

Revolutionizing Epigenetics with Direct Multiomic Mapping of DNA Methylation and Chromatin Proteins

Keith E. Maier¹, 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

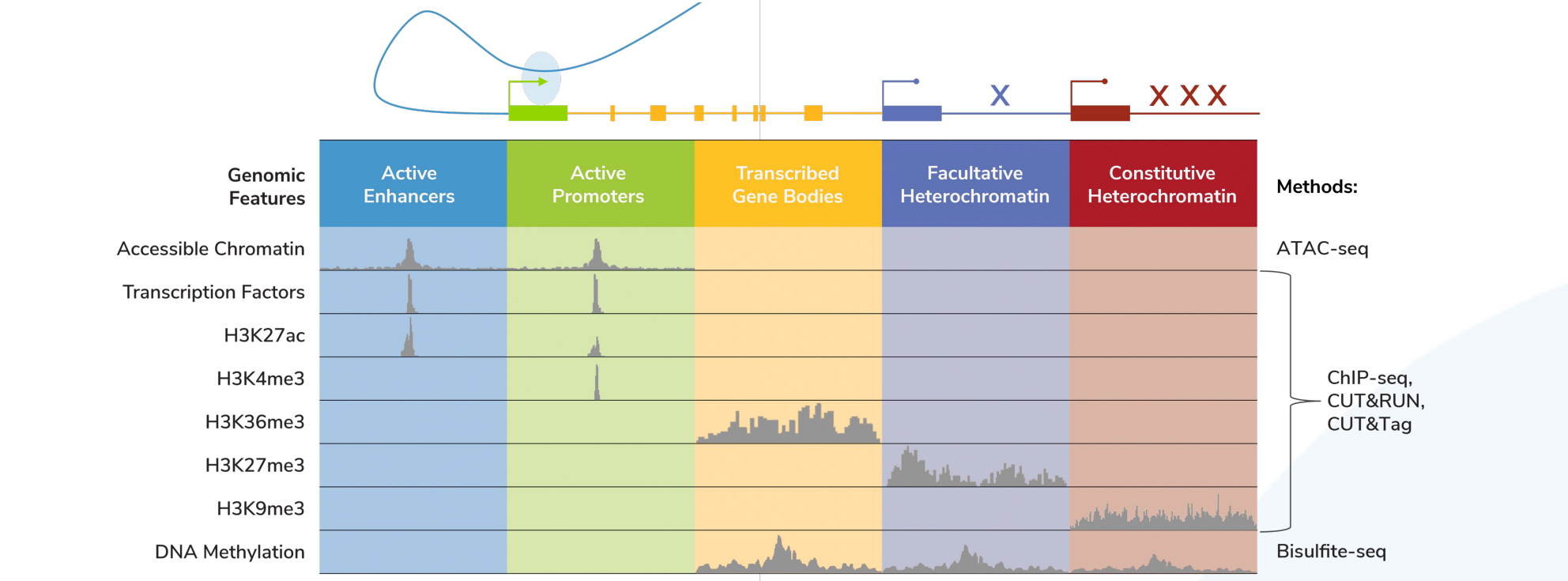


Figure 1. Chromatin proteins and DNAm define genomic features and reveal important regulatory mechanisms governing gene expression; however, currently available assays are correlative.

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

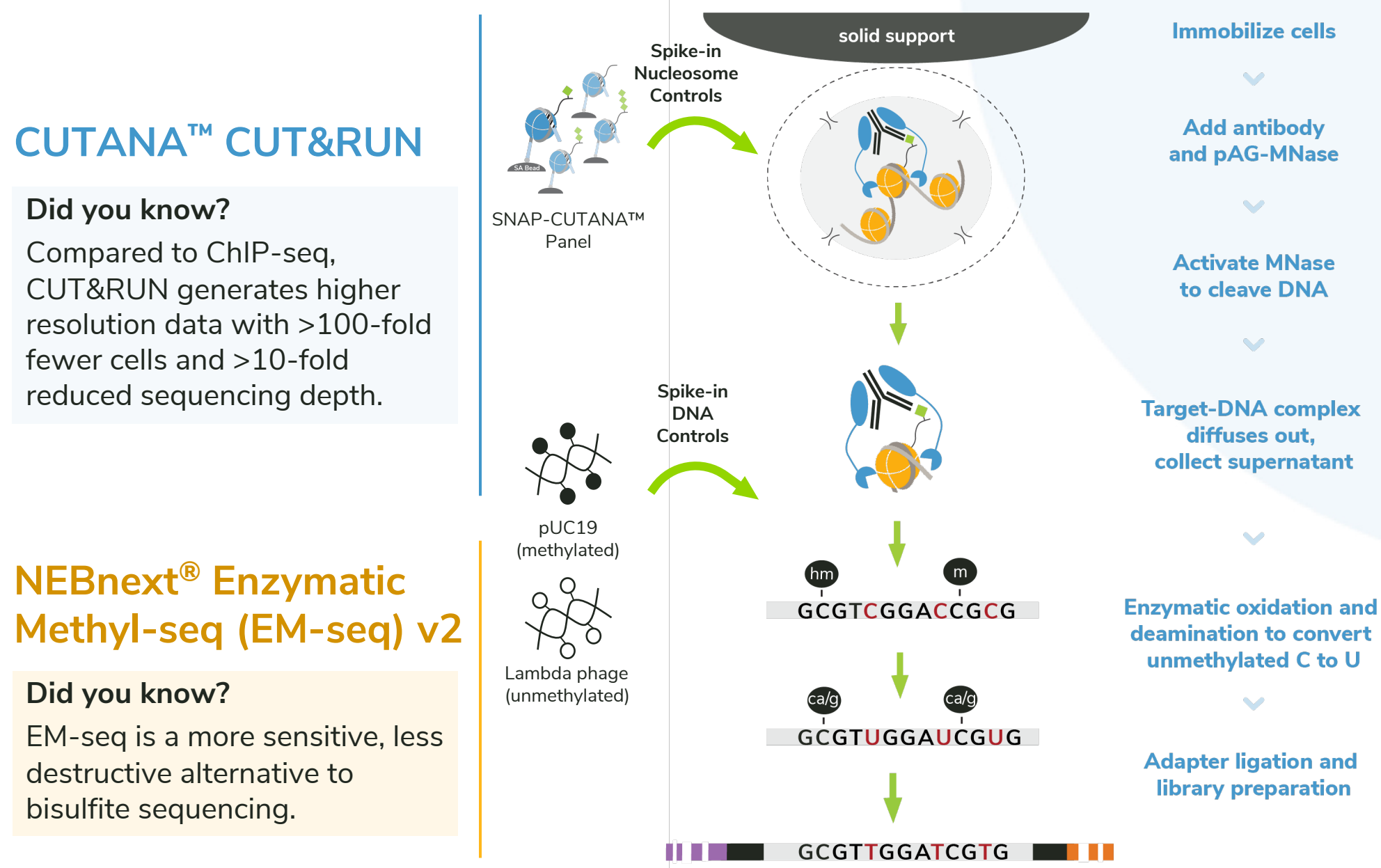


Figure 2. CUT&RUN-EM integrates EpiCypher CUT&RUN with NEB EM-seq v2. CUT&RUN isolates DNA associated with chromatin proteins of interest. EM-seq is then used to enzymatically convert unmethylated cytosines to uracils. Sequencing the resulting libraries enables CpG resolution of DNAm co-occurring with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

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

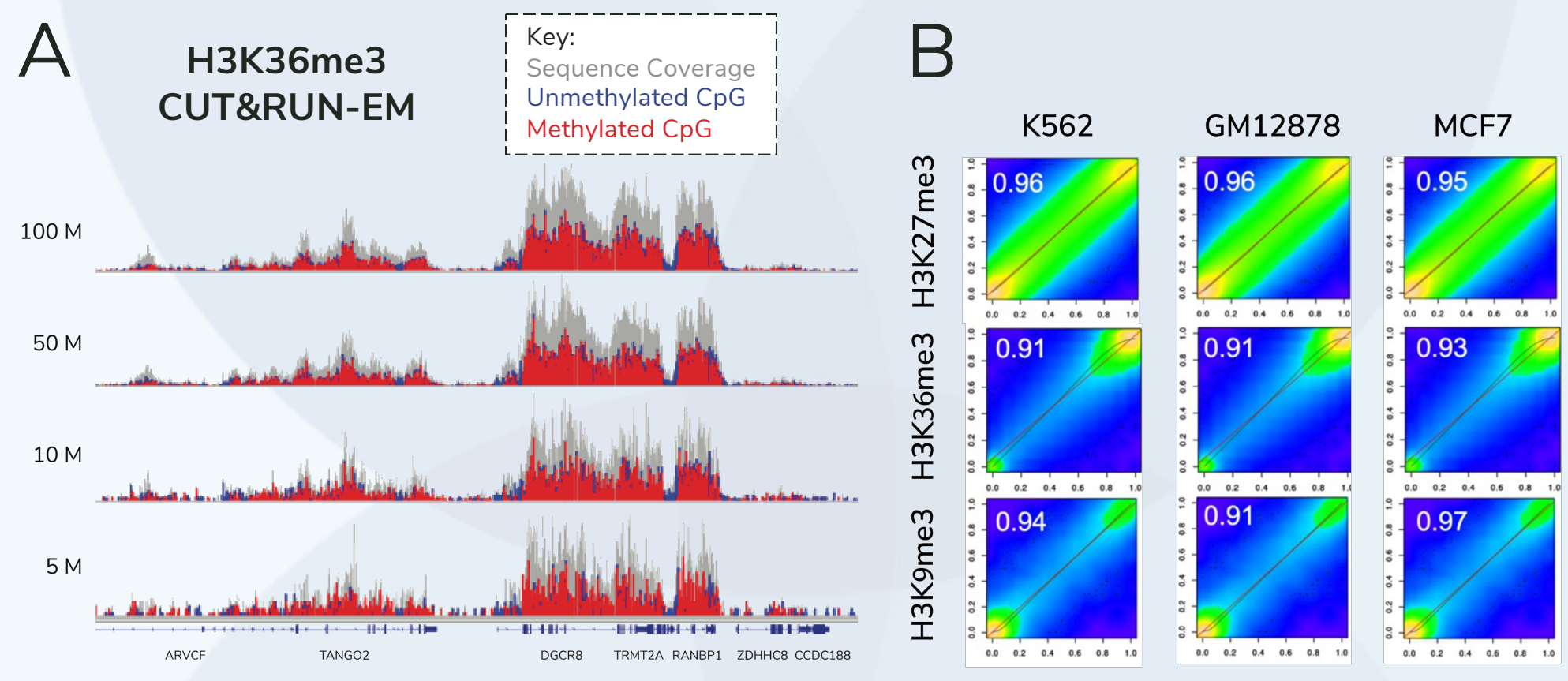
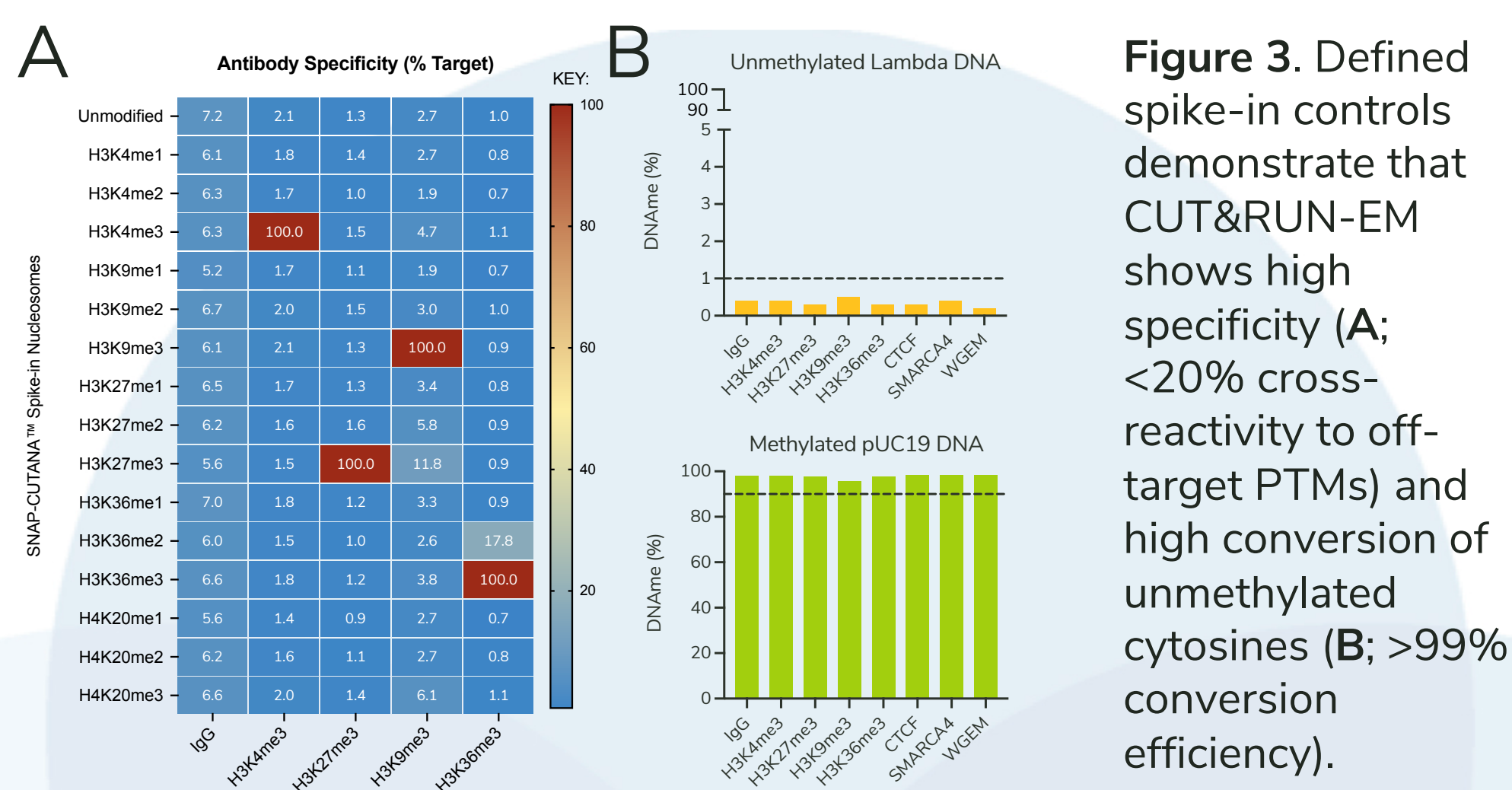


Figure 4. CUT&RUN-EM assays are highly reproducible across various sequencing depths (A; downscaled genome browser tracks) and replicates (B; Pearson correlation coefficients).

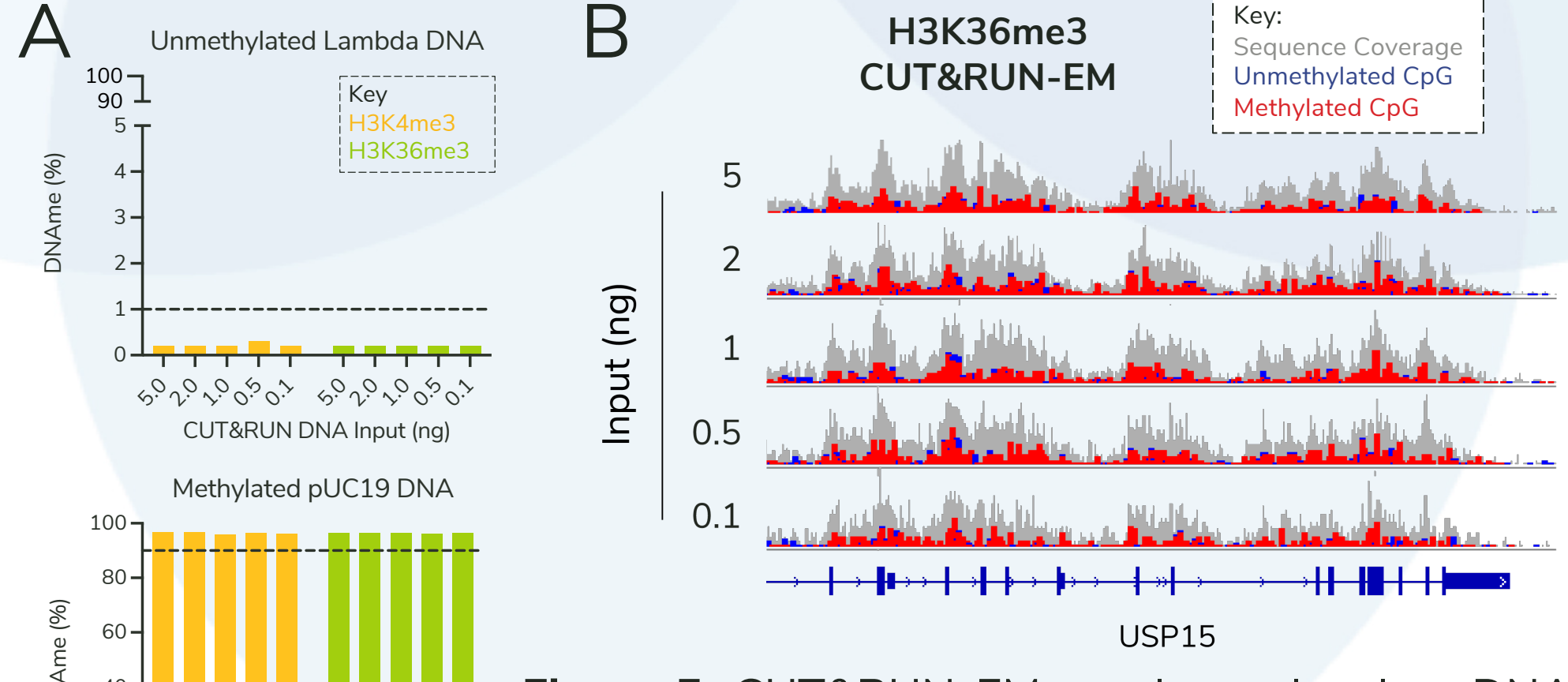


Figure 5. CUT&RUN-EM excels at ultra-low DNA inputs; demonstrating consistent conversion efficiency (A) and genomic enrichment (B) down to 0.1 ng input using the EM-seq v2 kit.

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

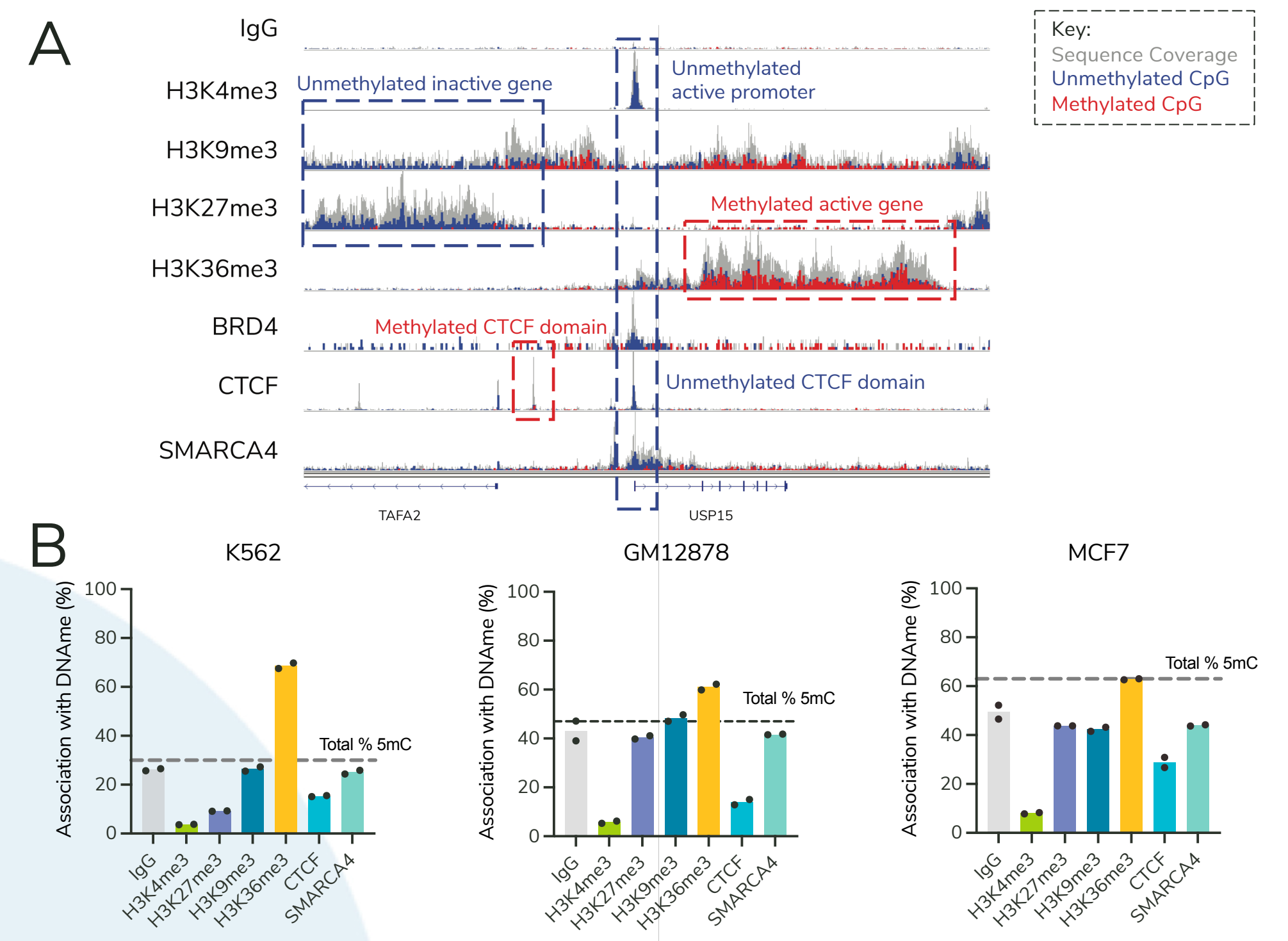


Figure 6. Representative genome browser tracks (A) demonstrate that CUT&RUN-EM reflects known biological functions of chromatin proteins (e.g. unmethylated H3K4me3 promoters, methylated H3K36me3 gene bodies) when overlaid with DNAm. The association of chromatin proteins with DNAm varies across cell lines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

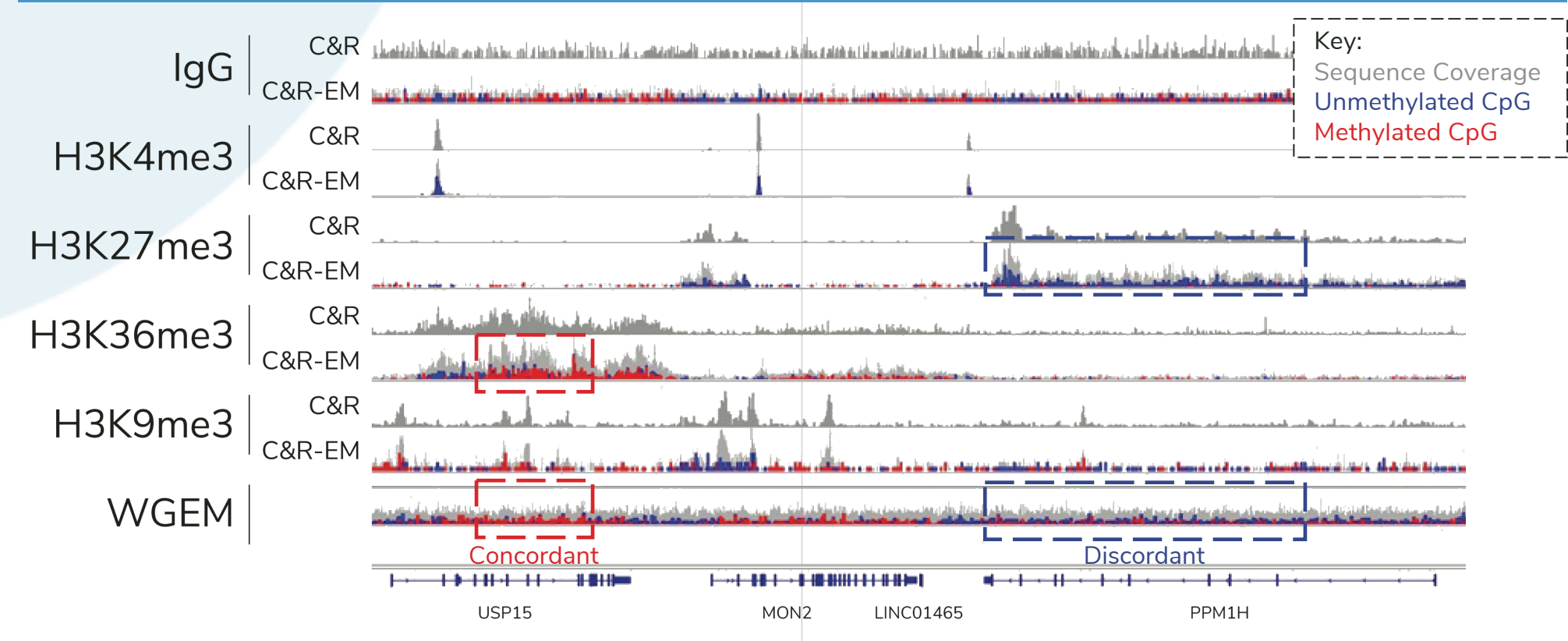


Figure 7. CUT&RUN-EM stratifies subpopulations of cells with unique PTM-DNAm crosstalk signatures that are obscured in whole genome EM-seq (WGEM). IGV genome browser comparison of CUT&RUN (C&R), CUT&RUN-EM (C&R-EM), and WGEM in K562 cells. C&R-EM loci with concordant results (red box) recapitulate the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

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

Modified CUT&RUN excises methylDNA to reduce seq depth

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

MeCP2 captures high-density 5mC with 34x less sequencing than WGEM

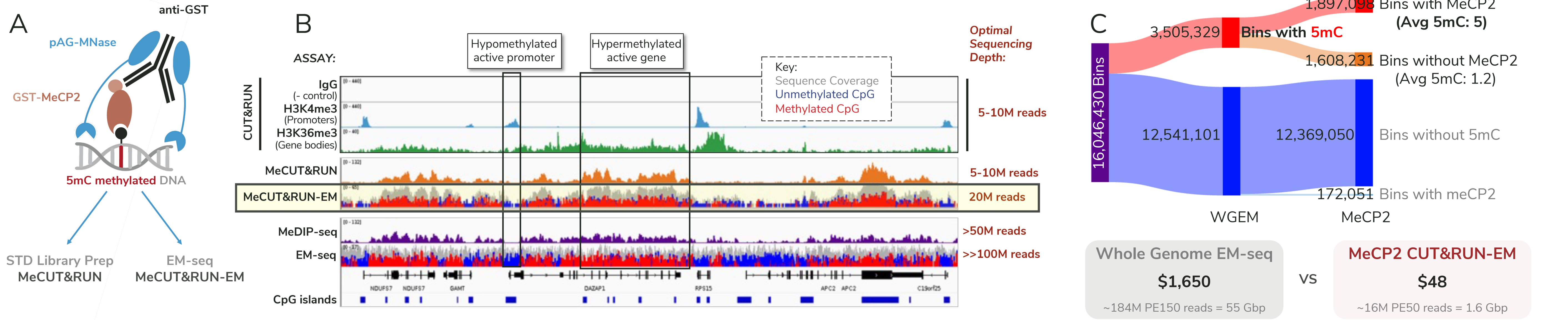


Figure 8. (A) Schematic representation of a GST-tagged MeCP2 methyl binding domain used in lieu of an antibody in traditional CUT&RUN (MeCUT&RUN) to enrich methylated DNA. Sequencing libraries can be prepared via standard library prep to provide a snapshot of genome-wide DNAm at 150 bp resolution, or unmethylated cytosines can be converted with EM-seq to achieve CpG resolution. (B) MeCUT&RUN (orange) with 500k K562 cells shows high concordance with Methyl-DNA Immunoprecipitation sequencing (MeDIP-seq; purple) at >2.5-fold reduced sequencing depth with 10-fold fewer cells. MeCP2 CUT&RUN-EM (yellow boxed track) generates similar DNA methylation profiles compared to whole genome EM-seq (bottom track). (C) Sankey plot of the binned human genome (200 bp) comparing the presence of 5mC in WGEM and MeCUT&RUN-EM. MeCP2 enriches for regions of high 5mC concentration with 34x less sequencing than WGEM, greatly reducing sequencing costs.

Conclusions

- CUT&RUN-EM enables a direct interrogation of DNAm and chromatin protein crosstalk
- MeCUT&RUN provides fast and efficient mapping of global DNA methylation levels with ~150 bp resolution
- MeCUT&RUN-EM resolves base pair resolution of DNA methylation with 34x less sequencing than WGEM

See EpiCypher assays in action

- AML therapy selection San José-Enériz, Nat Commun 2024 (PMID: [38956053](#))
- Epigenomics of stress Torres-Berrio, Neuron 2024 (PMID: [38959894](#))
- SCA1 disease mechanism Coffin, Neuron 2023 (PMID: [36577402](#))
- Aging & chromatin opening Patrick, Cell Metab 2024 (PMID: [38959897](#))
- iPSC differentiation by BRD4 Padmanabhan, Nat Cardio Res 2024 (PMID: [39196112](#))
- H3.3K36M disrupts DNAm Sinha, Mol Cell 2024 (PMID: [39368466](#))

Download Poster

