

# CUTANA™ CUT&Tag : Powerful Platform for Streamlined, Ultra-Sensitive Epigenomics



James Anderson<sup>1</sup>, Laylo Mukhsinova<sup>1</sup>, Allison R Hickman<sup>1</sup>, Vishnu U. Sunitha Kumary<sup>1</sup>, Katherine Novitzky<sup>1</sup>, John Pettingill<sup>1</sup>, Eva Brill<sup>1</sup>, Zu-Wen Sun<sup>1</sup>, Martis W. Cowles<sup>1</sup>, Bryan J. Venters<sup>1</sup>, & Michael-Christopher Keogh<sup>1</sup>

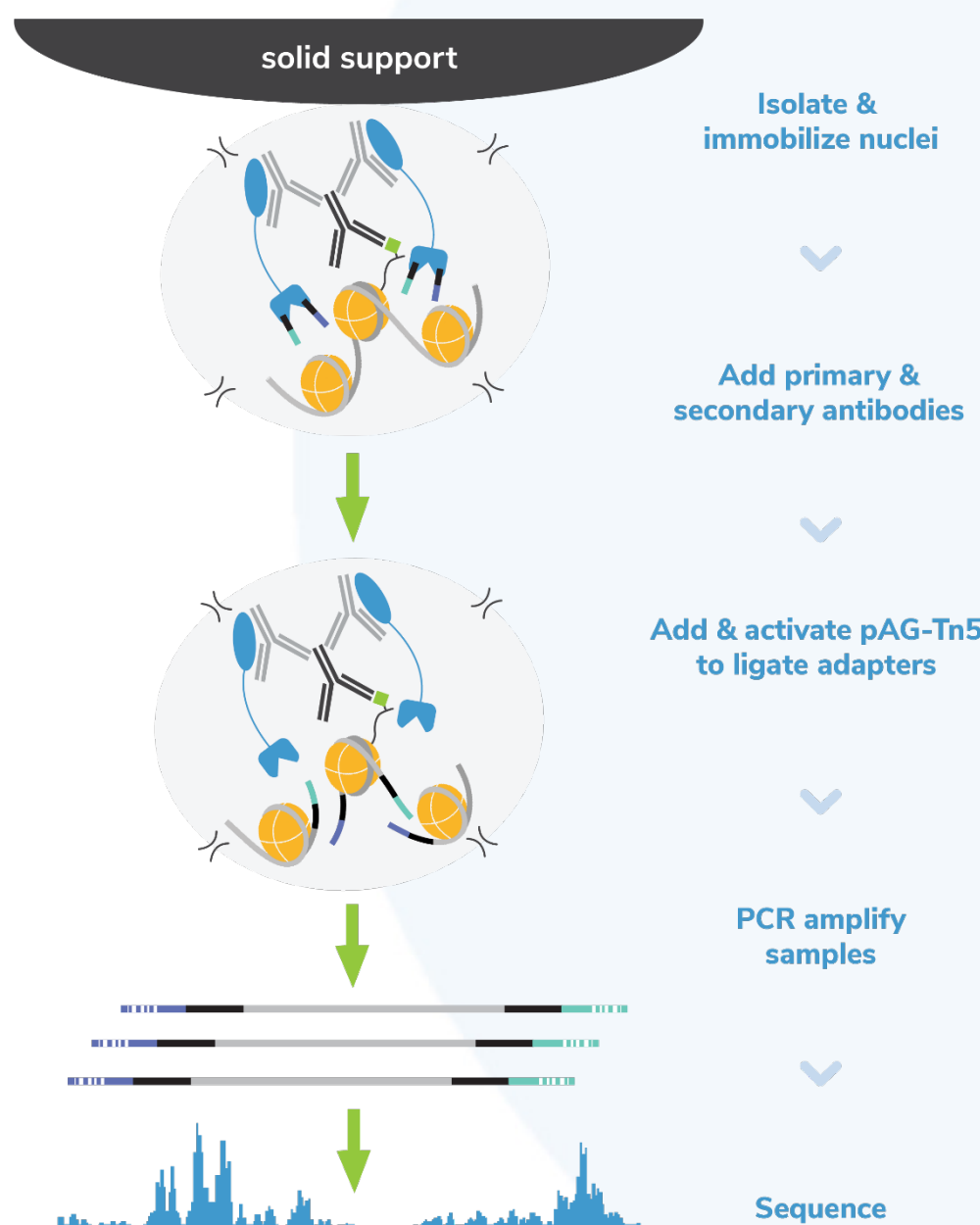
<sup>1</sup> EpiCypher Inc, Durham, NC 27709

## 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<sup>1,2</sup>
  - ✗ Requires large cell numbers – unsuitable for clinical or rare cell samples
  - ✗ Poor data quality – low signal to noise ratio, poor reproducibility
  - ✗ Lacks defined controls – crucial for reliable, quantitative results
- Compared to ChIP-seq, CUT&Tag<sup>3</sup> provides higher quality sequencing data with improved sensitivity and dramatically reduced background.

## Advantages of CUTANA™ CUT&Tag over ChIP-seq

### (A) CUT&Tag Workflow



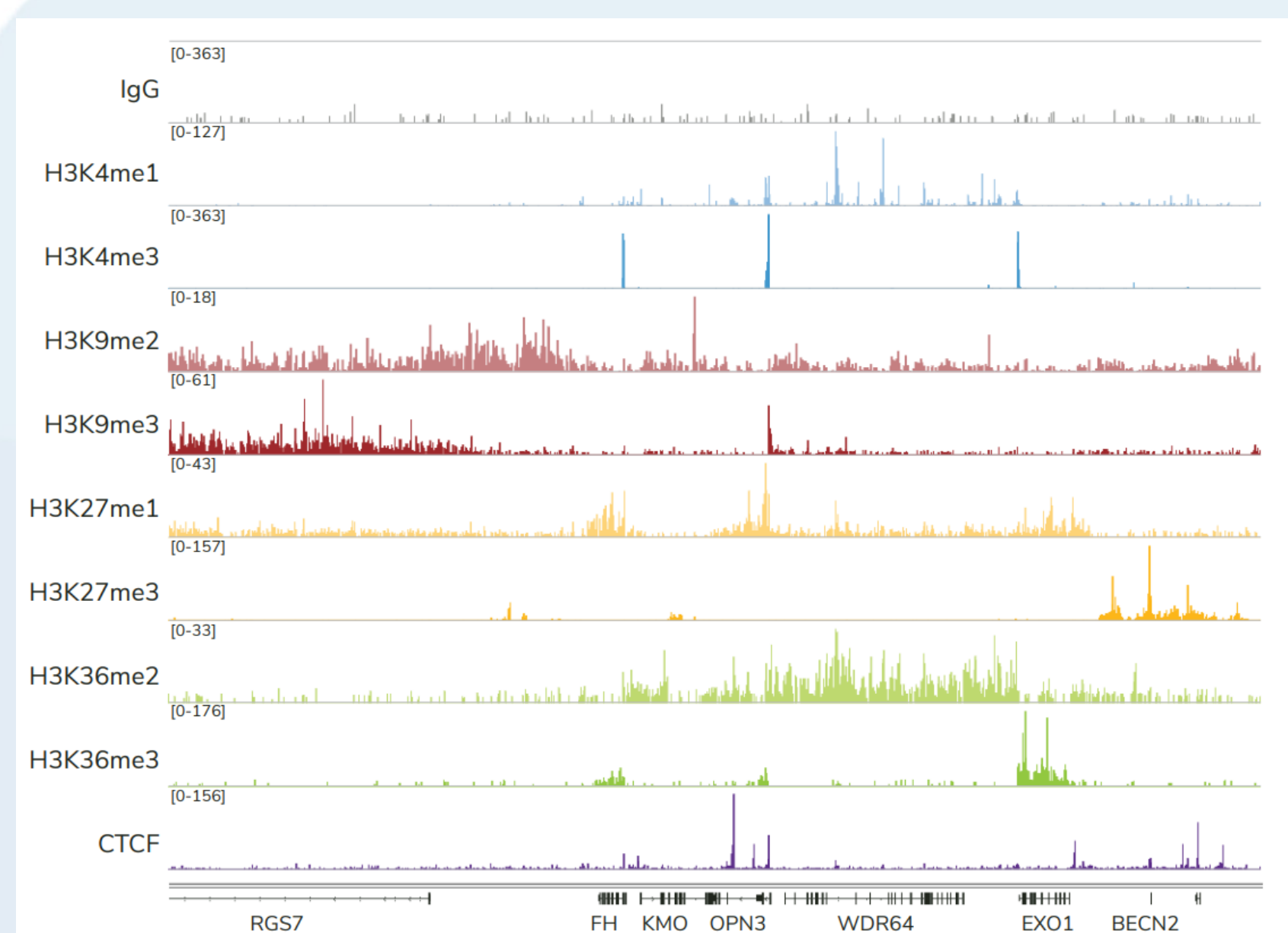
### (B) CUTANA™ vs. ChIP-seq

Platform Comparison	ChIP-seq	CUTANA™ CUT&RUN	CUTANA™ CUT&Tag
Required Cells	>1 million	5,000-500,000	10,000-100,000
Ideal for Profiling	Histone PTMs, TFs	Histone PTMs, TFs & remodelers	Histone PTMs
Sequencing Depth (Reads)	>30 million	3-8 million	3-8 million
Experimental Throughput	Low	High	High
Signal-to-Noise	Low	High	High
Library Prep	Standard	Standard	Streamlined Direct-to-PCR

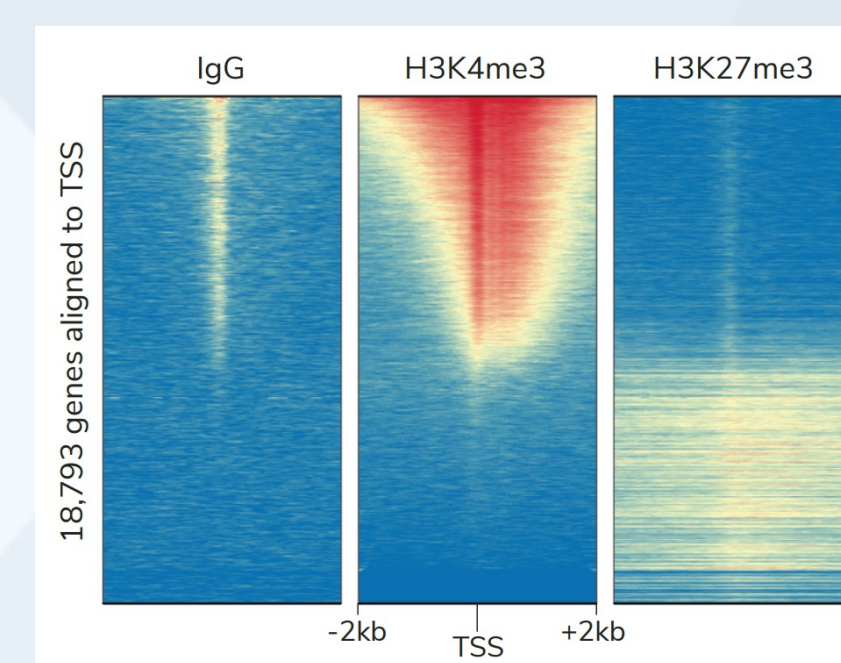
**Figure 1.** Overview of the CUTANA™ CUT&Tag workflow and advantages compared to ChIP-seq. Because CUTANA™ technologies release antibody bound fragments into solution (A), it has improved signal-to-noise even with significantly reduced cell numbers and sequencing depth (B). In CUT&Tag, pAG-Tn5 inserts sequencing adapters at antibody bound chromatin in intact nuclei, streamlining library preparation.

## CUTANA™ CUT&Tag is ideal for mapping histone PTMs

### (A) CUT&Tag delivers robust data across diverse PTM targets



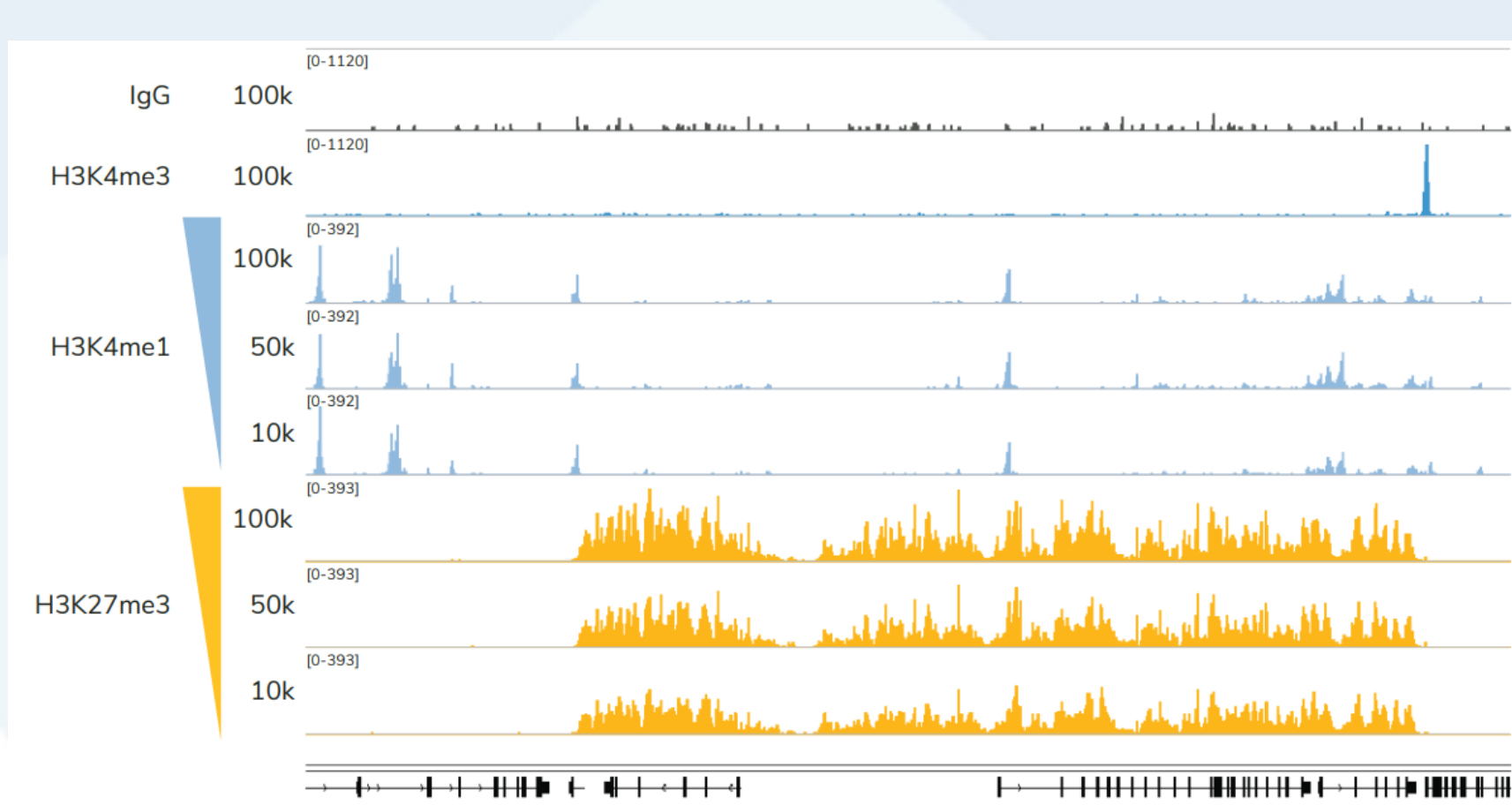
### (B) TSS Enrichment



**Figure 2.** (A) CUTANA CUT&Tag generates high-quality profiles for targets in both active and repressed chromatin regions and select proteins. Rabbit IgG is shown as a negative control. (B) Expected results from CUTANA CUT&Tag assays. Genes are aligned across targets and ranked by H3K4me3 intensity at transcription start sites (TSSs) from top (high signal, red) to bottom (low signal, yellow).

## CUT&Tag excels at epigenetic profiling from low cell numbers

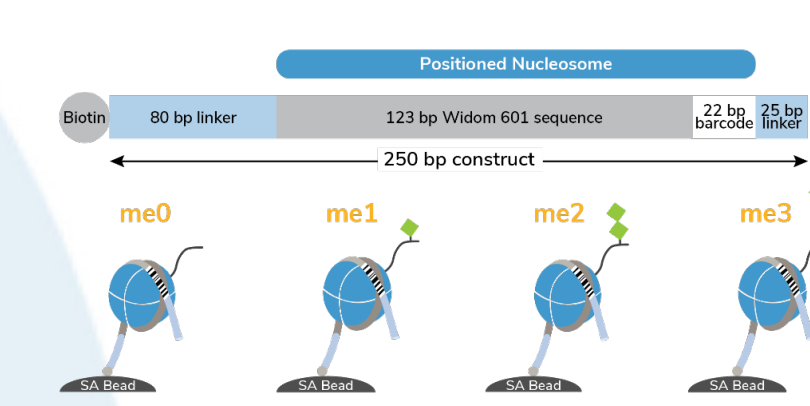
### (A) CUT&Tag generates reliable profiles down to 10K cells



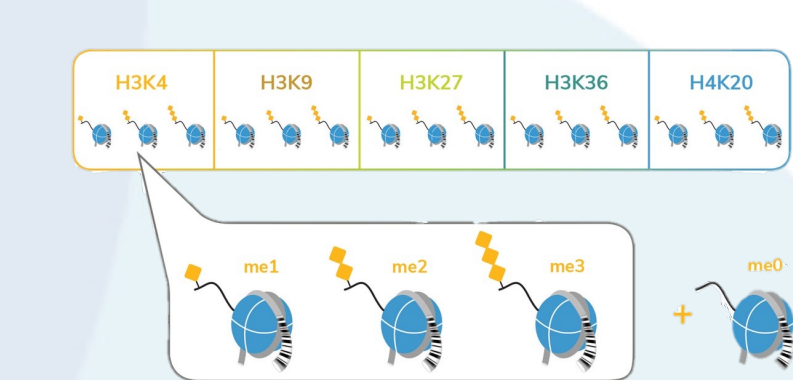
**Figure 3.** (A) CUTANA CUT&Tag enables reliable chromatin profiling from low cell numbers. CUT&Tag was used to map H3K4me1 and using decreasing numbers of K562 cells. Data quality at 10,000 cells is comparable to standard inputs of 100,000 cells.

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

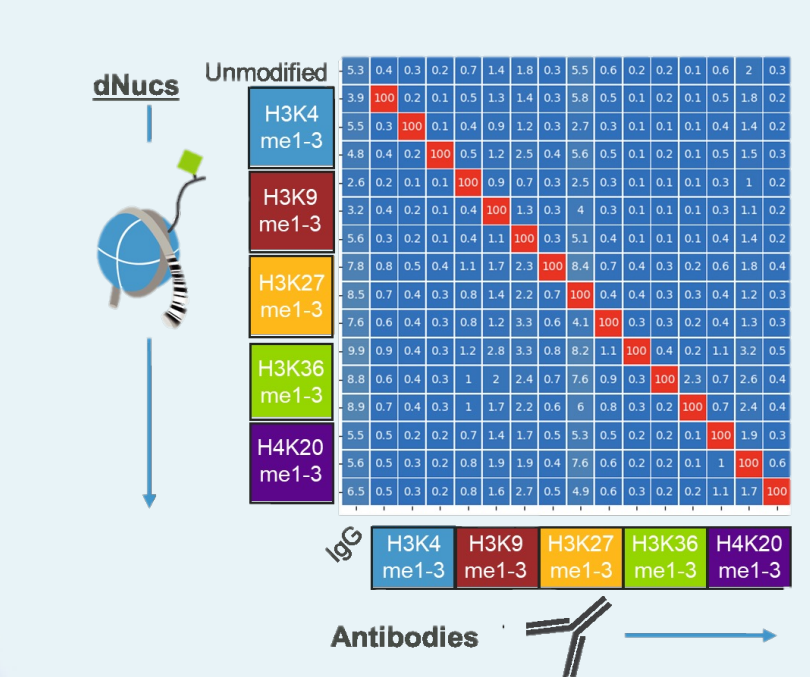
### (A) SNAP Spike-ins for CUTANA Technologies



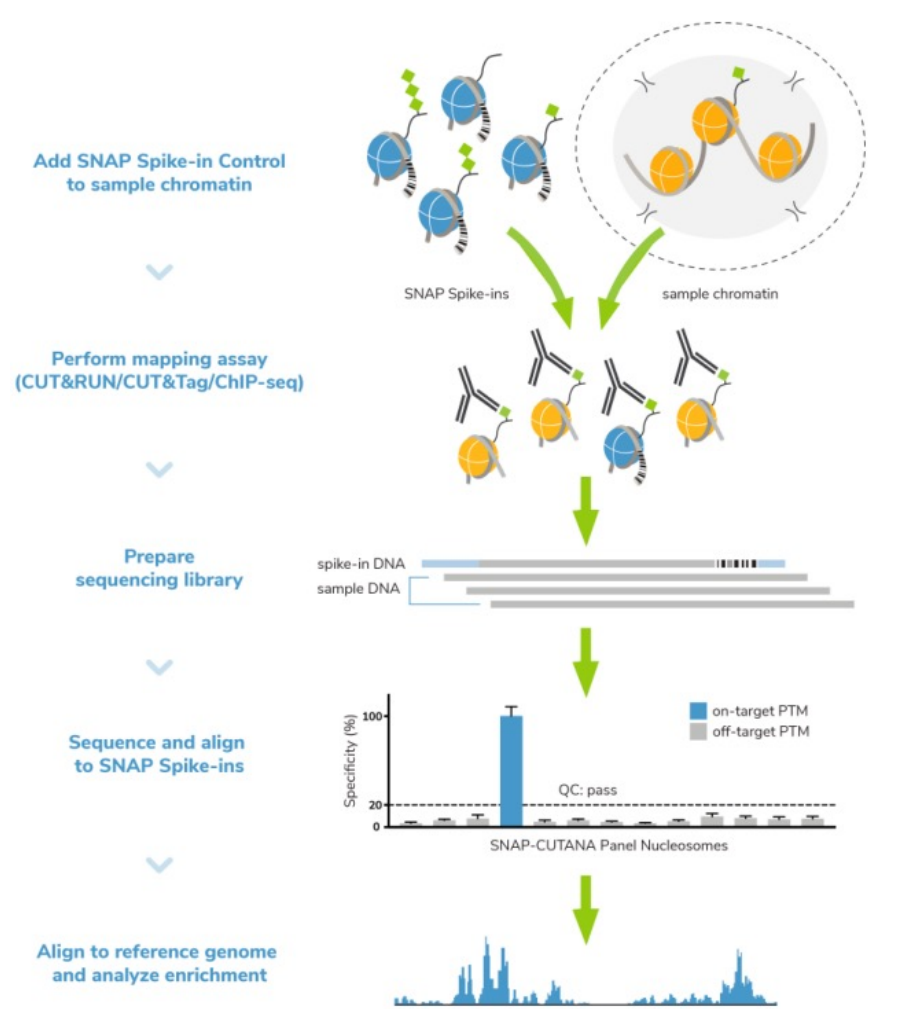
### (B) SNAP testing identifies specific antibodies



### (D) SNAP-CUTANA Specificity Heatmap



### (C) Spike-In Workflow



**Figure 4.** SNAP CUTANA controls can be used for Sample Normalization & Antibody Profiling in CUTANA assays. (A) SNAP-CUTANA™ panels consist of defined, DNA-barcoded nucleosomes spiked-in to sample chromatin. (B) KMetStat panel that would be processed alongside sample as an ideal internal control. (C) Panel is spiked into CUT&RUN workflows just prior to antibody addition. (D) They provide a quantitative readout of on- vs. off-target recovery that predicts non-specific peaks in genomic data.

## Conclusions

- CUTANA™ technologies (CUT&RUN and CUT&Tag) are poised to rapidly replace ChIP-seq.
- CUT&Tag uses a streamlined protocol to produce high quality data for diverse targets with low cell number requirements and reduced sequencing costs.
- SNAP Spike-in controls inform antibody specificity and monitor assay success.
- CUTAC recapitulates published ATAC-seq datasets but with significantly improved sensitivity and signal-to-noise.

## References

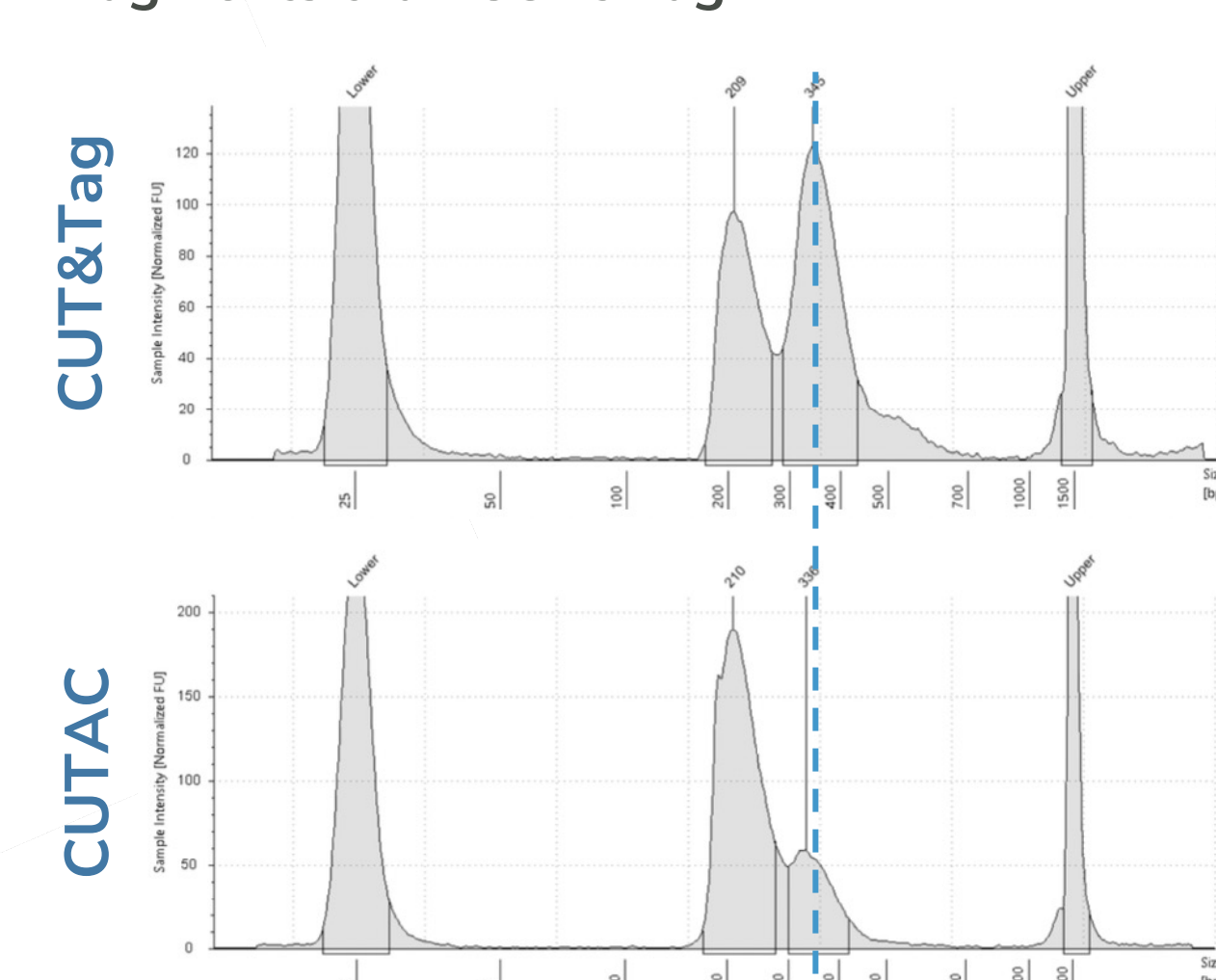
- Shah et al. Examining the roles of H3K4 methylation states with systematically characterized antibodies. *Mol. Cell* 72, 162-177 (2018).
- Small et al. Chromatin Immunoprecipitation (ChIP) to study DNA-Protein Interactions. *Methods Mol. Biol.* 2261, 323-343 (2021).
- Kaya-Okur HS et al. CUT&Tag for efficient epigenomic profiling of small samples and single cells. *Nat Commun* 10, 1930 (2019).
- Steven Henikoff/Lojza G Henikoff/Hatice S Kaya-Okur/Kami Ahmad (2020) Efficient chromatin accessibility mapping in situ by nucleosome-tethered tagmentation. *eLife* 9:e63274.

# CUTAC : Identifying Open Chromatin with Modified CUT&Tag

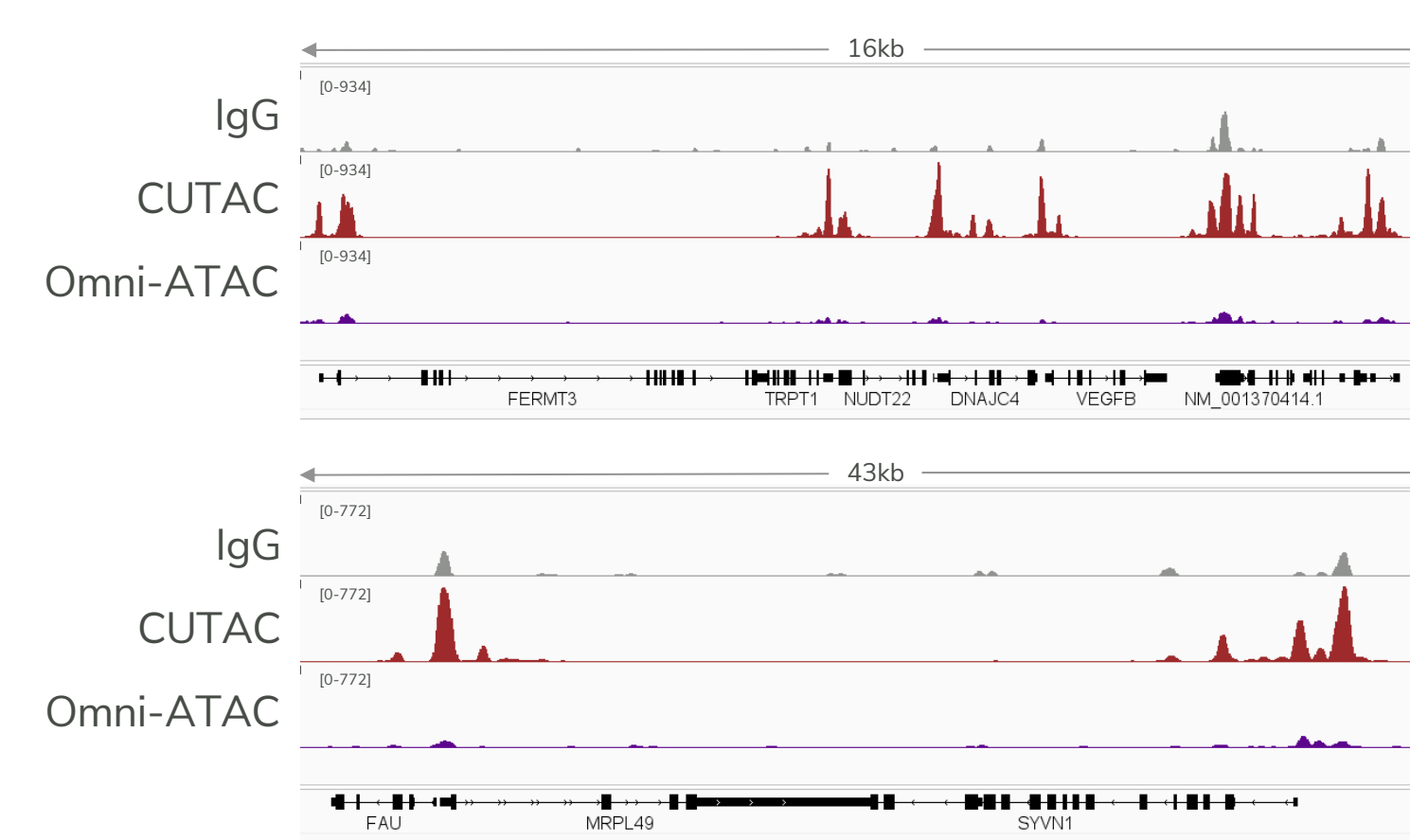
### (A) CUTAC conditions differ from CUT&Tag

Platform Comparison	CUT&Tag	CUTAC
Feature Enrichment	Histone PTMs	Chromatin Accessibility
Detection Reagent	PTM-specific Antibody	H3K4me2 Antibody
Tagmentation Salt Conditions	300mM NaCl, 10mM MgCl <sub>2</sub>	0mM NaCl, 5mM MgCl <sub>2</sub>
Tagmentation Time	1hr	20min
Tagmentation Temperature	37C	37C

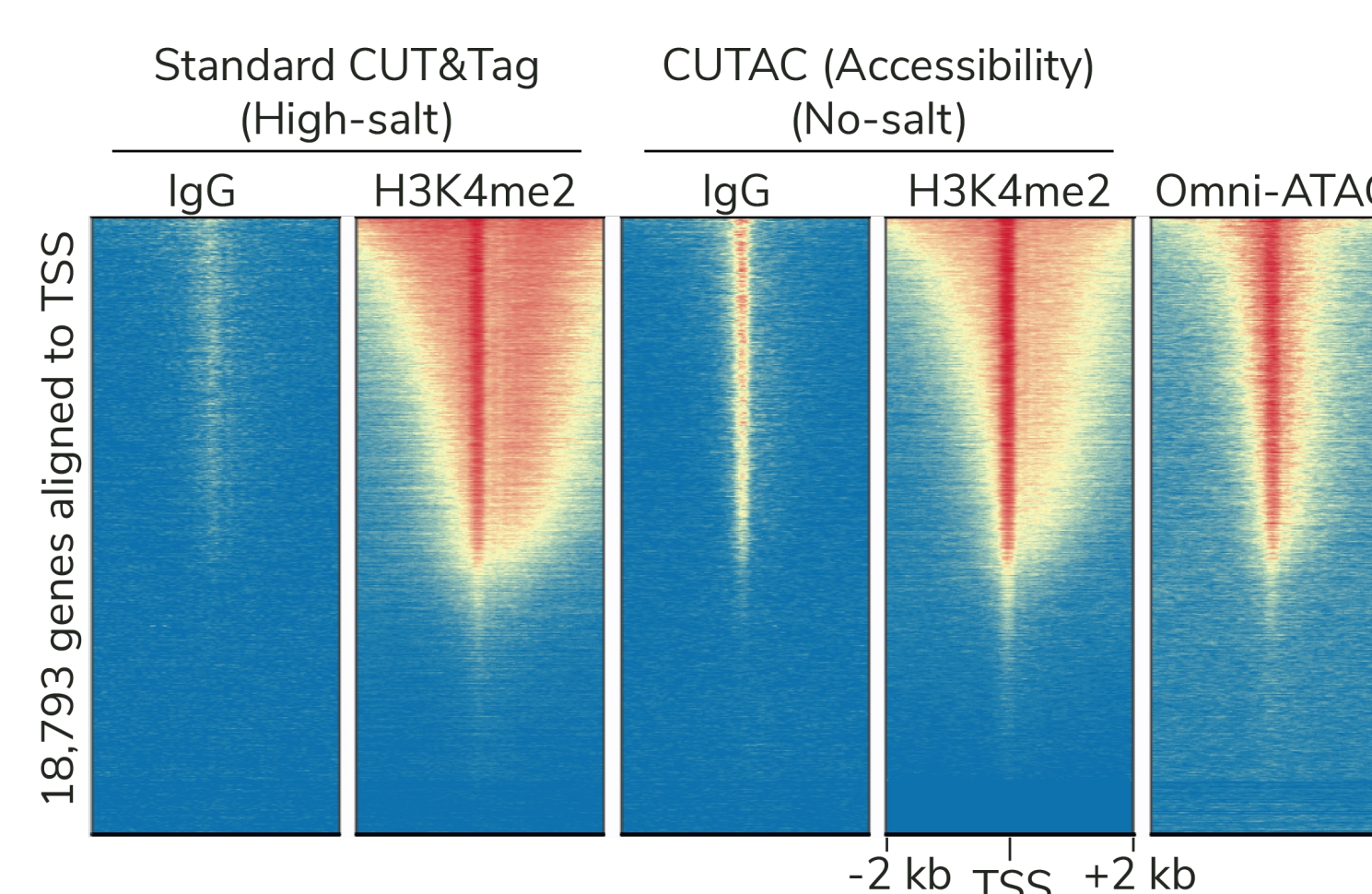
### (B) CUTAC releases smaller genomic fragments than CUT&Tag



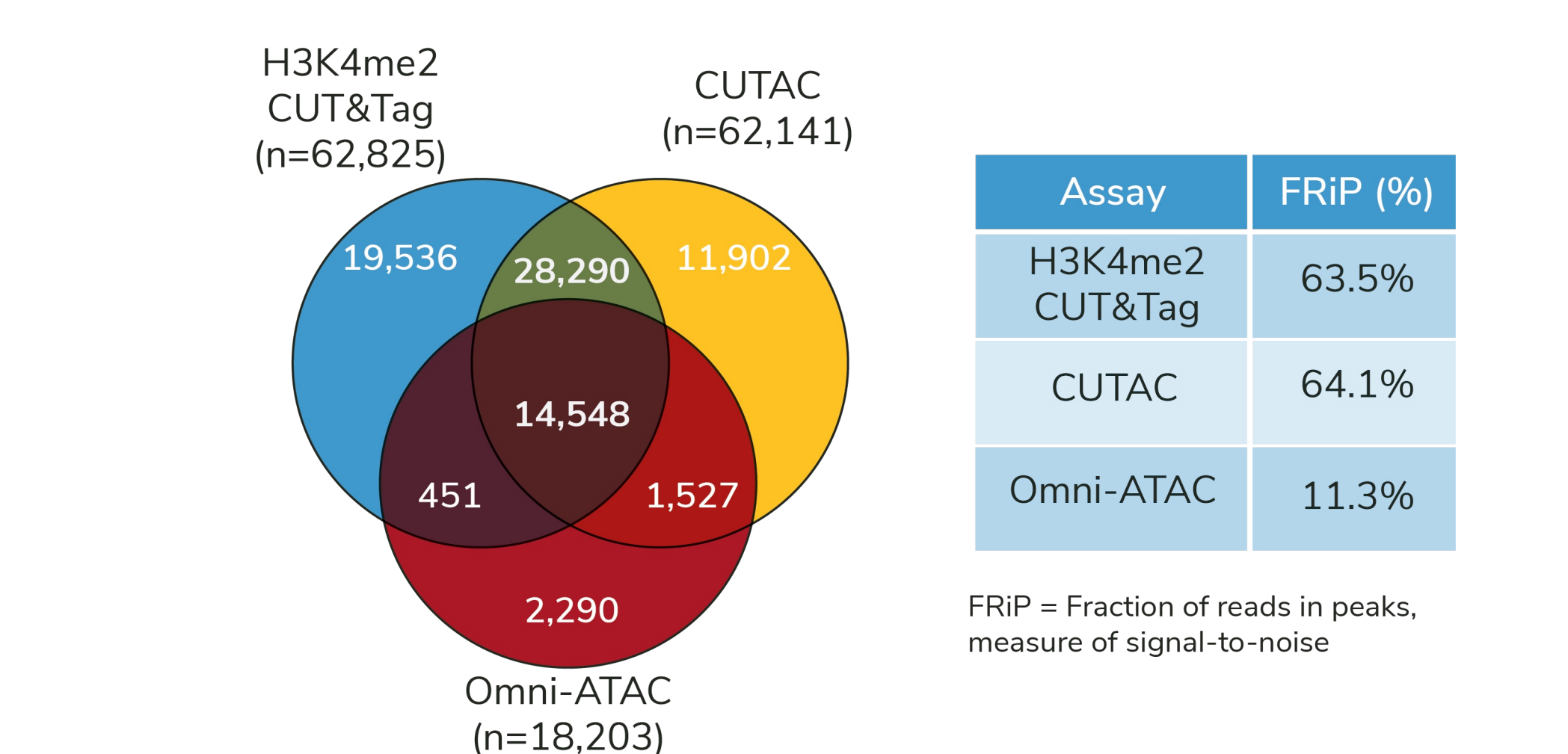
### (C) CUTAC shows stronger signal than Omni-ATAC



### (D) CUTAC, CUT&Tag, and Omni-ATAC enrichment relative to TSS



### (E) Peak calling comparisons across CUTAC, CUT&Tag, and Omni-ATAC



**Figure 5.** CUTAC<sup>4</sup> is an exciting new application of CUT&Tag that identifies open chromatin regions by restricting tagmentation to accessible DNA at nearby transcriptionally active chromatin. (A) CUTAC experiments use less salt and a shorter tagmentation time compared to CUT&Tag. (B) TapeStation traces from CUTANA CUT&Tag and CUTAC libraries prepared using H3K4me2 antibodies. While CUT&Tag libraries predominantly enrich for mononucleosome-sized fragments (~300bp fragments), CUTAC libraries enrich for smaller fragments (~200bp). (C) IGV screenshots of CUTAC and Omni-ATAC data compared to non-specific IgG. (D) Heatmaps show signal relative to TSS in each assay. Gene rows are aligned across conditions. (E) Venn diagram showing peak overlap between H3K4me2 CUT&Tag, CUTAC, and Omni-ATAC-seq. FRiP scores show that CUT&Tag and CUTAC data have higher signal-to-noise vs. Omni-ATAC-seq.

### Other Applications of CUT&Tag

- SWI/SNF in SARS-CoV-2 Infection (PMID: 36894709)
- CD4+ T cell calibration (PMID: 37188942)
- 3D genome in AML (PMID: 36289338)

Scan to download this poster now:

