

Exploring Antibodies and Rationally Designed Readers for Genomic Mapping of Histone Monoubiquitin



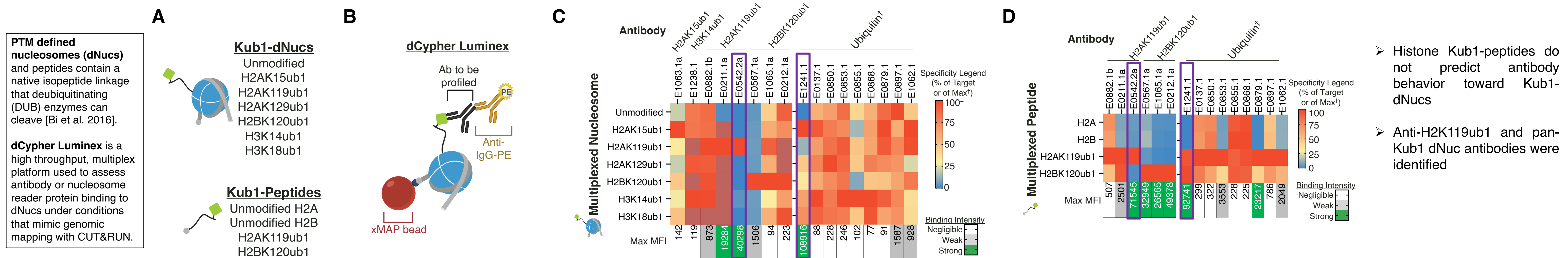
Susan L. Gloor¹, Tessa M. Firestone¹, Allison R. Hickman¹, Catherine E. Smith¹, Matthew R. Marunde¹, Nathan Hall¹, Laiba Khan¹, Matthew J. Meiners¹, Zachary B. Gillespie¹, Anup Vaidya¹, Sabrina R. Hunt¹, Ryan J. Ezell¹, Jonathan M. Burg¹, Hailey F. Taylor¹, Ugochi C. Onuoha¹, Marcus A. Cheek¹, Emily F. Patteson¹, Hannah E. Willis¹, Danielle N. Maryanski¹, Chitra Mohan³, Carolina dos Santos Passos¹, Jessica K. Tyler³, Robert E. Cohen², Tingting Yao², Martis Cowles¹, Zu-Wen Sun¹ & Michael-Christopher Keogh¹

¹ EpiCypher, Inc., Research Triangle Park, NC 27709, USA. ² Colorado State University, Fort Collins, CO 80523, USA. ³ Weill Cornell Medicine, New York, NY 10065, USA.

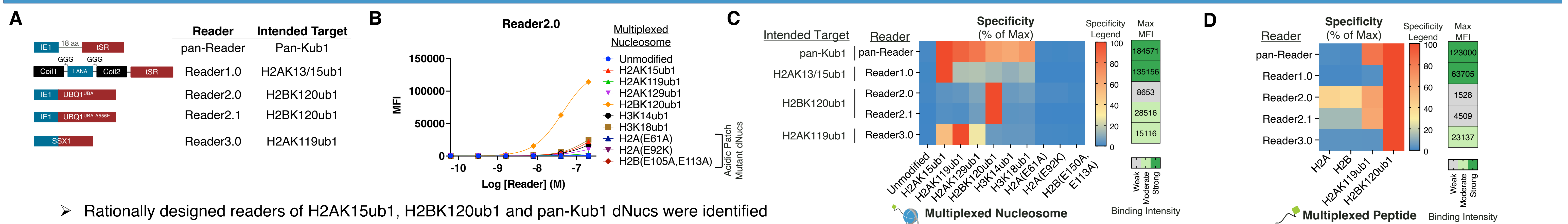
Background

Chromatin function is regulated by reversible histone post-translational modifications (PTMs) including lysine monoubiquitination (Kub1). Multiple Kub1 sites are involved in gene silencing (e.g., H2AK119) or activation (e.g., H2BK120), and marking DNA double-strand breaks (e.g., H2AK13/15). Misregulation of histone Kub1 has been linked to the pathogenesis of diverse diseases. However, the PTM class is a challenging study target due to a lack of physiologically relevant substrates and reliable, well-characterized detection reagents. Here we describe the characterization of diverse tools to overcome these issues and enable the accurate genomic mapping of histone Kub1. We set out to identify truly-capable genomic mapping reagents from two classes: 1) ubiquitin and ubiquitinated-histone antibodies; and 2) recombinant Kub1-nucleosome readers from a rational structure-guided approach. We used a high throughput, multiplexed platform (dCypher™ Luminex) to assess antibody / nucleosome reader binding to Kub1-peptides or PTM defined nucleosomes (Kub1-dNucs™) under conditions that mimic genomic mapping with CUT&RUN (Cleavage Under Targets and Release Using Nuclease). Most antibodies failed to bind Kub1-dNucs or lacked specificity for their purported target; exceptions being an **anti-H2AK119ub1** and a **pan-Kub1**. However, we also identified rationally designed Kub1-dNuc readers of **H2AK15ub1**, **H2BK120ub1** and **pan-Kub1**. Unsurprisingly, **histone Kub1-peptides did not predict antibody or Kub1-nucleosome reader behavior towards Kub1-dNuc substrates**. We next explored using antibodies / nucleosome readers to study the genomic enrichment of Kub1 PTMs by CUT&RUN, including a spike-in panel of fully defined dNuc standards to provide an *in situ* metric of reagent capability and assay performance. We showed that reagent behavior with dCypher Luminex was recapitulated by CUT&RUN and are able to effectively map Kub1 associated with gene repression (H2AK119ub1), gene activation (H2BK120ub1) and DNA DSB repair (H2AK15ub1, pan-Ub1). Such findings demonstrate the power of PTM-defined physiological substrates to characterize histone-PTM antibodies / readers and perform truly insightful genomic studies.

Survey of Ub and ubiquitinated-histone antibodies identifies Kub1-nucleosome specific antibodies



Novel Kub1-nucleosome readers engage the nucleosome acidic patch and monoubiquitin



➤ Rationally designed readers of H2AK15ub1, H2BK120ub1 and pan-Kub1 dNucs were identified

Figure 2. Binding strength and specificities of rationally designed Kub1-nucleosome readers were determined using dCypher Luminex. A) Schematic representation of Kub1-nucleosome readers and their intended targets. Acidic patch Anchor domains: blue, Linkers: black, ubiquitin binding domain (UBD): red. B) The binding strength and specificity of all readers was tested by dCypher Luminex (Figure 1B) including three additional acidic patch mutant dNucs. Characterization of Reader2.0 shown under conditions that mimic genomic mapping with CUT&RUN. Binding to the unmodified nucleosome is disrupted by acid-patch mutations (as would be expected given the mode of engagement). C) The binding strength and specificity of each Kub1-nucleosome reader compared under optimized conditions as determined in (B): pan-Reader, 200 nM; Reader1.0, 8 nM; Reader2.0 / 2.1, 40 nM; Reader3.0, 200 nM. MFI were normalized to max MFI to determine percent specificity. D) The binding strength and specificity of each Kub1-nucleosome reader toward Kub1-histone peptides was similarly tested as in (C): (pan-Reader; 40 nM, Reader1.0; 8 nM, Reader2.0 / 2.1 / 3.0; 200 nM). Excluding the pan-Reader, Kub1-peptide binding profiles lack a clear link to their target nucleosome, underscoring the importance of both the acidic patch anchor and the UBD for Kub1-dNuc readers' binding to the Kub1 nucleosomal target.

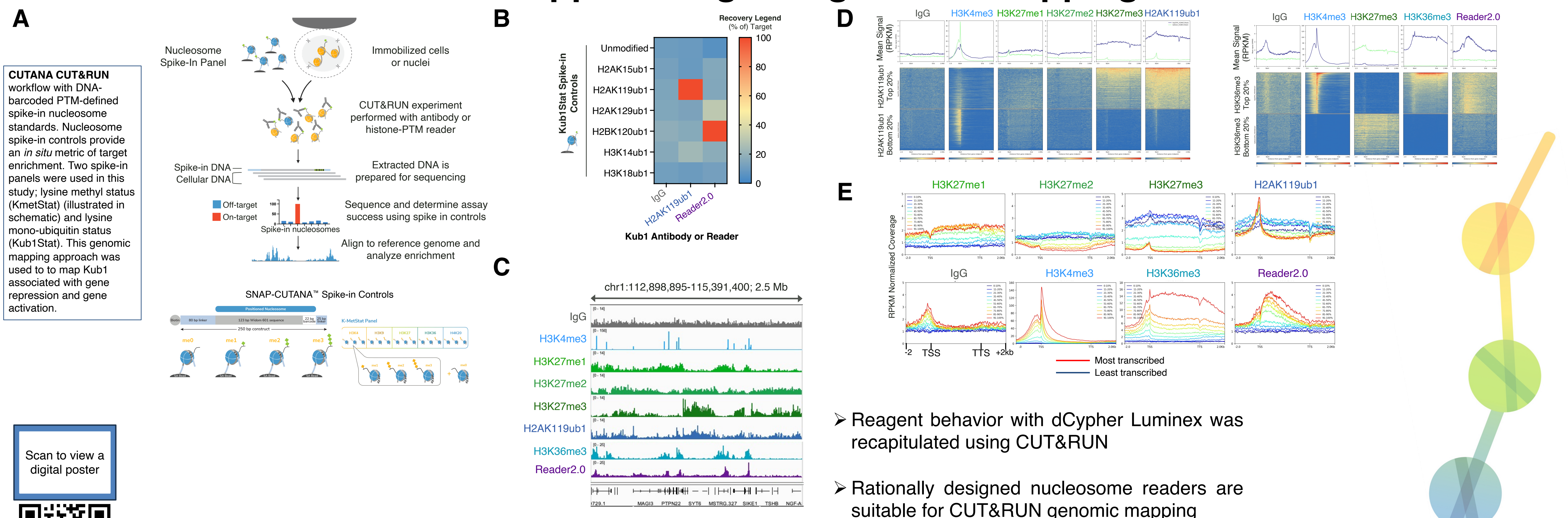
Conclusions

- Kub1-dNucs are an invaluable tool for studying histone lysine mono-ubiquitin.
- A survey of Ub and histone Kub1 antibodies identified an H2AK119ub1 and pan-Kub1 dNuc antibody while the majority failed to bind to Kub1-dNucs or were not specific for their target *in vitro*.
- Rationally designed readers of H2AK15ub1, H2BK120ub1, and pan-Kub1 dNucs were identified that can enable CUT&RUN studies where antibodies are unavailable.
- Kub1-histone peptides do not predict antibody or reader behavior toward Kub1-dNucs.
- CUTANA CUT&RUN genomic mapping with *in situ* SNAP-CUTANA spike-in controls using Luminex triaged reagents provides a powerful platform for studying Kub1-histone-PTMs.

References

- 1 Bi et al (2016) Semisynthetic UbH2A reveals different activities of deubiquitinases and inhibitory effects of H2A K119 ubiquitination on H3K36 methylation in mononucleosomes. *Organic & Biomolecular Chemistry* **14**, 835–839.
- 2 Dawson and Kouzarides (2012) Cancer epigenetics: from mechanism to therapy. *Cell* **150**:12–27.
- 3 Dos Santos Passos et al (2021) Design of genetically encoded sensors to detect nucleosome ubiquitination in live cells. *J Cell Biol* **220(4)**:e201911130.
- 4 ENCODE Project Consortium (2012) An integrated encyclopedia of DNA elements in the human genome. *Nature* **489(7414)**:57–7.
- 5 Sergey et al (2022) The complete sequence of a human genome. *Science* **376**:44–53.
- 6 Skene and Henikoff (2017) An efficient targeted nuclease strategy for high-resolution mapping of DNA binding sites. *eLife* **6**:e21856.
- 7 Skrajna et al (2020) Comprehensive nucleosome interactome screen establishes fundamental principles of nucleosome binding. *Nucleic Acids Res* **48(17)**:25.
- 8 Vaughn et al (2021) Chromatin regulation through ubiquitin and ubiquitin-like histone modifications. *Trends in Biochem Sci* **46(4)**:258.

Kub1-dNuc standards support insightful genomic mapping of Kub1 PTMs



➤ Reagent behavior with dCypher Luminex was recapitulated using CUT&RUN

➤ Rationally designed nucleosome readers are suitable for CUT&RUN genomic mapping

Scan to view a digital poster

