

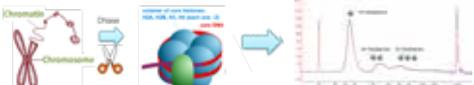


Correlation between mutations found in FFPE tumor tissue and paired cfDNA samples

Brittany Niccum PhD, Cindy Heath, Lauren Saunders PhD, Antonia Hur, and Asmita Patel
Beckman Coulter Life Sciences, Indianapolis, IN

Introduction

Liquid biopsies represent a promising area of facilitating cancer research as blood collection is less invasive than tumor biopsies. Cell free DNA (cfDNA) consists of small (150 – 500 bp) DNA fragments that circulate in the blood. cfDNA levels tend to be low in healthy, non-pregnant patients, and increase in patients with cancer, pregnancy, or extensive damage to tissue. cfDNA is believed to be derived mostly from apoptotic cells for which biomarkers for a variety of diseases have been found in cfDNA.



Apoptotic or necrotic cell death results in near-complete digestion of native chromatin and can be recovered from peripheral blood plasma as cfDNA.

Each 160-175 bp DNA is wrapped ~1.67 times around one nucleosome. These protein-bound DNA fragments preferentially survive digestion and are released into the circulation, and can be recovered from peripheral blood plasma as cfDNA.

Typical cfDNA peaks characterized by Agilent 2100 Bioanalyzer, with a main peak at 175 bp, second and third peaks at 350 and 525 bp.

FFPE tissue is often used to look for cancer-associated mutations despite invasiveness; however it does not always correlate with the mutations seen in cfDNA. In this poster we present a comparison of matched FFPE and plasma samples to determine how many mutations are seen in both tissues. We also look at where the mutational mismatches appear in the chromosome. Different chromosomal regions can have different mismatch rates, and we use this to draw conclusions about the best chromosomal locations for biomarkers. We automated from extraction through sequencing in collaboration with Swift Biosciences.

As cfDNA is extracted from blood, it is a non-invasive way to detect disease. However, there is some concern that cfDNA does not contain the same biomarkers as tumor tissue. Tumor tissue is typically removed and stored as formalin-fixed, paraffin-embedded tissue, a process that preserves the morphological structures well but chemically modifies and degrades the nucleic acids.

Methods

Sample Preparation

DNA was extracted from FFPE tissues using *FormaPure XL* Total automated on a Biomek i5 multichannel workstation. DNA concentrations for 4 10 µl curts were estimated using the Quant-iT PicoGreen assay; varied between blocks. Some blocks had very low yields, most likely due to tissue distribution in the blocks; for these samples the extraction was repeated with 7 10 µl curts. DNA from FFPE was sheared on a Covaris S220 following the 200bp shear protocol. cfDNA was extracted from 1ml of plasma using the *MinElute* High Efficiency cfDNA kit. cfDNA yield varied between samples as expected. Concentration was estimated using the Kapa Hg-Quant kit.

Library Construction

cfDNA and sheared FFPE DNA was processed with *Swift Accer-NGS 25 Hyd DNA Library Kit*. DNA input was 100ng. Following the library construction, the *Swift Pan-Cancer Hyb Panel* was used for gene enrichment. The hybridized libraries were sequenced on the Illumina NextSeq 550.

Analysis

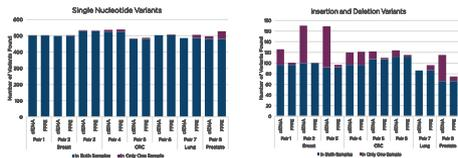
The sequencing data was analyzed using the BWA-enrichment tool on Illumina BaseSpace mapped to the genes in the panel.

The two extractions, library construction and hybridization panel have all been automated on a Biomek Workstation. The approximate time for an automated workflow is compared to manual workflows in the table below.

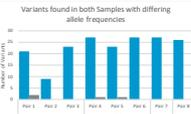
	Throughput per run	Manual Timings		Automated Timings	
		Hands on Time	Total Time	Hands on Time	Total Time
• <i>Adaptive High Efficiency cfDNA Kit</i>	24 Manual/ 96 Automation	45 minutes	75 minutes	30 minutes	5.3 hours
• <i>FormaPure XL Total</i>	24 Manual/ 96 Automation	3.5 hours	6.5 hours	30 minutes	8.5 hours
• <i>Swift Accer-NGS 25 Hyd DNA Library Kit</i>	96	2.7 hours	4.6 hours	30 minutes	6 hours
• <i>Swift Pan-Cancer Hyb Panel</i>	16	2.5 hours	8.5 hours	30 minutes	8.8 hours

Majority of Variants are identified in both sample types

We identified both single nucleotide variants (SNV) and insertion and deletion (indel) events in the sequencing results from all the samples. First we compared the SNVs found in both sample types. The majority of the SNVs and indels were identified in both sequencing methods and all tissues (Figure 1 and Figure 2). In all of the samples unique SNVs and indels were identified in both cfDNA and FFPE DNA. More SNVs were identified when sequencing DNA from FFPE than cfDNA, however this difference wasn't statistically significant. Conversely, more indels were identified when sequencing cfDNA and this difference is statistically significant.

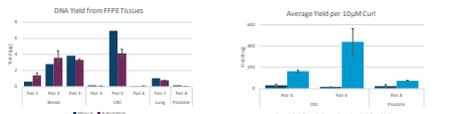


We also wanted to verify that the variants that were found in both sample types were found at equal frequency. This was determined by comparing the allele frequency identified by sequencing cfDNA and DNA from FFPE tissue. We compared the ratio of the two allele frequencies. For all 8 pairs the ratio of cfDNA allele frequency to FFPE allele frequency was 1; the allele frequency for variants identified by cfDNA and FFPE were not significantly different than each other. Interestingly though, when looking at variants with allele frequency less than 2 standard deviations from the mean, cfDNA showed allele frequencies higher than FFPE for more variants.



Methods Continued: Extraction Results and Sequencing Coverage

DNA extracted from FFPE was done both manually and on a Biomek i5 multichannel workstation. Both the manual user and the liquid handler extracted similar amounts of DNA (Figure 1); they are not significantly different as determined by a student t-test (P=0.79). Three of the blocks had very low yields and were extracted in the manual. The initial low yield were most likely due to in block tissue distribution. In Figure 2, the first 12 curts had much lower yield of DNA per curts than the second 14 curts did.



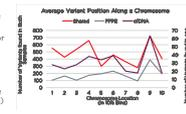
Below is a table of the sequencing coverage for all the samples. All samples had sequencing coverage of 50x at least 95% of the bases sequenced.

Tumor Tissue	Sample Material	Mean Region Coverage	Uniformity of Coverage	Target Coverage at			
				1X	10X	20X	50X
Breast	Pair 1	cfDNA 1946.5	99.50%	100.00%	100.00%	100.00%	100.00%
	FFPE 1322.4	98.00%	100.00%	100.00%	100.00%	99.50%	
	Pair 2	cfDNA 1677.2	89.50%	100.00%	100.00%	99.50%	99.50%
	FFPE 1306.5	99.00%	100.00%	100.00%	100.00%	100.00%	
	Pair 3	cfDNA 1051.4	90.50%	100.00%	99.50%	99.80%	99.20%
	FFPE 1931	98.50%	100.00%	100.00%	100.00%	100.00%	
	Pair 4	cfDNA 1973.2	99.30%	100.00%	100.00%	100.00%	100.00%
	FFPE 1239.8	89.80%	100.00%	100.00%	99.80%	99.50%	
CRC	Pair 5	cfDNA 2068.1	89.40%	100.00%	100.00%	100.00%	100.00%
	FFPE 1480.1	98.70%	100.00%	100.00%	100.00%	100.00%	
Lung	Pair 6	cfDNA 1772.4	98.00%	100.00%	100.00%	100.00%	100.00%
	FFPE 2656	99.60%	100.00%	100.00%	100.00%	100.00%	
Prostate	Pair 7	cfDNA 1775.8	99.50%	100.00%	100.00%	100.00%	100.00%
	FFPE 1359	87.90%	100.00%	99.60%	98.60%	92.30%	
Prostate	Pair 8	cfDNA 1436.4	99.00%	100.00%	100.00%	100.00%	99.90%
	FFPE 598.6	91.70%	99.90%	99.80%	99.60%	98.60%	

Variant location

To determine if mutations were evenly distributed throughout chromosomes, each chromosome was divided into 10 bins; the first 10% of each of the 23 chromosomes were treated as one bin. The mutations were pooled into the 10 bins for each chromosome. This was done separately for mutations found in only DNA from only FFPE or cfDNA and found in both FFPE DNA and cfDNA.

The graph to the right shows how the mutations mapped across the 10 bins. A Pearson's correlation was done to test for how differently the mutations mapped across the 10 bins. Mutations found only in FFPE tissue did not correlate (P=0.3) as well as mutations found only in cfDNA (P=0.8) to the position of mutations found in both FFPE DNA and cfDNA.



Variants Found at ClinVar Pathogenic Locations

We wanted to determine if we found any pathogenically relevant variants. To do this we compared against a list of variants from the NCBI ClinVar database. We only used variants that had the clinical significance of conflicting interpretations, uncertain significance, likely pathogenic, pathogenic and risk factor. We did this for all four of the tumor types. We were able to identify 2 variants in the ClinVar database only found using the DNA from FFPE tissue.

Tissue	Pair	cfDNA, FFPE, or Both	Gene	Chromosome	Position	Mutation	ClinVar Mutation	Conditions associated
CRC	Pair 4	FFPE	APC	chr5	112E-08	C>T	C>T	Familial adenomatous polyposis
CRC	Pair 6	FFPE	TP53	chr17	757840G	C>T	C>T	Carcinoma of colorectal/ovarian/colonic carcinoma, hereditary/Fraser's syndrome

In most of the samples however, we were unable to find any variants identified in the ClinVar database. We were however able to find variants at the same position as variants in the ClinVar database. These could also alter the encoded protein and would be good candidates for further study into pathology. Interestingly all of these variants were found in the FFPE DNA sequences; one of them was only found in FFPE DNA sequences; higher sequencing coverage of the cfDNA could help to identify these variants.

Tissue	Pair	cfDNA, FFPE, or Both	Gene	Chromosome	Position	Mutation	ClinVar mutation
Breast	Pair 1	Both	BRC1A1	chr17	4323394	T>C	GCC insertion
		Both	BRC1A1	chr17	4324436	G>A	G Duplication or TT insertion
		Both	CHEK2	chr22	2913048	T>C	T Deletion
		Both	TP53	chr17	7579472	G>C	G>T
	Pair 2	Both	BRC1A1	chr17	4323394	T>C	TA A
		Both	BRC1A1	chr17	4324400	T>C	TA A
		Both	BRC1A1	chr17	4324435	T>C	TA A
		Both	STK11	chr19	122022	G>C	G>A or G>T
Pair 3	Both	BRC1A1	chr17	4323394	T>C	T Deletion	
	Both	BRC1A1	chr17	4323994	T>C	TA A	
	Both	BRC1A1	chr17	4324435	T>C	TA A	
	Both	TP53	chr17	7579472	G>C	G>T	
Lung	Pair 7	FFPE	TP53	chr17	7579471	G>C	G Duplication
	Both	BRC1A1	chr17	4323994	T>C	GCC insertion	
Prostate	Pair 8	Both	BRC1A1	chr17	4324436	G>A	G Duplication or TT insertion
	Both	CHEK2	chr22	2913048	T>C	T Deletion	

Conclusions

Here we show:

- Sequencing of cfDNA captures the majority of variants that were found in sequenced FFPE DNA
- More indels are identified using cfDNA than with FFPE DNA, especially with breast tissue
- Distribution of variants across the genome differs when sequencing FFPE DNA
- More previously identified clinically relevant variants, as identified by the ClinVar database were found when sequencing FFPE DNA

This study is small and further work should be done using larger data sets to gain more conclusive information.



Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to these protocols, 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. These protocols are for demonstration only, and are not validated by Beckman Coulter.

© 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. AAG-6156PST11.9