

Captify™ Screening of BPTF Reader Domains Show Nucleosomes, Not Peptides, Dictate the Histone Code

EpiCypher®

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 Captify 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 Captify 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.

