

CUT&RUN-EM: An Ultra-Sensitive Multiomic Method that Directly Links Chromatin Features to DNA Methylation

Keith E. Maier¹, Vishnu U. Sunitha Kumary¹, Bryan J. Venters¹, Allison Hickman¹, Anup Vaidya¹, Ryan Ezell¹, Jonathan M. Burg¹, Louise Williams², Chaithanya Ponnaluri², Hang Geong Chin², Pierre Esteve², Isaac Meek², Sriharsa Pradhan², Zu-Wen Sun¹, Martis W. Cowles¹, Michael-Christopher Keogh¹

¹EpiCypher Inc., Durham, NC, USA; ²New England Biolabs, Ipswich, MA, USA



Gene expression is controlled by complex molecular crosstalk between DNA methylation and chromatin proteins

- Precise regulation of DNA methylation (DNAm) and chromatin proteins underlies many gene expression programs
- DNAm and chromatin proteins are associated with specific genomic features (Figure 1)
- Current methods are low-resolution or rely on correlating parallel assays
- Improved technologies are needed to directly resolve DNAm - chromatin protein crosstalk**

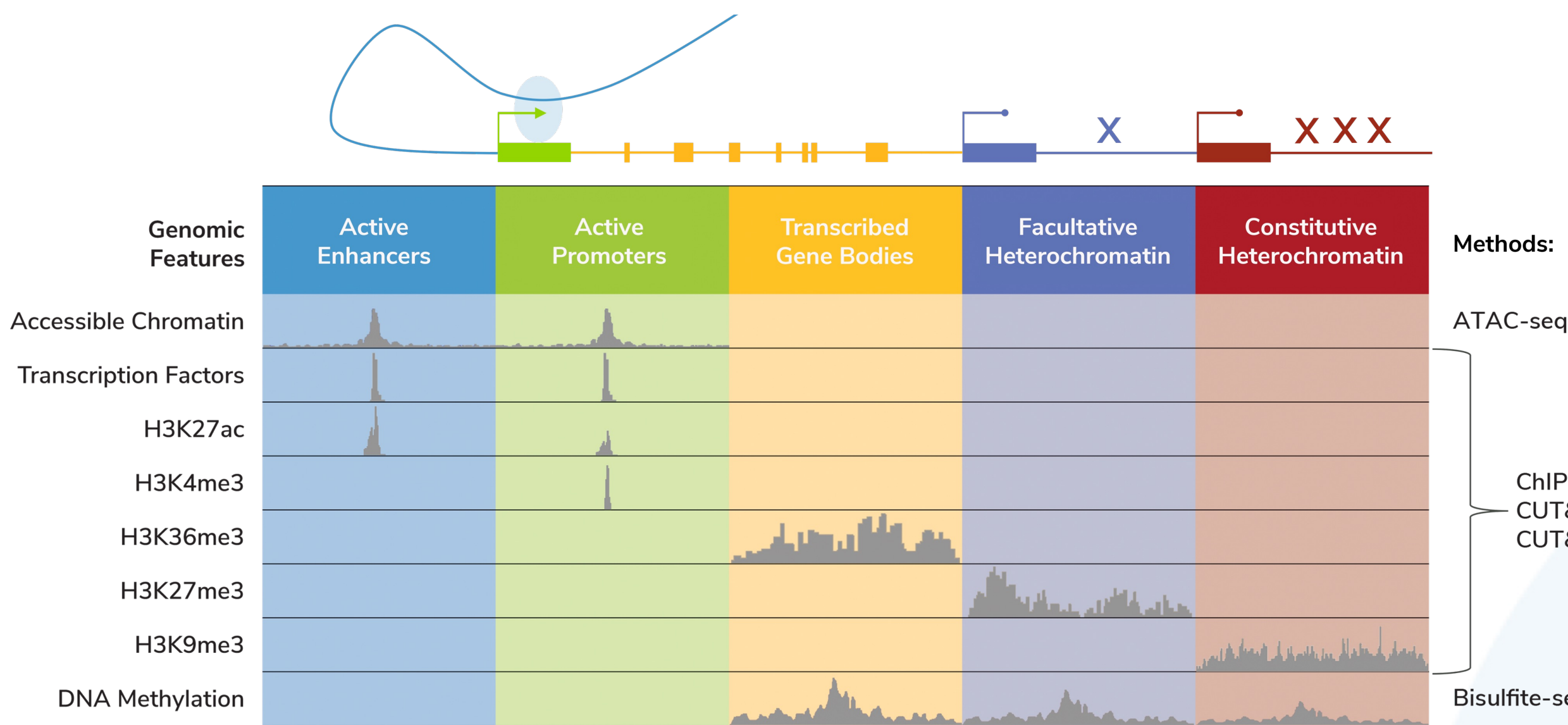


Figure 1. Chromatin proteins and DNAm define genomic features and reveal important regulatory mechanisms governing gene expression; however, currently available assays are correlative.

CUT&RUN-EM is a powerful multiomic assay to simultaneously profile DNAm and chromatin proteins

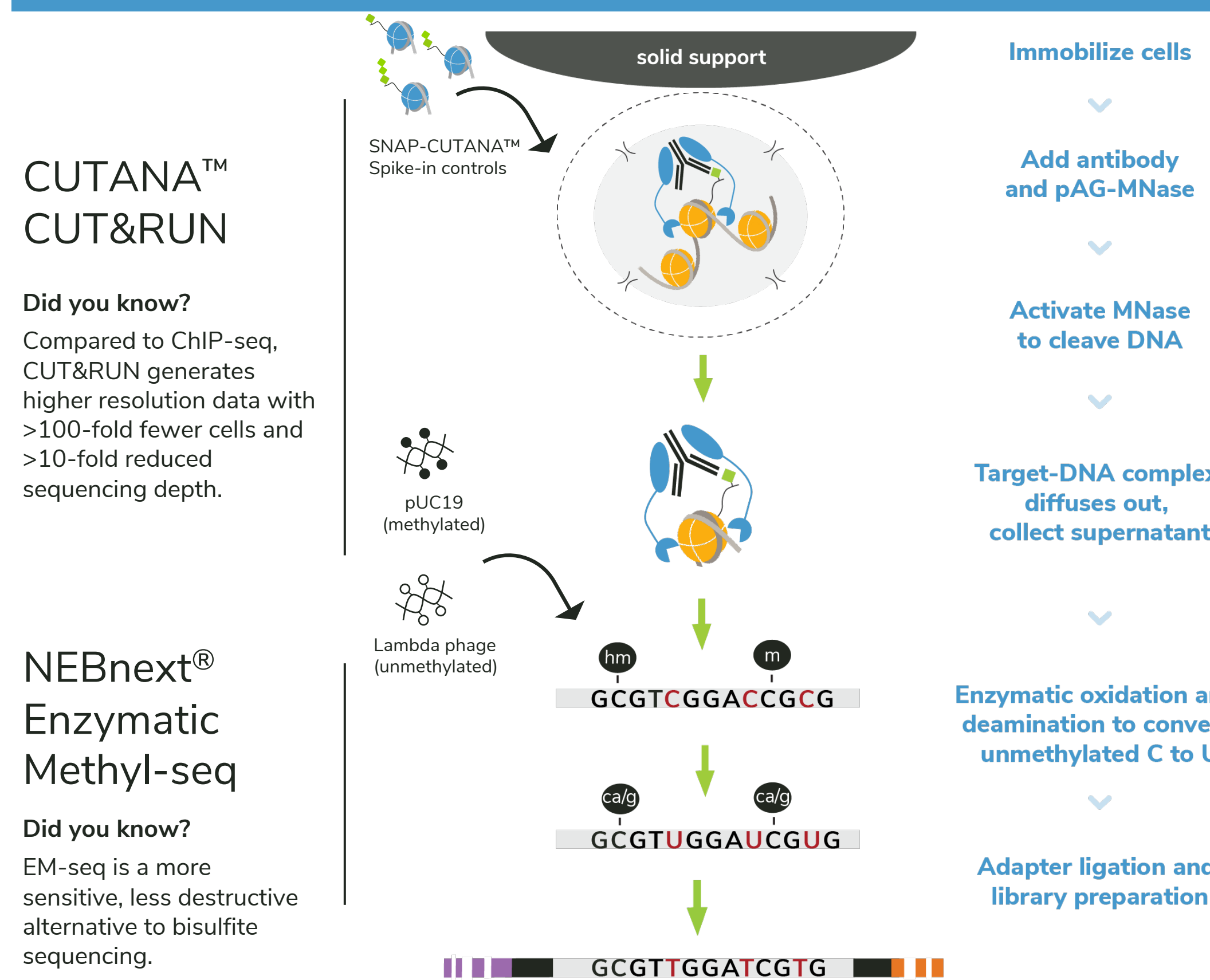


Figure 2. CUT&RUN-EM integrates EpiCypher CUT&RUN with NEB EM-seq to deliver a direct multiomic assay enabling the simultaneous profiling of DNAm and chromatin proteins. CUT&RUN has significantly improved signal to noise compared to ChIP-seq, resulting in reduced cellular input and sequencing requirements. Unmethylated cytosines from the isolated DNA is enzymatically converted, enabling CpG resolution of DNA methylation that co-occurs with the chromatin protein of interest.

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

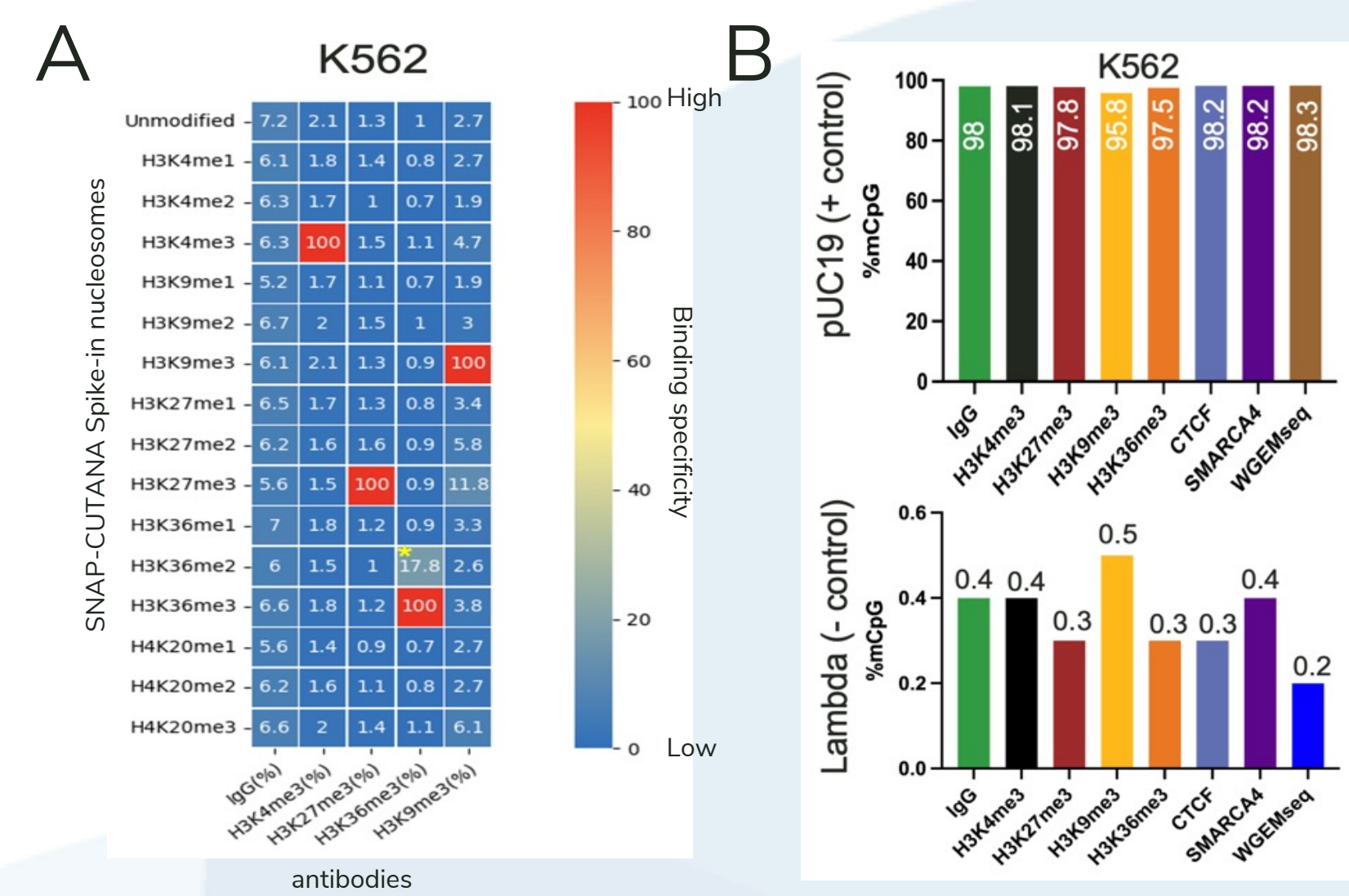


Figure 3. CUT&RUN-EM shows high specificity and percent conversion for a variety of chromatin proteins as determined using defined spike-in controls. (A) Heatmap of SNAP-CUTANA™ Spike-in nucleosomes showing the percent enrichment of each antibody tested in CUT&RUN. Red and blue indicate high and low binding specificity, respectively. (B) Methylated and Unmethylated EM-seq spike-in DNAs were used to monitor enzymatic conversion of unmethylated cytosines. As expected, pUC19 was found to be >95% methylated and Lambda DNA methylation was <0.5%, consistent with the standard for whole genome EM-seq (WGEMseq).

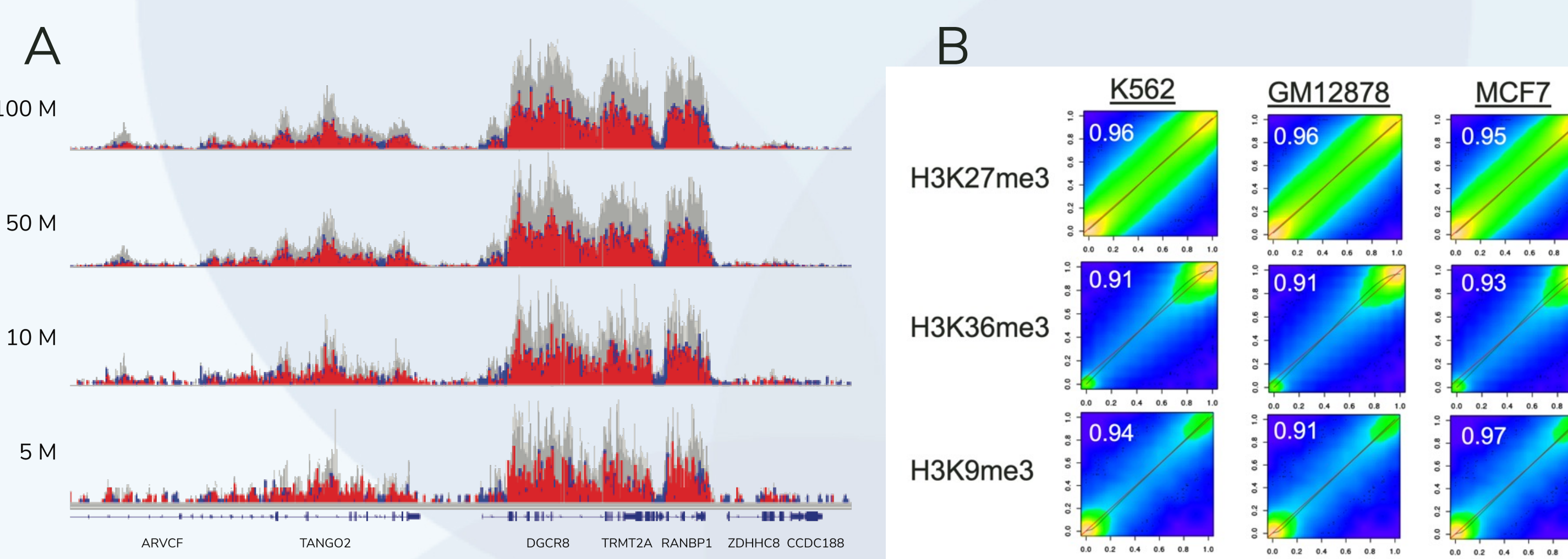


Figure 4. CUT&RUN-EM assays are highly reproducible across sequencing depths and replicates. (A) Representative genome browser tracks from H3K36me3 CUT&RUN-EM data, bioinformatically downsampled, shows qualitatively similar DNAm distribution using as few as 5M reads. (B) Biological replicates in K562, GM12878, and MCF7 cells shows that CUT&RUN-EM assays are highly reproducible ($r > 0.91$).

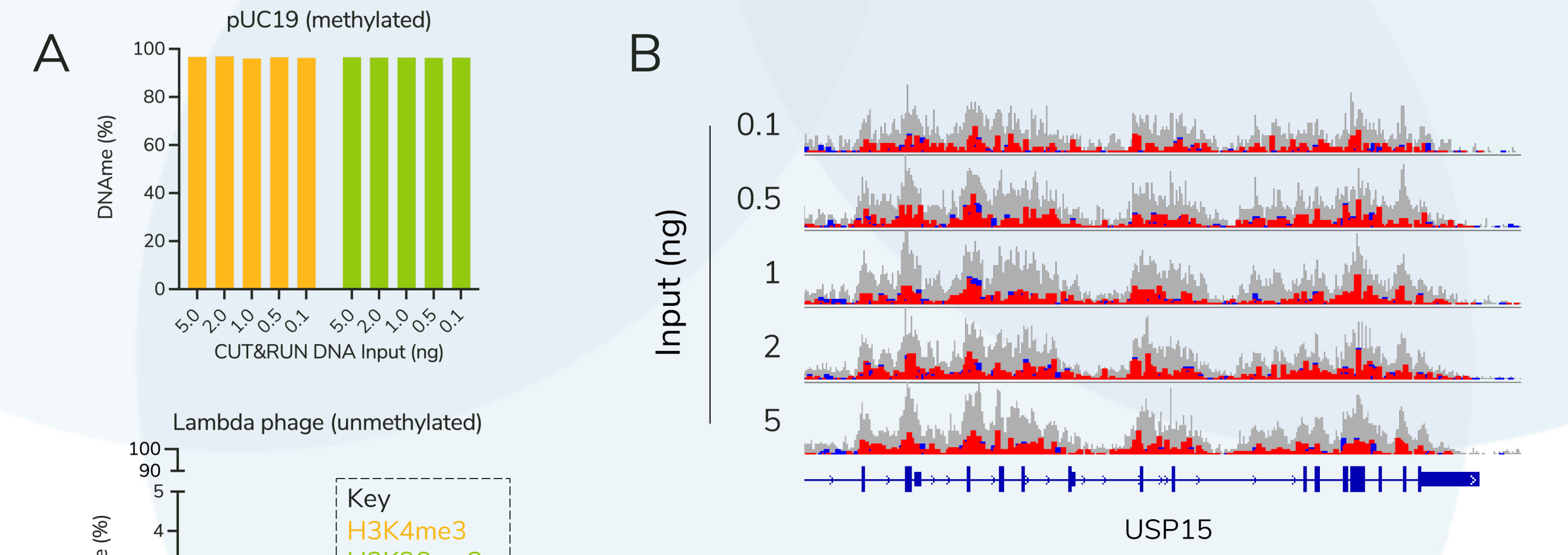
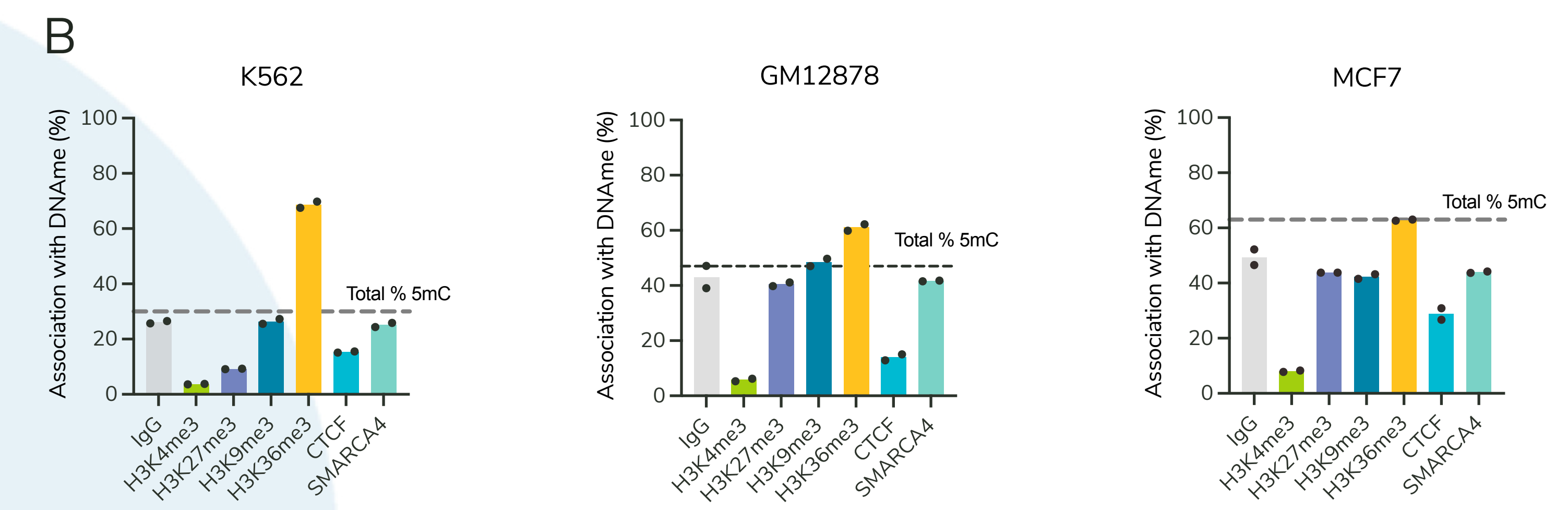
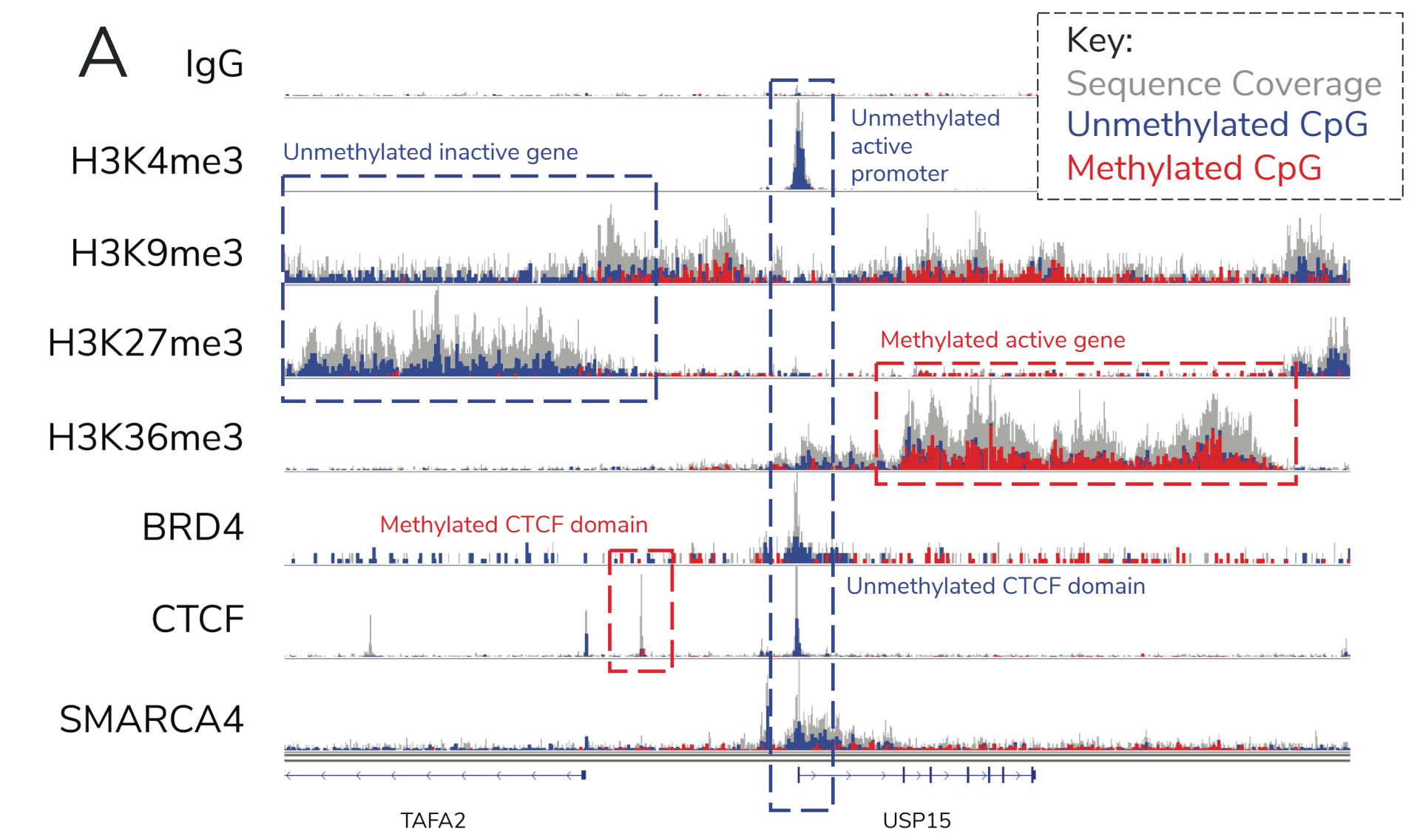


Figure 5. Input titrations reveal CUT&RUN-EM excels at ultra-low DNA inputs. (A) Conversion efficiency of EM-seq controls are unaffected across a range of inputs for H3K4me3 and H3K36me3 CUT&RUN outputs. (B) H3K36me3 genome browser tracks are qualitatively equivalent down to 0.1 ng of EM-seq input.

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

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



CUT&RUN-EM deconvolutes mechanisms of chromatin protein-DNAm crosstalk that would be masked in correlative assays

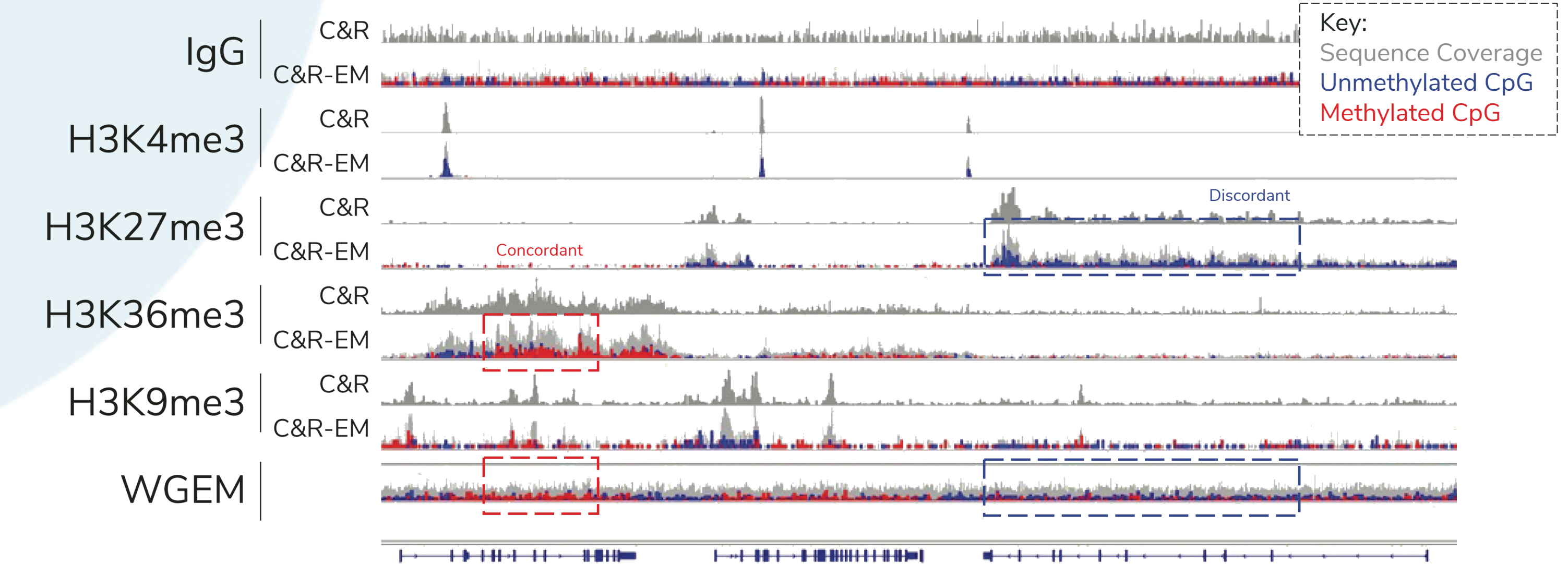


Figure 7. CUT&RUN-EM stratifies specific subpopulations of cells with unique PTM-DNAm crosstalk signatures that are obscured in the global whole genome EM-seq (WGEM) profile. IGV genome browser comparison of CUT&RUN (C&R), CUT&RUN-EM (C&R-EM), and WGEM in K562 cells. C&R-EM loci with concordant results (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 a low cost

Higher quality data with reduced cells & sequencing vs. MeDIP

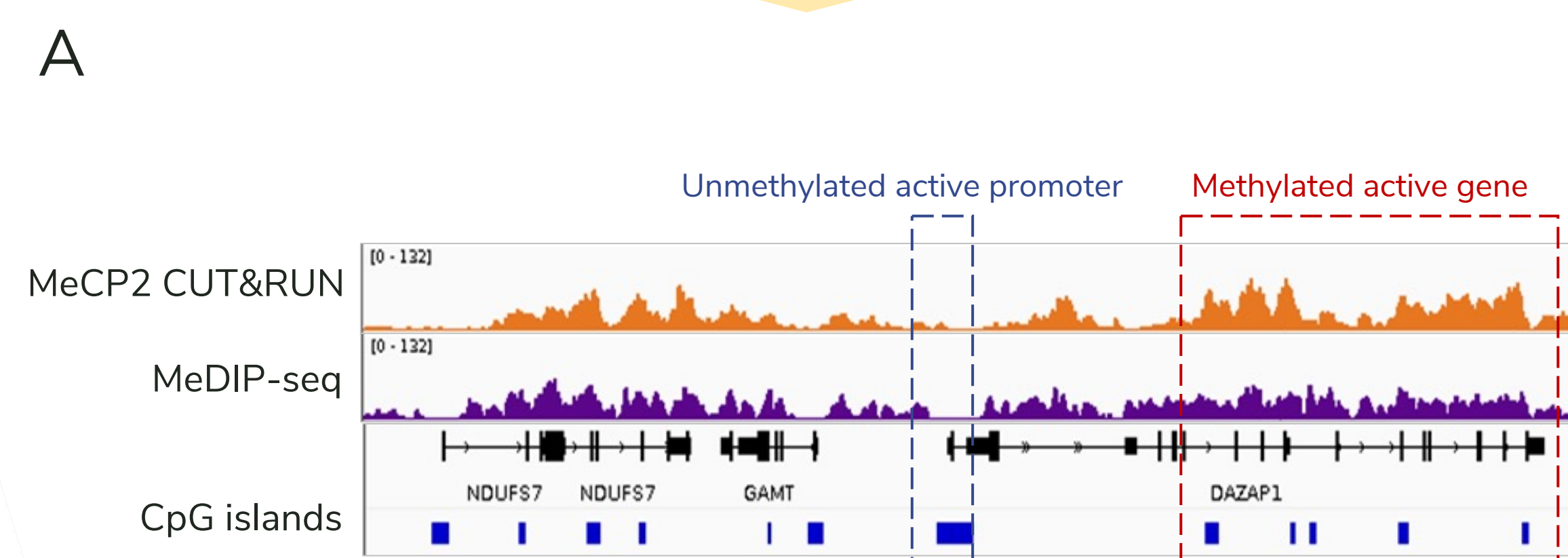
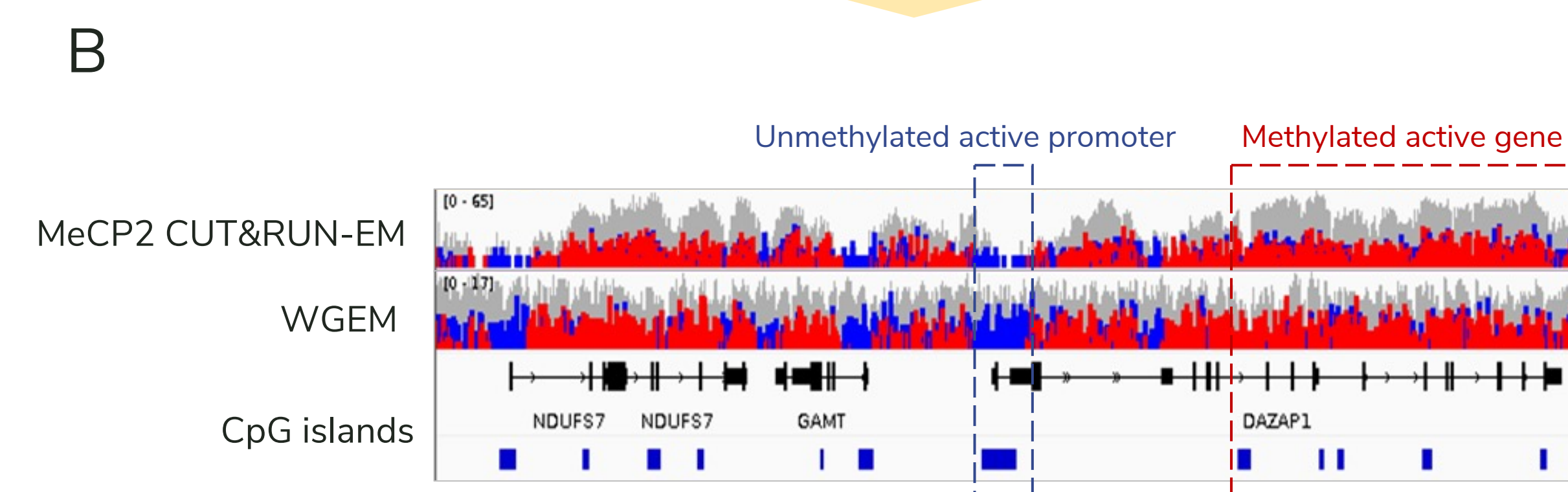


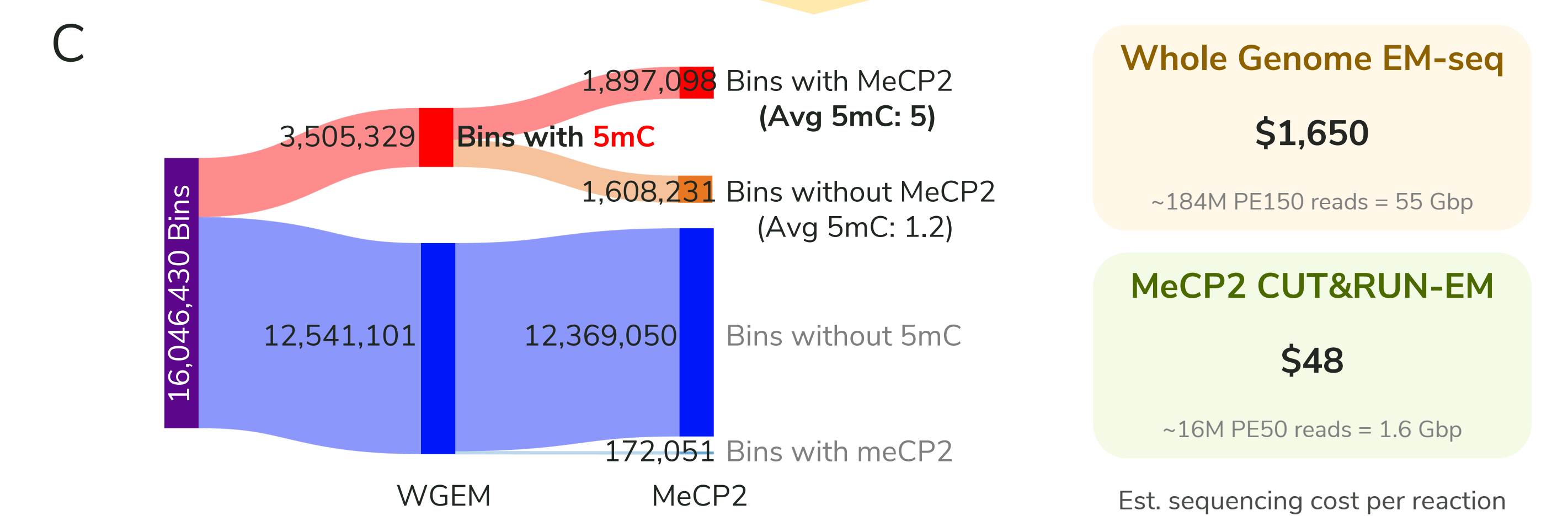
Figure 7. (A) GST-tagged MeCP2 methyl binding domain used in lieu of an antibody in traditional CUT&RUN with 500k K562 cells (orange) provides ~150bp resolution of genome-wide DNAm. MeCP2 CUT&RUN showed high concordance with Methyl-DNA Immunoprecipitation sequencing (MeDIP-seq; purple) at >5-fold reduced sequencing depth. (B) To achieve CpG resolution of DNAm, MeCP2 was used in CUT&RUN-EM. MeCP2 CUT&RUN-EM generates similar DNA methylation profiles compared to WGEM. (C) Sankey plot of the binned human genome (200 bp) comparing the presence of 5mC in WGEM and MeCP2 CUT&RUN. MeCP2 enriches for regions with higher concentrations of 5mC, identifying 83% of 5mCs found in WGEM with 34x less sequencing, greatly reducing sequencing costs.

Combine with EM-seq for base-pair resolution



- ### 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.

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



See EpiCypher assays in action

- AML therapy selection PMID: [38956053](#)
- Epigenomics of stress PMID: [38959894](#)
- SCA1 disease mechanism PMID: [36577402](#)
- Aging & chromatin opening PMID: [38959897](#)
- iPSC differentiation by BRD4 PMID: [39196112](#)
- H3.3K36M disrupts DNAm PMID: [39368466](#)

Download poster

