

Multiplex Exome Enrichment from Pooled Barcoded Libraries Yields Efficient SNP and Indel Detection on the SOLiD™ System

Gavin D Meredith, Christopher Adams, Gary Bee, Loni Pickle, Miroslav Dudas, Jennifer Kilzer, Adam Harris, Marie Callahan, George Marnellos, Yongming Sun, Vrunda Sheth, Clarence Lee, Tanya Sokolsky, Chris Clouser, Kim Mather, Maryam Shenasa, Andrew Hutchinson, Jeffrey Ichikawa, Timothy Harkins, and Rob Bennett, Life Technologies Corp., 5791 Van Allen Way, Carlsbad, CA, 92008

ABSTRACT

The identification of genetic variation associated with human disease requires the development of a robust and cost-effective approach for systematic resequencing of candidate regions in the human genome. Even though the cost of sequencing a human genome continues to drop, the demand for increased sample throughput continues to increase. Higher sample throughput is considered necessary to enable larger patient cohort studies which hold the key to identifying rare disease-related alleles. Thus, scalable and automatable workflows for target enrichment and sequencing are needed to facilitate cancer and other genetic disease research. Described here is a targeted resequencing workflow that employs pooled barcoded fragment libraries, multiplexed exome enrichment, and multiplexed sequencing on the Applied Biosystems™ SOLiD™ System. To validate the performance of this multiplexed workflow, barcoded fragment libraries were made from HuRef DNA using the new 5500 SOLiD™ fragment library protocol. Resulting libraries were then pooled in multiples of 4 for exome capture with the Agilent SureSelect™ Human All Exon 50 Mb Kit. The 4-plex data obtained from 2 quads of SOLiD™ 4 fragment sequencing yielded an average depth of coverage over the targets of 23.1x. The 8-plex data from a full slide yielded average depth of 29.6x. Overall, good barcode balance, similar mapping efficiencies and similar SNP/indel calls were observed for 4-plex and 8-plex exome capture samples. The percentage of on-target reads varied from 71.2% to 74.0% which is comparable to numbers reported by others. For the 8-plex samples, the concordance of SNP calls (average of 31,474 SNPs) to dbSNP was 98.7% (sd=0.1%) for homozygous and 90.3% (sd=0.3%) for heterozygous variants and the concordance of small indels (average of 1560 indels) was 55.9% (sd=0.9%). Of particular note, sequencing a single (38 Mb) exome on a single lane of a 5500x SOLiD™ System flow-cell yielded an average coverage of 66.9x (76.9% of target bases covered at >= 20x depth) with only 5.1% of target bases left uncovered. The combination of multiplexed exome enrichment and multiplexed on the SOLiD™ System provides an efficient and economical solution for the high-throughput detection of genetic variation in multiple human genomes.

INTRODUCTION

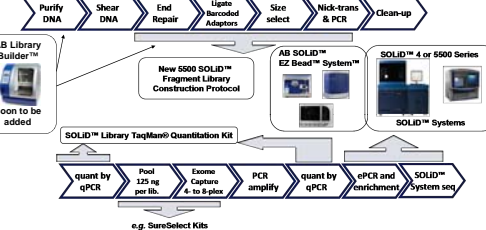
Next-generation sequencing technology has brought high throughput sample processing to genome sequencing, but an accompanying solution for high throughput target enrichment is lacking. Target enrichment is a term used to describe the ability to selectively enrich specific regions of a genome. The method employed by the Agilent SureSelect Human All Exon 50 Mb Kit extracts target regions from genomic libraries by hybridization to in-solution biotinylated cDNA probes, or "baits". Post-enrichment material is amplified and used directly for downstream steps, including emulsion PCR (ePCR) and sequencing on the SOLiD™ System (Figure 1). The inherent scalability and flexibility for automation of the SureSelect in-solution enrichment system coupled with the ultra-high throughput of the SOLiD™ sequencing platform provides an integrated approach to targeted resequencing. The new Agilent SureSelect Human All Exon 50 Mb Kit builds upon previous exon products with additional validated novel content developed by the Wellcome Trust Sanger Institute. The new design encompasses coding exons annotated by the GENCODE project and also includes all exons annotated in the consensus CCDS (March 2009) database. In addition, the content contains small non-coding RNAs from miRBase (v13) and Rfam.

MATERIALS AND METHODS

HuRef genomic DNA, purchased from the Coriell Institute for Medical Research, was fragmented to a mean length of ~200 bp with the Covaris S2 System, then 3 µg amounts were used for library construction using a new protocol that included the use of 5500 SOLiD™ System compatible barcoded adapters. After nick translation, libraries were PCR amplified for 6 cycles, quantified and pooled in 4-plex (BC1-BC4) or 8-plex (BC1-BC8) using 125 ng of each library, based on Bioanalyzer estimates of average size and qPCR determinations of amplifiable molecules. The pooled 500 ng (for 4-plex) or 1 µg (for 8-plex) of library DNA was mixed with adaptor barcodes, dried-down, and handled as described in the Agilent protocol, using 1X capture probes for 4-plex and 2X probes for 8-plex. After hybridization, capture, elution, and clean-up, the enriched libraries were amplified by 10 more cycles of PCR. Standard steps were taken thereafter to create enriched, templated beads for SOLiD™ System sequencing. The beads were sequenced as 50-color fragment tags (FS) and the data was progressively mapped in color-space and target enrichment and variant calling statistics were generated with the Targeted Resequencing pipeline in SOLiD™ BioScope™ 1.3 software.

RESULTS

Figure 1. Workflow for multiplexed exome capture with SOLiD™ System sequencing



An exome enrichment workflow that permits pre-capture pooling of barcoded libraries and incorporates many of the most recent SOLiD™ System innovations has been developed. A working protocol has been established and may be available upon request.

Table 1. 4-plex library pooling prior to dry-down

Barcode	Number of mapped reads	Percent on target	Fold Enrichment	Percent target bases not covered	coverage >=1x	coverage >=5x	coverage >=10x	coverage >=20x	average coverage depth
BC1	34,421,202	74.4%	44.7	15.6%	84.3%	71.6%	60.8%	42.2%	21.3
BC2	39,463,644	74.9%	45.0	14.9%	85.1%	73.5%	63.9%	46.9%	24.7
BC3	33,740,286	75.9%	45.6	15.9%	84.1%	71.3%	60.5%	42.0%	21.3
BC4	40,992,057	73.2%	44.0	15.4%	84.6%	73.5%	63.0%	46.9%	25.0
Total	148,617,394								
average	37,154,239	74.6%	44.8	15.4%	84.5%	72.2%	62.1%	44.5%	23.1
std dev	3,614,173	1.1%	0.7	0.4%	0.4%	1.0%	1.6%	2.8%	2.0

Figure 2. Exome-enriched 4-plex library

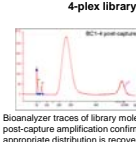


Figure 3. 4-plex barcode representation

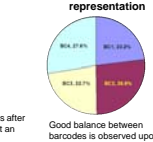


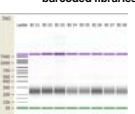
Table 2. 4-plex mapping and enrichment stats (SureSelect 50 Mb exome)

Barcode	Number of mapped reads	Percent on target	Fold Enrichment	Percent target bases not covered	coverage >=1x	coverage >=5x	coverage >=10x	coverage >=20x	average coverage depth
BC1	34,421,202	74.4%	44.7	15.6%	84.3%	71.6%	60.8%	42.2%	21.3
BC2	39,463,644	74.9%	45.0	14.9%	85.1%	73.5%	63.9%	46.9%	24.7
BC3	33,740,286	75.9%	45.6	15.9%	84.1%	71.3%	60.5%	42.0%	21.3
BC4	40,992,057	73.2%	44.0	15.4%	84.6%	73.5%	63.0%	46.9%	25.0
Total	148,617,394								
average	37,154,239	74.6%	44.8	15.4%	84.5%	72.2%	62.1%	44.5%	23.1
std dev	3,614,173	1.1%	0.7	0.4%	0.4%	1.0%	1.6%	2.8%	2.0

Table 3. 4-plex variant calls and dbSNP132 concordance (SureSelect 50 Mb exome)

Barcode	total SNPs	homo dbSNP	heter dbSNP	het SNPs	total indels	homo dbSNP	heter dbSNP	het indels	total concord
BC1	34364	14997	98.7%	20467	96.5%	1175	552	70.8%	1023
BC2	37839	16127	98.4%	21712	96.5%	1647	587	70.4%	1060
BC3	35819	15758	98.6%	20060	91.1%	1540	545	70.6%	995
BC4	37160	15849	98.6%	21311	91.1%	1715	582	71.0%	1133
average	36796	15908	98.6%	20913	96.9%	1619	567	70.7%	1053
std dev	887	137	0.3%	722	0.3%	78	21	0.3%	60

Figure 4. Pre-pooled 8-plex barcoded libraries



The protocol yields libraries of consistent size and concentration. Average library molecule size based on Bioanalyzer™ traces and molar concentration by qPCR are used to determine the "qPCR ng/µl"; this value is used to apportion 125 ng of each library into the pool.

Table 5. 8-plex mapping and enrichment stats (SureSelect 50 Mb exome)

Barcode	Number of mapped reads	Percent on target	Fold Enrichment	Percent target bases not covered	coverage >=1x	coverage >=5x	coverage >=10x	coverage >=20x	average coverage depth
BC1	48,326,122	72.4%	43.6	22.4%	77.6%	62.4%	54.4%	44.1%	28.8
BC2	53,597,196	72.5%	43.5	21.1%	78.5%	63.6%	55.6%	45.9%	31.9
BC3	48,066,332	74.0%	44.4	23.1%	78.5%	61.8%	53.9%	43.3%	27.3
BC4	53,581,754	71.2%	42.7	22.4%	77.6%	62.3%	54.4%	44.7%	33.4
BC5	54,065,056	72.4%	43.5	21.3%	78.7%	64.0%	56.2%	46.3%	32.1
BC6	44,246,062	72.5%	43.4	23.6%	78.4%	60.3%	51.9%	45.9%	26.2
BC7	52,462,469	71.8%	43.1	21.9%	78.1%	62.9%	55.0%	45.1%	30.8
BC8	49,969,589	72.3%	43.4	22.2%	78.8%	62.7%	54.6%	44.4%	29.7
Total	399,433,878								
average	49,929,197	72.4%	43.5	22.5%	77.7%	62.8%	54.9%	44.3%	29.6
std dev	4,365,958	0.8%	0.5	0.8%	0.8%	1.1%	1.3%	1.7%	2.4

Very similar enrichment statistics were obtained from a full slide (half a run) of the SOLiD™ 4 System sequencing on the 8-plex simultaneous exome enrichment sample as compared to those obtained from 2 quads (~40% of a slide) of sequencing on a 4-plex reaction (compare to Table 2). There may have been a minor degree of complexity loss upon scaling to 8-plex based on a slightly lower "on-target" rate and a larger percentage of "target bases not covered". Nonetheless, both samples have nearly identical numbers of bases (~44.4%) that are covered at 20x depth or greater.

Table 6. 8-plex variant calls and dbSNP132 concordance

Barcode	total SNPs	homo dbSNP	heter dbSNP	het SNPs	total indels	homo dbSNP	heter dbSNP	het indels	total concord
BC1	31,437	12,964	98.7%	18,473	90.3%	1509	547%	58.4%	1060
BC2	32,138	13,090	98.7%	19,048	90.3%	1600	566%	58.6%	1060
BC3	30,787	12,789	98.6%	17,998	90.3%	1508	566%	58.6%	1060
BC4	31,218	12,831	99.0%	18,387	90.9%	1552	564%	58.6%	1060
BC5	32,462	13,253	98.7%	19,209	90.2%	1628	566%	58.6%	1060
BC6	30,866	12,844	98.7%	17,712	89.8%	1488	566%	58.6%	1060
BC7	31,724	12,938	98.9%	18,786	90.4%	1514	577%	58.6%	1060
BC8	31,464	13,017	98.7%	18,447	90.6%	1589	565%	58.6%	1060
average	31,474	12,862	98.7%	18,612	90.3%	1589	565%	58.6%	1060
std dev	633	139	0.1%	611	0.3%	52.3	0.9%	0.9%	60

Again, the variant calls of this 8-plex exome capture compare well to the 4-plex capture overall (see Table 3), particularly in degree of concordance with dbSNP and number of indels called (96.3% as many); however, there are somewhat fewer total SNPs called (81.5% as many homozygous SNPs and 88.5% as many heterozygous SNPs). Taken along with the differences in enrichment statistics, this suggests, not surprisingly, that there is a trade-off made between the degree of multiplexing one can do and the degree to which one confidently identifies variants with exome enrichment.

Table 4. 4-plex library pooling prior to dry-down

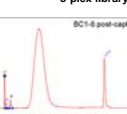


Figure 5. Exome-enriched 8-plex library

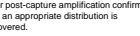


Table 7. SOLiD™ 4 and 5500x1 System enrichment on a single exome (SureSelect 38 Mb)

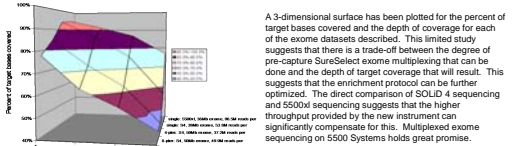
	Total Reads	Percent mapped	Percent on target	fold enrichment	Percent target bases not covered	coverage >=1x	coverage >=5x	coverage >=10x	coverage >=20x	average coverage
SOLID 4 (1 quad)	83,098,814	64.8%	72.0%	57.9	6.0%	94.0%	86.2%	76.0%	54.9%	33.2
5500x1 (1 lane)	123,464,339	73.9%	73.5%	59.2	5.1%	95.0%	91.1%	86.0%	76.9%	66.9

Table 8. SOLiD™ 4 and 5500x1 System variant calls on a single exome (SureSelect 38 Mb)

	total SNPs	homo dbSNP	heter dbSNP	het SNPs	total indels	homo dbSNP	heter dbSNP	het indels	total concord
SOLID 4 (1 quad)	26,681	12,164	97.9%	14,520	89.3%	606	215	89.8%	385
5500x1 (1 lane)	32,619	12,433	99.1%	20,186	89.9%	1,406	421	87.7%	639

The increased throughput of the 5500 Series SOLiD™ Sequencers permits more exome data to be obtained per run. As an example, one lane of 5500x1 sequencing (there are 6 lanes per flow-cell) yields an average coverage >=66x for a 38Mb version of the exome) that is approximately twice that of a quad of SOLiD™ 4 sequencing. This added depth permits more SNPs (~4,000 and indels ~80) to be called as well.

Figure 7. Landscape of exome target base coverage for all data shown



CONCLUSIONS

- Barcoded SOLiD™ System libraries can be pooled prior to exome enrichment with the Agilent SureSelect Human All Exon 50 Mb Kit. 4-plex and 8-plex simultaneous capture is possible.
- Multiplex capture yields reproducible results. Good barcode balance is observed and SOLiD™ System barcodes 1-8 do not bias performance.
- The 5500 Series SOLiD™ System yields the largest amount of high-quality data observed for single exome sequencing. Sequencing 2 exomes per 5500 flow-cell lane can yield >30X average coverage for the SureSelect 38 Mb exome. This is a throughput of ~24 exomes per 5500x run.
- Multiplex exome capture leads to gains in throughput that are balanced against depth of coverage. A full slide of SOLiD™ 4 System sequencing will yield an average coverage of ~30X for 8-plex and ~50X for 4-plex for the SureSelect 38 Mb exome.

REFERENCES

- http://www.appliedbiosystems.com/aba/us/en/home/applications-technologies/solid-next-generation-sequencing/targeted-resequencing.html
- Nosichenko A, et al. Nat. Genet. 2010;42(6):483-85. Epub 2010 May 2.
- Teer JK and Mulligan JC. Hum Mol Biol. 2010;19(14):151-151. Epub 2010 Aug 12.
- Beil C, et al. Sci Transl Med. 2011;3(65):114
- Blankenship MK, et al. Genome Biol. 2011;11(6):R82. Epub 2010 Jun 17.

TRADEMARKS/LICENSING

For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.
© 2011 Life Technologies Corporation. All rights reserved.
The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.