

Evaluation and Comparison of Liquid Biopsy Reference Materials from Commercial Sources using Oncomine™ Pan-Cancer Cell-free Assay

Ru Cao, Kris Lea, Jeff Schageman, Yanchun Li, Jian Gu, Priyanka Kshatriya, Varun Bagai, Khalid Hanif, Kelli S. Bramlett
Thermo Fisher Scientific, 2130 Woodward Street, Austin, Texas 78744

INSTRODUCTION

Over the last few years, analysis of liquid biopsies has grown rapidly to interrogate cell-free nucleic acid and characterize tumor mutation profiles from plasma. To date, a number of liquid biopsy research assays utilizing next generation sequencing technologies have been made commercially available to detect various types of variants at low allele frequency with minimal sample input. To ensure each assay is performed consistently and adopted successfully by researchers, reference materials are essential in assay validation and can be used as standards to compare technologies and standardize protocols across different laboratories.

MATERIALS AND METHODS

Reference material

Multiplex cfDNA reference standard set: this standard set (Horizon Discovery, Cat# HD780) contains 4 contrived DNA samples which mimic cfDNA type of material. Each sample covers 8 engineered single nucleotide variants (SNVs/SNPs) with allele frequency at 5%, 1%, 0.1%, and 0% (100% wild type), respectively.

Structural Multiplex cfDNA Reference Standard: this standard (Horizon Discovery, Cat# HD786) contains RET and ROS1 fusion variants, MYC and MET amplifications.

Seraseq® cfDNA complete™ mutation mix: four ctDNA mutation mix samples were provided by Seracare including wild type (Cat# 0710-0533), 1% (Cat# 0710-0530), 0.5% (Cat# 0710-0531), and 0.1% (Cat# 0710-0532). Each sample contains 25 unique multiplexed variants in 16 genes, covering 12 SNVs, 7 INDELs, 3 CNVs, and 3 SVs at different allele frequencies.

Seraseq® fusion RNA mix: this contrived fusion standard (Cat# 0710-0431) contains 16 clinically relevant RNA fusions mixed in total RNA from GM24385 cell line as background. In this study, it was further diluted to 10%, 5%, and 2% using cfDNA from healthy donor as background to generate a serial of standards with different copy numbers of fusion variants for limit of detection (LOD) evaluation.

Library preparation and sequencing:

Targeted libraries were prepared using Oncomine™ Pan-Cancer Cell-free assay reagents (Thermo Fisher Scientific, Cat# A37664).

For Horizon multiplex cfDNA reference standard, 25 ng was used as input in library preparation. In addition, input amount for 0.1% standard was increased to 50 ng to mimic 6000 copies of DNA allowing for improved sensitivity at lower LOD.

For Seraseq® cfDNA complete™ mutation mix, 25 ng was used as input in library preparation. For the serial dilutions from Seraseq® fusion RNA mix in cfDNA background, 20 ng was used as input in library preparation.

The Ion 540™ Kit-Chef was used for template preparation on Ion Chef™ (Thermo Fisher Scientific, Cat# 4484177) and followed by sequencing on Ion GeneStudio S5 XL system (Thermo Fisher Scientific, Cat# A27214) using the Ion 540™ Chip (Thermo Fisher Scientific, Cat# A27766).

Data analysis:

Data analysis was performed using Torrent Suite™ Software v5.6 and Ion Reporter™ v5.6 for simultaneous SNV/Indel, CNV, and fusion variant calls .

Table 1. Oncomine™ Pan-Cancer Cell-free Assay Content

Assay	Configuration	Unique Genes	DNA	RNA
Pan Cancer	TNA (DNA + RNA)	52	50	12

Hotspot Genes	Tumor Suppressor Genes	Copy Number Genes	Gene Fusions
AKT1 ALK AR ARAF BRAF CDKN2A CTNNB1 DDR2 EGFR ERBB2 ERBB3 ESR1 FGFR1 FGFR2 FGFR3 FGFR4 FLT3 GNA11 GNAQ GNAS	APC FBXW7 IDH1 IDH2 KIT KRAS MAPK1 MAPK2 MET MTOR NRAS NTRK1 NTRK3 PDGFRA PIK3CA RAF1 RET ROS1 SF3B1 SMAD4 SMO	CCND1 CCND2 CCND3 CDK4 CDK6 EGFR ERBB2 FGFR1 FGFR2 FGFR3 GATA1 GATA2 GATA3 GATA4 GATA5 GATA6 GATA7 GATA8 GATA9 GATA10 GATA11 GATA12 GATA13 GATA14 GATA15 GATA16 GATA17 GATA18 GATA19 GATA20 GATA21 GATA22 GATA23 GATA24 GATA25 GATA26 GATA27 GATA28 GATA29 GATA30 GATA31 GATA32 GATA33 GATA34 GATA35 GATA36 GATA37 GATA38 GATA39 GATA40 GATA41 GATA42 GATA43 GATA44 GATA45 GATA46 GATA47 GATA48 GATA49 GATA50 GATA51 GATA52 GATA53 GATA54 GATA55 GATA56 GATA57 GATA58 GATA59 GATA60 GATA61 GATA62 GATA63 GATA64 GATA65 GATA66 GATA67 GATA68 GATA69 GATA70 GATA71 GATA72 GATA73 GATA74 GATA75 GATA76 GATA77 GATA78 GATA79 GATA80 GATA81 GATA82 GATA83 GATA84 GATA85 GATA86 GATA87 GATA88 GATA89 GATA90 GATA91 GATA92 GATA93 GATA94 GATA95 GATA96 GATA97 GATA98 GATA99 GATA100	ALK BRAF ERG ETV1 FGFR1 FGFR2 FGFR3 MET NTRK1 NTRK2 NTRK3 RET ROS1

Table 1. Oncomine™ Pan-Cancer cell-free assay provides broad coverage across 52 unique genes that detects DNA and RNA variants across multiple cancer types. The workflow also offers multiplexing flexibility utilizing different chip configurations available to GeneStudio S5 systems to accommodate the needs for desired throughput..

RESULTS

Figure 1. Oncomine™ Pan-Cancer Cell-free Assay Workflow

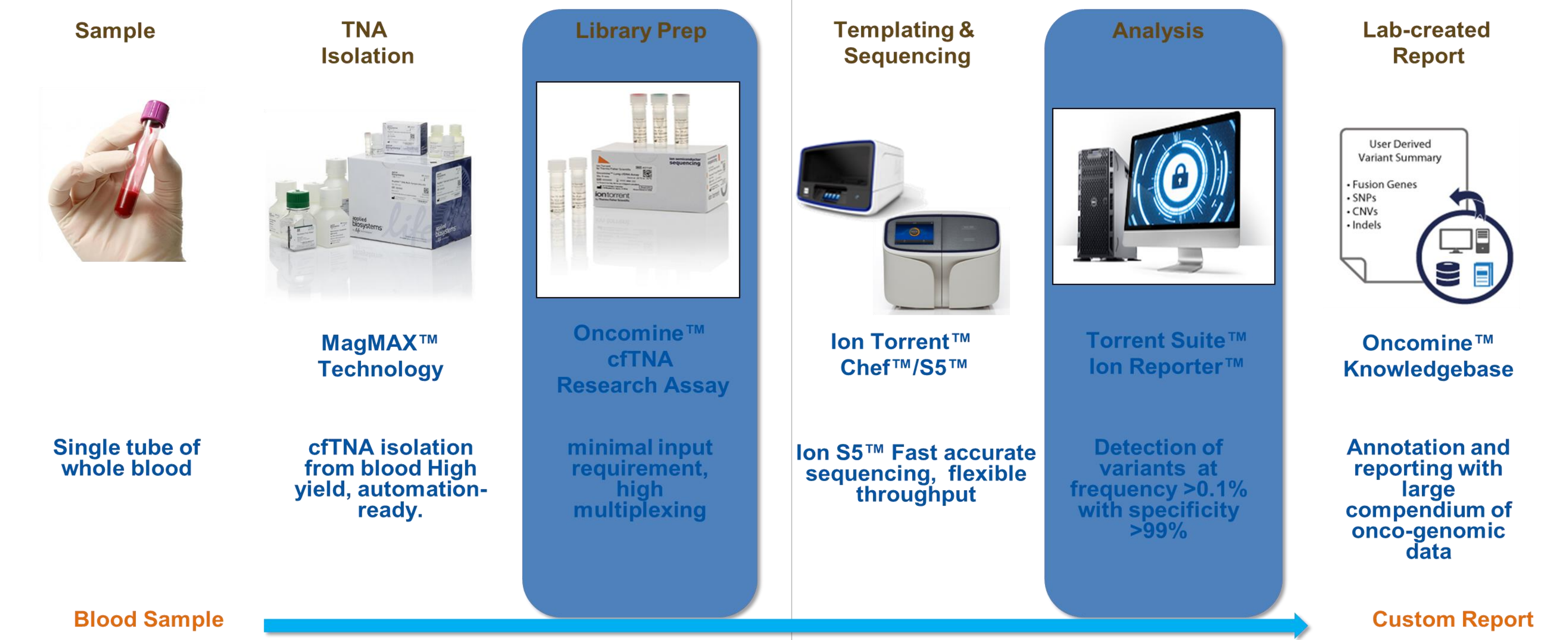


Figure 1. The Oncomine™ Pan-Cancer cell-free assay enables a fast 3-day workflow that starts with cell-free nucleic acid purification from a single tube of blood and library preparation to interrogate both DNA and RNA in a single reaction. The amplified library will undergo high-throughput semiconductor sequencing, with results analyzed and reported through an integrated bioinformatics solution.

Table 2. Result Summary for Horizon Multiplex cfDNA Reference Standards

Sample	Gene	AA Chg	SNV / Indel	AF % (R1)	AF % (R2)	AF % (R3)	AF % (R4)
Horizon cfDNA RS WT	CTNNB1	p.S33Y	24.0	24.3	26.7	24.2	
	PIK3CA	p.H1047R	24.9	24.0	26.5	24.7	
	EGFR	p.G719S	20.6	23.5	22.9	22.2	
	BRAF	p.V600E	29.9	29.3	30.7	33.5	
	MAP2K1	p.Q56P	22.4	23.7	23.7	22.4	
Sample	Gene	AA Chg	SNV / Indel	AF % (R1)	AF % (R2)	AF % (R3)	AF % (R4)
Horizon cfDNA RS 5% (25ng)	NRAS	p.Q61K	7.37	6.86	6.46	6.22	
	NRAS	p.A59T	7.20	7.22	6.34	6.96	
	PIK3CA	p.E545K	6.94	6.55	6.33	5.96	
	EGFR	p.Q746_A750del	5.48	6.98	6.70	5.80	
	EGFR	p.M766_A767insASV	3.60	3.78	3.37	3.12	
Horizon cfDNA RS 0.1% (50ng)	EGFR	p.T790M	3.72	3.53	3.01	3.98	
	EGFR	p.L858R	6.05	4.74	4.20	5.09	
	KRAS	p.G12D	6.33	5.28	6.22	6.27	
Sample	Gene	AA Chg	SNV / Indel	AF % (R1)	AF % (R2)	AF % (R3)	AF % (R4)
Horizon cfDNA RS 1.0% (25ng)	NRAS	p.Q61K	1.12	1.47	1.15	1.46	
	NRAS	p.A59T	1.46	0.78	1.77	1.19	
	PIK3CA	p.E545K	1.10	1.13	1.52	1.12	
	EGFR	p.Q746_A750del	1.12	1.53	1.45	1.71	
	EGFR	p.M766_A767insASV	0.74	0.62	0.69	0.68	
Horizon cfDNA RS 0.1% (50ng)	EGFR	p.T790M	0.81	0.80	0.69	0.58	
	EGFR	p.L858R	0.57	0.86	0.80	0.94	
	KRAS	p.G12D	0.90	1.43	1.24	1.49	
Sample	Gene	AA Chg	SNV / Indel	AF % (R1)	AF % (R2)	AF % (R3)	AF % (R4)
Horizon cfDNA RS 0.1% (50ng)	NRAS	p.Q61K	ND	0.21	0.17	0.13	
	NRAS	p.A59T	ND	0.13	0.17	0.13	
	PIK3CA	p.E545K	ND	0.11	0.19	0.15	
	EGFR	p.Q746_A750del	0.13	0.08	ND	ND	
	EGFR	p.M766_A767insASV	0.11	ND	0.08	0.21	
Horizon cfDNA RS 0.1% (50ng)	EGFR	p.T790M	0.20	0.16	0.11	0.15	
	EGFR	p.L858R	ND	ND	ND	ND	
	KRAS	p.G12D	ND	ND	0.09	ND	
Sample	Gene	AA Chg	SNV / Indel	AF % (R1)	AF % (R2)	AF % (R3)	AF % (R4)
Horizon cfDNA RS 0.1% (50ng)	NRAS	p.Q61K	ND	0.098			
	NRAS	p.A59T	0.096	0.117			
	PIK3CA	p.E545K	0.079	0.212			
	EGFR	p.Q746_A750del	0.101	0.120			
	EGFR	p.M766_A767insASV	0.073	ND			
Horizon cfDNA RS 0.1% (50ng)	EGFR	p.T790M	0.193	0.117			
	EGFR	p.L858R	0.136	0.117			
	KRAS	p.G12D	ND	ND			

Table 2. All 13 variants were successfully detected in Horizon Multiplex I cfDNA RS standards above 1% AF (25 ng input) using Pan-Cancer assay, including 8 SNV/Indels at several important gene loci and 5 additional SNVs from parental cell lines as shown in RS wild-type. For 0.1% RS, high DNA input at 50 ng that is equivalent 6000 copies of DNA molecules significantly improved variant detection sensitivity from ~73% to ~85% comparing to 25 ng input. No FPs were observed in any samples.

Table 3. Result Summary for Seraseq® cfDNA Complete™ Mutation Mix

Sample	Gene	AA Chg	SNV / Indel	R1 (%AF)	R2 (%AF)	R3 (%AF)
Seraseq cfDNA complete mix WT	FP		1	2	2	
	Gene	AA Chg	R1 (%AF)	R2 (%AF)	R3 (%AF)	
	NRAS	p.Q61R	1.09	1.33	0.92	
	ALK	p.G1202R	1.14	1.20	1.34	
	ALK	p.F1174L	0.74	1.26	1.05	
Seraseq cfDNA complete mix 1.0%	PIK3CA	p.H1047R	1.18	1.50	1.13	
	KIT	p.D816V	1.45	0.88	1.26	
	EGFR	p.E746_A750del	1.31	1.48	0.95	
	EGFR	p.T790M	1.29	1.43	0.91	
	EGFR	p.L858R	1.02	1.29	0.77	
Seraseq cfDNA complete mix 0.5%	BRAF	p.V600E	0.70	0.47	1.08	
	KRAS	p.Q61H	1.02	1.04	1.04	
	KRAS	p.G12D	0.93	0.81	1.06	
	KRAS	p.G12C	1.00	1.07	1.03	
	EGFR	p.L747_P753>S	1.77	1.24	1.68	
Seraseq cfDNA complete mix 0.1%	EGFR	p.S752_I759del	1.16	1.03	1.11	
	ERBB2	p.E770_A771insAVVM	0.94	1.00	0.87	
	AKT1	p.E17K	0.75	0.76	0.69	
	TP53	p.G245D	0.37	0.66	0.44	
	FP		1	2	2	
Seraseq cfDNA complete mix 0.1%	Gene	AA Chg	R1 (%AF)	R2 (%AF)	R3 (%AF)	
	NRAS	p.Q61R	0.70	0.68	0.54	
	ALK	p.G1202R	0.50	0.79	0.50	
	ALK	p.F1174L	0.50	0.60	0.46	
	PIK3CA	p.H1047R	0.51	0.58	0.52	
Seraseq cfDNA complete mix 0.1%	KIT	p.D816V	0.88	0.82	0.91	
	EGFR	p.E746_A750del	0.56	0.59	0.68	
	EGFR	p.T790M	0.20	0.42	0.45	
	EGFR	p.L858R	0.55	0.64	0.72	
	BRAF	p.V600E	0.28	0.64	0.55	
Seraseq cfDNA complete mix 0.1%	KRAS	p.Q61H	0.63	0.55	0.77	
	KRAS	p.G12D	0.36	0.59	0.50	
	KRAS	p.G12C	0.72	0.59	0.61	
	EGFR	p.L747_P753>S	0.88	0.59	0.74	
	EGFR	p.S752_I759del	0.88	0.59	0.59	
Seraseq cfDNA complete mix 0.1%	ERBB2	p.E770_A771insAVVM	0.18	0.28	0.20	
	AKT1	p.E17K	0.53	0.65	0.52	
	TP53	p.G245D	0.41	0.39	0.33	
	FP		1	1	2	
	Gene	AA Chg	R1 (%AF)	R2 (%AF)	R3 (%AF)	
Seraseq cfDNA complete mix 0.1%	NRAS	p.Q61R	0.24	0.13	0.21	
	ALK	p.G1202R	0.10	0.15	0.12	
	ALK	p.F1174L	0.09	ND	ND	
	PIK3CA	p.H1047R	ND	ND	0.24	
	KIT	p.D816V	0.09	0.09	ND	
Seraseq cfDNA complete mix 0.1%	EGFR	p.E746_A750del	ND	0.31	0.12	
	EGFR	p.T790M	ND	0.15	ND	
	EGFR	p.L858R	0.21	0.13	ND	
	BRAF	p.V600E	0.21	ND	ND	
	KRAS	p.Q61H	0.43	ND	ND	
Seraseq cfDNA complete mix 0.1%	KRAS	p.G12D	ND	ND	0.14	
	KRAS	p.G12C	ND	0.08	0.20	
	EGFR	p.L747_P753>S	0.09	0.15	ND	
	EGFR	p.S752_I759del	0.09	0.15	ND	
	ERBB2	p.E770_A771insAVVM	0.13	ND	ND	
Seraseq cfDNA complete mix 0.1%	AKT1	p.E17K	0.15	ND	0.14	
	TP53	p.G245D	0.65	0.29	0.41	
	FP		3	1	3	

Table 3. All 16 variants were successfully detected in Seraseq™ cfDNA Complete Mutation Mix with higher allele frequency ≥ 0.5%. As observed previously (data not shown) additional FP calls were made in these libraries.

Table 4. Result Summary for Seraseq® Fusion RNA Dilutions

Fusion Variants	Copies (20%)*		Molecule Count ^Δ			
	ddPCR_SeraCare	20%	10%	5%	2%	
TPM3-NTRK1	95	77	66	22	11	
LMNA-NTRK1	97	123	30	16	15	
SLC45A3-BRAF	132	110	35	6	2	
EM14-ALK	184	78	36	15	19	
FGFR3-BAIAP2L1	259	38	16	16	3	
FGFR3-TACC3	147	35	20	14	3	
SLC34A2-ROS1	101	24	4	3	ND	
CD74-ROS1	153	40	18	4	8	
MET Exon14 Skipping	166	68	45	15	9	
KIF5B-RET	83	76	45	15	5	
NCOA4-RET	144	127	30	13	14	
ETV6-NTRK3	128	87	51	15	8	
TMPS22-ERG	83	8	ND	ND	ND	
PAX8-PPARG1	259					
EGFR-SEPT14	133					
EGFR variant III	124					

* Estimated copy number in 20% dilution based on ddPCR measurement from SeraCare
Δ Measured copy number using Pan-Cancer assay

Table 4. All 13 fusion variants covered by Pan-Cancer assay were detected in Seraseq fusion standards diluted down to 5% in cfDNA background. Based on the molecule counts, LODs for each fusion target were able to achieve at 2% standard for Pan-Cancer assay.

Table 5. Result Summary for Horizon Structural Multiplex cfDNA Reference Standard

Sample	SNV / Indel		CNV		Fusion	
	Gene	AA Chg	AF % (R1)	Gene	CNV Ratio	Molecule Counts
Structural Multiplex cfDNA Reference Standard	CTNNB1	p.S33Y	5.4	FGFR3*	1.7	50
	PIK3CA	p.E545K	5.8	CDK6*	1.4	6
	PIK3CA	p.H1047R	15.7	MET	1.8	26
	EGFR	p.G719S	5.8	MYC	3.1	
	EGFR	p.Q746_A750del	5.7			
Structural Multiplex cfDNA Reference Standard	EGFR	p.M766_A767insASV	4.9			
	BRAF	p.V600E	16.4			
	KRAS	p.G13D	4.7			
	AKT1	p.E17K	5.0			
	MAP2K1	p.Q56P	4.7			
Structural Multiplex cfDNA Reference Standard	TP53	p.S241F	6.9			
	TP53	p.R175C	0.2			
	GNA11	p.Q209L	5.8			