

From Genome to Phenome: Revolutionizing Agricultural Epigenetics with Direct Multiomic Mapping of DNA Methylation and Chromatin Proteins

Andrea L Johnstone¹, Vishnu U Kumary¹, Allison Hickman¹, Anup Vaidya¹, Sriharsa Pradhan², Michael-Christopher Keogh¹ and Bryan J. Venters¹
¹EpiCypher Inc., Durham, NC, USA; ²New England Biolabs, Ipswich, MA, USA



Gene expression is controlled by crosstalk via DNA methylation and chromatin proteins

- ▶ DNA methylation (DNAm) and chromatin proteins (e.g. histone PTMs and transcription factors) oversee gene expression
- ▶ These factors are associated with specific genomic features (Figure 1)
- ▶ Current methods are low-resolution or rely on indirect parallel assays
- ▶ Improved technologies are needed to directly resolve DNAm / chromatin protein crosstalk

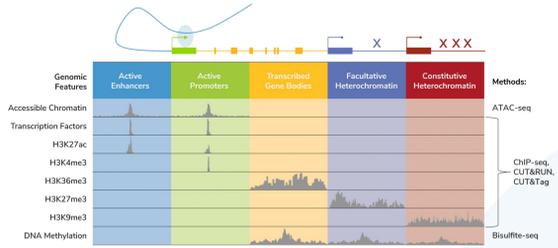


Figure 1. Chromatin proteins and DNAm define genomic features and reveal important regulatory mechanisms governing gene expression; however, currently available assays are correlative. Enrichment profiles shown are characteristic for mammals; some differences exist in plants.

CUT&RUN-EM simultaneously profiles direct DNAm / chromatin protein interactions

CUTANA™ CUT&RUN

Did you know?
 Compared to ChIP-seq, CUT&RUN generates higher resolution data with >100-fold fewer cells and >10-fold reduced sequencing depth.

NEBnext® Enzymatic Methyl-seq (EM-seq) v2

Did you know?
 EM-seq is a more sensitive, less destructive alternative to bisulfite sequencing.

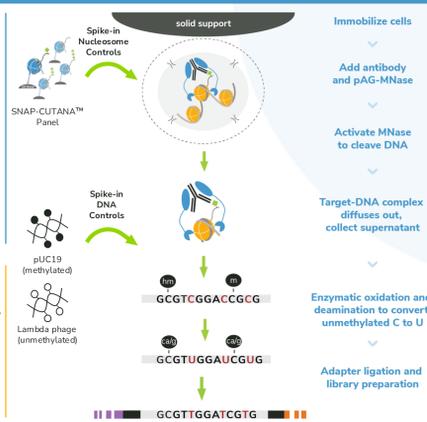


Figure 2. CUT&RUN-EM integrates EpiCypher CUT&RUN with NEB EM-seq v2 kit. CUT&RUN isolates DNA associated with chromatin proteins of interest. EM-seq is then used to enzymatically convert unmethylated cytosines to uracils. Sequencing the resulting libraries enables CpG resolution of DNAm co-occurring with the chromatin mark of interest.

CUT&RUN-EM is highly reproducible across targets, sequencing depths, and DNA inputs

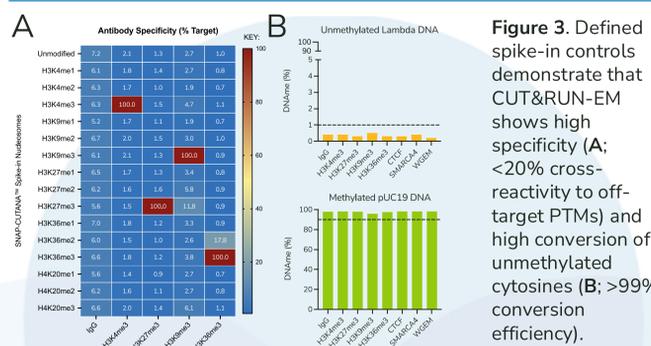


Figure 3. Defined spike-in controls demonstrate that CUT&RUN-EM shows high specificity (A; <20% cross-reactivity to off-target PTMs) and high conversion of unmethylated cytosines (B; >99% conversion efficiency).

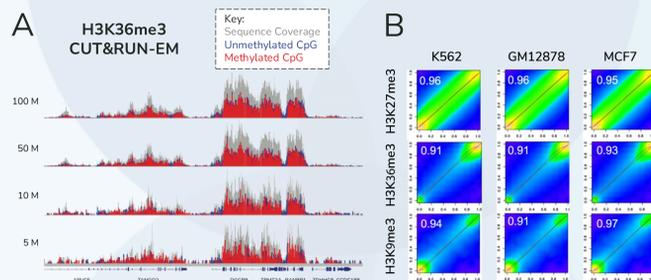


Figure 4. CUT&RUN-EM assays are highly reproducible across various sequencing depths (A; downsampled genome browser tracks) and replicates (B; Pearson correlation coefficients).

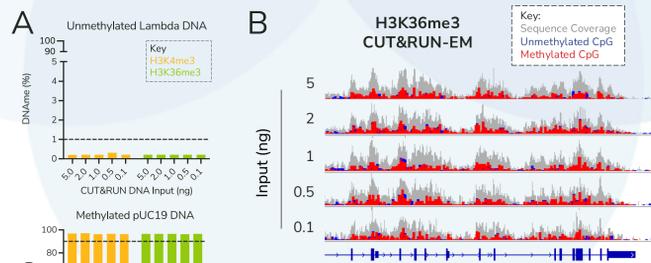


Figure 5. CUT&RUN-EM excels at ultra-low DNA inputs; demonstrating consistent conversion efficiency (A) and genomic enrichment (B) down to 0.1 ng input using the EM-seq v2 kit.

CUT&RUN-EM reveals distinct DNAm profiles at defined genomic features

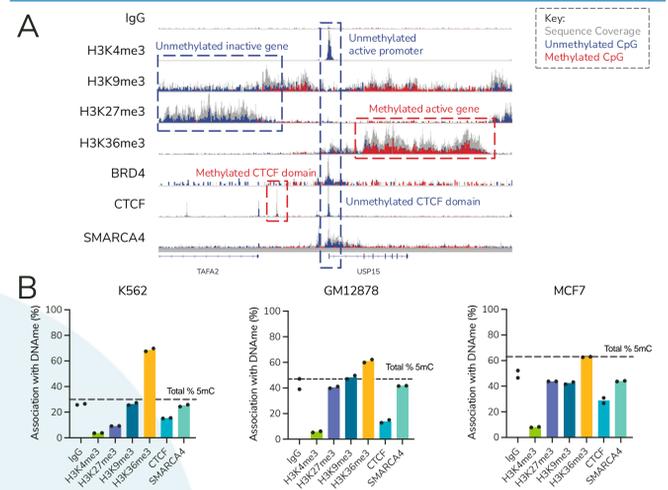


Figure 6. Representative genome browser tracks (A) demonstrate that CUT&RUN-EM reflects known biological functions of chromatin proteins (e.g. unmethylated H3K4me3 promoters, methylated H3K36me3 gene bodies) when overlaid with DNAm. The association of chromatin proteins with DNAm varies across cell lines (B), highlighting the utility of CUT&RUN-EM to provide deep gene regulatory insights.

CUT&RUN-EM deconvolutes epigenetic crosstalk that is masked in correlative assays

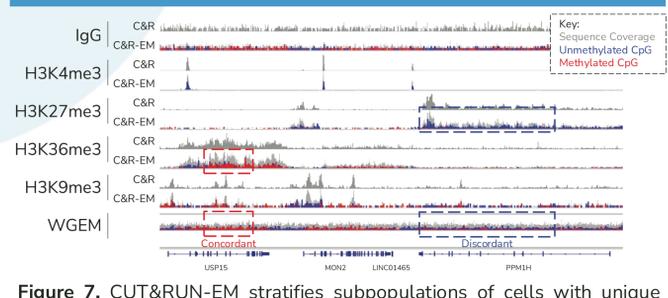


Figure 7. CUT&RUN-EM stratifies subpopulations of cells with unique PTM-DNAm crosstalk signatures that are obscured in whole genome EM-seq (WGEM). IG genome browser comparison of CUT&RUN-EM (C&R), CUT&RUN-EM (C&R-EM), and WGEM in K562 cells. C&R-EM loci with concordant results (red box) recapitulate the findings of WGEM, while discordant loci (blue box) resolve PTM-DNAm specific signatures.

MeCP2-targeted CUT&RUN-EM generates ultra-sensitive, global DNA methylation profiles at low cost

Higher quality, bp resolution data using reduced cells & sequencing vs. MeDIP

MeCP2 captures 83% of 5mC with 34x less sequencing vs. WGEM

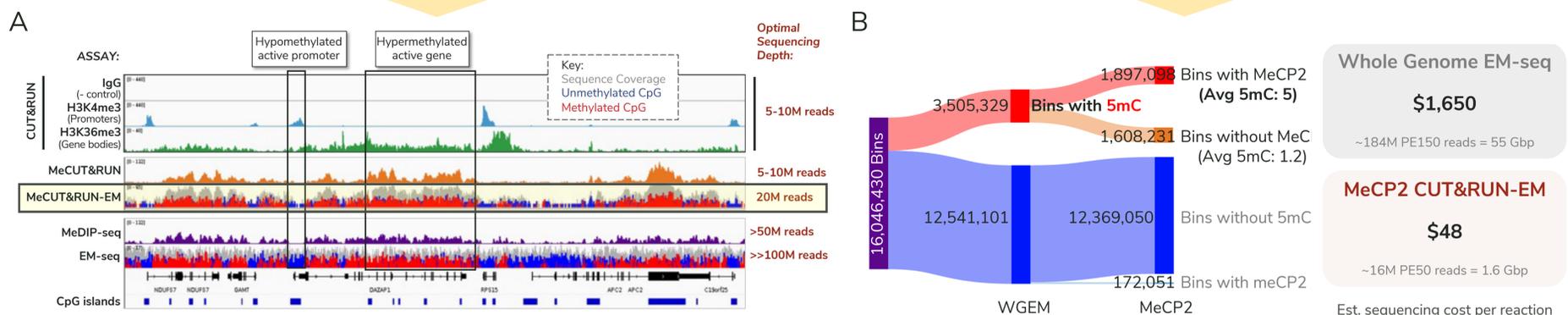


Figure 8. (A) GST-tagged MeCP2 methyl binding domain used in lieu of an antibody in traditional CUT&RUN ("MeCUT&RUN"; orange) with 500k K562 cells provides ~150bp resolution of genome-wide DNAm. MeCUT&RUN shows high concordance with Methyl-DNA Immunoprecipitation sequencing (MeDIP-seq; purple) at >2.5-fold reduced sequencing depth. To achieve CpG resolution of DNAm, MeCP2 CUT&RUN was followed by EM-seq ("MeCP2 CUT&RUN-EM"; yellow boxed track). MeCP2 CUT&RUN-EM generates similar DNA methylation profiles compared to whole genome EM-seq (bottom track). (B) Sankey plot of the binned human genome (200 bp) comparing the presence of 5mC in WGEM and MeCUT&RUN-EM. MeCP2 enriches for regions with higher concentrations of 5mC, identifying 83% of 5mCs found in WGEM with 34x less sequencing, greatly reducing sequencing costs.

Conclusions

- ▶ CUT&RUN-EM reveals the direct association of DNAm and chromatin proteins in a single workflow
- ▶ CUT&RUN-EM can deconvolute bulk cell heterogeneity through multiomic filtering
- ▶ MeCP2 CUT&RUN-EM provides base pair resolution of global DNA methylation using 34x less sequencing than WGEM

See CUT&RUN/Tag in Ag Research

- Arabidopsis imprinting León, *Nucleic Acids Res* 2024 (PMID: [38967011](#))
- Porcine development Innis, *Epigen Chrom* 2024 (PMID: [38773546](#))
- Maize heat tolerance Li, *Plant Cell* 2024 (PMID: [38573521](#))
- Arabidopsis development Zhang, *Mol Plant* 2024 (PMID: [38825830](#))
- Rice fungal infection Xue, *Plant Cell Rep* 2024 (PMID: [38393406](#))
- Wheat regeneration Liu, *Nat Plants* 2023 (PMID: [37142750](#))
- Chicken development Sun, *Poult Sci* 2023 (PMID: [36764138](#))
- Cotton methods Tao, *Plant Methods* 2020 (PMID: [32884577](#))

Download Poster

