

# Nucleosome spike-in controls enable reliable next-generation epigenomic mapping

Andrea L. Johnstone<sup>1</sup>, Danielle N. Maryanski<sup>1</sup>, Keli L. Rodriguez<sup>1</sup>, Ellen Weinzapfel<sup>1</sup>, Bryan J. Venters<sup>1</sup>, Katherine Novitzky<sup>1</sup>, Matthew R. Marunde<sup>1</sup>, Carolina P. Lin<sup>1</sup>, Zachary B. Gillespie<sup>1</sup>, Augustus Adeleke<sup>1</sup>, Saarang Gopinath<sup>1</sup>, Leslie Lewis<sup>1</sup>, Marcus A. Cheek<sup>1</sup>, Matthew J. Meiners<sup>1</sup>, Zu-Wen Sun<sup>1</sup>, Michael-Christopher Keogh<sup>1</sup>

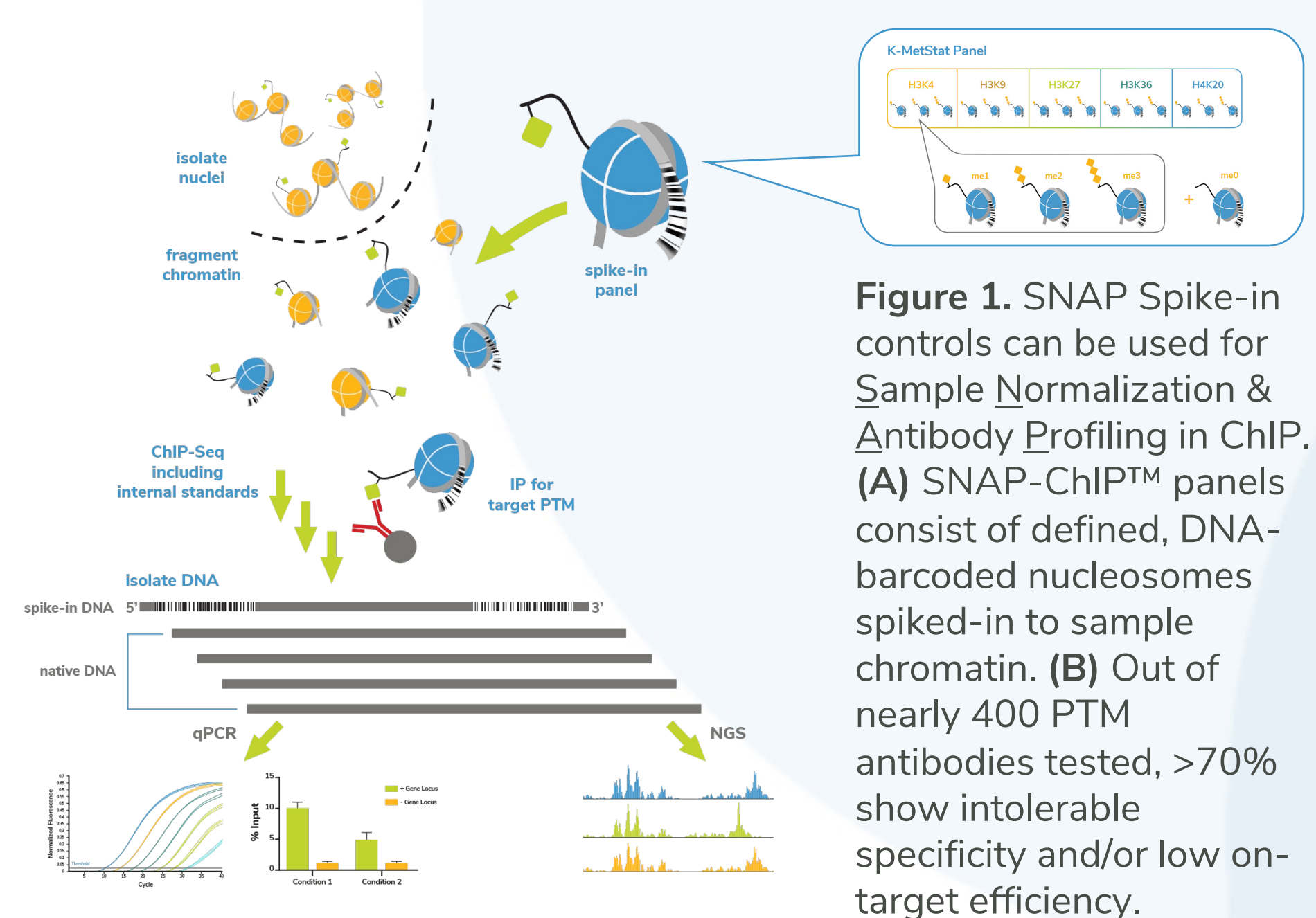
<sup>1</sup>Epicpyher Inc., Research Triangle Park, Durham NC 27709, USA

## The Problem: Improved epigenomics assays are needed to advance cancer research

- Epigenomic mapping for histone post-translational modifications (PTMs) can be used to discover gene regulatory mechanisms in cancer
- ChIP-seq is the most widely used epigenomic mapping assay, but has major limitations:
  - ✗ Depends on PTM antibodies – which are notoriously cross-reactive<sup>1,2</sup>
  - ✗ Requires large cell numbers – unsuitable for clinical or rare cell samples
  - ✗ Poor data quality – low S:N, 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 cancer research

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

### (A) Fully defined ChIP controls mimic physiological targets

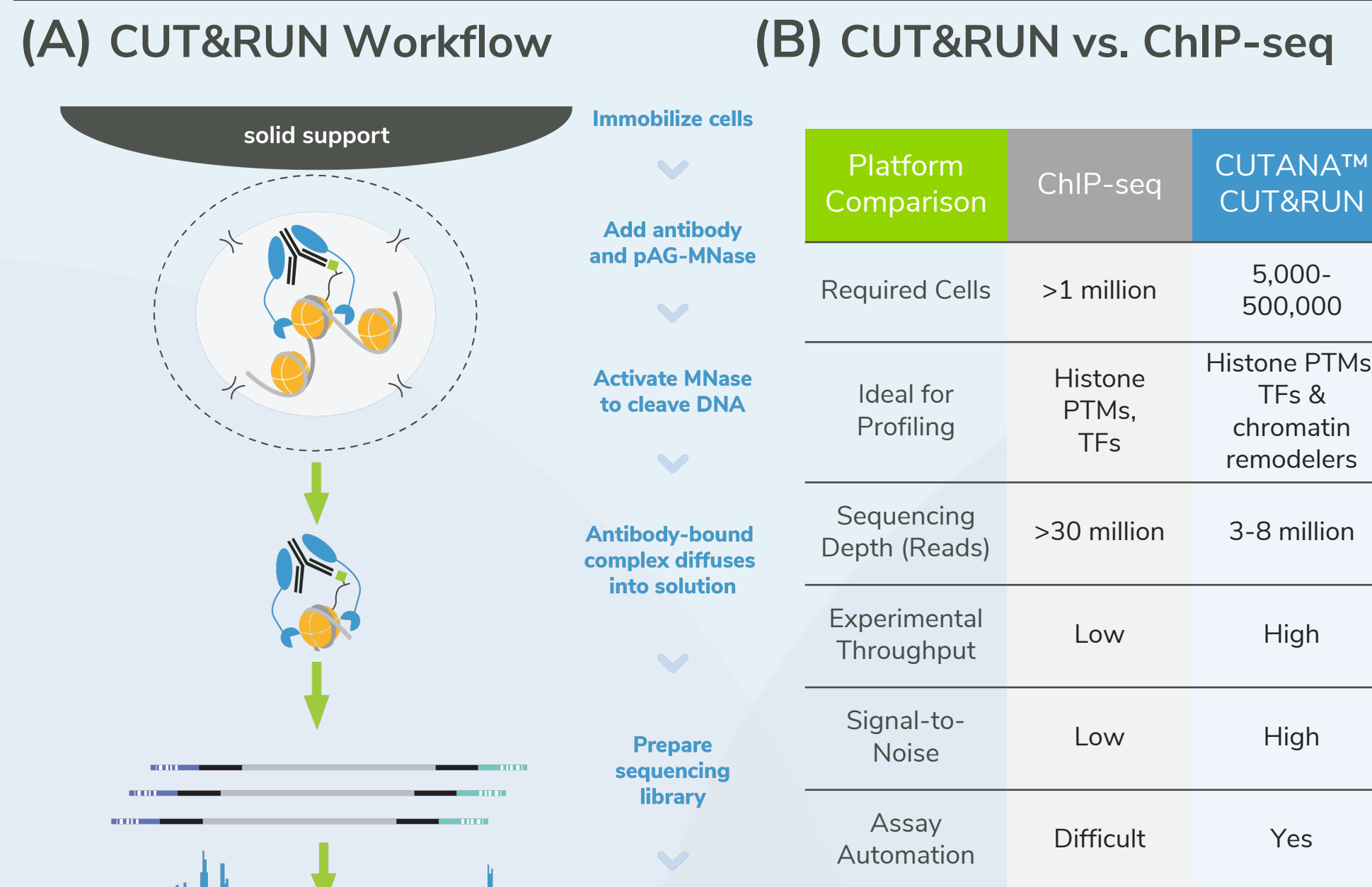


### (B) Most commercial antibodies are unfit for epigenomics

Antibodies Tested by SNAP-ChIP™	Lysine Methylation	Lysine Acylation	Total
Total Tested	263	129	392
Failure Rate	74.5%	64.5%	<b>71.2%</b>
\$ 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%	<b>79.2%</b>

Data available at: [www.ChromatinAntibodies.com](http://www.ChromatinAntibodies.com)<sup>3</sup>

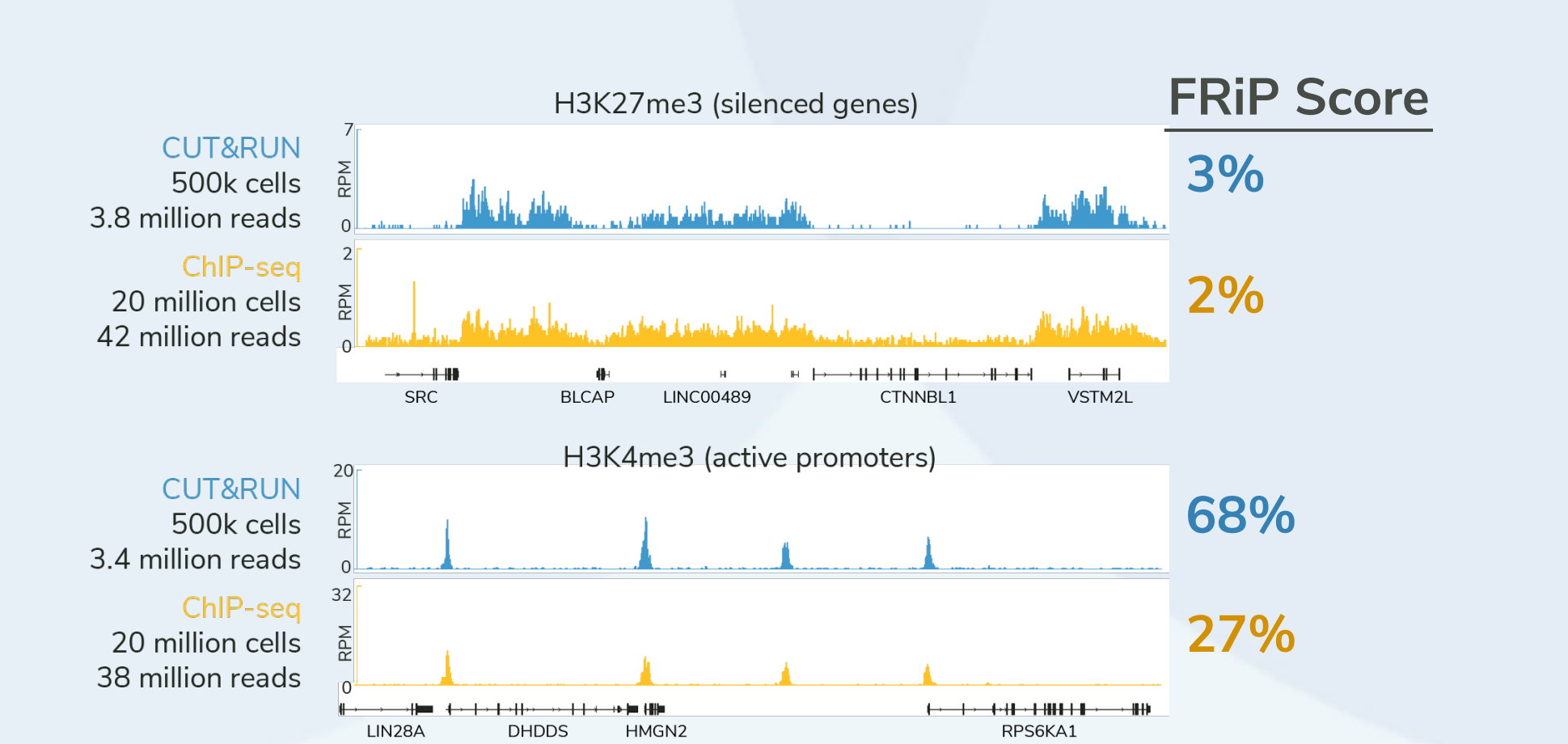
## CUTANA™ CUT&RUN has major advantages over ChIP



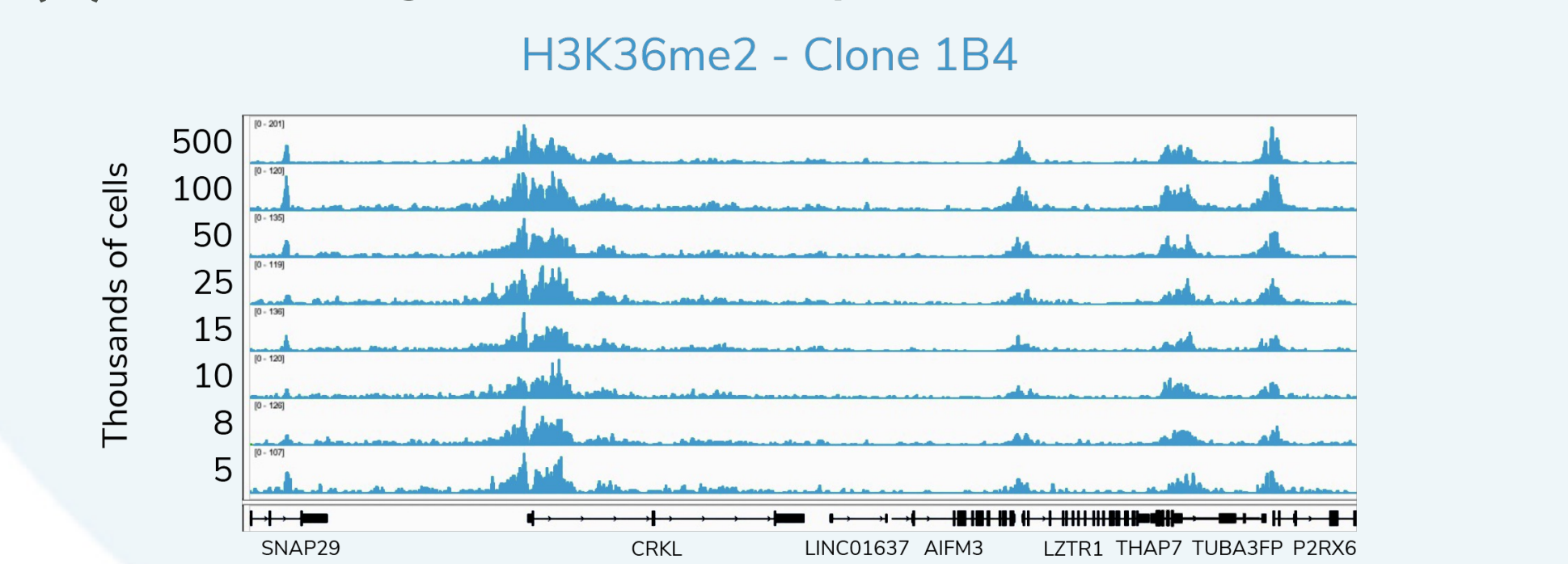
**Figure 2.** Overview of the CUTANA™ CUT&RUN workflow and advantages compared to ChIP-seq. Because CUT&RUN releases antibody-bound fragments into solution (A), it has improved signal-to-noise even with significantly reduced cell numbers and sequencing depth (B).

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

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



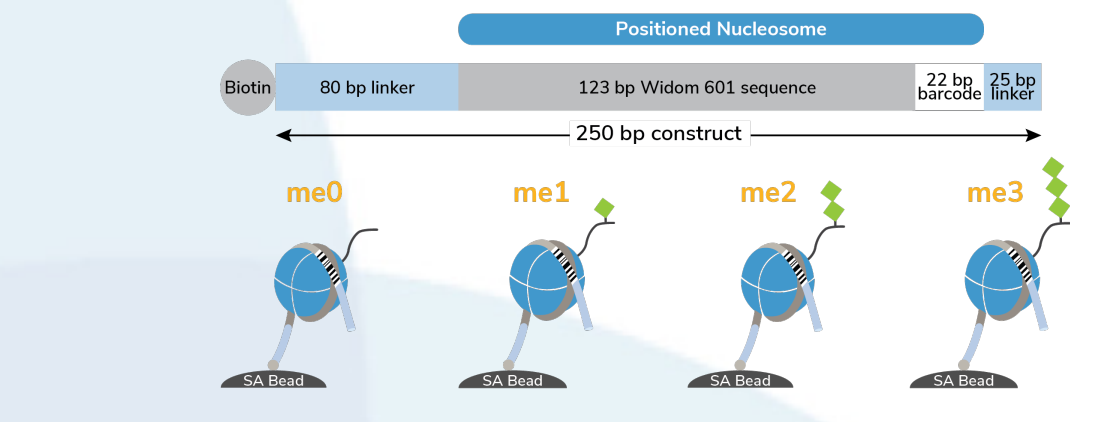
### (B) CUT&RUN generates reliable profiles down to 5,000 cells



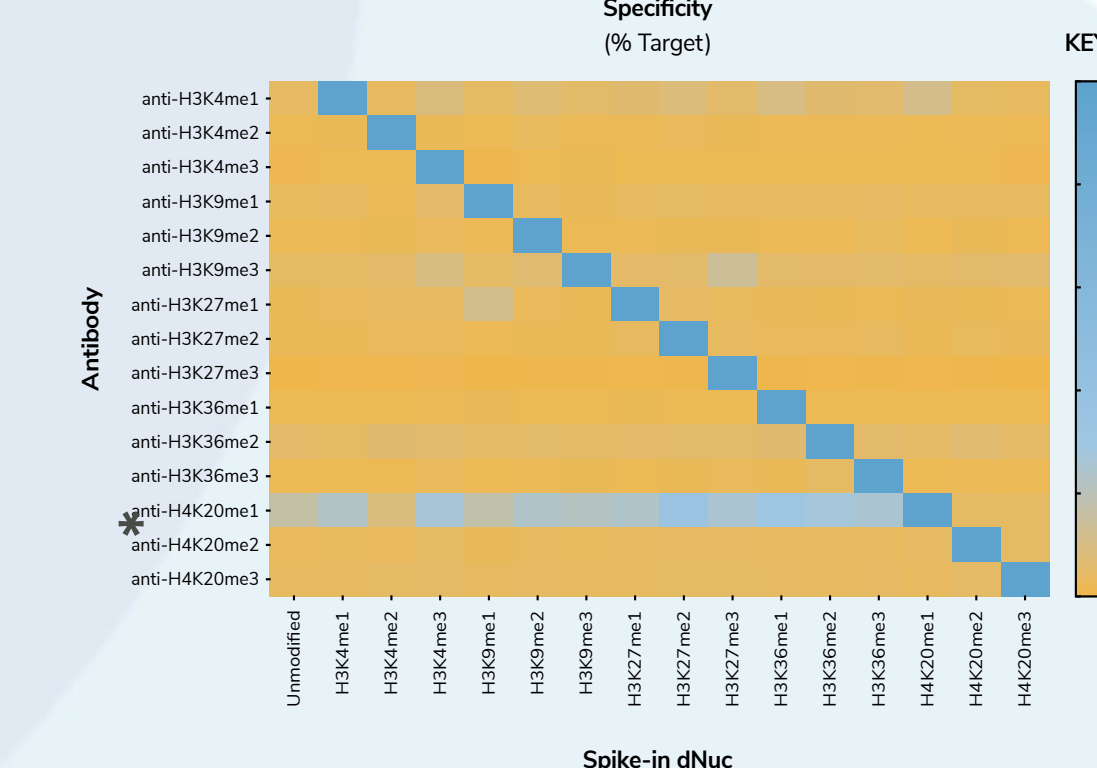
**Figure 3.** (A) The Fraction of Reads on Peaks (FRiP) score, a measure of signal-to-noise (S:N), is higher in CUT&RUN vs. ChIP-seq, despite using fewer cells and reduced sequencing depths. (B) Enrichment profiles are conserved when inputs are titrated from 500k to 5k cells.

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

### (A) SNAP Spike-ins for CUT&RUN

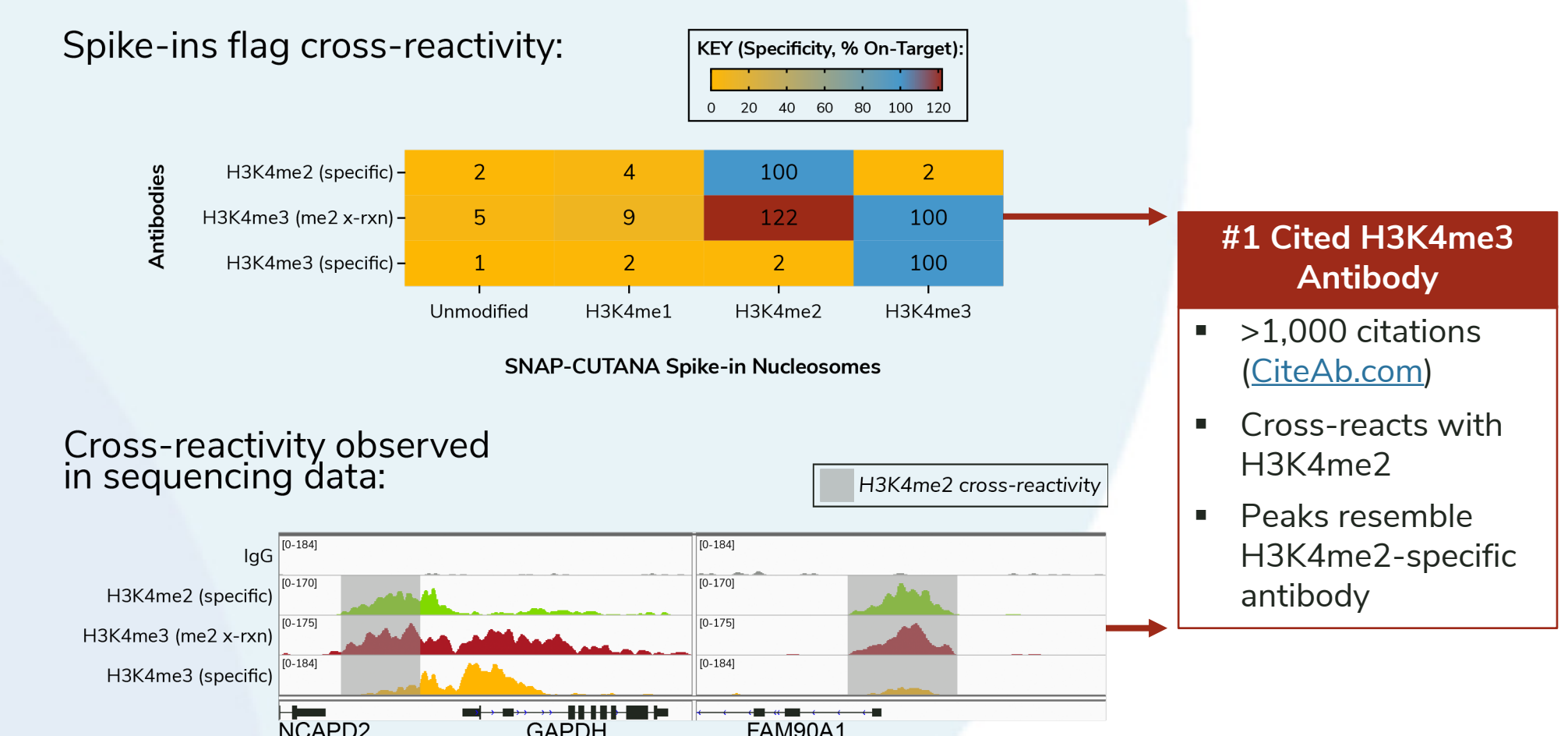


### (B) SNAP testing identifies specific antibodies for CUT&RUN



**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](mailto:info@epicypher.com) for up-to-date information.

### (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 cancer epigenomics, 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!)

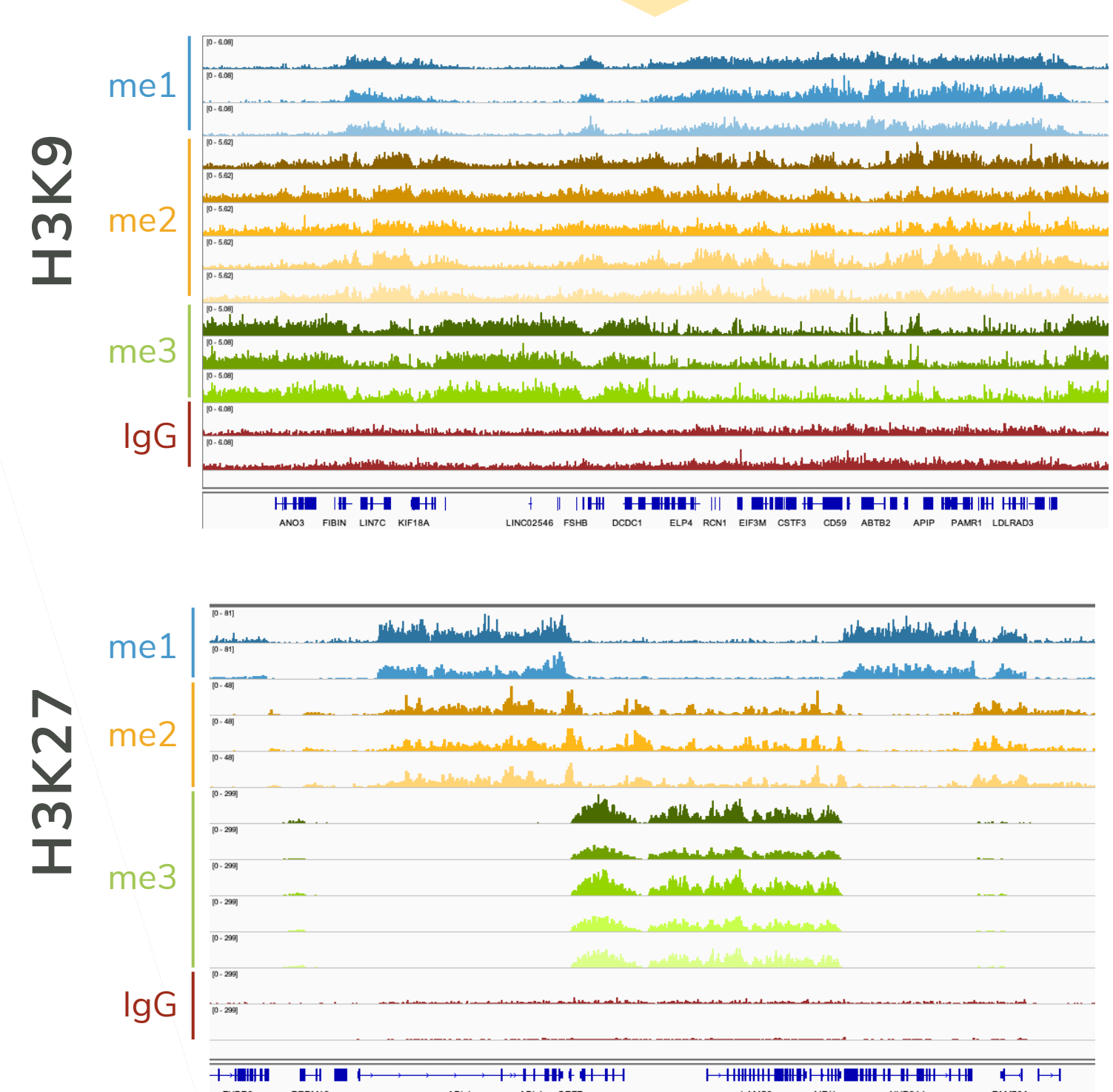
## References

- 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](http://www.ChromatinAntibodies.com)

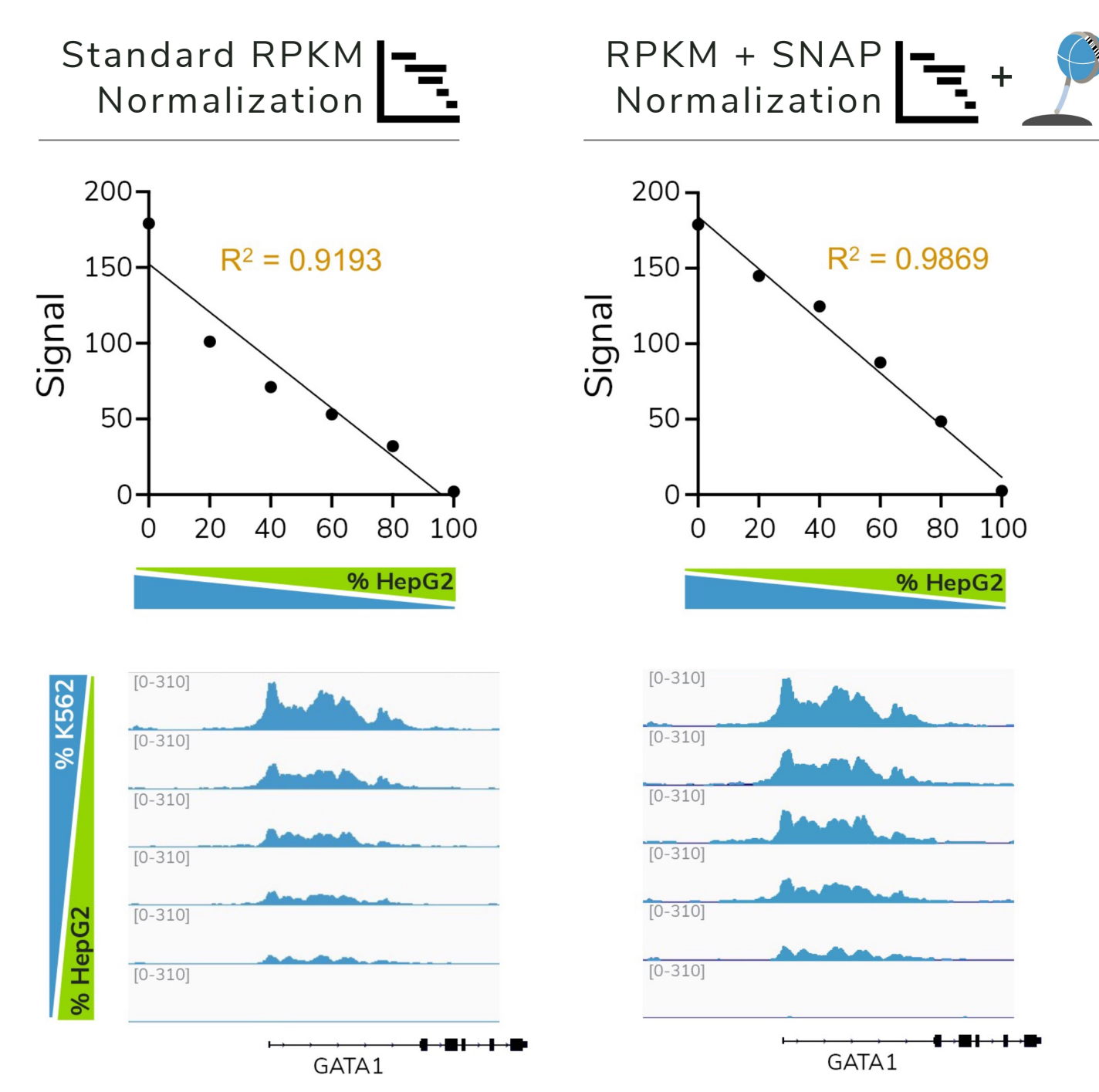
# Ultra-sensitive epigenomics advances cancer research

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)

SCAN TO DOWNLOAD THIS POSTER NOW

