

Method Biography for Illumina TruSight™ Oncology 500 DNA Only on Biomek NGenius System

Presented by Zach Smith | 1/26/2024

This method has been demonstrated for use on the Biomek NGenius System but has not been validated by Beckman Coulter for use in the diagnosis of disease or other clinical conditions. All products are for Research Use Only. Not for use in Diagnostic Procedures.

Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this App Template, including but not limited to warranties of fitness for a particular purpose or merchantability or that the App Template is non-infringing. All warranties are expressly disclaimed. Your use of the App Template is solely at your own risk, without recourse to Beckman Coulter. This App Template has been demonstrated for use on the Biomek NGenius system for the chemistry kit version and the release date shown at the time when the App is selected and created but has not been validated by Beckman Coulter for use in the diagnosis of disease or other clinical conditions.

©2024 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein, including Biomek and Biomek NGenius, 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.

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

Scoping

Illumina TruSight Oncology 500 DNA Only Kit

Kit Highlights:

The kit is optimized to provide high sensitivity and specificity for low-frequency somatic variants across 523 genes.

DNA biomarkers include the following:

- Single nucleotide variants (SNVs)
- Insertions
- Deletions
- Gene amplifications
- Multinucleotide variants (MNVs)

TruSight Oncology 500 also detects immunotherapy biomarkers for tumor mutational burden (TMB) and microsatellite instability (MSI) in DNA.

Figure 1 TruSight Oncology 500 Kit DNA Only Workflow (Part 1)

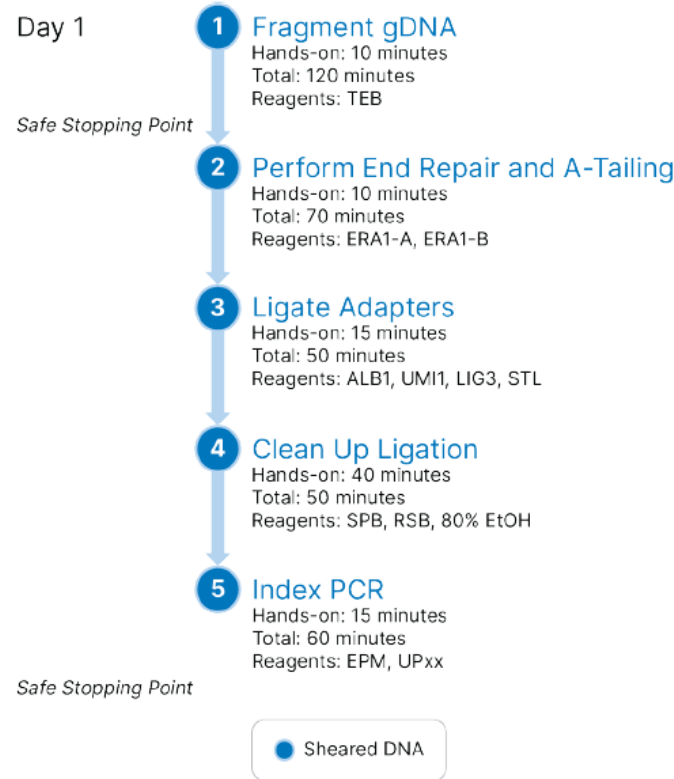
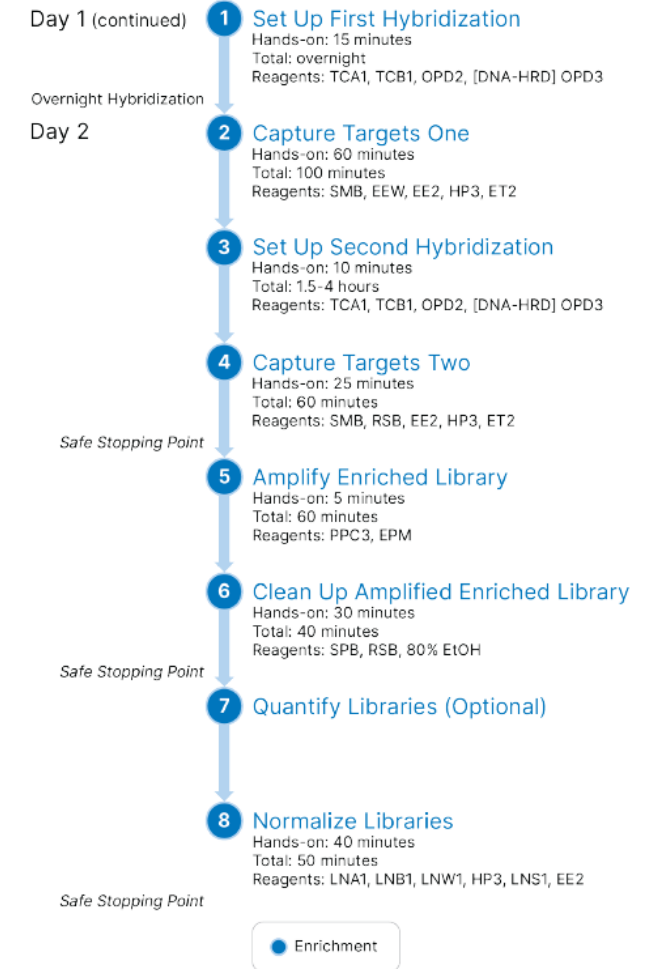


Figure 2 TruSight Oncology 500 Kit DNA Only Workflow (Part 2)



Above: Illumina TSO 500 DNA Only Workflow (adapted from TruSight Oncology 500 Reference Guide)

App Template Description

The Illumina* TruSight* Oncology 500 DNA Automation Kit App Template allows for the creation of Illumina TruSight Oncology 500 DNA Only libraries compatible with Illumina sequencing platforms. The App Template allows the user to produce between 4 to 16 libraries in a single batch utilizing pre-sheared DNA samples. Refer to the Illumina TruSight Oncology 500 Reference Guide for details concerning DNA input mass, sample quality, and shearing parameters. Combined DNA/RNA workflow is not supported by the App Template. Supplementary HRD enrichment workflow is not supported by the App Template. 80% ethanol wash volumes have been reduced to 50 μ L from 200 μ L to reduce tip consumption. First and Second Hybridization times have been limited to the minimum time listed in the manual protocol (8 hours for First Hybridization) to reduce application run time and cannot be changed. The App Template was designed using the Illumina TruSight Oncology 500 Reference Guide (Document# 1000000067621 v10). The App Template utilizes the TruSight Oncology 500 DNA Automation Kit (16 indexes, 64 Samples) (Illumina Part Number 20045504).

*Illumina and TruSight are registered trademarks of Illumina, Inc.

Document ID 2023-GBL-EN-101637-v1

TSO500 DNA Only Workflow on Biomek NGeniusS System

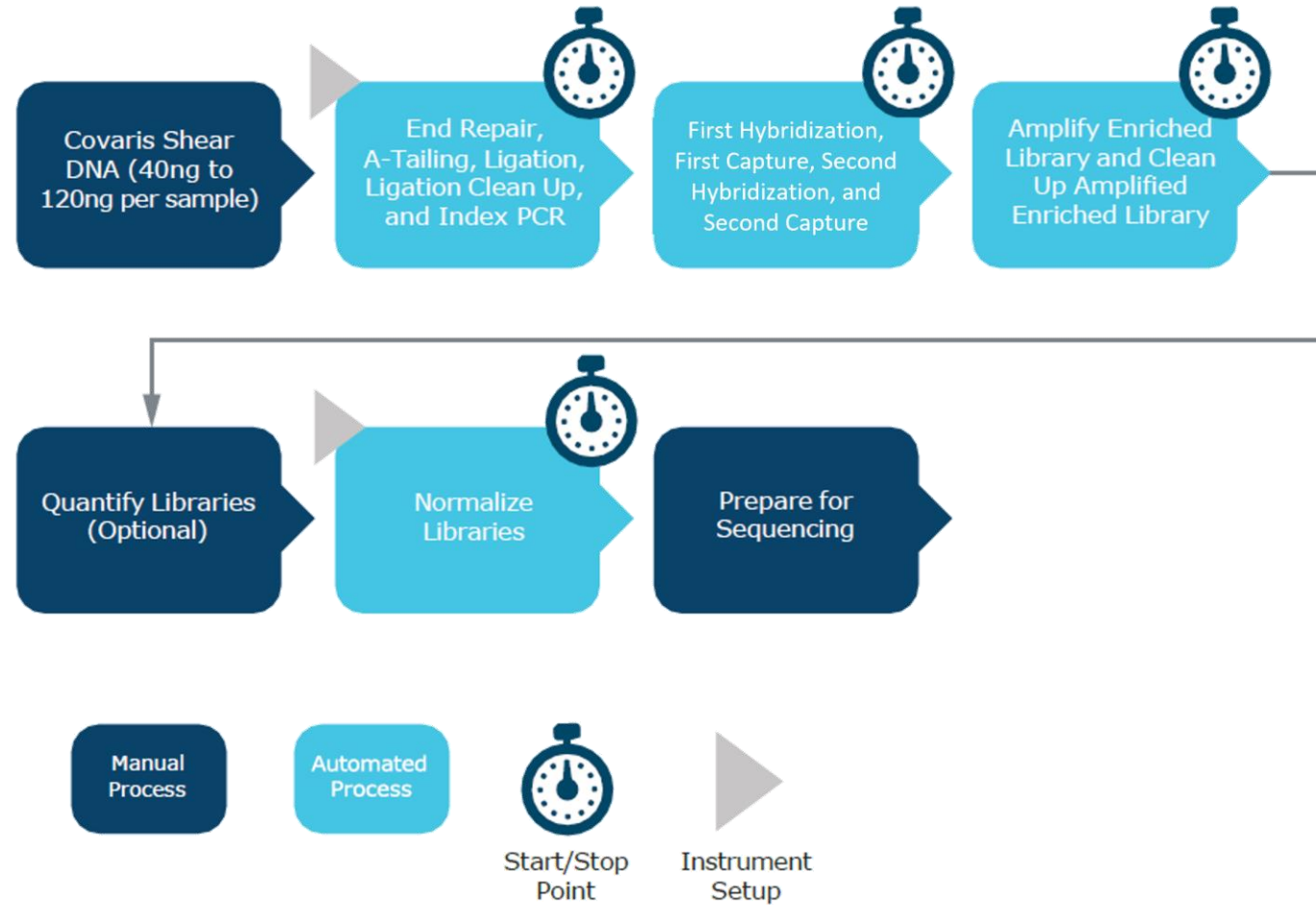
Features of the TSO500 DNA Only App

- Process any number of samples from 4 to 16, including odd numbers of samples.
- Can process a 16-sample batch with samples on sequencer by the end of Day 2.
- Built for the TruSight Oncology 500 DNA Automation Kit (16 indexes, 64 Samples) (Illumina Part Number 20045504).
- Can process up to four 16-sample batches from a single kit.

Not Included in the TSO500 DNA Only App

- Support for processing RNA and DNA samples in the same batch.
- Support for the supplemental HRD workflow.

TSO500 DNA Only Workflow on Biomek NGeniusS System



Expectation (From Reference Guide and Other Materials)

Metric Location	Metric Source	Results Metrics
Pre-Sequencing	Illumina TruSight Oncology 500 Reference Guide (1000000067621 v10)	Pre-library normalization yield should be 3ng/ul or higher as determined by Qubit.
Post Sequencing	TSO500 DRAGEN Evaluation App MetricsOutput.tsv	Sequencing run percent pass filter reads should be 80% or higher
Post Sequencing	TSO500 DRAGEN Evaluation App MetricsOutput.tsv	Post Sequencing: Contamination Score should be less than 1457
Post Sequencing	TSO500 DRAGEN Evaluation App MetricsOutput.tsv	Median Insert Size should be greater than or equal to 70bp
Post Sequencing	TSO500 DRAGEN Evaluation App MetricsOutput.tsv	Median Exon Coverage (Count) should be greater than or equal to 150X
Post Sequencing	TSO500 DRAGEN Evaluation App MetricsOutput.tsv	Percent Exon Coverage 50% should be greater than or equal to 90%
Post Sequencing	TSO500 DRAGEN Evaluation App MetricsOutput.tsv	Usable MSI Sites (Count) should be greater than or equal to 40
Post Sequencing	TSO500 DRAGEN Evaluation App MetricsOutput.tsv	Gene_Scaled_MAD (Count) should be less than 0.134
Post Sequencing	TSO500 DRAGEN Evaluation App MetricsOutput.tsv	Median Bin Count CNV Target (Count) should be greater than or equal to 1
Post Sequencing	TruSight Oncology 500 and TruSight Oncology 500 High-Throughput Data Sheet (M-GL-00173 v5.0)	5% Variant Allele Frequency (VAF) for small variants
Post Sequencing	TruSight Oncology 500 and TruSight Oncology 500 High-Throughput Data Sheet (M-GL-00173 v5.0)	2.2X fold-change for Copy Number Variants (CNV)
Post Sequencing	TruSight Oncology 500 and TruSight Oncology 500 High-Throughput Data Sheet (M-GL-00173 v5.0)	Analytical sensitivity greater than 96% for all variant types at 5% VAF or higher

Operating the Illumina TruSight Oncology 500 DNA Only App Template

TSO500 DNA Only Workflow on Biomek NGeniusS System

Batches can be executed with a minimum of three user interactions

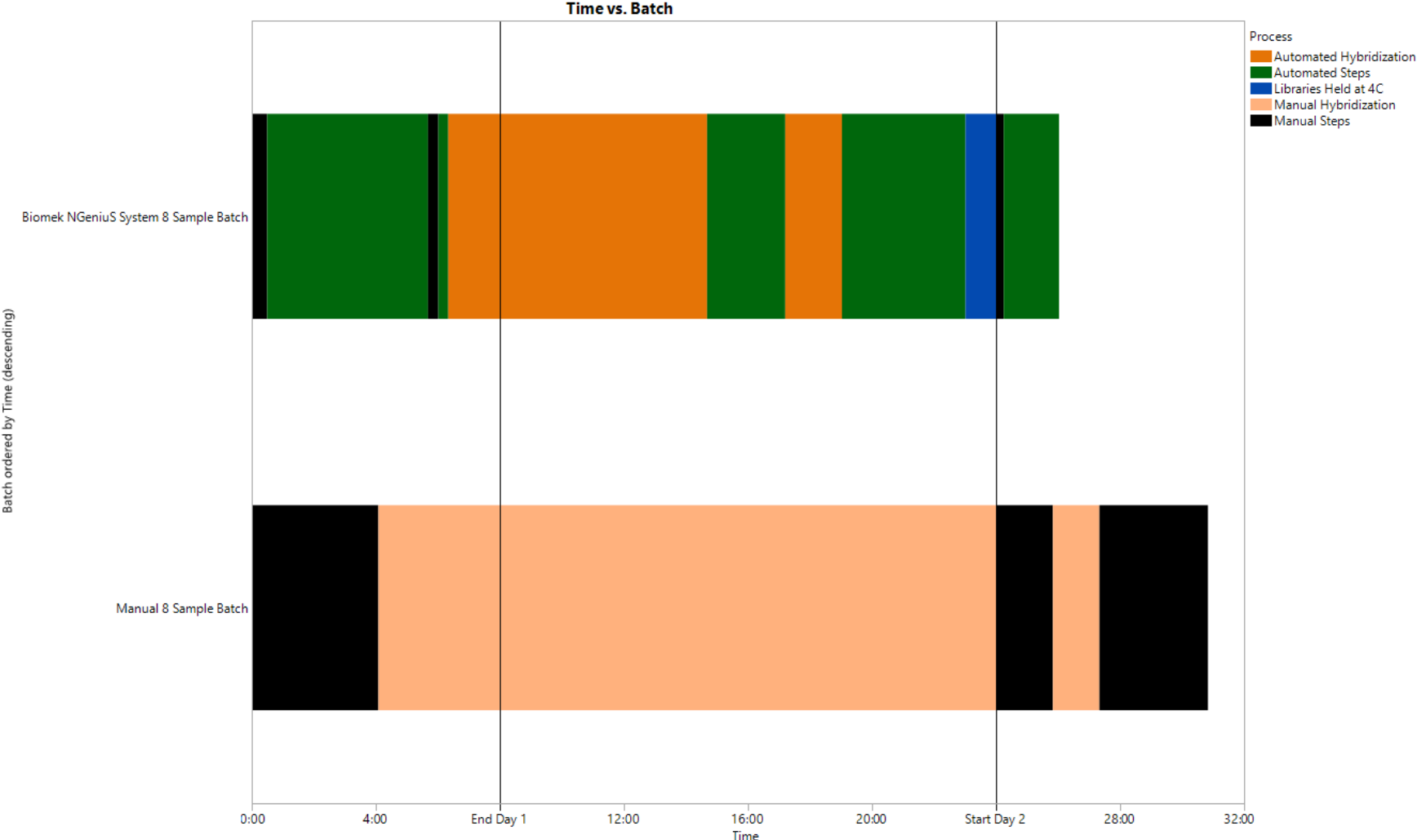
Sections		
Start at section		
1. End Repair, A-Tailing, Ligate Adapters, Clean Up Ligati...		
#	Section	Status
1	End Repair, A-Tailing, Ligate Adapters, Clean Up Ligation, and Index PCR	—
2	Retrieve ALS Product From Cold Storage 3	Off System
3	First Hybridization, Capture Targets One, Second Hybridization, and Capture Targets Two	—
4	Amplify Enriched Libraries and Clean Up Amplified Enriched Library	—
5	Quantify Libraries Off System	Off System
6	Normalize Libraries	—

Sections		
Start at section		
3. First Hybridization, Capture Targets One, Second Hyb...		
#	Section	Status
1	End Repair, A-Tailing, Ligate Adapters, Clean Up Ligation, and Index PCR	Skipped
2	Retrieve ALS Product From Cold Storage 3	Off System
3	First Hybridization, Capture Targets One, Second Hybridization, and Capture Targets Two	—
4	Amplify Enriched Libraries and Clean Up Amplified Enriched Library	—
5	Quantify Libraries Off System	Off System
6	Normalize Libraries	—

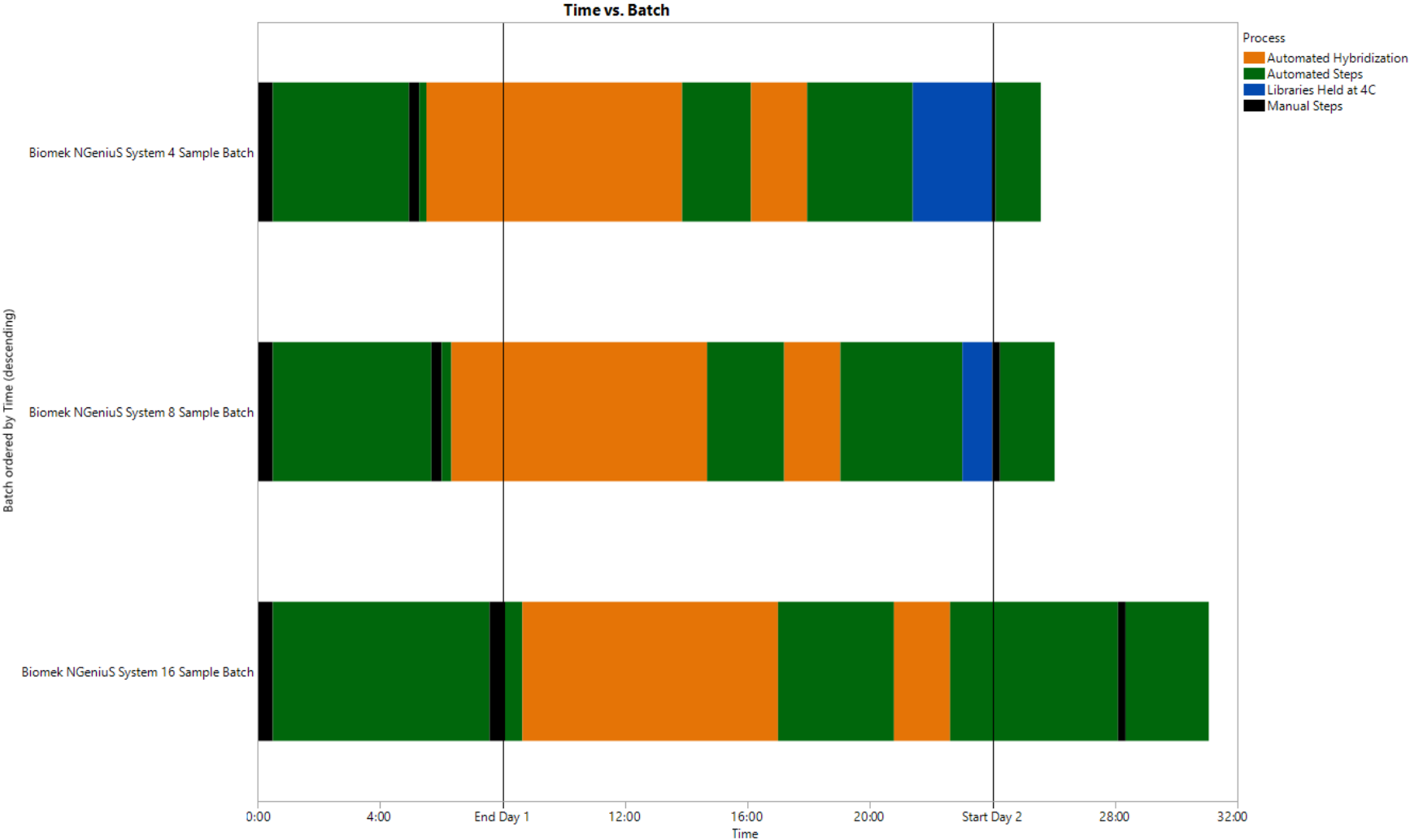
Sections		
Start at section		
6. Normalize Libraries		
#	Section	Status
1	End Repair, A-Tailing, Ligate Adapters, Clean Up Ligation, and Index PCR	Skipped
2	Retrieve ALS Product From Cold Storage 3	Off System
3	First Hybridization, Capture Targets One, Second Hybridization, and Capture Targets Two	Skipped
4	Amplify Enriched Libraries and Clean Up Amplified Enriched Library	Skipped
5	Quantify Libraries Off System	Off System
6	Normalize Libraries	—

There are two off-system sections that require a user interaction. In the first one, the user is shown in the App Template Instructions for Use how to retrieve the ALS product that isn't being used for the first hybridization. The second off-system section is for the quantification of the libraries prior to bead-based library normalization.

TSO500 DNA Only on Biomek NGeniusS System Run Time and Hands-On Time



TSO500 DNA Only on Biomek NGeniusS System Run Time and Hands-On Time



TSO500 DNA Only on Biomek NGenius System Consumables Required By Batch Size

Part Number	Name	Number Required	Number Required	Number Required	Quantity of Part Number to Purchase
		by App Template (tip boxes, reaction vessels, etc.) for 4 Samples	by App Template (tip boxes, reaction vessels, etc.) for 8 Samples	by App Template (tip boxes, reaction vessels, etc.) for 16 Samples	
C62705	Biomek Reaction Vessel, 24 well	19	19	19	1
C62706	Biomek Lid, 24 well	15	15	15	1
C62707	Biomek Bulk Reservoir, 25 mL/section	5	5	5	1
C59585	Case, Tips, 1025µL, Conductive, Filtered, 480 Tips	1	3	6	1
C62712	Case, Tips, 70 µL, Conductive, Filtered, 3840 Tips	2	2	4	1
C70672	Case, Biomek 1025 µL Tip Box, Empty, 5 Racks	1	1	1	1
C70673	Case, Biomek 70 µL Tip Box, Empty, 10 Racks	1	1	1	1
C70665*	Biomek Seal Plate, 24 well	3	3	3	1

* For the updated seal plate information please see Notification letter – Biomek Seal Plate Revision

TSO500 DNA Only on Biomek NGeniusS System Cost per Sample

Manual Processing	Per Sample*
Library Prep Kit	\$389.96
Plastic Consumables	\$46.34
Total Consumables	\$436.30

Biomek NGeniusS System Processing	Per Sample*
Library Prep Kit	\$389.96
Plastic Consumables	\$32.33
Total Consumables	\$422.29



Per Sample	
NGeniusS	\$422.29
Manual	\$436.30
Difference	-\$14.01



* Prices calculated for a 16-sample batch. Prices calculated as of Jan 2024; subject to change; for reference only

Batch runs per kit

Batch Size	4	5	6	7	8	9	10	11	12	13	14	15	16
Number of Batches per Kit	6	6	6	6	6	6	5	5	5	4	4	4	4
Samples	24	30	36	42	48	54	50	55	60	52	56	60	64
Largest Batch with Leftover Volume	16	16	16	11	5	0	7	0	0	9	5	0	0
Total Samples from Kit	40	46	52	53	53	54	57	55	60	61	61	60	64

The *Batch Size* can be run *Batches per kit* times, leaving enough reagent volume to do one additional batch with *Largest batch with leftover volume* samples.

Run combinations calculated based on published reagent vial volumes.

Demonstration Data

TSO500 DNA Only Demonstration Runs

TSO500 DNA Only 16-Sample Demo Run

Nine 40 ng Sera Care DNA shear replicates, two HD798 replicates, two HD799 replicates, two HD803 replicates, and one negative control (water) were used for the run.

TSO500 DNA Only 9-Sample Demo Run

Five 80 ng Sera Care DNA shear replicates, one 80 ng HD798 replicate, one 80 ng HD799 replicate, one 80 ng HD803 replicate, and one negative control (water) were used for the run.

TSO500 DNA Only 4-Sample Demo Run

Three 120 ng Sera Care DNA shear replicates and one negative control (water) were used for the run.

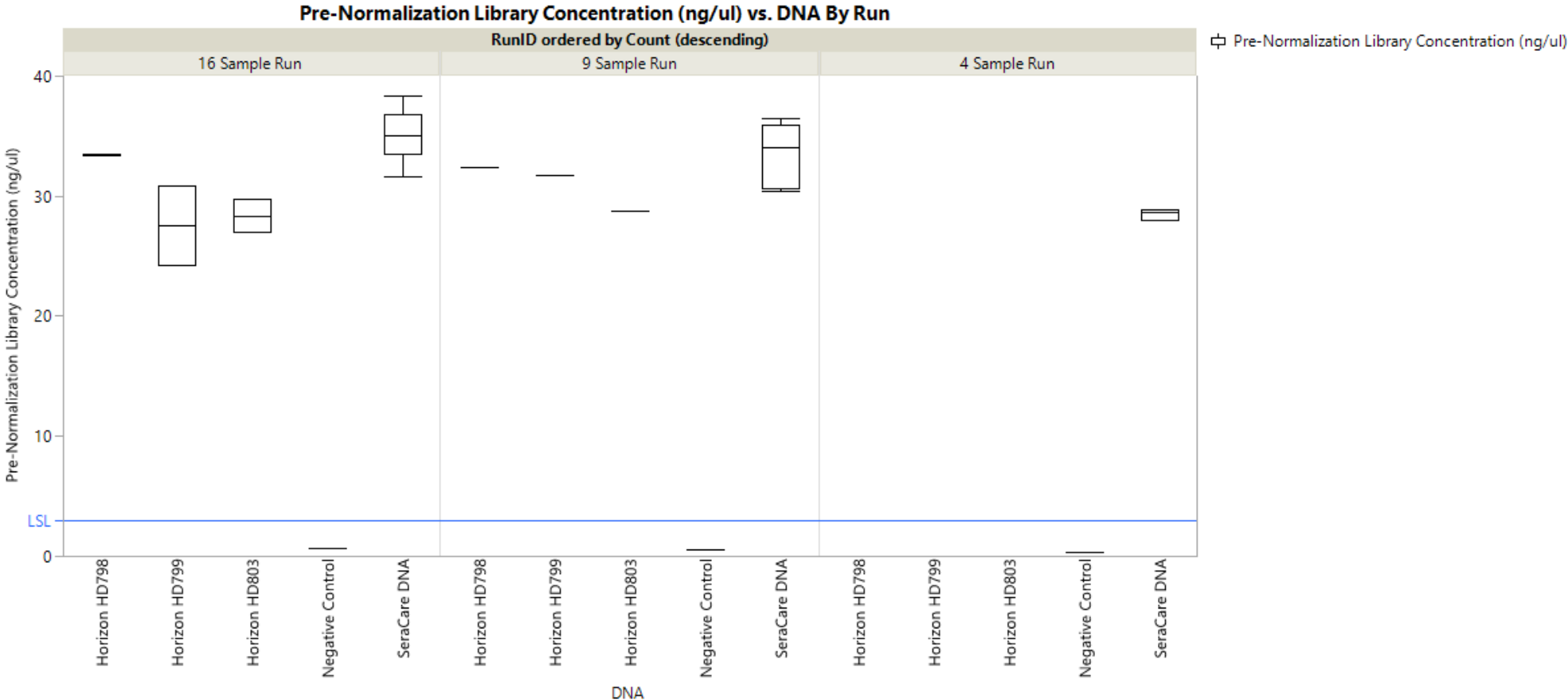
<https://www.seracare.com/Seraseq-TriLevel-Tumor-Mutation-DNA-Mix-v2-HC-0710-0097/>

<https://horizondiscovery.com/en/reference-standards/products/quantitative-multiplex-reference-standard-fcdna-mild?catalognumber=HD798>

<https://horizondiscovery.com/en/reference-standards/products/quantitative-multiplex-reference-standard-fcdna-moderate?catalognumber=HD799>

<https://horizondiscovery.com/en/reference-standards/products/quantitative-multiplex-reference-standard-fcdna-severe>

TSO500 DNA Only Demonstration Data: Pre-Normalization Library Concentration



Sequencing Runs

TSO500 DNA Only 16-Sample Demo Run

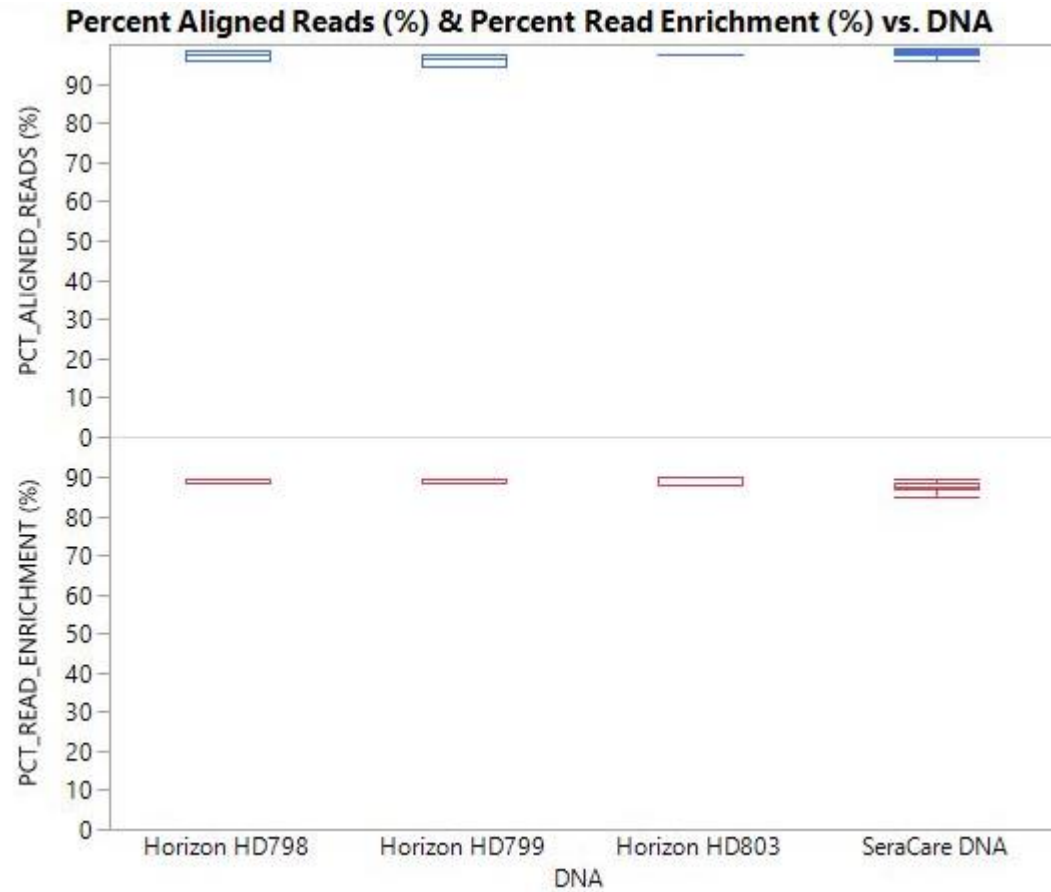
The sequencing run generated 205,866,836 pass filter reads (92.66% of total reads), of which 85% of reads were successfully identified. 94.35% of all bases were Q30 or higher. Total sequencing output was 88.25 Gb. Sequencing results should have 80% of reads pass filter with 80% of bases Q30 or higher.

TSO500 DNA Only 9-Sample and 4-Sample Demo Runs

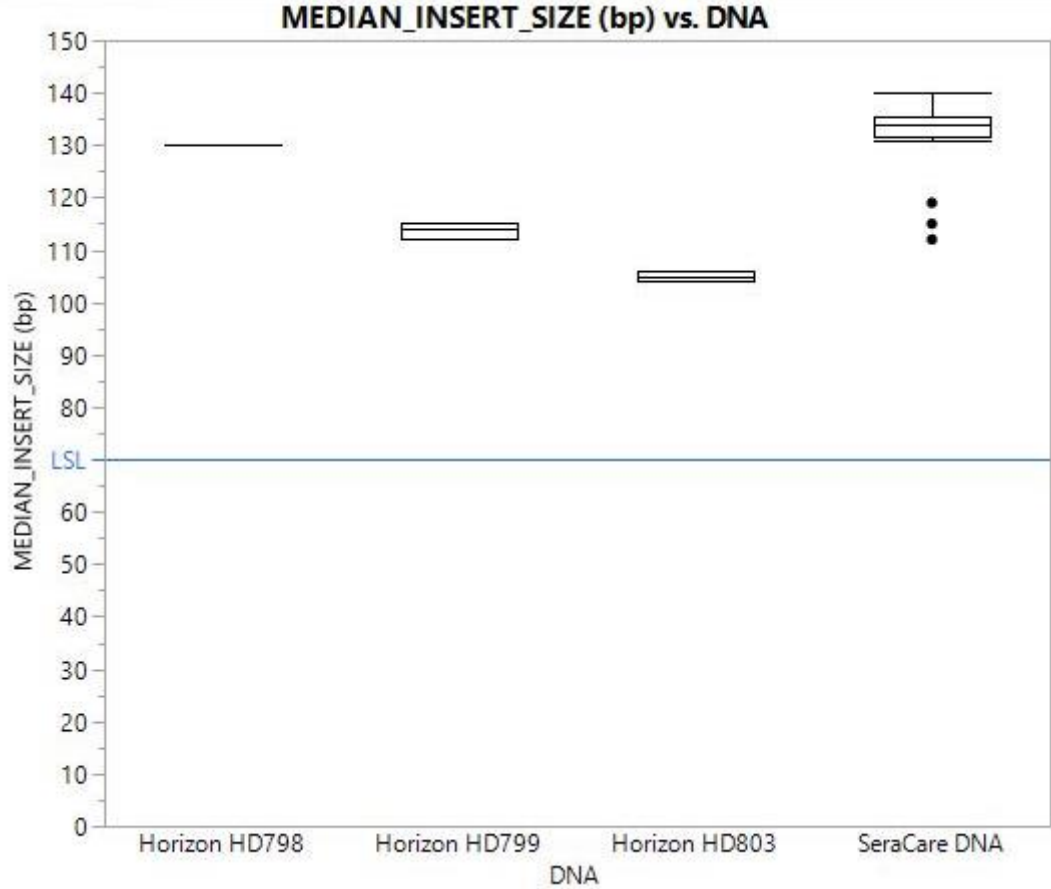
The sequencing run generated 271,579,046 pass filter reads (86.54% of total reads), of which 87.29% of reads were successfully identified. 91.37% of all bases were Q30 or higher. Total sequencing output was 115.45 Gb. Sequencing results should have 80% of reads pass filter with 80% of bases Q30 or higher.

Sequencing results were analyzed on BaseSpace using the TSO500 DRAGEN Evaluation App. The results were then entered into JMP and graphed along with Upper and Lower Spec Limits when appropriate.

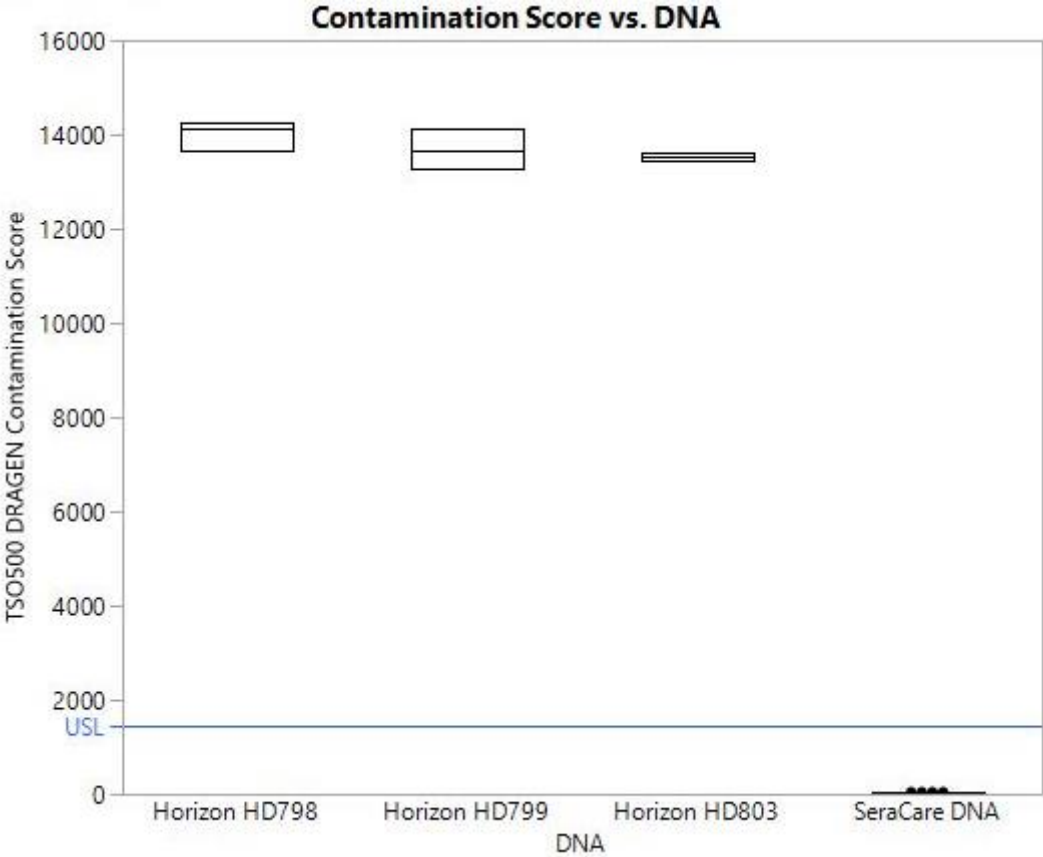
DNA Metrics



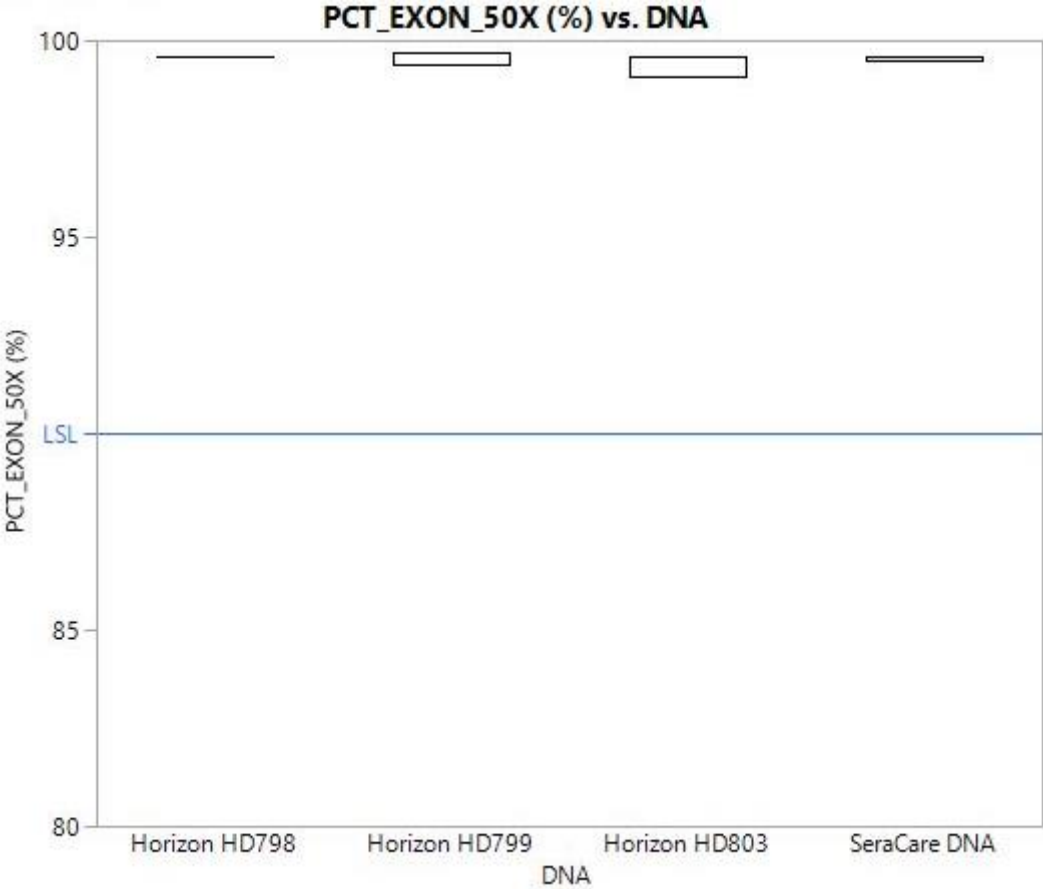
DNA Metrics



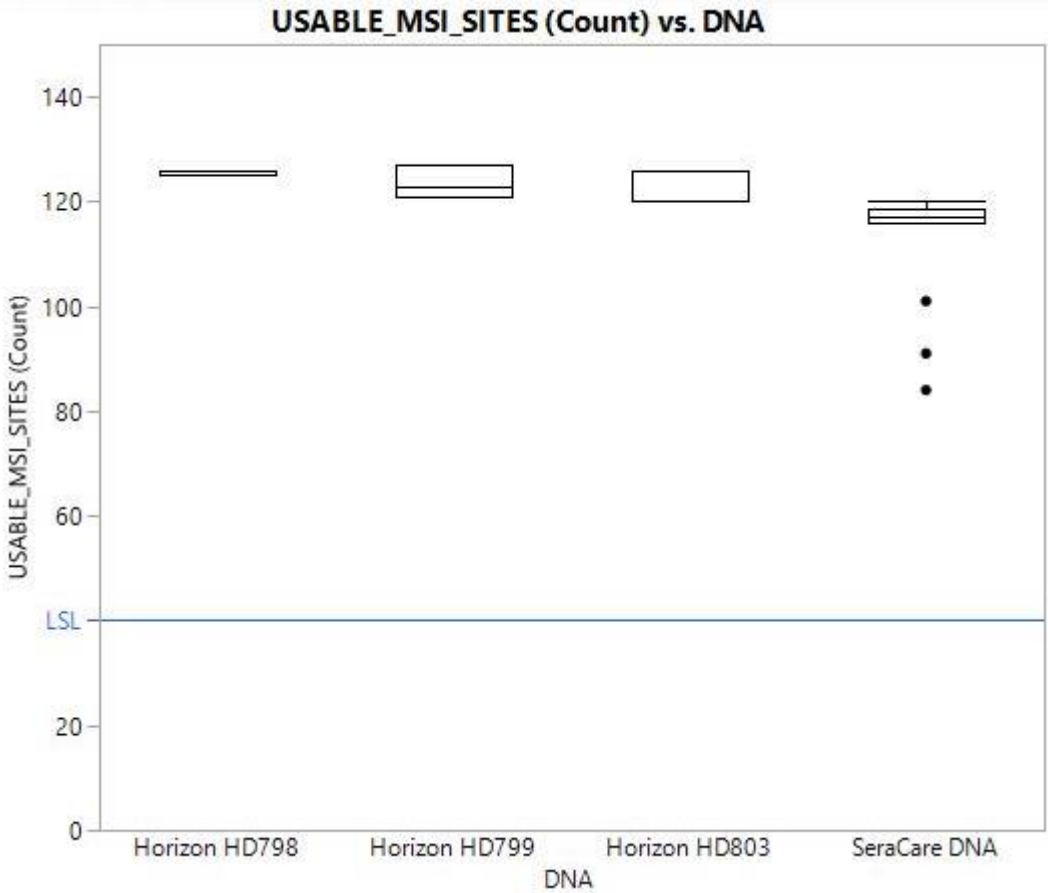
DNA Metrics



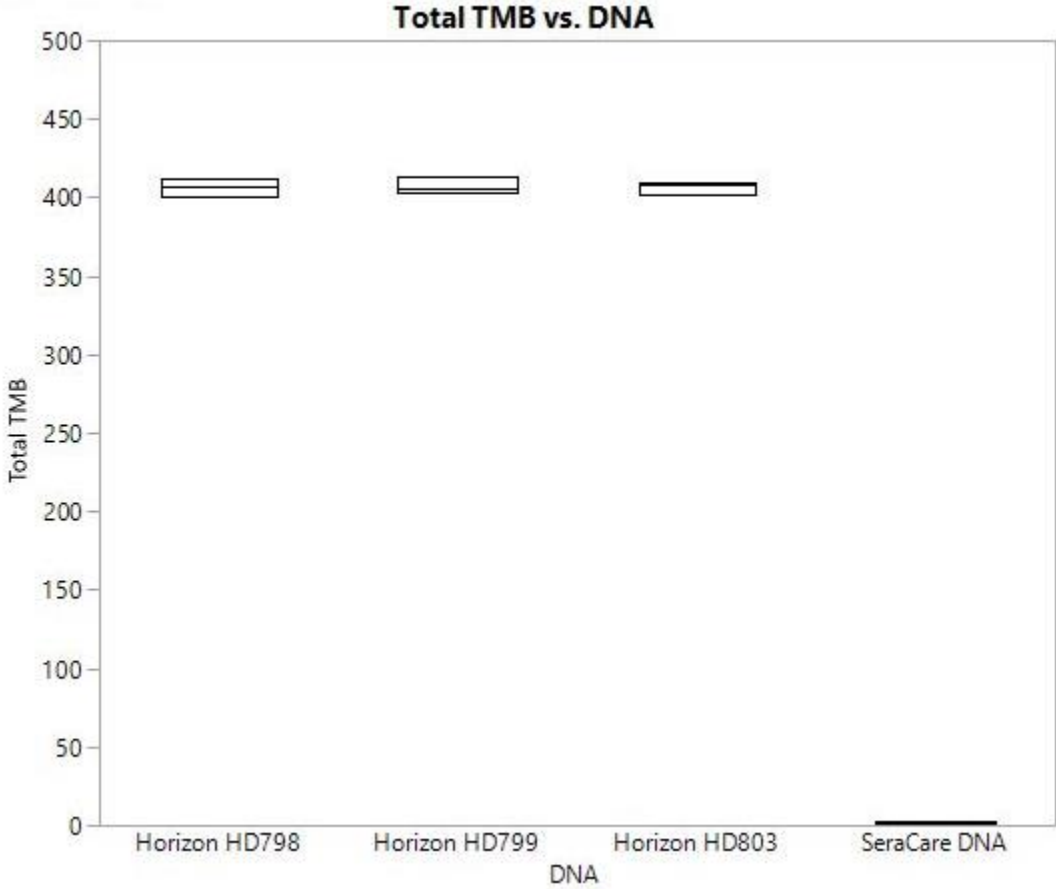
DNA Metrics



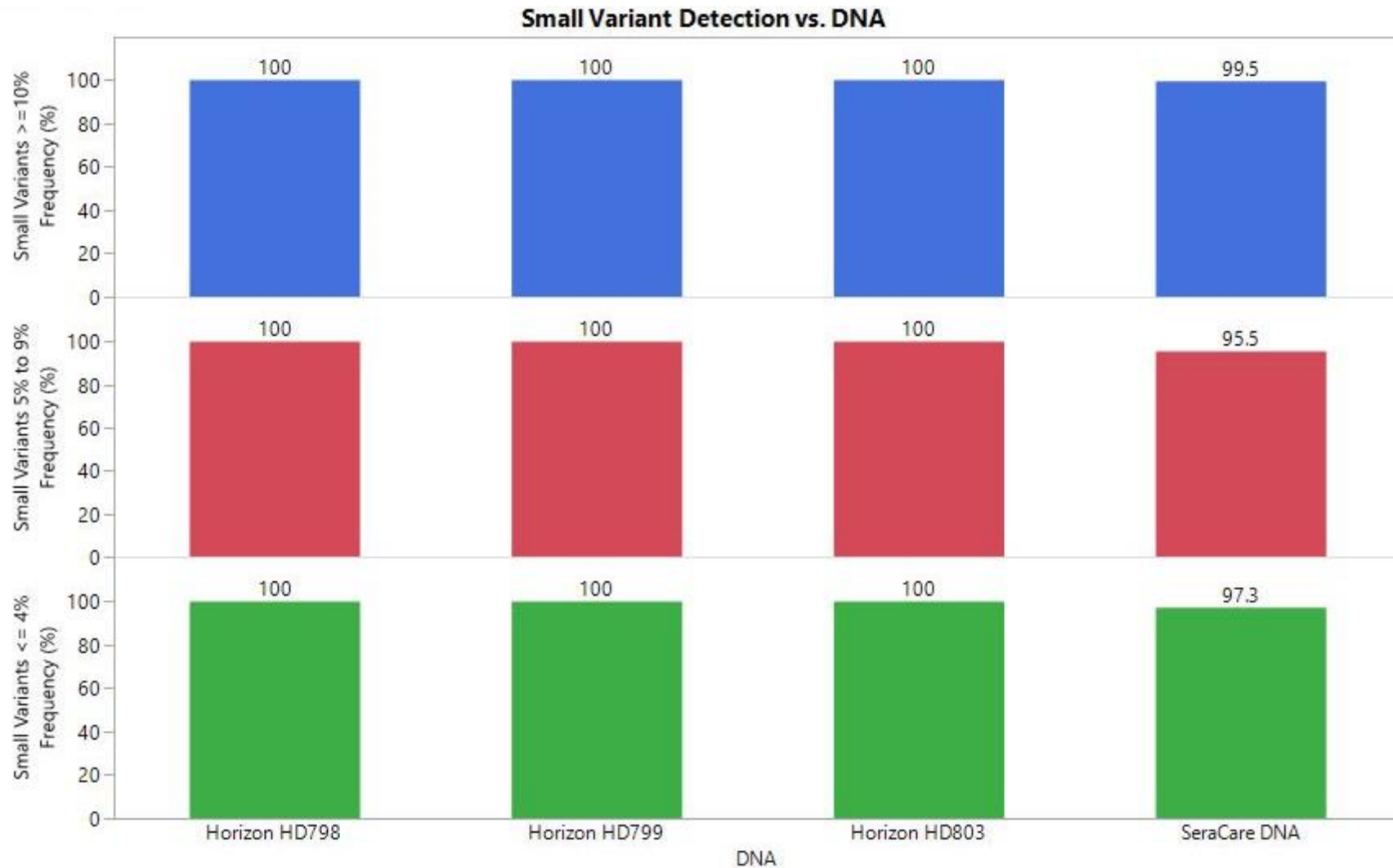
DNA Metrics



DNA Metrics



DNA Metrics: Sera Care and Horizon Discovery Small Variants



Metric Results

Results Metrics

DNA Libraries

Pre-Sequencing: Pre-library normalization yield should be 3ng/ul or higher as determined by Qubit.

Passed (26/26)

Post Sequencing: Sequencing run percent pass filter reads should be 80% or higher

Passed (85% and 87.2%)

Post Sequencing: Contamination Score should be less than 1457

Passed (26/26)*

Post Sequencing: Median Insert Size should be greater than or equal to 70bp

Passed (26/26)

Post Sequencing: Median Exon Coverage (Count) should be greater than or equal to 150X

Passed (26/26)

Post Sequencing: Percent Exon Coverage 50% should be greater than or equal to 90%

Passed (26/26)

Post Sequencing: Usable MSI Sites (Count) should be greater than or equal to 40

Passed (26/26)

Post Sequencing: Gene_Scaled_MAD (Count) should be less than 0.134

Passed (26/26)

Post Sequencing: Median Bin Count CNV Target (Count) should be greater than or equal to 1

Passed (26/26)

Post Sequencing: 5% Variant Allele Frequency (VAF) for small variants

Pass. Variants as low as 4% frequency detected

Post Sequencing: 2.2X fold-change for Copy Number Variants (CNV)

Passed (26/26)

Post Sequencing: Analytical sensitivity greater than 96% for all variant types at 5% VAF or higher

Pass (98.2% of all 10% and 7% variants identified)

*Horizon Discovery samples are known to have higher contamination scores than the 1457 upper limit due to the mix of cell lines used per Illumina.

Considerations

General automation considerations

- Do not use other sized kits as their reagents might come in an unexpected tube size
 - Chance for misread OCR token on vial
 - Chance for failed chemistry
 - Chance for damaged instrument
- Avoid bubbles in reagent tubes to ensure accurate liquid level sensing and aliquoting
 - Sample Purification Beads (SPB) and Streptavidin Magnetic Beads (SMB)
- Dead volume will be left behind in some storage wells
 - The nature of automation, tolerance stack-ups, and environment necessitates some overage
- Make sure bulk reagents wet the entire length of reservoir
 - Ensures accurate liquid volume sensing
- Prepare samples while Biomek NGenius System is aliquoting reagents
 - Avoids sample evaporation while Biomek NGenius System is preparing run

Conclusion

Illumina TruSight Oncology 500 DNA Only Libraries prepared on the Biomek NGeniusS Next Generation Library Prep System meets all specifications for successful sequencing.



This document contains proprietary information owned by Beckman Coulter, Inc.

Any disclosure, use, copying or distribution of this information without the express written consent of Beckman Coulter, Inc. is strictly prohibited.

© 2024 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.

