

Eva Brill¹, Allison Hickman¹, Bryan J. Venters¹, Andrea L. Johnstone¹, Katherine Novitzky¹, Danielle N. Maryanski¹, Keli L. Rodriguez¹, Ellen Weinzapfel¹, Matthew R. Marunde¹, Carolina P. Lin¹, Ashton Gladney¹, Rachel Watson¹, Zachary B. Gillespie¹, Augustus Adeleke¹, Saarang Gopinath¹, Leslie Lewis¹, Hannah Willis¹, Lu Sun¹, Vishnu Sunitha Kumary¹, Marcus A. Cheek¹, Matthew J. Meiners¹, James R. Bone¹, Martis W. Cowles¹, Zu-Wen Sun¹, Kami Ahmad² & Michael-Christopher Keogh¹

¹EpiCypher Inc, ²Fred Hutchinson Cancer Center

Improved assays and reagents are needed to advance epigenetic research

- Epigenomic mapping for histone post-translational modifications (PTMs) is essential for driving biological discovery
- ChIP-seq is the most widely used epigenomic mapping assay, but has major limitations:
 - Depends on PTM antibodies – which are notoriously cross-reactive^{1,2}
 - Requires large cell numbers – unsuitable for clinical or rare cell samples
 - Poor data quality – low signal to noise ratio, poor reproducibility
 - Expensive and laborious – making it difficult to scale
 - Lacks defined controls – crucial for reliable, quantitative results
- Defined controls and improved methods are needed to fully leverage epigenomics in innovative research

Nucleosome spike-in controls show that most histone PTM antibodies fail in ChIP-seq

(A) Fully defined ChIP controls mimic physiological targets

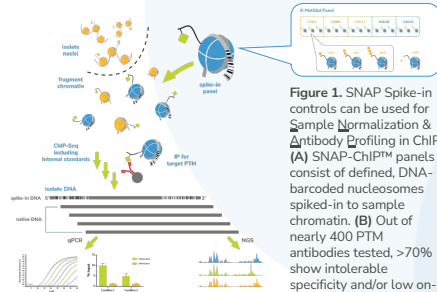


Figure 1. SNAP Spike-in controls can be used for Sample Normalization & Antibody Profiling in ChIP. (A) SNAP-ChIP panels consist of defined, DNA-barcode nucleosomes spiked-in to sample chromatin. (B) Out of nearly 400 PTM antibodies tested, >70% show intolerable specificity and/or low on-target efficiency.

(B) Most commercial antibodies are unfit for epigenomics

Antibodies Tested by SNAP-ChIP TM	Lysine Methylation	Lysine Acylation	Total
Total Tested	263	129	392
Failure Rate	74.5%	64.5%	71.2%
\$ Spent	\$105,200	\$51,600	\$156,800
\$ Wasted	\$78,374	\$33,282	\$111,656
Top Cited Citations w/ChIP [®]	3,720 (N=18)	1,031 (N=6)	4,751 (N=24)
Top Cited Failure Rate	74.5%	64.5%	79.2%

Data available at: www.ChromatinAntibodies.com³

CUTANA[™] CUT&RUN has major advantages over ChIP

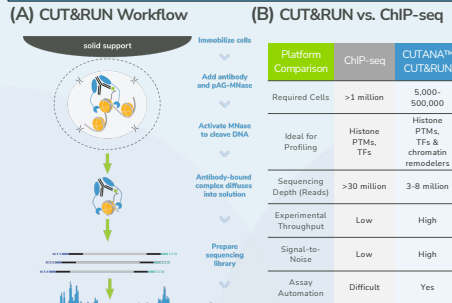
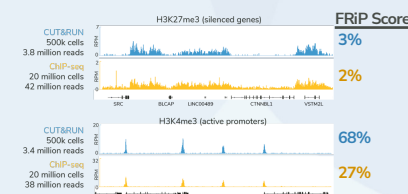


Figure 2. Overview of the CUTANA[™] CUT&RUN workflow and advantages compared to ChIP-seq. Because CUT&RUN releases antibody-bound fragments into solution (A), it has improved signal-to-noise even with significantly reduced cell numbers and sequencing depth (B).

CUT&RUN enables ultra-sensitive epigenomic profiling from low cell numbers

(A) CUT&RUN vs. ChIP-seq: better data, low sequencing costs



(B) CUT&RUN generates reliable profiles down to 5,000 cells

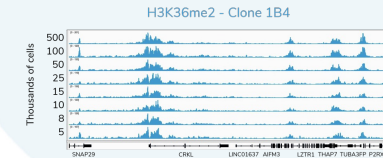


Figure 3. (A) The fraction of reads in peaks (FRIP) score, a measure of signal-to-noise (S/N), is higher in CUT&RUN vs. ChIP-seq, despite using fewer cells and reduced sequencing depths. (B) Enrichment profiles are conserved when inputs are titrated from 500k to 5k cells.

SNAP-CUTANA[™] Spike-ins are critical controls for reliable chromatin mapping

(A) SNAP Spike-ins for CUT&RUN

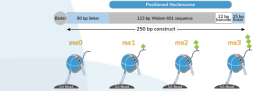
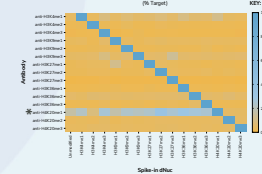
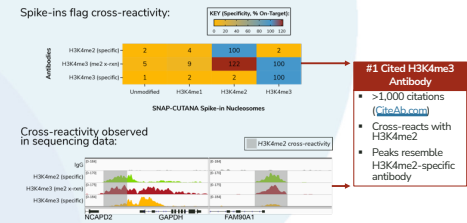


Figure 4. SNAP-CUTANA spike-in controls optimized for use in CUT&RUN (A) can be used to identify highly specific antibodies to widely studied histone PTMs (B). By spiking them into CUT&RUN workflows just prior to antibody addition, they provide a quantitative readout of on-vs. off-target recovery that predicts non-specific peaks in genomic data (C). *Ongoing testing for H4K20me1. Inquire at info@epicypher.com for up-to-date information.

(B) SNAP testing identifies specific antibodies for CUT&RUN



(C) SNAP Spike-ins predict non-specific recovery in CUT&RUN



Conclusions

- CUTANA[™] CUT&RUN is poised to rapidly replace ChIP-seq
- Our data illustrate how CUT&RUN could be applied for epigenomics research, particularly for low sample inputs and/or high-throughput applications
- SNAP Spike-in controls address pervasive antibody specificity problems while enabling a direct readout of assay success and quantitative normalization (see below)

References

- Shah et al. Examining the roles of H3K4 methylation states with systematically characterized antibodies. *Mol. Cell* 72, 162-177 (2018).
- Small et al. Chromatin Immunoprecipitation (ChIP) to study DNA-Protein Interactions. *Methods Mol. Biol.* 2261, 323-343 (2021).
- Maryanski et al. Novel nucleosome-based methods for rapid screening and identification of best-in-class antibodies: a community resource to improve genomic mapping. In preparation. www.ChromatinAntibodies.com

Ultra-sensitive epigenomics drives biological discovery

Uncover novel biology with highly specific antibodies

Quantitative readout enables drug & clinical development applications

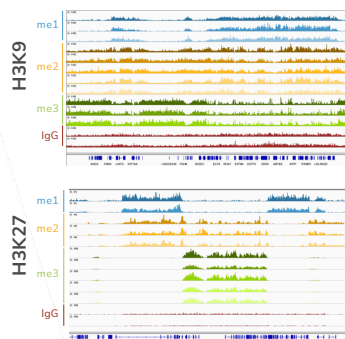


Figure 5. GENOMICALLY DISTINCT = FUNCTIONALLY DISTINCT. The use of highly specific histone PTM antibodies in CUT&RUN (Figure 4B) enables novel insights into the histone code. Distinct genomic profiles observed by mono-, di- and tri-methyl states imply distinct biological functions not previously appreciated.

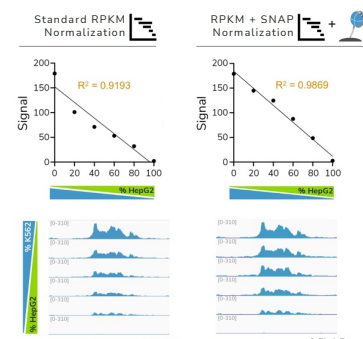


Figure 6. CUT&RUN linearity is improved by SNAP normalization. Cells with distinct lineages were mixed at defined intervals. Linear regression analysis of an H3K4me3 peak shows that SNAP normalization (right) shows improved linearity of a K562 cell-specific peak over standard RPKM normalization (left).

● K562 cells (mesodermal) ● HepG2 cells (ectodermal)

