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



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

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

Molecular crosstalk between DNA methylation and other chromatin features

DNA methylation (DNAme) is an epigenetic mark that includes the modification of cytosine resides (5mC) within CpG islands. In addition to well characterized roles regulating gene expression, imprinting and silencing parasitic DNA elements, the misregulation of DNAme is implicated in multiple diseases. Evidence is emerging that DNAme 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 conversation of DNA, which introduces DNA breaks and systemic biases.

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



TEM-seq allows sensitive profiling to relate DNAme 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 gray. Blue = unmethylated CpG; Red = methylated CpG. TEM-seq generates high quality genomic maps for both histone PTMs and ChAPs and also reveals differential DNAme at various genomic regions, (e.g. hypermethylation at gene bodies / hypomethylation at promoters).

TEM-seq excels at sub-nanogram amounts of input DNA



In vitro characterization of Rett syndrome **MeCP2** mutants with Luminex assay

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



High reproducibility and sensitivity of TEM-seq 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 DNAme distribution observed using as few as 5M reads

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





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 AT-like hooks (DNA-binding elements). Adapted from (8).

Box: Rett syndrome X-linked disorder neurodegenerative resulting from mutation of the DNAme reader MeCP2 (8), a transcriptional repressor that binds methylated CpG residues via its Methyl Binding domain (MBD).

Α

where DNA

labels a

DNA

EM-sea

box

and

locus

400000 ¬

300000

100000

-12

₩ 200000-

MeCP2 titration on Luminex

MeCP2 MBD WT binding to DNA

Antibody to Luminex be profiled Assay Anti-IgG - P xMAP Bead - dNuc

MeCP2 MBD WT binding to nucleosomes



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



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

TEM-seq reveals distinct DNAme 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 PTM 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 spike-ins provide quantitative metrics for rigorous assay monitoring



Figure 7: (A) dNuc spike-ins are compatible with TEM-seq assays. Heatmap showing the percent enrichment of each antibody tested (H3K4me3, H3K27me3, H3K36me3, or H3K9me3). Red and blue colors indicate high and low binding specificity, respectively. (B) DNA spike-ins are compatible with TEM-seq assays. Spike-in DNAs were used to monitor enzymatic conversion across different targets. As expected, we observed a >95% conversion of our methylated DNA control and <0.5% conversion of the unmethylated DNA control. EM-seq datasets Benchmarking with WGEMseq) shows that these results are consistent with conventional EM-seg reactions

Figure 10. Luminex 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 in a buffer containing 200mM NaCl (right panel). (B) Impact of Rett syndrome mutations on binding of MeCP2 MBD to free DNA (left) under different NaCl concentrations. Arrow indicates conditions chosen for reader TEM-seg 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



(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-methyl, hemi-methyl and symmetric-methyl Nucs [22-N-22; [0x (un) - 3x (hemi) - 6x (symmetric) 5meC in each linker region] as spike in controls. MeCP2 showed preferential enrichment of the symmetric-methyl spike-in compared to non-methyl (135-fold) and hemi-methyl (19-fold) Nucs. Anti-H3K4me3 and SNAP-CUTANA K-MetStat mini-panel (Figure 2) positive control showed 81-fold enrichment of H3K4me3 over unmodified Nuc. (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 methyl-CpG (red) over non methyl CpG (blue).

Conclusions

> TEM-seq is a novel multi-omic approach that directly links DNA methylation to chromatin features

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



- \succ TEM-seq delivers high quality data at low sequence depths.
- > Spike-in controls are important to monitor TEM-seq performance.
- Cypher 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., MDB of MeCP2)

References

1. Jones (2012) Functions of DNA methylation: islands, start sites, gene bodies and beyond. Nature Rev. Genet. 13:484.

2. Cedar & Bergman (2009) Linking DNA methylation and histone modification: patterns and paradigms. Nature Rev. Genet. 10:295.

3. Shah *et al* (2018) Examining the Roles of H3K4 Methylation States with Systematically Characterized Antibodies. Mol. Cell. 72:162.

4. Xu & Pozzo-Miller (2013) A novel DNA-binding feature of MeCP2 contributes to Rett syndrome. Front Cell Neurosci. 7:64.

5. Marunde et al (2022) The dCypher Approach to Interrogate Chromatin Reader Activity Against Posttranslational Modification-Defined Histone Peptides and Nucleosomes. *Methods* Mol. Biol. 2458:231.

Acknowledgement: NIH-NHGRI for financial support

