

Keith E. Maier<sup>1</sup>, Andrea L. Johnstone<sup>1</sup>, Matthew R. Marunde<sup>1</sup>, Irina K. Popova<sup>1</sup>, Jonathan M. Burg<sup>1</sup>, Nathan Hall<sup>1</sup>, Katherine Novitzky<sup>1</sup>, Eileen McAnarney<sup>1</sup>, Harrison A. Fuchs<sup>2,3</sup>, Matthew J. Meiners<sup>1</sup>, Zachary Gillespie<sup>1</sup>, Marcus A. Cheek<sup>1</sup>, Sarah A. Howard<sup>1</sup>, Catherine Musselman<sup>2,3</sup>, Zu-Wen Sun<sup>1</sup>, Bryan J. Venters<sup>1</sup> & Michael-Christopher Keogh<sup>1</sup>

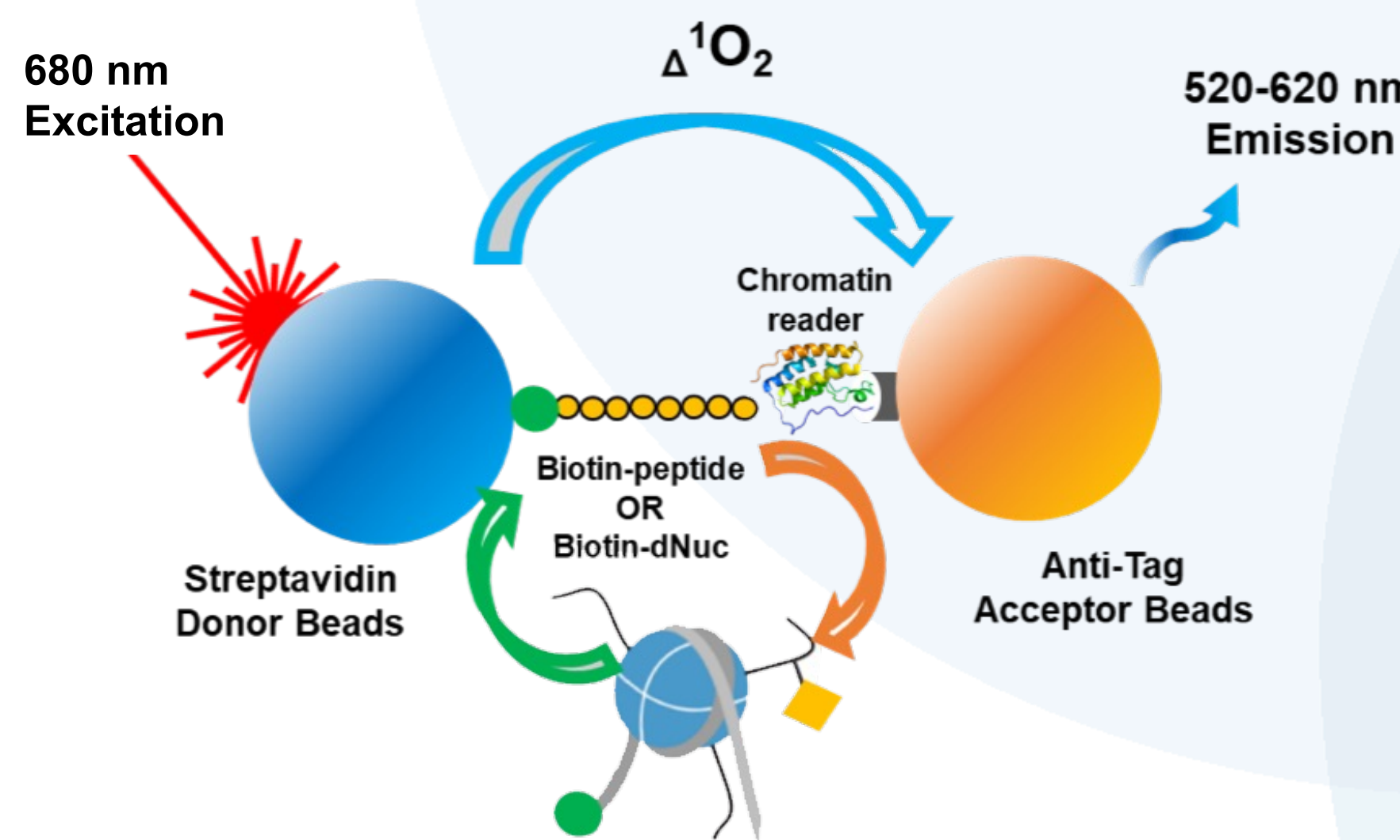
<sup>1</sup>EpiCypher, Inc., Research Triangle Park, NC 27713, USA. <sup>2</sup>University of Iowa, Iowa City, IA 52246. <sup>3</sup>University of Colorado, Aurora CO 80045

## BACKGROUND

Epigenetic regulators recognize histone post-translational modifications (PTMs) through evolutionarily conserved binding domains (aka. readers), thereby recruiting nuclear complexes to specific genomic loci. Modified histone peptides are frequently used to model chromatin and enable characterization of [reader - histone PTM] specificity, and in doing so, decipher the histone code; however, peptides are structurally very different from chromatin and their use assumes PTM specificity is unaffected by higher order factors. To assess the limitations and shortcomings of this reductive approach, EpiCypher® has developed the dCypher assay, a high-throughput discovery platform for the rapid screening & detailed interrogation of chromatin interactors (readers, enzymes, and antibodies) against comprehensive libraries of modified histone peptides and designer nucleosomes. Here, we used dCypher assays to show PTMs presented on peptides engage individually with a dual PHD-BD domain of BPTF (subunit of the Nucleosome Remodeling Factor [NURF] complex); however, when the same modifications are presented on nucleosomes, combinatorial engagement drives a highly specific interaction. Further, genomic mapping with CUT&RUN recapitulates the specificity of the nucleosome-reader interaction, where reader-mediated enrichment is only seen when both methyl and acetyl marks are present. These findings demonstrate the critical importance of using nucleosome substrates to garner accurate insights into *in vivo* binding mechanisms.

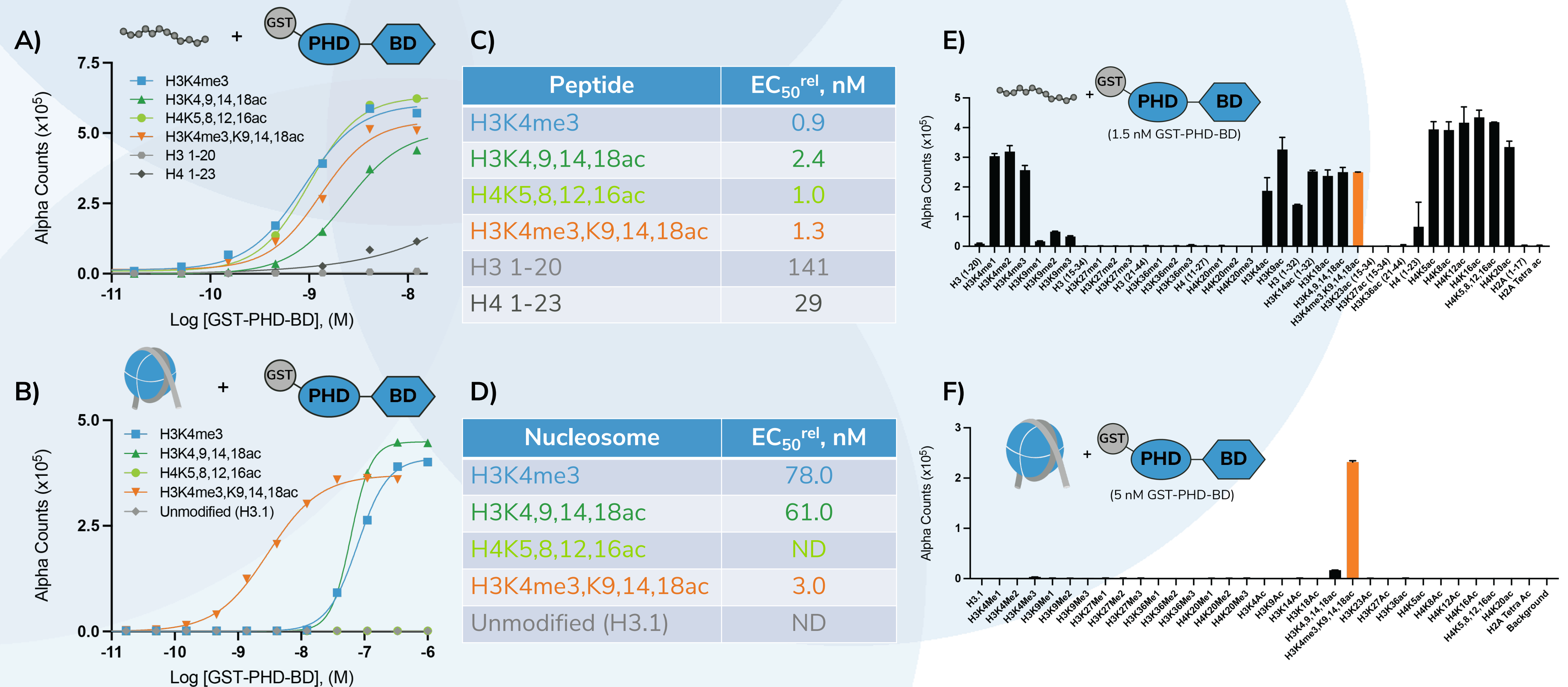
### dCypher Binding Assay

- Robust signal / background
  - Identifies hits missed by peptide arrays
- Highly sensitive no-wash format
  - Lower protein input requirements vs peptide arrays
- Accommodates full length & multidomain proteins
  - Assess cooperative binding
- Access comprehensive libraries
  - >287 modified histone peptides
  - >100 recombinant nucleosomes
- Unprecedented epigenetic diversity
  - PTMs: Kme, Rme, Kac, S/T/Y-phos
  - Other: DNAm, histone variants, onco-mutations
- Custom services by expert scientists



**Figure 1.** dCypher combines the benefits of peptides, nucleosomes and AlphaScreen® (Perkin Elmer) in one powerful assay. Above, a graphical depiction of dCypher: a no-wash liquid phase assay compatible with any biotinylated histone peptide or nucleosome. The approach has been optimized for use with GST-, 6HIS-, or FLAG-tagged query proteins.

### BPTF PHD-BD Dual Domain Binds Cooperatively to Nucleosomes

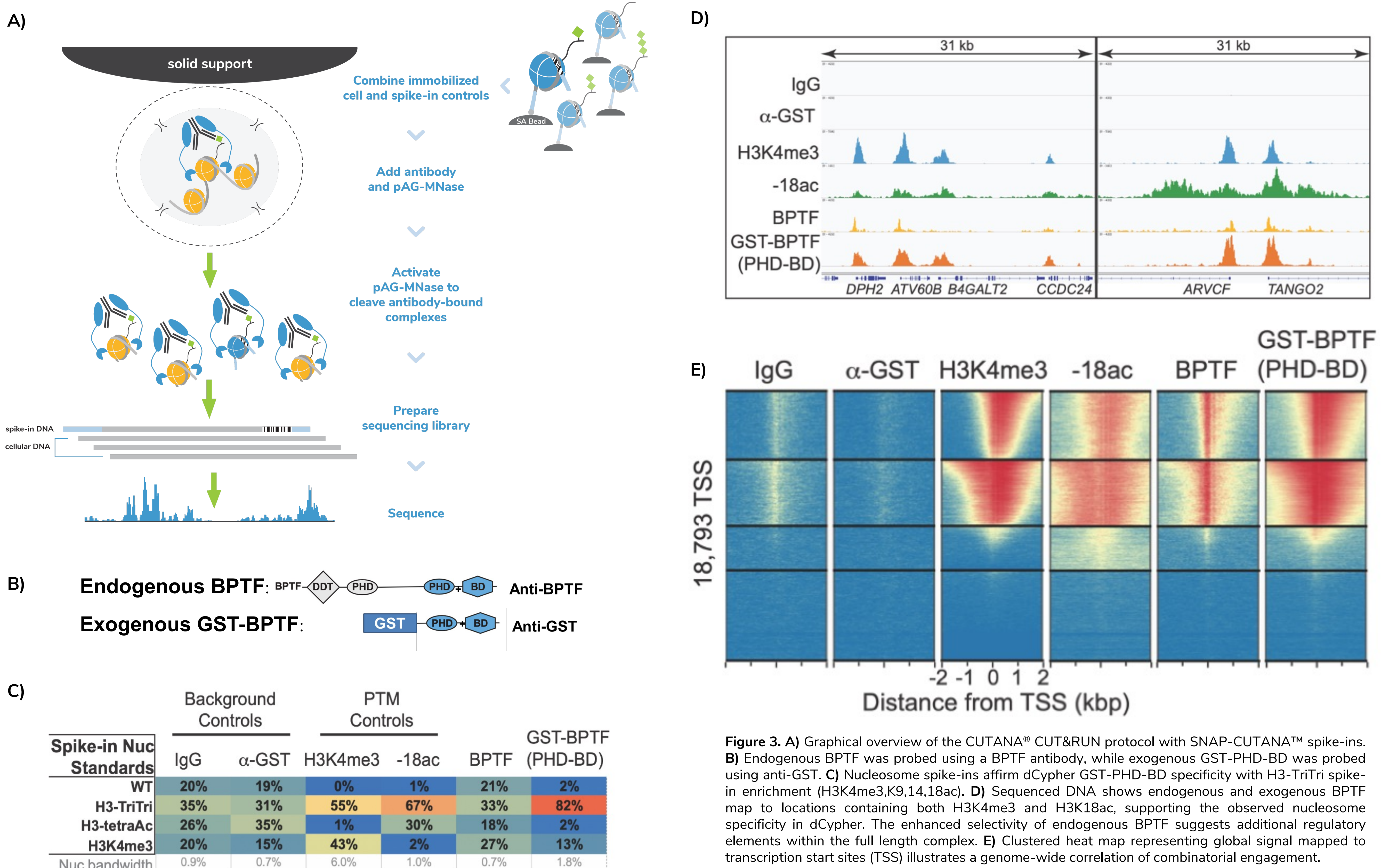


**Figure 2.** BPTF PHD-BD binds H3K4me3,K9,14,18ac cooperatively on nucleosomes. **A)** Titration of GST-BPTF PHD-BD against H3 and H4 histone peptides. Robust binding observed against methyl and acetyl peptides, consistent with expected domain function, though no cooperative binding was observed. **B)** Titration of GST-BPTF PHD-BD against H3 and H4 modified nucleosomes. Strong combinatorial engagement is observed on the H3K4me3,K9,14,18ac substrate (orange). **C,D)** Relative EC<sub>50</sub> values from A & B determine optimal query protein concentration for E & F. **E)** Representative data from a 287-member dCypher discovery screen shows broad binding to H3K4methyl and H3/H4 acetyl peptides. **F)** Representative data from a 65-member nucleosome screen shows highly selective engagement with H3K4me3,K9,14,18ac.

## CONCLUSIONS

- dCypher™ has major benefits compared to histone peptide arrays
- Nucleosome context is critical to decipher the histone code
- BPTF PHD-BD cooperatively engages me & ac PTMs on nucleosomes but not peptides
- QUESTION: Do nucleosomes model *in vivo* specificity?

## Reader-CUT&RUN Confirms PHD-BD Combinatorial Engagement *In Vivo*



**Figure 3.** **A)** Graphical overview of the CUTANA® CUT&RUN protocol with SNAP-CUTANATM spike-ins. **B)** Endogenous BPTF was probed using a BPTF antibody, while exogenous GST-PHD-BD was probed using anti-GST. **C)** Nucleosome spike-ins affirm dCypher GST-PHD-BD specificity with H3-TriTri spike-in enrichment (H3K4me3,K9,14,18ac). **D)** Sequenced DNA shows endogenous and exogenous BPTF map to locations containing both H3K4me3 and H3K18ac, supporting the observed nucleosome specificity in dCypher. The enhanced selectivity of endogenous BPTF suggests additional regulatory elements within the full length complex. **E)** Clustered heat map representing global signal mapped to transcription start sites (TSS) illustrates a genome-wide correlation of combinatorial engagement.



Download PDF