Method Overview

Below is a summary of the steps for the *WGA* method. For complete program details, the automation program can be downloaded from www.sigmaldrich.com/automation

1. 5× Fragmentation Buffer (2 µl) is dispensed into each well of the amplification plate containing DNA samples and mixed 4 times.
2. The user is prompted to transfer the amplification plate to a thermal cycler for incubation at 95 °C for 4 minutes.
3. Return the amplification plate back to the deck of the FX and click the “OK” button on the dialog box to continue the method.
4. Library Preparation and Stabilization Solution mixture (2 µl) is dispensed into each well of the amplification plate and mixed 4 times.
5. The user is prompted to transfer the amplification plate to a thermal cycler for incubation at 95 °C for 2 minutes.
6. Return the amplification plate back to the deck of the FX and click the “OK” button on the dialog box to continue the method.
7. Library Preparation Enzyme and Buffer mixture (2 µl) is dispensed into each well of the amplification plate and mixed 4 times.
8. The user is prompted to transfer the amplification plate to a thermal cycler for amplifications.
9. Return the amplification plate back to the deck of the FX and click the “OK” button on the dialog box to continue the method.
10. Amplification Master Mix (60 µl) is dispensed into each well of the amplification plate and mixed 4 times.
11. The user is prompted transfer the amplification plate to a thermal cycler for amplification.
12. Click the “OK” button on the dialog box to finish the program.

X. Recommended Parameters for PCR Amplification:

Step	Temperature	Time	Cycles
Initial Denaturation	95 °C	3 minutes	1
Denaturation	94 °C	15 seconds	14
Annealing/Extension	65 °C	5 minutes	
Hold	4 °C	Indefinitely	

XI. Method Customization

A. Use of a different PCR plate for Amplification

The automated method was created using half-skirt 96-well PCR amplification plates ABgene. Other PCR plates may be used in this method, but may require the creation of a new labware in the Biomek software.

B. Use of a different PCR plate or Reservoir for the Reagents

The automated method was created using full-skirt 96-well PCR plates from Sigma. Other PCR plates or reservoirs may be used in this method, but may require the creation of new labware in the Biomek software.

C. Use of a different Reservoir for the Amplification Master Mix

The automated method was created using a 12-column reservoir for holding the Amplification Master Mix. Other reservoirs may be used in this method, but may require the creation of new labware in the Biomek software.

XII. Performance Characteristics

Whole Genome Amplification Formalin-fixed, Paraffin-embedded Tissue

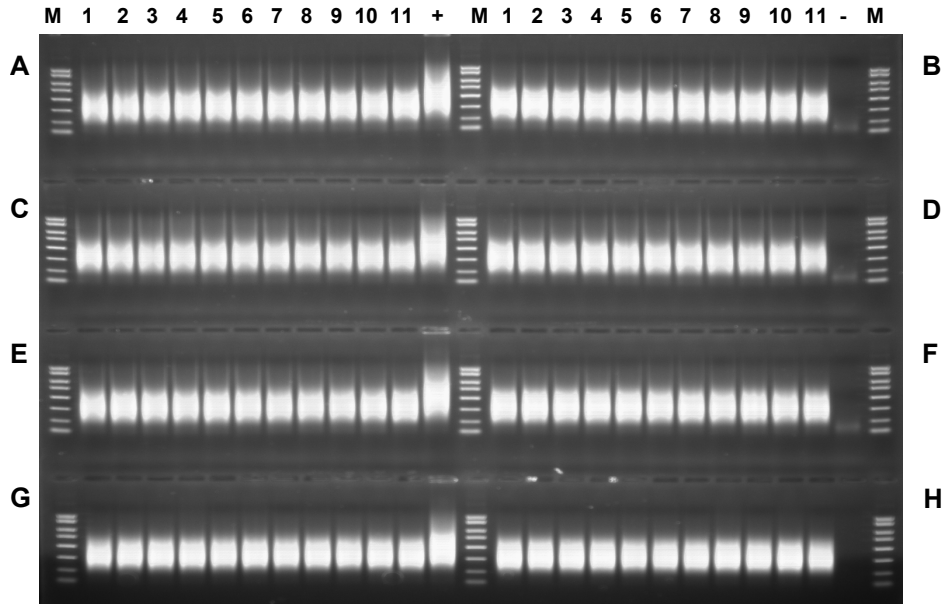


Figure 1. 10 ng of DNA isolated from 88 samples of formalin-fixed, paraffin-embedded (FFPE) rat liver tissue were amplified using the automated WGA procedure on the Biomek FX workstation. 6 μ l of each amplified product were analyzed on a 1% agarose gel. M: PCR marker. (+): Human genomic DNA control. (-): No DNA template control.

Whole Genome Amplification Different Sources

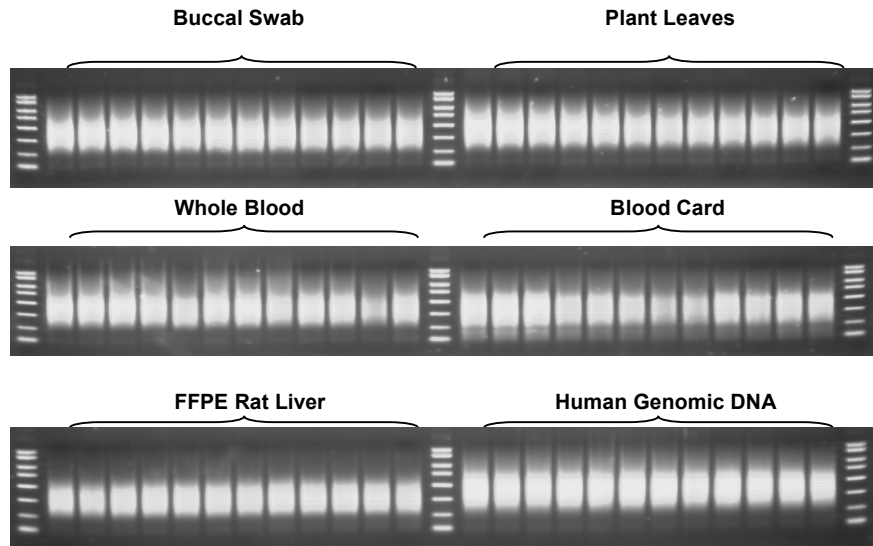


Figure 2. DNA isolated from buccal swab, tomato leaves, whole blood, blood card, and FFPE rat liver tissue was amplified using the automated WGA procedure on the Biomek FX workstation. 6 μ l of each amplified product were analyzed on a 1% agarose gel.

Cross-contamination Analysis

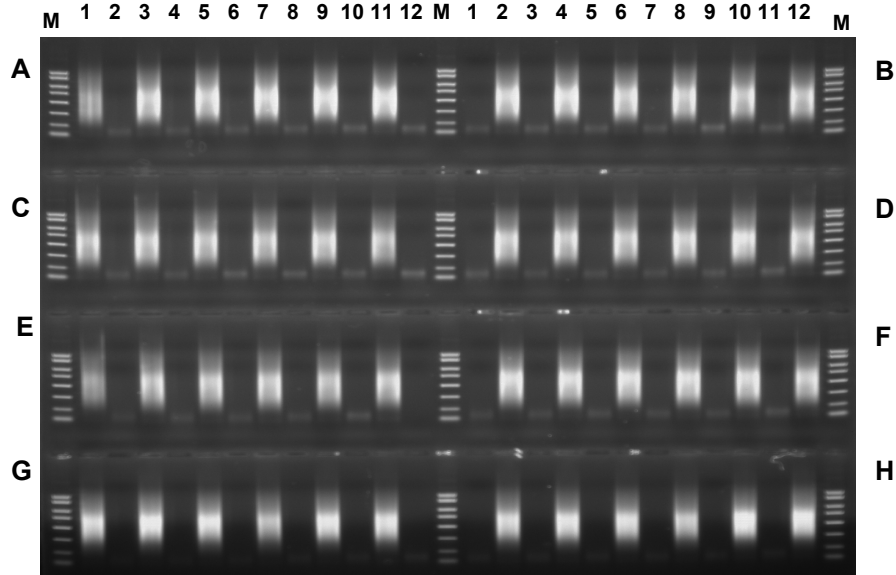


Figure 3. 10 ng of human genomic DNA samples or 8 μ l of water were placed in alternating wells of an amplification plate. The plate was processed using the automated WGA procedure on the Biomek FX workstation. All samples were then subjected to amplification and 6 μ l of the resultant products were electrophoresed on a 1% agarose gel. Amplified products were not observed in the wells containing water.

XIII. Troubleshooting

Problem	Cause	Solution
Low yield after cycling	Sample contains PCR inhibitors, or high buffer salts.	Dialyze samples in a suitable microdialysis unit to remove the inhibiting components. Requantitate the dialyzed product as loss of DNA may occur in the dialysis process.
	Input DNA is degraded or less than 10 ng of DNA is present in the sample.	Amplification of insufficient DNA quantities often result in poor yield or poor representation in the final product. Some templates can be rendered amplifiable by increasing input DNA. Successful WGA amplification has been performed with degraded samples by increasing starting template from 25–100 ng.
	More enzyme is required.	WGA yield suffers when limiting amounts of DNA polymerase are used. We recommend a minimum of 12 units hot start DNA polymerase per 75 µl reaction. This is preferable to adding cycles as the resulting DNA may suffer from amplification bias.
	Liquid handler is not dispensing low volumes accurately.	Decrease the tip height from the bottom of the well in the dispensing steps so that the reagents are added into the solution. Check the performance of liquid handler at low volume range.
	The mixing of reagents is not sufficient due to inefficient mixing by the liquid handler.	Increase the aspiration and dispensing speed and/or cycle times in the mixing steps. Decrease the tip height from the bottom of the well in the mixing steps.
	Others	Refer to the Technical Bulletin of GenomePlex WGA Kit.
Negative control shows an amplified product.	Reagents are contaminated.	Use new labware and new batch of reagents. Test a reagent blank without DNA template to determine if the reagents are contaminated.

XIV. Contact Information

Technical Service
(800) 325-5832
email: techserv@sial.com

Customer Service
(800) 325-3010
(800) 588-9160
www.sigma-aldrich.com/order

GenomePlex and OmniPlex are registered trademarks of Rubicon Genomics, Inc.
Thermowell is a registered trademark of Corning, Inc.
PicoGreen is a registered trademark of Invitrogen, Inc.

JV,KTA 06/06-1

Sigma brand products are sold through Sigma-Aldrich, Inc.
Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications.
Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply.
Please see reverse side of the invoice or packing slip.