

Development of a Targeted NGS Oncology Assay for Detection and Reporting Comprehensive Genomic Profiling

Paul Williams, Vinay Mittal, Jennifer Kilzer, Gary Bee, Santhoshi Bandla, Dinesh Cyanam, Sameh El-Difrawy, Aren Ewing, Nickolay Kahanov, Anelia Kraltcheva, Denis Kaznadzey, Cristina Van Loy, Scott Myrand, Warren Tom, Yu-Ting Tseng, James Veitch, Chenchen Yang, Janice Au-Young, Zheng Zhang, Fiona Hyland, Elaine Wong-Ho, Seth Sadis Thermo Fisher Scientific, Ann Arbor, MI, USA; South San Francisco, CA; Austin, TX, Carlsbad, CA.

INTRODUCTION

Next-generation sequencing (NGS) is a research tool that allows for the detection of cancer-associated somatic mutations, focal copy number aberrations, and gene fusions.<sup>1</sup> Current scientific and clinical research indicates these genomic variants may be associated with favorable or unfavorable responses to specific targeted therapeutic regimens.<sup>2,3,4</sup> The recent emergence of immunotherapeutic agents that enable the immune system to destroy cancer cells has driven research to identify additional biomarkers that may be associated with therapeutic response to these agents. Speculation that immune cells are recruited to the tumor by the presentation of mutant antigens<sup>5</sup> has led to the investigation of several biomarkers associated with increased somatic mutation rates. Comprehensive sequencing of the genes involved in DNA repair pathways may identify variants that lead to deficiencies in these pathways. These deficiencies may manifest themselves as microsatellite instability<sup>6</sup> (MSI), the extension of small repetitive sequence elements within the genome, or as an increase in overall tumor mutational burden (TMB)<sup>7</sup>, an estimate of the total number of mutations observed in a tumor sample, expressed in mutations/Megabase of DNA.

We developed an NGS assay for FFPE tissue samples that can detect a variety of DNA variants. Variants detected by the assay have been associated with response to targeted therapies and immune checkpoint inhibitors.<sup>2,3,6,7</sup> This assay covers over 500 genes, including several target genes that have been associated with oncology clinical research, and requires low amounts of input FFPE material. This assay detects important biomarkers associated with oncology research and is part of an automated sample-to-report workflow that allows streamlined utilization for clinical research.

MATERIALS AND METHODS

Gene content was prioritized based on variant prevalence in solid tumors. To support robust TMB estimation, additional genomic regions were added in order to bring the coding sequence footprint to >1MB. To enable MSI status assessment, coverage of a diverse set of microsatellites throughout the genome was added. The MSI analysis solution makes use of in-sample standards incorporated as internal references.

The assay uses Ion AmpliSeq™ technology with automated templating on the Ion Chef™ System and sequencing on the Ion Torrent GeneStudio™ S5 sequencing platform. An automated variant calling and TMB reporting workflow is provided within Ion Reporter™ Software.

RESULTS

Over 500 genes with DNA based alterations and over 50 RNA fusion drivers are included in the assay. Of the genes measured for DNA alterations, 169 cover important cancer hotspots, 333 cover copy number variants (CNV), and 227 genes have full coding sequence (CDS) coverage for detection of truncating variants. The total genomic coverage of the assay is 1.50MB with 1.1MB of exonic sequence, to support high-confidence TMB estimation. A diverse set of microsatellite markers targeting MSI locations comprising of mono- and di-nucleotide repeats ranging from 7 to 34 bp are included for MSI status assessment. RNA content for the assay is more fully described in poster ST091 “Development of a comprehensive next-generation sequencing assay for gene-fusions detection in solid tumors”.

FFPE tumor samples from a variety of tissue types were sequenced using the assay. The assay displays high (>95%) uniformity and consistent read depth (>2200x) to support robust variant calling at low allele frequency. An excellent variant calling (SNV/INDEL) performance was observed, with sensitivity ranging from 98%-100% and PPV ranging from 88%-100%, depending upon the variant type. CNV detection sensitivity and specificity were 89%-99% and 100%, respectively. In-silico TMB assessment using publicly available whole-exome cancer sequencing data resulted in a correlation of R<sup>2</sup> > 0.90 (0-40 mut/mb) in pan-cancer and specific cancer types including lung, colorectal and melanoma. MSI test performance was assessed using 192 samples with known MSI status. An overall performance of 96% sensitivity and 99% specificity was observed.

Figure 1. Summary of the Assay Content

CATEGORIZED BY SOMATIC ALTERATION TYPE

500+ GENES

169 Hotspot genes

333 Focal CNV gains and CNV loss

227 Full CDS genes

>1M Exonic bases

76 Microsatellite genes

50 Fusion drivers

CATEGORIZED BY RELEVANCE

15 Genes on 66 labels

22 Genes in 26 guidelines

180 Genes are used in 1,034 global clinical trials

Table 1. Oncomine Comprehensive Assay Plus Cancer Gene Table

CNV gain genes										Full CDS coverage																																																																			
CNV gain only										CNV loss																																																																			
CNV gain and hotspot					Hotspot only					CDS only																																																																			
ABCB1	ABL1	FGFR4	PIK3C2B	ACVR1	MYO10	ABRAXAS1	CDKN1A	FANCA	MAP3K1	POLE	SOX9	CALR	CTNND2	ABL2	FLT3	PIK3CA	ATP11A	NSD2	ACVR1B	CDKN1B	FANCC	MAP3K4	POT1	SPEN	GATA	CYP2D28																																																			
DDR1	AKT1	FLT4	PIK3CB	AURKB	NTSG2	ACVR2A	CDKN2A	FANCD2	MAPK8	PPM1D	STAG2	CYP2D28	EMSY	AKT2	FOXA1	PIK3R2	BCR	NTRK2	ADAMTS12	CDKN2B	FANCE	MEN1	PPP2R2A	STK11	ERCC25	FGF3	AKT3	GATA2	PIK1	BMP5	NUP205	ADAMTS2	CDKN2C	FANCF	MOA	PRDM1	STU1	FAS	ID3																																						
FGF4	ALK	GNAS	PLCG1	BTX	PAS5	AMER1	CDKN2A	FANCG	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK2	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9																										
FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3																										
FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK2	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3													
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC	CHEK1	FANCI	MLH3	PRKARIA	TAP2	KULH13	FGF19	ARAF	HSP3B	PPP2C	CD79B	PIK3CG	ARHGAP35	CIC	FANCL	MRE11	PTCH1	TBX3	MTUS2	FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9
FGF23	AURKA	IDH2	PRKACA	CSF1R	PTPRD	ARID1A	CSMD3	FAT1	PTEN	TCF7L2	PSMB8	PSMB9	GLI3	AXL	ILTR	PXN1L	CUL1	RHOA	ARID2	CTCF	FBXW7	MSH6	RAD50	TGFB2	PSMB10	IGF1R	BCL2	RAC1	CTSL	CYSLTR2	SLK1	ASXL1	CDKN2A	FANCD1	MLH1	PRDM9	TAP1	ID3	FGF9	ARL	HSP3A	PPP2R1A	CACNA1D	PIK3CD	APC																																