

Sequencing of Methylated Human DNA Enabled through MBD-Affinity is a Cost-Effective Alternative to Whole Genome Bisulfite Conversion



Gavin D. Meredith, Mirosław Dudas, Elizabeth Levandowsky, Tamara Gilbert, Daniel Krissinger, George Marnellos, Vrunda Sheth, Christopher Adams, Clarence Lee, Timothy Harkins and Rob Bennett, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, CA, USA, 92008

ABSTRACT

Aberrant DNA methylation is characteristic of many cancers and other complex diseases and differences in methylation have been observed in a wide variety of genomic contexts; for example, both within "classic" promoter-associated CpG islands and also in distal, non-CpG island regions including CpG island shores [1, 2]. High-throughput sequencing of DNA fragment libraries from the well-characterized and epigenetically profiled IMR-90 fibroblast cell-line [3] proves that methyl-CpG binding domain (MBD) proteins bind methylated DNA fragments sensitively (binding fragments with as few as 2 methyl-CpGs per MBD) and selectively (~50-fold over non-methylated fragments). Furthermore, when MBD-captured DNA is sub-fractionated by step-wise salt elution, genomic DNA fragments separate based on their average methyl-CpG content. Peak analysis of the mapped reads permitted accurate identification of thousands of regions of methylation in different genomic locus classes. The lowest salt fraction (eluted with 350 mM NaCl) is depleted of CpG islands but otherwise samples the genome (introns, promoters, shores, repeats, and intergenic regions) evenly, intermediate salt fractions (eluted successively with 450 mM and 600 mM NaCl) show moderate degrees of enrichment (2- to 3-fold) for exons and CpG island shores, and the highest salt fraction (eluted with 2M NaCl) is greatly enriched (14-fold) for CpG islands and moderately enriched (4- to 10-fold) for shores and exons. Methylation at a large number of these positions was confirmed by whole-genome bisulfite sequencing on the SOLiD™ System along with comparison to the published methylation. Bisulfite-sequencing on the SOLiD™ System yields high-quality data on the exact positions of methylation as evidenced by high concordance (~94%) of our results with the published dataset; this high similarity between biological samples cultured and analyzed independently indicates a remarkably stable methylation pattern for this differentiated cell-line. MBD-affinity enrichment coupled with high-throughput sequencing enables efficient surveying of genomic DNA methylation and yields 5- to 10-fold improvement in throughput and cost.

INTRODUCTION

DNA methylation plays a critical role in gene regulation that influences normal organism development and many diseases including cancer. Profiling the DNA methylation patterns of higher organisms is challenging because methylation patterns vary between tissues and with developmental state, hence there are far more methylomes to be analyzed than genomes. Furthermore, in order to map methylation positions with high precision and accuracy, greater depth of sequencing is required than for normal genome sequencing [4]. Affinity-based enrichment of methylated DNA sequences prior to high-throughput sequencing, as with the SOLiD™ System, provides an avenue to pursue this kind of genome-wide information in a minimally biased and cost-efficient manner. The workflow described here using MethylMiner™ enrichment with stepwise salt gradient elution enables the partitioning of the genome into low, moderate, and high density of methylation. This permits blind discovery of methylated regions and permits detection of differentially methylated regions (DMRs) between samples and across genomic feature subsets that harbor differing degrees of methylation density.

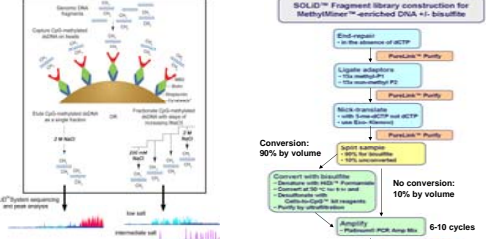
MATERIALS AND METHODS

Methylated DNA enrichment and SOLiD™ System sequencing. Genomic DNA from cultured IMR-90 cells was purified with PureLink® columns. Purified HuRef genomic DNA was purchased from the Coriell Institute for Medical Research. Genomic DNA was fragmented to 50-400 bp (mean ~250 bp) with a Covaris® S2 System (Woburn, MA). Methyl-CpG binding-domain protein affinity capture was with MethylMiner™ Methylated DNA Enrichment kits (Invitrogen, Carlsbad, CA) following the manufacturer's protocol. For salt-gradient elution of IMR-90 DNA, successive fractions were obtained by elution using buffer containing the following NaCl concentrations: 200 mM, 350 mM, 450 mM, 600 mM, and finally 2 M NaCl. For HuRef DNA [5], successive elutions were done with buffer containing 450 mM, 600mM, and 2M NaCl. Each elution step consisted of 2-3 serial incubations of the MethylMiner™ beads at each salt concentration.

MATERIALS AND METHODS (continued)

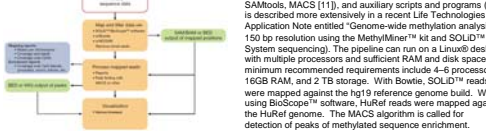
Whole-genome SOLiD™ bisulfite sequencing. Genomic DNA from cultured IMR-90 cells was purified with PureLink® columns. Genomic DNA was fragmented to 50-400 bp (mean ~250 bp) with a Covaris® S2 System (Woburn, MA). The workflow is depicted in Figure 2. Five (5) micrograms of genomic DNA fragments was end-repaired using reagents from a SOLiD™ Fragment Library Construction kit with a dNTP-mix lacking dCTP, then PureLink® bisulfite-S2-seq on methylation-enriched DNA is described in the Methods section of Ondov et al and is based on a workflow described by Borman Chung et al [6, 7] and a protocol is available on-line [8].

Figure 1. MethylMiner™ Kit Fractionation of CpG-methylated DNA for SOLiD™ System Sequencing



Fragmented double-stranded CpG-methylated genomic DNA is directly and specifically captured on MethylMiner™ MBD-coated magnetic beads then eluted all-at-once with buffer containing 2M NaCl or separated into complementary fractions by step-wise elution with buffers containing progressively increasing concentrations of NaCl up to 2 M. Sequencing after single-step elution using 2M NaCl shows greatest enrichment for densely methylated regions of the sample (lower left). Elution using step-wise salt gradient buffers yields subsets of the methylome with differing degrees of methylation density: sparsely methylated fragments (light blue) elute with low salt, more densely methylated fragments (purple) elute with higher salt, and heavily methylated fragments (red) elute at maximal salt. Selective enrichment prior to sequencing permits clearer identification of these subclasses of methylated genomic DNA fragments.

Figure 3. SOLiD™ System DNA Methylation Analysis Tool



RESULTS

Figure 3. MethylMiner™ Elution profile from 25 µg of human cell-line IMR-90 genomic DNA (and 500 ng, inset)

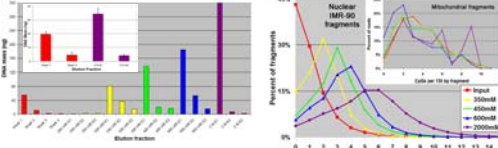
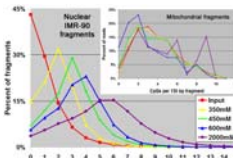
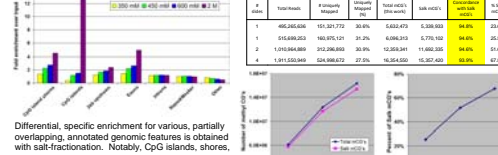


Figure 4. SOLiD™ System DNA Methylation Analysis Tool



Mass of human DNA recovered using a step-wise salt gradient (350mM - 2M NaCl) elution. The majority of input DNA is not captured on the MethylMiner™ beads because only about 0.5-1% of all bases are 5-methyl-C in CpG dinucleotides. This fact, and the observation that diminishing amounts of DNA are recovered in successive wash and step-wise elutions (e.g., fractions 350a, 350b, and 350c in Fig. 2A-B), indicate that the methylated DNA-enrichment protocol worked properly. Importantly, as shown in the inset of Figure 2, MethylMiner™ bead enrichment from as little as 500 ng of fragmented genomic DNA can yield enough methylated DNA for SOLiD™ library construction.

Figure 5. Enrichment trends for human genome features



Differential, specific enrichment for various, partially overlapping, annotated genomic features is obtained with salt-fractionation. Notably, CpG islands, shores, and exonic sequences increase in relative representation with increasing ionic strength. Single-fraction elution with high-salt (2M NaCl) yields an average behavior in this respect (not shown).

Figure 7. MethylMiner™ Elution profiles from 25 µg of human genomic DNA

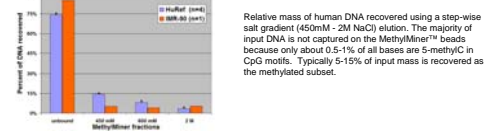


Table 2. Paired-end mapping stats for HuRef libraries with and without ECC

sample	pairings	P1	P2	non-redundant		reads per start		reads per end	
				pairings	point #3	point #3	point #3	point #3	point #3
IMR90_B5_S1P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1
IMR90_B5_S2P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1
IMR90_B5_S3P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1
IMR90_B5_S4P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1
IMR90_B5_S5P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1
IMR90_B5_S6P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1
IMR90_B5_S7P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1
IMR90_B5_S8P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1
IMR90_B5_S9P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1
IMR90_B5_S10P	53,577,406 (50.70%)	49,493,054 (46.82%)	49,493,054 (46.82%)	2,397,712	1.15	1.1	1.1	1.1	1.1

Bisulfite converted

Non-converted

Paired-end sequencing (75x35) shows very high rates of non-redundant pairings for unmethylated MethylMiner™ bead-enriched libraries. ECC consistently permits higher yields of unique mapping too. For bisulfite-converted libraries, there is a reduction in pairing efficiency at higher ionic strengths, presumably due to the reduced sequence complexity both of the selected fragments and caused by bisulfite treatment.

Table 3 and Figure 8. Methyl-Cytosines in HuRef detected using BioScope™ Software SNP-calls

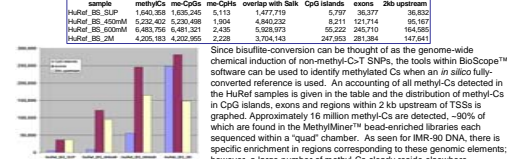


Figure 9 Example: Apparent Differential Methylation in the 5' region of the ISOC2 gene



CONCLUSIONS

- A) Bisulfite-sequencing of human methylomes is tractable on the SOLiD™ A system; it requires 3-4 full runs per methylome. The results are highly concordant with published data.
- B) MBD-based enrichment of methylated sequences with the MethylMiner™ kit is an efficient means to focus of SOLiD™ System sequencing on genomic feature subsets.
- C) MBD-based enrichment permits sub-fractionation (by varying ionic strength) of the genome based on the density of methylated CpGs.
- D) Methylated DNA enrichment can be coupled with bisulfite-sequencing permitting single-nucleotide resolution validation of specific positions of methylation and 4-fold reduction in sequencing cost.

TRADEMARKS/LICENSING

For Research Use Only. Not for release or further distribution or diagnostic use.
© 2011 Life Technologies Corporation. All rights reserved.
The trademarks shown here are the property of Life Technologies Corporation or their respective owners.