

Biomek Automated NGS Solutions Accelerate Genomic Research

Biomek i-Series¹ Automated Illumina TruSeq² Exome Library Prep Kit

Introduction

The Illumina TruSeq Exome Library Prep kit is a workflow to generate indexed paired-end libraries and is specially designed to work with Formalin Fixed Paraffin Embedded (FFPE) samples. Solid Phase Reversible Immobilization (SPRI) bead chemistry is utilized throughout the protocol for size selection and sample cleanup steps. The process can be laborious and error-prone and is therefore ideal for Biomek automation. Choose from low to high throughput Biomek liquid handling workstations to automate all steps of the workflow providing the option for a complete walk-away solution. The methods are proven to deliver sequence-ready libraries through demonstrated performance using scientifically relevant samples. These ready-to-implement methods are compatible with Illumina TruSeq Exome Library Preparation kits (Illumina P/N FC-150-001, FC-150-002 or FC-150-003).

This application is automated on multiple Biomek workstations and provides:

- Standardized workflow for improved results
- Reduction in costly errors
- Reduced hands-on-time and increased throughput
- Quick implementation with ready-to-implement methods available delivered by knowledgeable support teams

Spotlight: Biomek i7 Hybrid Genomics Workstation

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

- 300 uL or 1200 uL Multichannel head with 1-250 uL and 1-1000 uL pipetting range
- Span-8 with 0.5-5000uL 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. 45 position deck capacity maximizes efficiency and increases walk away time.



Demonstrated Method Interface (DMI):

Three simple modules that provide the user full instructions to better ensure error-free method setup and provides users maximum flexibility for scheduling their day

1. **Biomek Method Launcher (BML)** — secure interface for selecting methods without affecting method integrity and manual control



Figure 2. Biomek Method Launcher provided an easy interface to start the method



Figure 3. Manual Control can be run through the launcher interface

2. **Method Options Selector (MOS)** — Select run-time options and maximize flexibility in daily scheduling and method execution

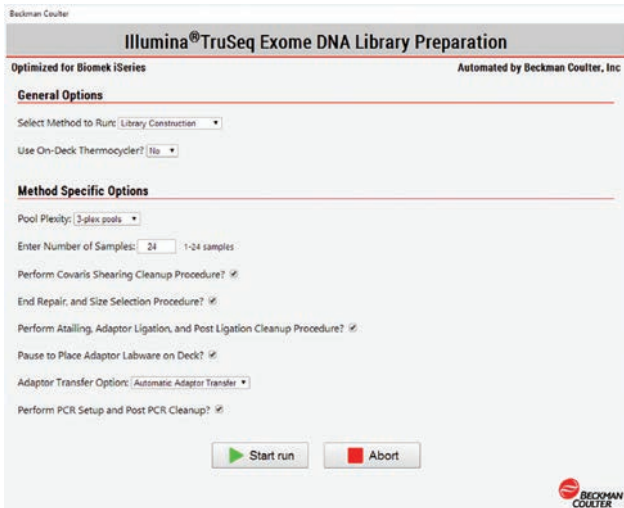


Figure 4. Illumina TruSeq Exome Method Options Selector showing the different features and run options. The method is broken into modules based on Illumina's recommended stop points, so the user has the flexibility to run specific modules or full method.

3. **Guided Labware Setup (GLS)** — Generated based on options selected in the MOS, and provides the user specific text and graphical setup instructions with reagent calculation

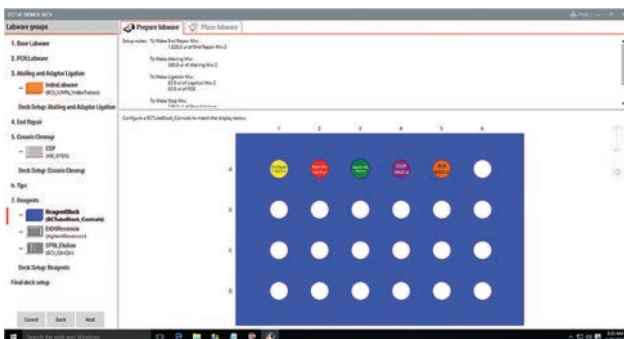


Figure 5. Guided Labware Setup showing reagent volumes and setup notes and guides the user for correct deck setup

Modular Design:

Built-in flexibility for scalable throughput and scheduling.

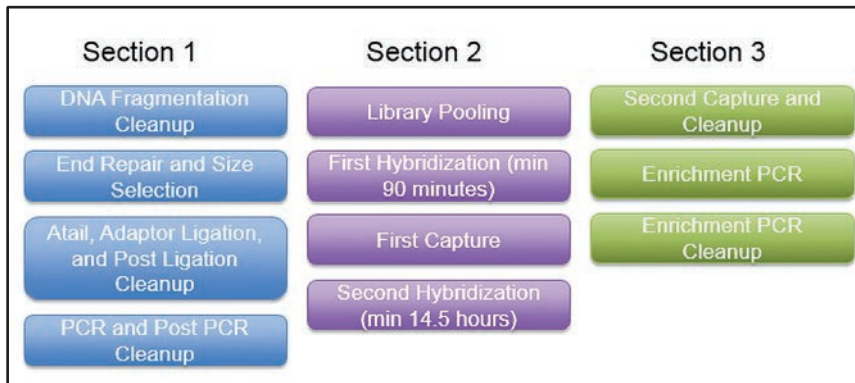


Figure 6. Illumina TruSeq Exome workflow.

Major Process Description for 96 samples/8 pools*	Section 1	Section 2	Section 3
TruSeq Exome	6 hrs 23 mins	2 hrs 16 min	3 hrs 28 mins
Hands On Time	30 mins	15 mins	15 mins
**Timing does not include thawing of reagents, hybridization or Library QC.			

Table 1. Illumina TruSeq Exome turnaround time on i7 Hybrid Genomics Workstation. The method is modular design based on Illumina's recommended stop points. Users can choose to run between 1-96 samples and 1 to 8 pools.

Experimental Design and Results³

DNA was extracted from FFPE samples using FormaPure DNA Extraction Kit from Beckman Coulter. Horizon Quantitative Multiplex Reference Standard for FFPE (HD200), breast (FFPE 76 and FFPE 96), lung (FFPE 110 and FFPE 121) and liver (FFPE 107) blocks were used to generate libraries using Illumina's TruSeq Exome Kit and sequenced with NextSeq. Library quality was assessed using the Agilent TapeStation High Sensitivity D5000 kit.

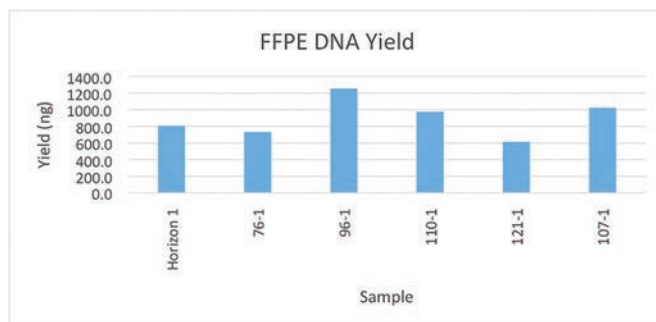


Figure 7. Library Yield in ng as measured by Quant-IT Picogreen assay.

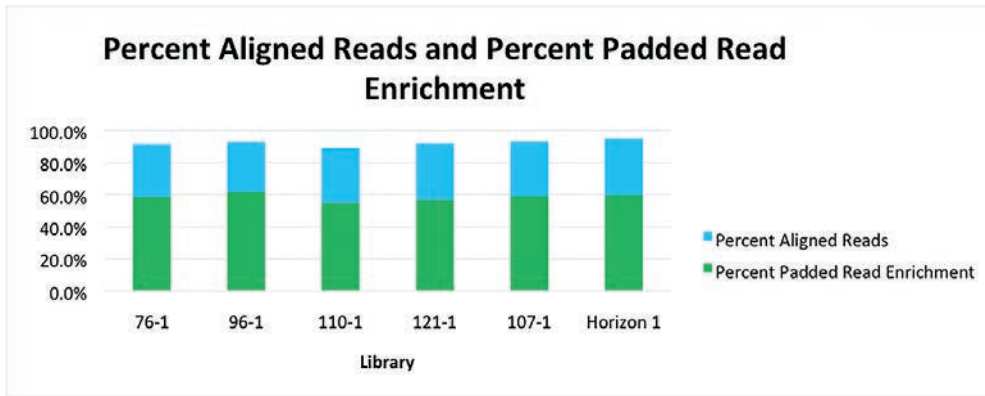


Figure 8. FFPE TruSeq Exome Sequence Alignment Results

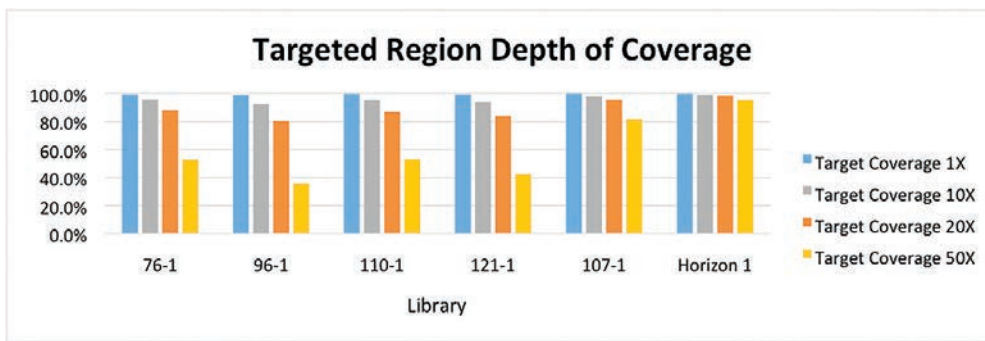


Figure 9. FFPE TruSeq Exome Depth of Targeted Region Analysis

Summary

The sequencing run generated 112.9M passed filter reads with 96% of those reads identified. 80% of targeted regions were covered to the a depth of at least 20X for the FFPE block samples, while over 95% of targeted regions for the Horizon sample were covered at a depth of 20X. In conclusion, Biomek i-Series liquid handlers offer a flexible solution for complex NGS workflows with challenging sample types.

1. Product in development
2. All trademarks are properties of their respective owners
3. Data obtained during development

For research use only, not for use in diagnostic procedures.

©2016 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the Stylized Logo, and Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.



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

AAG-2356FLY02.17