

CUTANA™ CUT&RUN and CUT&Tag epigenomic mapping assays for agricultural research

Kelsey E. Noll¹, Katherine Novitzky¹, Matthew R. Marunde¹, Danielle N. Maryanski¹, Catherine Smith¹, Allison Hickman¹, Kateřina Adamusová², Karina Chalenko², Swetha Chenna², Quentin Thomas², Jinghan Liu², Andrea L. Johnstone¹, Keith E. Maier¹, Ellen Weinzapfel¹, Irina K. Popova¹, Bryan J. Venters¹, Keli L. Rodriguez¹, Carolina P. Lin¹, Ashton Gladney¹, Rachel Watson¹, Zachary B. Gillespie¹, Augustus Adeleke¹, Saarang Gopinath¹, Leslie Lewis¹, Hannah Willis¹, Lu Sun¹, Marcus A. Cheek¹, Matthew J. Meiners¹, James R. Bone¹, Martis W. Cowles¹, Zu-Wen Sun¹, Sebastian Marquardt², Michael-Christopher Keogh¹

¹EpiCypher Inc., Durham, NC, USA; ²Department of Plant and Environmental Sciences, Copenhagen Plant Science Centre, University of Copenhagen, Denmark



Epigenomic mapping provides key insights into gene regulation, but many commonly used approaches are not sufficient

- Many genomic strategies for cellular profiling focus on transcription; however, RNA-seq reveals the **outcomes** – not driving **mechanisms**
- **Epigenomics is the solution:** mapping the location of chromatin features provides mechanistic insights that are central to cell fate and function
- Existing epigenomic technologies (e.g., ChIP-seq) lack standardized controls, require high cell inputs, are low-throughput, and have high costs

CUTANA™ CUT&RUN and CUT&Tag chromatin mapping assays are highly enabling to advance epigenetic studies

Compared to ChIP-seq, CUT&RUN and CUT&Tag generate data with >100-fold fewer cells and >10-fold reduced sequencing depth, with >3-fold cost improvement per reaction

- CUT&RUN and CUT&Tag are next-generation immunotethering-based mapping approaches
- Nuclei extraction and genomic mapping can be performed across separate days and worksites

Defined nucleosome spike-in controls are necessary for reliable epigenomic mapping and data analysis

SNAP-CUTANA™ nucleosome spike-in controls are specifically designed for compatibility with CUTANA™ CUT&RUN and CUT&Tag assays, and are essential reagents to enable:

1. Antibody validation
2. Technical monitoring
3. Quantitative normalization

>70% of histone PTM antibodies are unfit for genomic mapping due to poor specificity and/or low efficiency
search your antibody of interest at www.chromatinantibodies.com

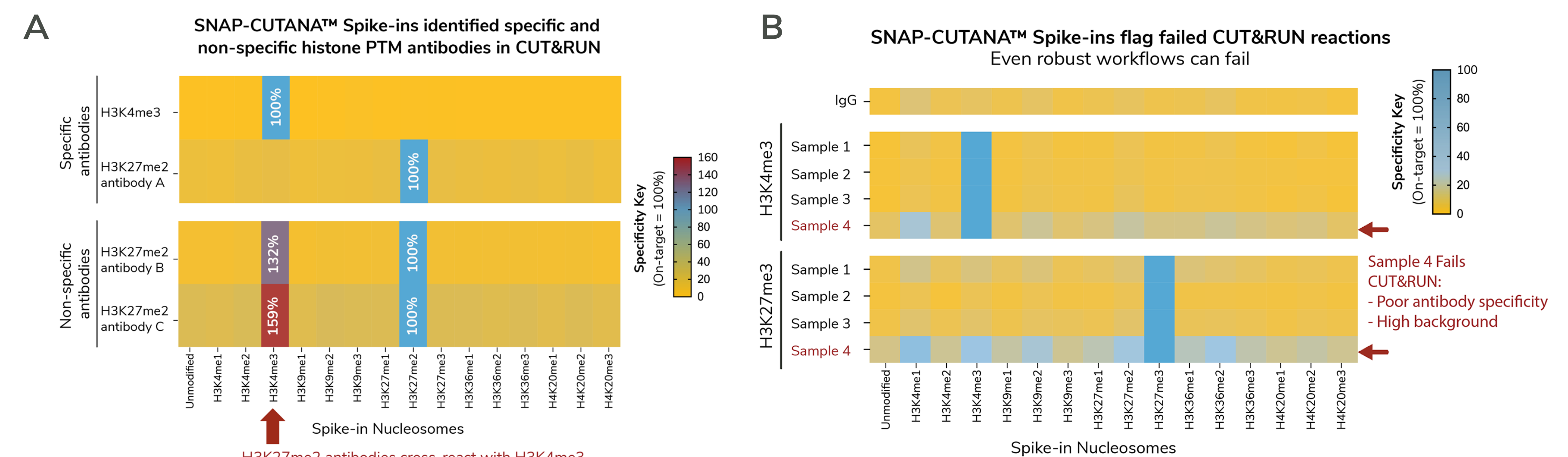


Figure 2: SNAP-CUTANA™ spike-in control heatmaps highlighting applications to (A) validate antibody specificity in-assay, and (B) monitor technical performance for sample quality control.

Automated CUTANA™ assays enable genomic mapping at scale

Platform comparison	ChIP-seq	CUTANA™ CUT&RUN and CUT&Tag	Implications for automation
Sample input	Fragmented chromatin	Intact cells or nuclei	Streamlined workflow
Required cells	>1 million	500k - 5k	More data from precious samples
Defined controls	Uncommon	SNAP-CUTANA spike-ins	Standardized protocols
Seq depth (reads)	>30 million (mammalian)	3-5 million (mammalian)	Greater multiplexing
Assay cost (per reaction)	~\$225	~\$72	70% cost savings
Background	High	Low	Better data quality
Experimental throughput	Low	High	Compatible with 96-well plates

Automation halves hands-on time and increases throughput >8-fold

- End-to-end optimization: sample prep to library QC
- Reduced variance and reaction volumes with 96-well liquid handling
- Standardized workflow for native and cross-linked cells and nuclei
- Buffer optimization for improved sample handling
- AutoCUT&RUN assays generate reliable signal down to 5,000 mammalian cells
- AutoCUT&Tag assays in active development

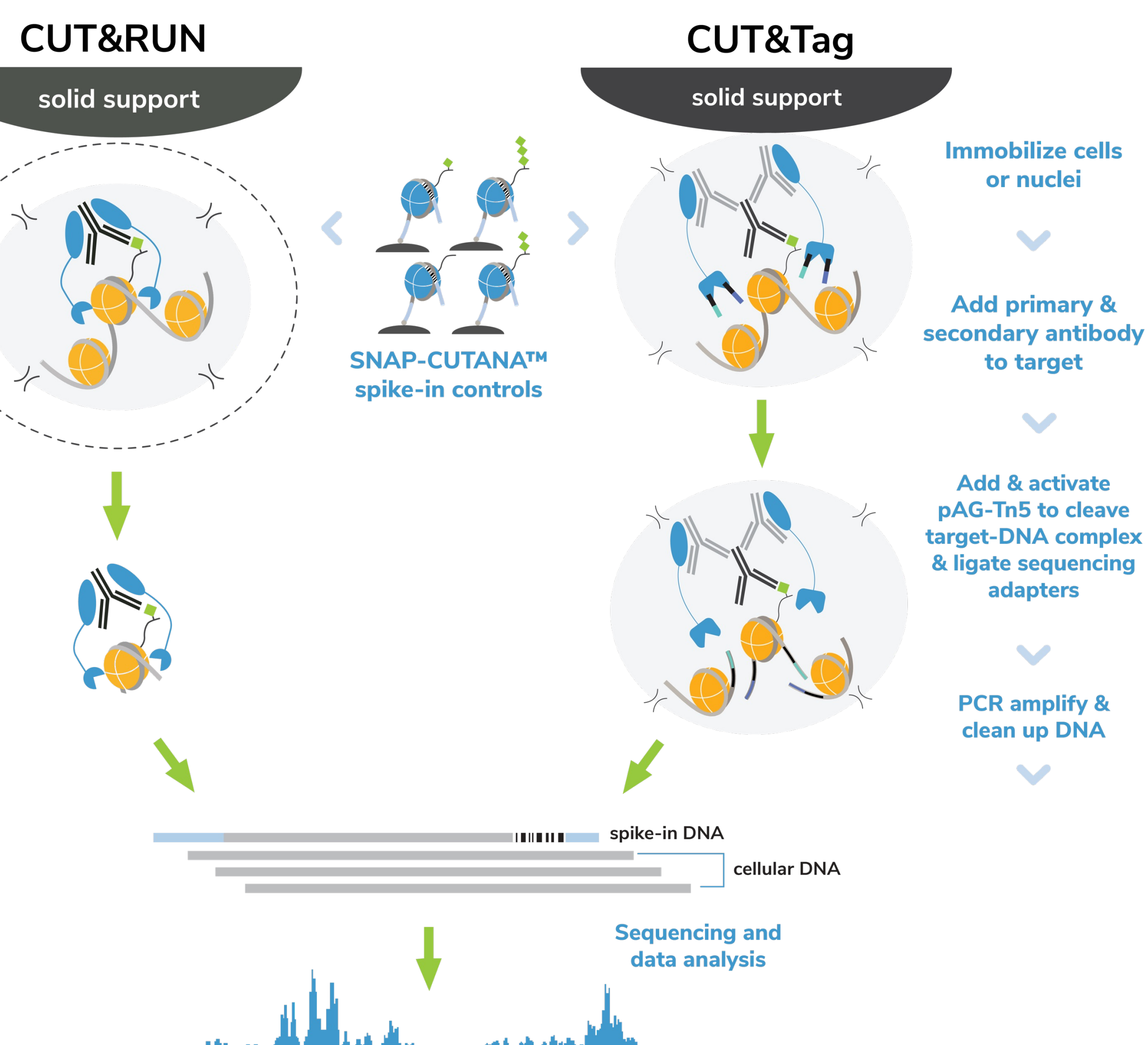


Figure 1: CUTANA™ CUT&RUN and CUT&Tag workflows for chromatin mapping

CUTANA™ CUT&RUN and CUT&Tag assays are enabling for next-generation agricultural research

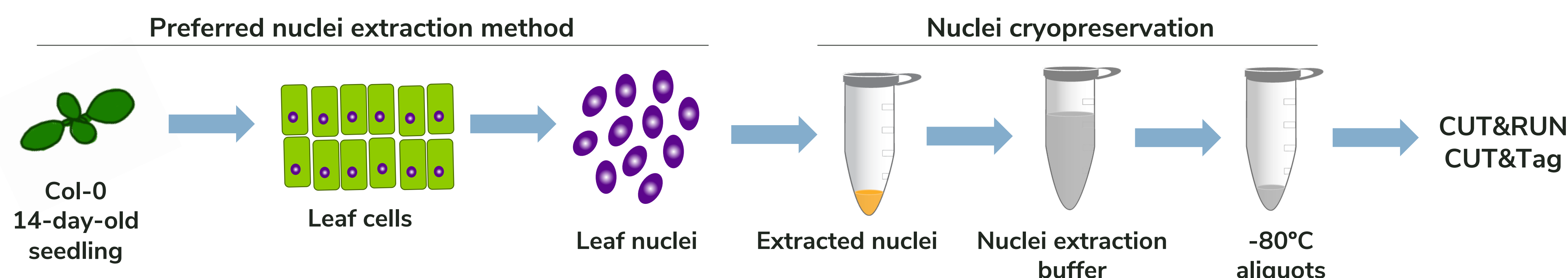


Figure 4: Workflow for sample processing of *Arabidopsis thaliana* nuclei upstream of CUTANA™ CUT&RUN and CUT&Tag assays. Cryopreservation in EpiCypher nuclei extraction buffer (Cat. No. 21-1026) is critical to preserve plant nuclei integrity and enables flexibility to perform the assay across multiple days and/or collaborative sites.

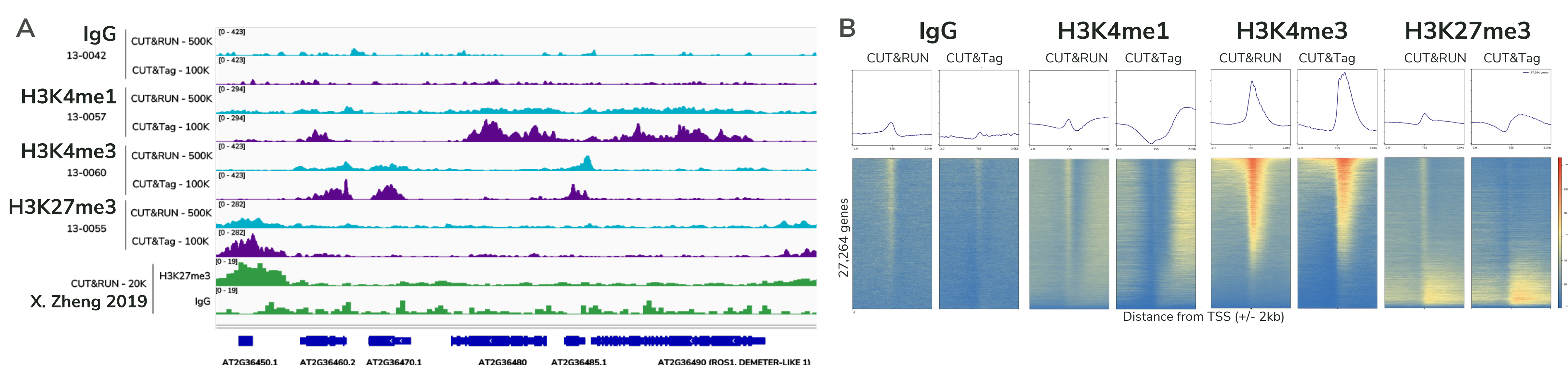


Figure 5: CUTANA™ CUT&RUN and CUT&Tag assays map epigenomic landscape of *Arabidopsis thaliana* seedlings. (A) IGV tracks of histone PTMs; H3K27me3 aligns with previously published tracks (Zheng, 2019). (B) Heatmaps show PTM signal relative to transcription start sites (TSS), genes ordered by H3K4me3 signal. (C) Metagene plots show mean signal for each PTM relative to annotated genomic features.

- Both CUT&RUN and CUT&Tag are compatible with plant cells
- Preliminary experiments indicate CUT&Tag may generate higher signal-to-noise

