



Biomek Automated Genomic Sample Prep Accelerates Research

Biomek i-Series Automation of the Beckman Coulter GenFind V3 Blood and Serum DNA Isolation Kit

Introduction

The GenFind V3 Blood and Serum DNA Isolation Kit from Beckman Coulter isolates high quality genomic DNA from fresh or frozen whole blood and serum containing Citrate, EDTA, or Heparin anticoagulants, as well as from cultured cells. The kit produces a high recovery of DNA for downstream applications such as agarose gel analysis, restriction enzyme digestion, PCR amplification, and next-generation sequencing library construction. GenFind V3 uses an improved cell lysis buffer and Proteinase K treatment to rupture cell membranes and digest proteins. DNA is then immobilized using Beckman Coulter's patented SPRI (Solid Phase Reverse Immobilization) paramagnetic bead technology by adding a magnetic bead reagent. This differential binding of DNA enables separating DNA with the magnetic field, followed by rinsing away contaminants with multiple simple wash procedures, leaving the genomic DNA ready for elution. SPRI also enables fast separation, easy manipulation and simple automation compared to traditional vacuum filtration and centrifugation technologies. In this technical note, we demonstrate the automated performance of the Beckman Coulter GenFind V3 kit on the Biomek i7 Hybrid Genomics Workstation.

When compared to manual operations, the Beckman Coulter GenFind V3 kit automated on Biomek platforms provide

- Reduced hands-on time and increased throughput
- Option to run the method end-to-end with only setup and tear-down touch points
- Reduction in pipetting errors
- Standardized workflow for improved results
- Quick implementation with demonstrated methods
- Knowledgeable support for reagents, automation and methods all from a single vendor

Spotlight

Biomek i7 Hybrid Genomics Workstations

System features provide the highest throughput and flexibility of all Biomek Genomics workstations for maximized efficiency and reliability to increase user confidence and walk-away time

- 300 μ L or 1200 μ L Multichannel head with 1-300 μ L and 1-1200 μ L pipetting capability
- Span-8 with 1-1000 μ L pipetting capability
- Enhanced Selective Tip pipetting to transfer custom array of samples
- Two Independent 360° rotating gripper with offset fingers
- 45 positions
- Orbital Shakers, Peltiers and Tip washing for controlling sample processing
- Optional Enclosure



Figure 1. Biomek i7 Hybrid Genomics Workstation with optional enclosure on a Biomek Cart. The deck has a maximum capacity of 45 positions to maximize efficiency and increase walk away time

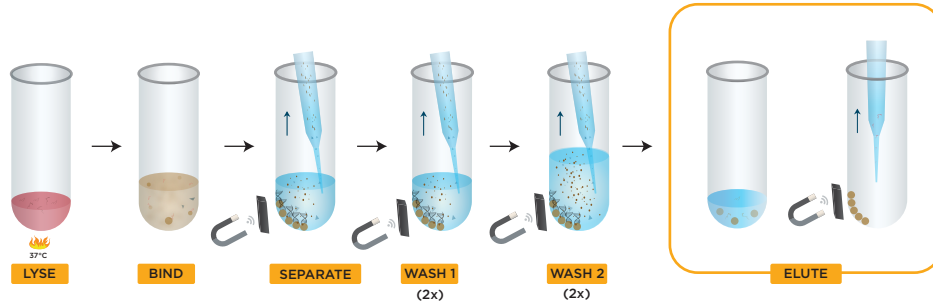


Figure 2. Beckman Coulter GenFind V3 kit protocol

Automated Method

We automated the Beckman Coulter GenFind V3 kit protocol on the Biomek i7 Hybrid, incorporating on-deck lysis. In general, the automated protocol enables DNA extraction from 1-96 blood and serum samples up to 400 μ L, in approximately 4 hours (Table 1).

The handling of blood and blood waste require special precautions and waste disposal procedures. Therefore, the automated GenFind V3 method isolates blood waste into waste plates, facilitating proper disposal of biohazardous waste.

Major Process Description	Automated/ Hands on Time	
	24 Samples	96 Samples
Sample Extraction - 400 μL input		
Prepare Reagents/ Set up Inst	30 min	30 min
Method Run Time	2 hr, 29 min	3 hr, 24 min
Total	2 hr, 59 min	3 hr, 54 min

Major Process Description	Automated/ Hands on Time	
	24 Samples	96 Samples
Sample Extraction - 200 μL input		
Prepare Reagents/ Set up Inst	30 min	30 min
Method Run Time	2 hr, 8 min	2 hr, 51 min
Total	2 hr, 38 min	3 hr, 21 min

* Timing does not include reagent and sample thawing, blood/serum collection, homogenization or dissection.

Table 1. . Estimated run times for Beckman Coulter GenFind V3 automated method on the Biomek i7 Hybrid.

Demonstrated Method Interface (DMI)

Three modules used to set up the GenFind V3 protocol provide the user with full instructions to set up the method and reduce set up errors.

1. Biomek Method Launcher (BML)

BML is a secure interface for selecting methods without affecting method integrity and manual control (Figure 3).

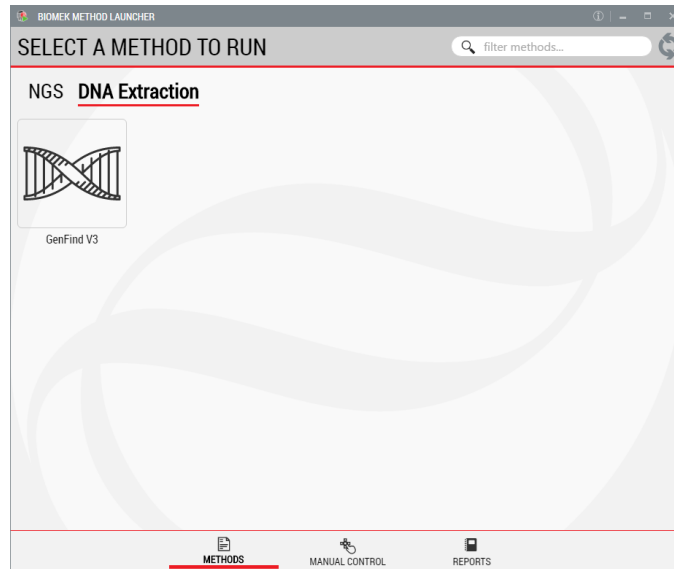


Figure 3. Biomek Method Launcher provides an easy interface to start the method

2. Method Options Selector (MOS)

MOS enables the selection of run-time options to maximize the flexibility in the set up of the method execution (Figure 4). For the GenFind V3 method, the automated protocol allows for 1-96 samples in the volume range of 40-400 μ L for processing.

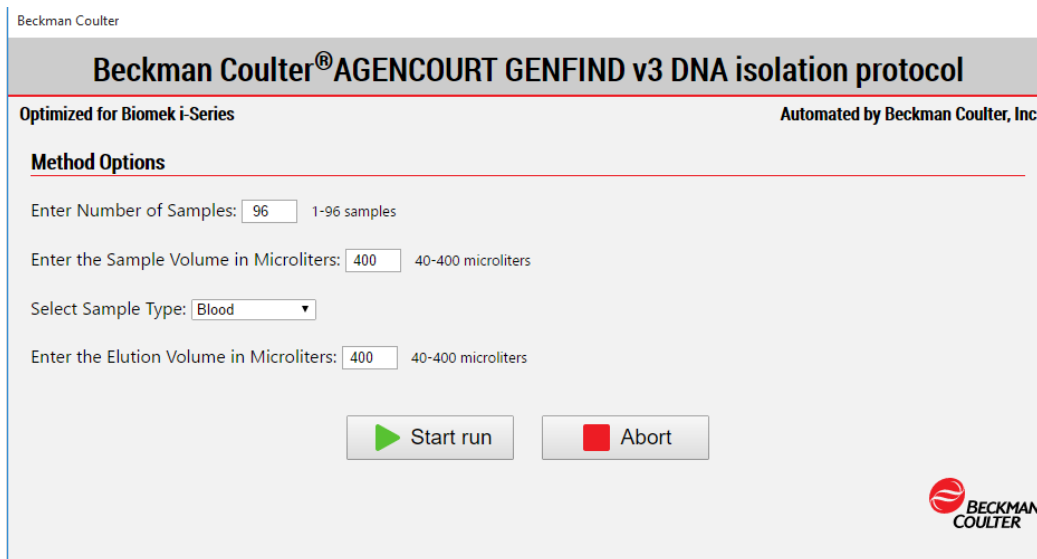


Figure 4. Biomek Method Options Selector indicates sample number and processing options

3. Guided Labware Setup (GLS)

GLS is generated based on user selected options in the MOS and provides the user specific text and step-by-step graphical setup instructions with reagent calculations (Figure 5).

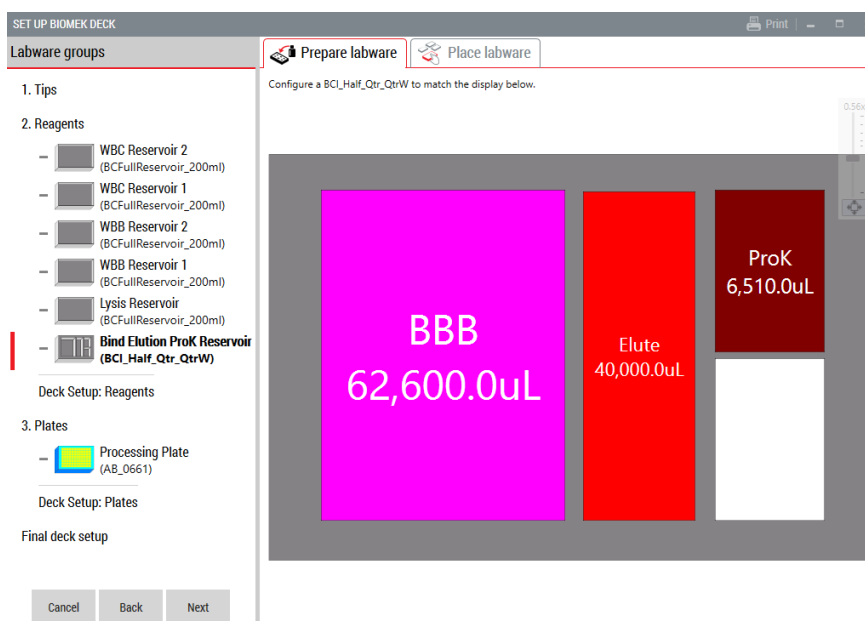


Figure 5. Guided Labware Setup indicates reagent volumes and guides the user for correct deck setup

Experimental Design

Whole blood was collected from four individual donors in blood collection tubes coated with EDTA, Citrate, and Heparin anticoagulants. Four technical replicates for each donor were used for a total of 48 samples of 400 μ L aliquots (16 EDTA, 16 Citrate, 16 Heparin). Genomic DNA was then extracted using the GenFind V3 Blood and Serum DNA Isolation automated method implemented on the Biomek i7 Hybrid platform. All extracted genomic DNA samples were analyzed using a NanoDrop™ 8000 (Thermo Fisher Scientific). 14 samples (four EDTA, five Citrate, and five Heparin) were randomly selected and analyzed using 2200 TapeStation using the Genomic DNA ScreenTape kit (Agilent Technologies). Those same 14 genomic DNA samples were then converted to NGS sequencing libraries using the Kapa HyperPlus Kit implemented on the Biomek i5 NGS Workstation. NGS library QC was performed using the 2100 Bioanalyzer with the High Sensitivity DNA kit (Agilent Technologies).

Results

The quantity as well as quality of DNA extracted from the automated protocol were comparable to those of the manual protocol (Figure 6). The DNA Integrity Number (DIN) for 14 selected samples were also calculated by TapeStation to further estimate DNA quality. Both methods indicate the extraction of high quality DNA, as indicated by their high DIN scores (Table 2).

To demonstrate the use of the extracted DNA in downstream applications, we constructed NGS libraries from the same 14 selected samples using the Kapa HyperPlus Kit. Successful library preparation was verified using a Bioanalyzer. Certain anticoagulants such as Heparin act as PCR inhibitors. The resulting traces indicate successful library construction and that the inhibitors were successfully removed through the GenFind V3 protocol (Figure 7).

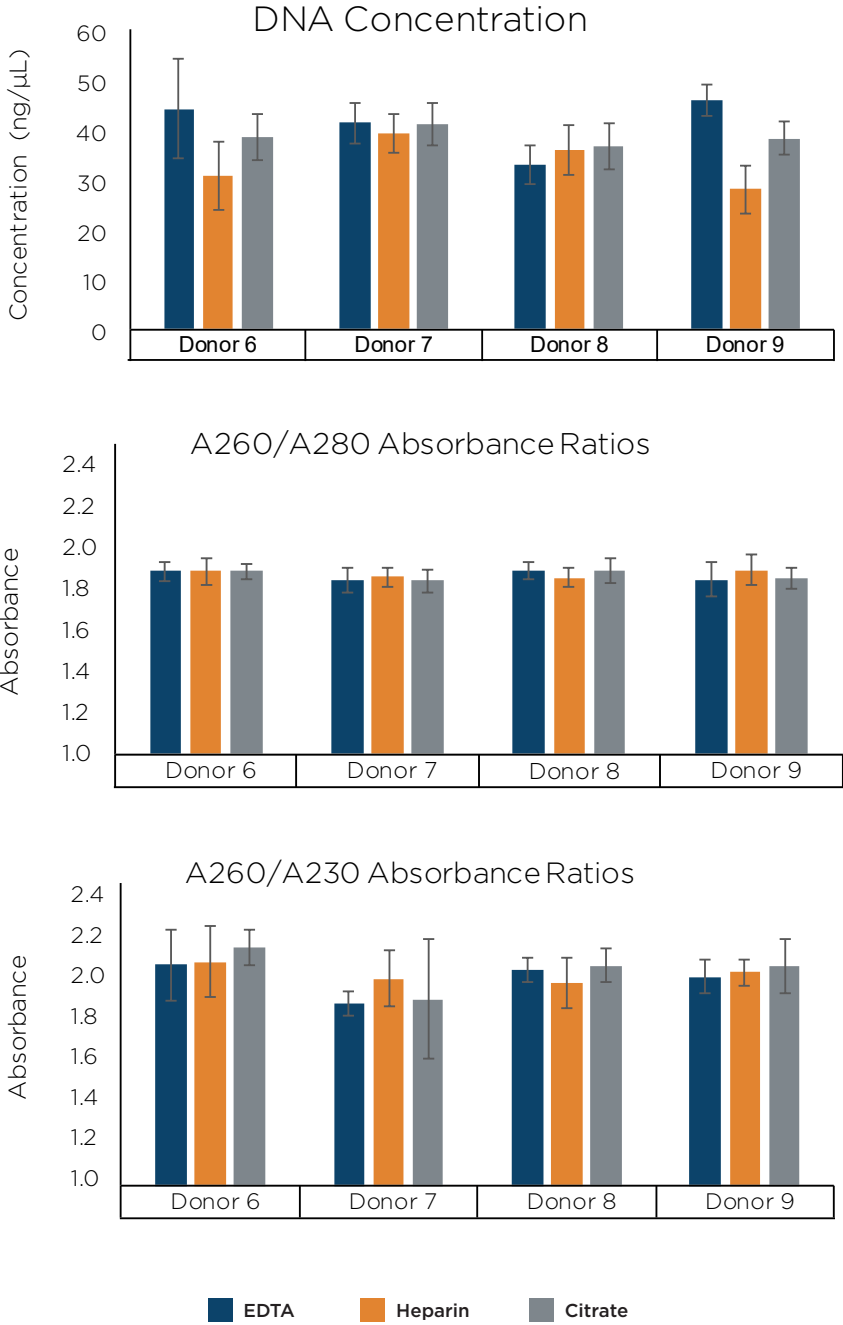


Figure 6. Average DNA concentration and purity ratios of replicate samples and blood tube types as indicated by NanoDrop 8000

Donor	Blood Tube Type	DIN
D6	EDTA	9.6
D6	Heparin	9.5
D6	Citrate	9.6
D7	EDTA	9.2
D7	Heparin	9.4
D7	Heparin	9.6
D7	Citrate	9.5
D8	EDTA	9.1
D8	Heparin	9.6
D8	Citrate	9.5
D8	Citrate	9.2
D9	EDTA	9.5
D9	Heparin	9.6
D9	Citrate	9.2

Table 2. DIN scores for randomly selected samples analyzed on Agilent TapeStation 2200

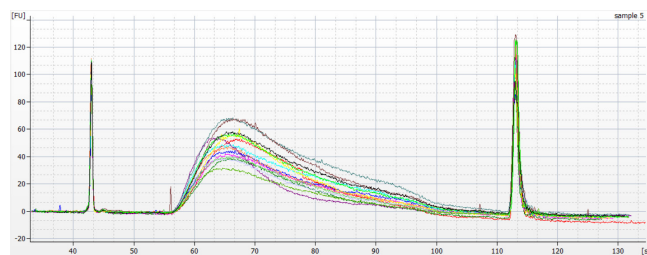


Figure 7. Bioanalyzer traces of NGS libraries constructed from samples in Table 2.

Summary

We have demonstrated automation of Beckman Coulter GenFind V3 on the Biomek i7 Hybrid Genomics Workstation results in high quality DNA and expected yields from whole blood. Furthermore, sequence-ready libraries can be generated when using these samples as input for NGS processes.

Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this protocol, including but not limited to warranties of fitness for a particular purpose or merchantability or that the protocol is non-infringing. All warranties are expressly disclaimed. Your use of the method is solely at your own risk, without recourse to Beckman Coulter. Not intended or validated for use in the diagnosis of disease or other conditions.

This protocol is for demonstration only, and is not validated by Beckman Coulter.

Biomek i-series are not labeled for IVD use and are not intended or validated for use in the diagnosis of disease or other conditions.

©2019 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

Content previously approved from AAG-5106PST03.19.



For Beckman Coulter worldwide office locations and phone numbers, please visit Contact Us at beckman.com

AAG-5653FLY07.19