

# Direct multi-omics for the masses: Linking DNA methylation to chromatin targets via TEM-seq

Keith E. Maier<sup>1</sup>, Bryan J. Venters<sup>1</sup>, Vishnu U. Sunitha Kumary<sup>1</sup>, Allison Hickman<sup>1</sup>, Anup Vaidya<sup>1</sup>, Ryan Ezell<sup>1</sup>, Jonathan M. Burg<sup>1</sup>, Louise Williams<sup>2</sup>, Chaitanya Ponnaluri<sup>2</sup>, Hang Geong Chin<sup>2</sup>, Pierre Esteve<sup>2</sup>, Isaac Meeke<sup>2</sup>, Sriharsa Pradhan<sup>2</sup>, Zu-Wen Sun<sup>1</sup>, Martis W. Cowles<sup>1</sup>, & Michael-Christopher Keogh<sup>1</sup>

<sup>1</sup>EpiCypher Inc., Research Triangle Park, Durham NC 27709, USA, <sup>2</sup>New England Biolabs, Ipswich, MA 01938

## Molecular crosstalk between DNA methylation and other chromatin features

DNA methylation (DNAm) is an epigenetic mark that includes the modification of cytosine residues (5mC) within CpG islands. In addition to well characterized roles regulating gene expression, imprinting and silencing parasitic DNA elements, the misregulation of DNAm is implicated in multiple diseases. Evidence is emerging that DNAm is not an independent epigenetic mark but rather closely linked to the post-translational modification (PTM) of histone proteins. However, examining the direct relationships between 5mC and PTMs are hampered by correlated analyses of separate assays that cannot establish a direct mechanistic linkage. Furthermore, the traditional approach to measure 5mC relies upon harsh bisulfite chemical conversion of DNA, which introduces DNA breaks and systemic biases.

To address these limitations, we developed a Targeted Enzymatic Methylation-sequencing (TEM-seq) approach, an ultra-sensitive multi-omic genomic mapping technology that delivers high resolution DNAm profiles at epitope-defined chromatin features. Importantly this assay examines the direct molecular link between 5mC and histone PTMs or chromatin associated proteins (ChAPs).

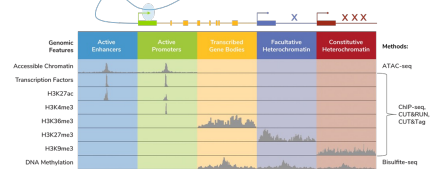


Figure 1: Gene expression is regulated by chromatin structure and accessibility. PTMs provide useful markers of chromatin elements, such as promoters, enhancers, and gene bodies. Adapted from (1).

## TEM-seq is a powerful multi-omic assay that directly links 5mC to PTMs and/or ChAPs

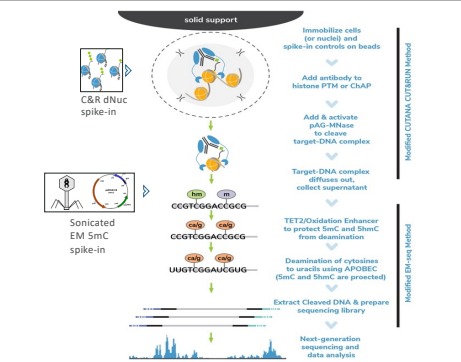


Figure 2: Workflow for TEM-seq (Targeted Enzymatic Methylation-sequencing).

## TEM-seq reveals distinct DNAm levels linked to different histone PTMs and ChAPs

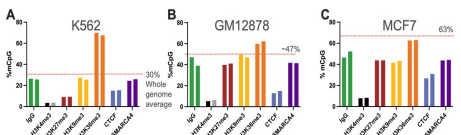


Figure 3: Percent methylation analysis from TEM-seq data sets are consistent with known biological functions of antibody targets. TEM-seq was performed using a set of antibodies to different histone PTMs or chromatin targets. For each data set, the percent of methylated CpGs was calculated to provide a global view of relative methylation directly linked to each antibody target.

## TEM-seq allows sensitive profiling to relate DNAm to Histone PTMs and ChAPs

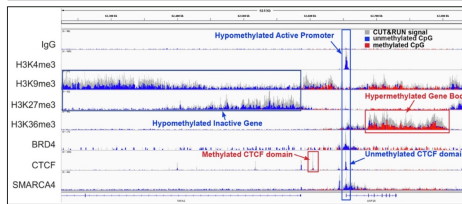


Figure 4: TEM-seq genome browser tracks using a range of histone PTM and ChAP (Chromatin Associated Protein) targets. Composite histone PTM or ChAP signal (i.e. CUT&RUN data) is in grey. Blue = unmethylated CpG; Red = methylated CpG. TEM-seq generates high quality genomic maps for both histone PTMs and ChAPs and also reveals differential DNAm at various genomic regions, (e.g. hypermethylation at gene bodies / hypomethylation at promoters).

## TEM-seq: Highly reproducible and sensitive at low seq depths

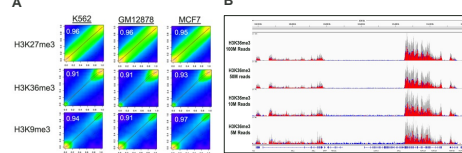
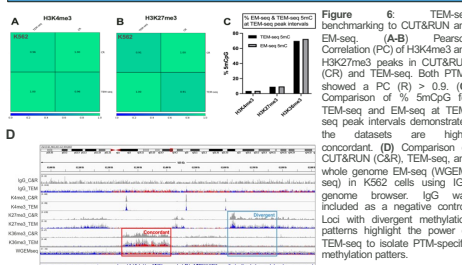
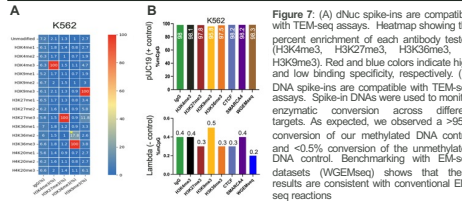


Figure 5: (A) TEM-seq assays are highly reproducible. Biological replicate analysis in K562, GM12878, and MCF7 cells shows that TEM-seq assays are highly reproducible ( $r = 0.91$ ). (B) TEM-seq genome browser tracks for H3K36me3 were randomly downsampled / analyzed at represented sequence depths and similar DNAm distribution observed using as few as 5M reads.

## Benchmarking TEM-seq to CUT&RUN and EM-seq

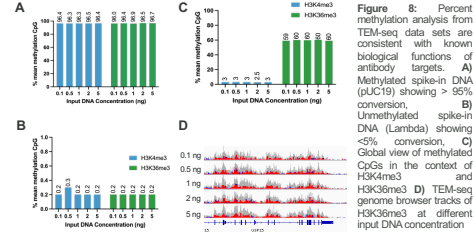


## TEM-seq spike-ins provide quantitative metrics for rigorous assay monitoring



## TEM-seq spike-ins provide quantitative metrics for rigorous assay monitoring

## TEM-seq excels at sub-nanomgram amounts of input DNA



## In vitro characterization of Rett syndrome MeCP2 mutants with Luminescence Assay

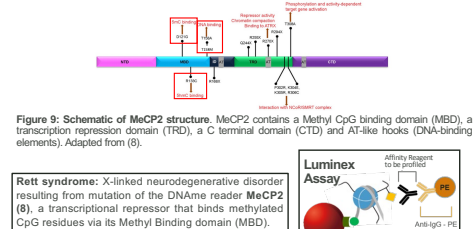
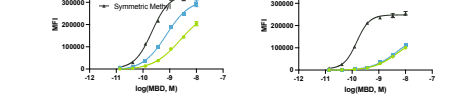


Figure 9: Schematic of MeCP2 structure. MeCP2 contains a Methyl CpG binding domain (MBD), a transcription repression domain (TRD), a C terminal domain (CTD) and Atf-like hooks (DNA-binding elements). Adapted from (8).

## MeCP2 titration on Luminescence Assay



## Specificity of MeCP2 (WT & Rett syndrome alleles) for symmetric 5mC (salt dependence)

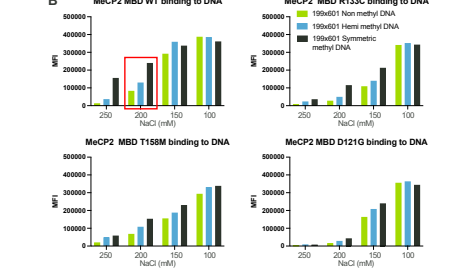
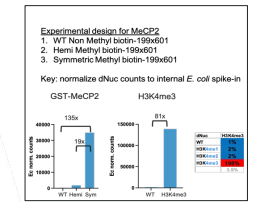


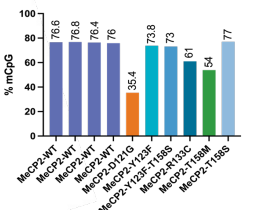
Figure 10: Luminescence characterization of MeCP2 and Rett syndrome mutations. (A) GST-MeCP2 MBD WT shows a strong preference for symmetrically methylated 5mC, most particularly in the nucleosome context (right panel) compared to free DNA (left panel). (B) Impact of Rett syndrome mutations on binding of MeCP2 MBD to free DNA under decreasing NaCl concentrations. The red box indicates conditions chosen for reader TEM-seq studies with GST-MeCP2 MBD WT (see Figure 11).

# TEM-seq provides multi-omic insights into Rett syndrome

## (A) EpiCypher dNucs can be used as spike-in controls to monitor specific enrichment in reader TEM-seq



## (B) TEM-seq reveals differences in DNAm levels linked to Rett MeCP2 mutants



## (C) Rett MeCP2 mutants display defective genome distribution.

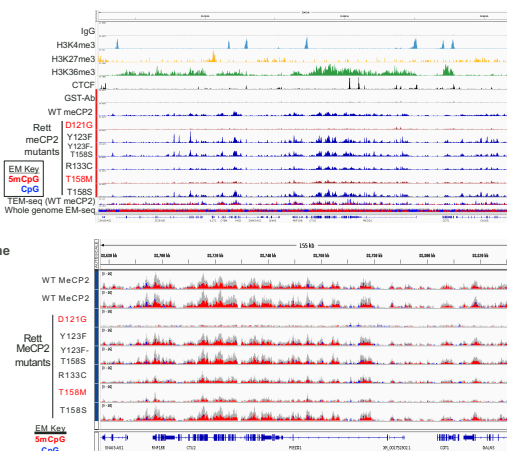


Figure 11. Reader TEM-seq using GST-MeCP2. (A) MeCP2 MBD WT was used in a reader TEM-seq experiment with non-methylated, hemi-methylated and symmetrically methylated DNA (pUC19) as spike-in controls. MeCP2 showed preferential enrichment of the symmetric-methylated spike-in compared to non-methylated (135-fold) and hemi-methylated (19-fold) DNAs. Anti-H3K4me3 and SNA-CUT&RUN, K-MeSati mini-panel (Figure 2) positive control showed 81-fold enrichment of H3K4me3 on unmethylated DNA. (B) Percent methylation analysis reveals the effect of Rett mutants in the MeCP2 binding (C) Genome browser tracks (CUT&RUN and TEM-seq) for GST-MeCP2 MBD using K562 cells show its localization to methylated CpG (red) over non-methylated CpG (blue).

## Conclusions

- TEM-seq is a novel multi-omic approach that directly links DNA methylation to chromatin features.
- TEM-seq delivers high quality data at low sequence depths.
- Spike-in controls are important to monitor TEM-seq performance.
- dCypher 3.0 optimizes conditions for reader TEM-seq experiments.
- TEM-seq can be leveraged to gain mechanistic insights into clinically relevant chromatin reader domains (e.g., MBD of MeCP2).

## References

- Jones (2012) Functions of DNA methylation: islands, start sites, gene bodies and beyond. *Nature Rev. Genet.* 13:484.
- Cedar & Bergman (2009) Linking DNA methylation and histone modification: patterns and paradigms. *Nature Rev. Genet.* 10:255.
- Shah et al (2018) Examining the Roles of H3K4 Methylation States with Systematically Characterized Antibodies. *Mol. Cell.* 72:162.
- Xu & Krutz-Miller (2013) A novel DNA-binding feature of MeCP2 contributes to Rett syndrome. *Front Cell Neurosci.* 7:84.
- Manunze et al (2022) The dCypher Approach to Interrogate Chromatin Reader Activity Against Posttranslational Modification-Defined Histone Peptides and Nucleosomes. *Methods Mol Biol.* 2488:231.

Acknowledgment: NIH-NHGR for financial support

