

Enzymatic Methyl-seq: Next Generation Methylomes

Fiona J. Stewart¹, V K Chaithanya Ponnaluri¹, Brittany S. Sexton¹, Louise Williams¹, Bradley W. Langhorst¹, Romualdas Vaisvila¹, Yvonne Helbert², Lei Zhang², Timothy Harkins², Kevin J. McKernan², Eileen T. Dimalanta¹, Theodore B. Davis¹

¹ New England Biolabs, Inc., Ipswich, MA 01938 ² Medicinal Genomics Corp., Woburn, MA 01801

Introduction

DNA methylation is important for gene regulation. The ability to accurately identify 5-methylcytosine (5mC) and 5-hydroxymethylcytosine (5hmC) gives us greater insight into potential gene regulatory mechanisms. Bisulfite sequencing (BS) is traditionally used to detect methylated Cs, however, BS does have its drawbacks. DNA is commonly damaged and degraded by the chemical bisulfite reaction resulting in libraries that demonstrate high GC bias and are enriched for methylated regions. To overcome these limitations, we developed an enzymatic approach, NEBNext[®] Enzymatic Methyl-seq (EM-seq[™]), for methylation detection that minimizes DNA damage, resulting in longer fragments and minimal GC bias.

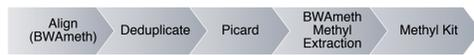
Illumina libraries were prepared using bisulfite and EM-seq methods with 50 ng DNA from *Arabidopsis thaliana* and *Cannabis sativa* DNA. Libraries were sequenced using Illumina's NextSeq 500 (2x75). Reads were aligned using BWA-Meth 0.2, and methylation information was extracted from the alignments using MethyDackel. Total 5mC levels were compared between the sequencing data from EM-seq and WGBS libraries and LCMS (*Liquid Chromatography Mass Spectrometry*). 5mC levels determined by EM-seq are close to those from LCMS, whereas, WGBS results in an overestimation of 5mC. Additionally, EM-seq libraries produce higher quality sequencing metrics such as longer inserts, lower duplication rates, a higher percentage of mapped reads and less GC bias compared to bisulfite converted libraries. We conclude that EM-seq is superior to WGBS and delivers higher library yields, more accurate methylation information, reduced DNA damage, increased sequencing length, and decreased GC-bias.

Sample Preparation

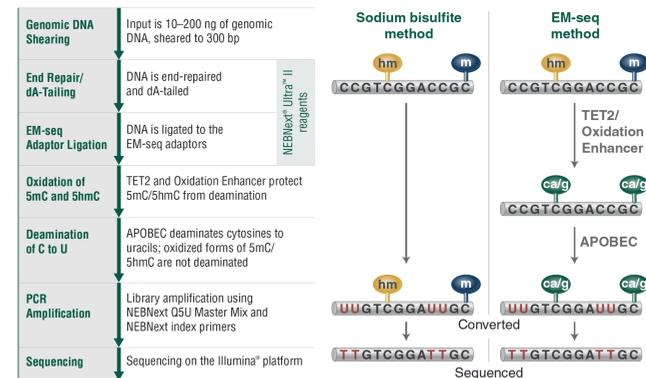
- Two plant DNAs were used to make EM-seq libraries
 - Cannabis sativa* genomic DNA (Jamaican Lion): female clones (leaf, seeded and unseeded flowers) & male sibling (flowers) plants
 - Arabidopsis thaliana*
- Libraries were made using 50 ng genomic DNA, spiked with control DNA (unmethylated lambda & CpG-methylated pUC19)
- Libraries were sequenced using an Illumina NextSeq 500, 2x76 base paired reads. 5caC is sequenced as C and deaminated C as T.
- Bisulfite conversion was performed using Zymo Research EZ DNA Methylation-Gold[™] kit

Data Analysis

- Reads were aligned to Jamaican Lion reference genome (August 2018 assembly) or the *Arabidopsis* reference genome (TAIR10) (for Jamaican Lion, four miscellaneous contigs were removed from methylation analysis)
- Data were analyzed using the tools in above flowchart.



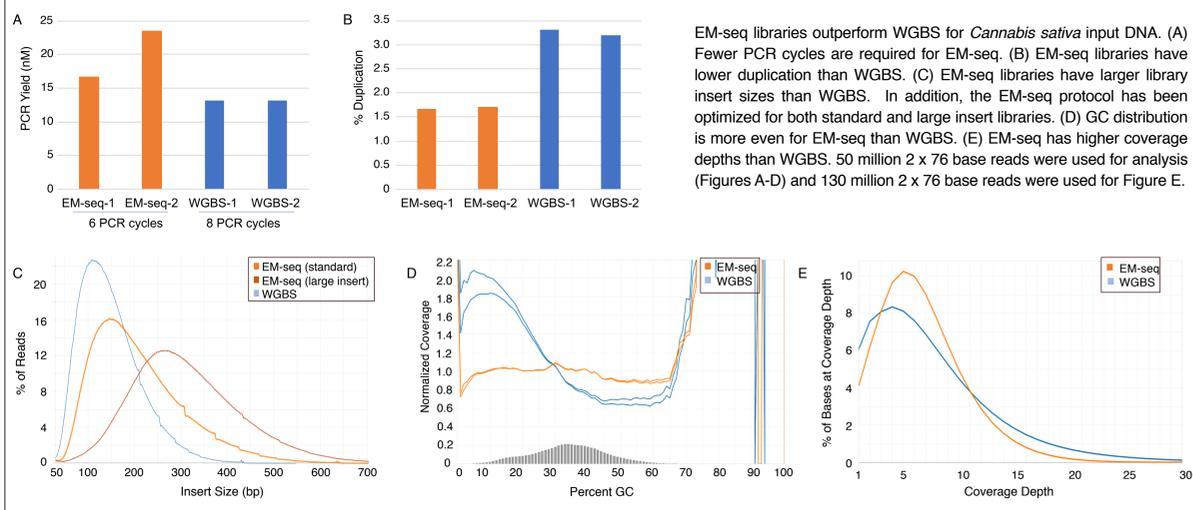
Methods



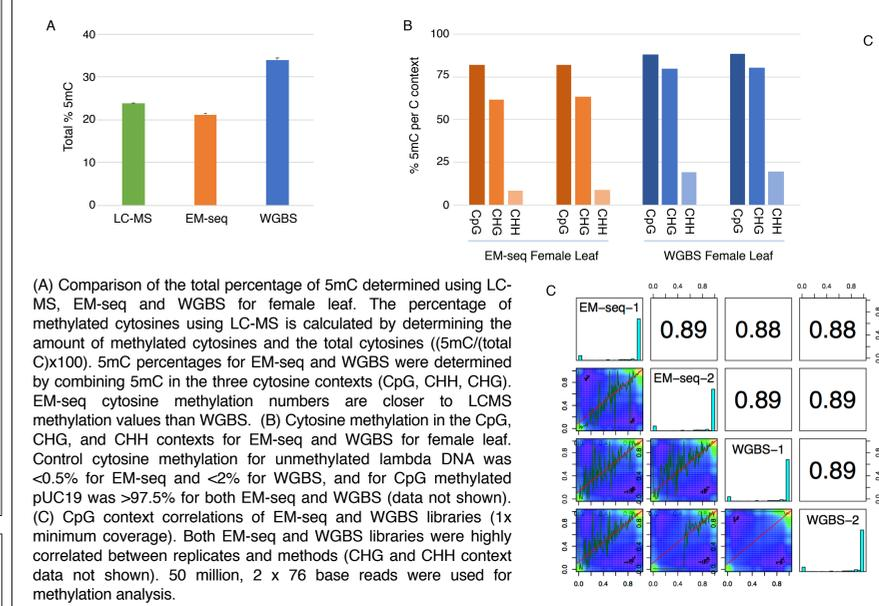
Methods

Cannabis sativa

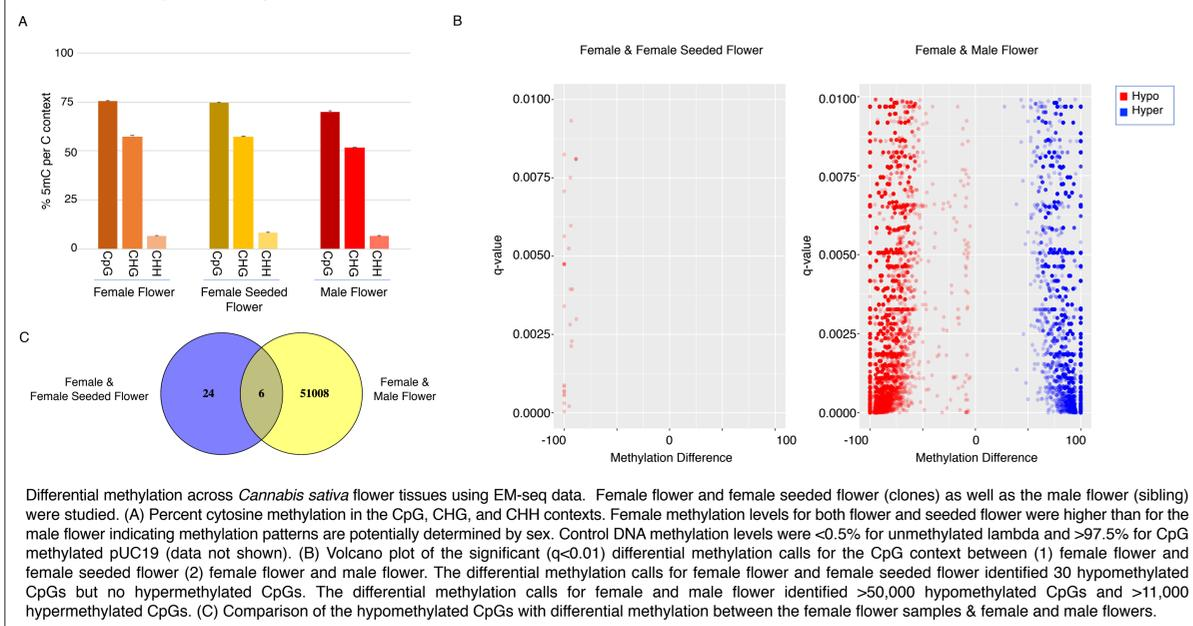
Cannabis sativa: Higher Quality Sequencing Metrics with EM-seq compared to WGBS



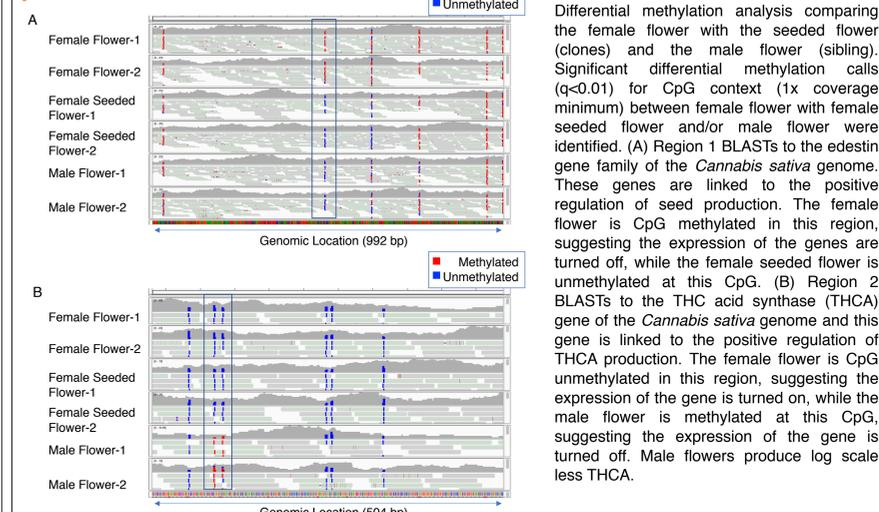
Cannabis sativa female leaf EM-seq libraries are superior to WGBS



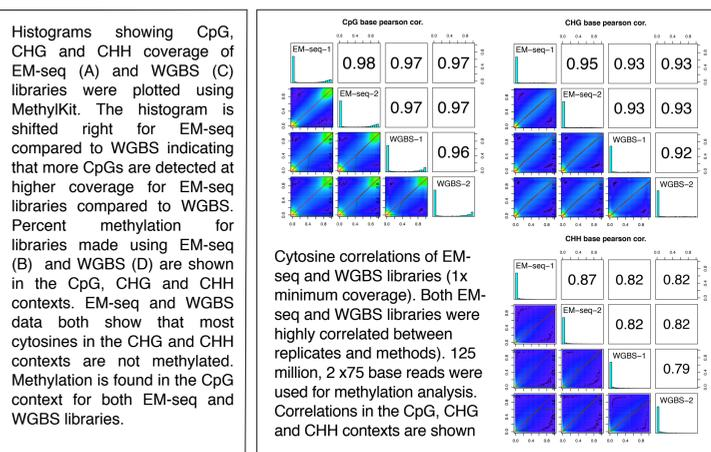
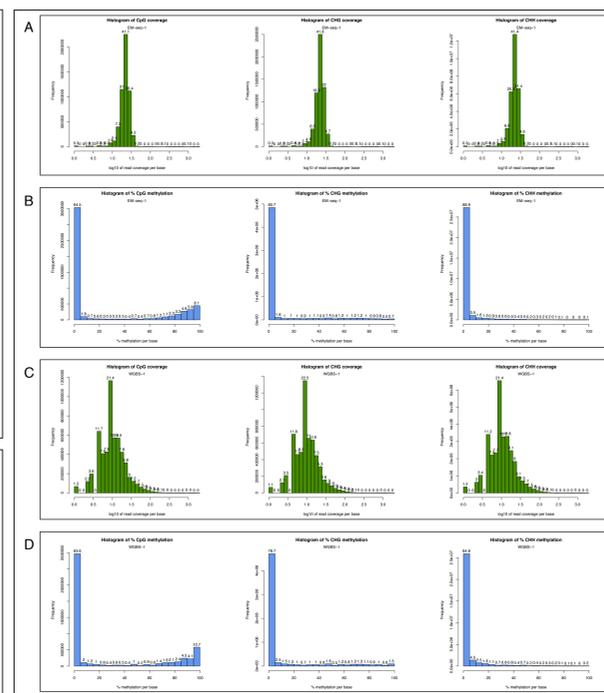
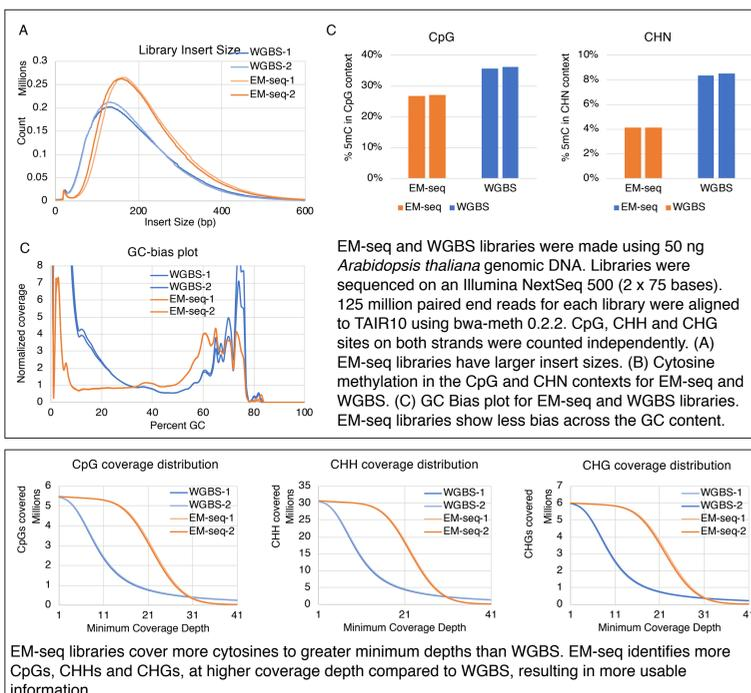
Differential CpG methylation identified between Cannabis flower tissues



Methylation profiles identified genes involved in seed & cannabinoid production



Arabidopsis thaliana



Conclusion

- EM-seq can be used to investigate plant genomic DNA
 - analysis of the *Cannabis sativa* methylome identified genes involved in seed and THC production
 - the *Arabidopsis thaliana* methylome was successfully probed

- EM-seq libraries compared to WGBS libraries had:
- Higher library yields with fewer PCR cycles
 - Lower percent duplication
 - More even base coverage
 - Larger library insert sizes
 - Less GC bias
 - Similar percentage methylation as LC-MS