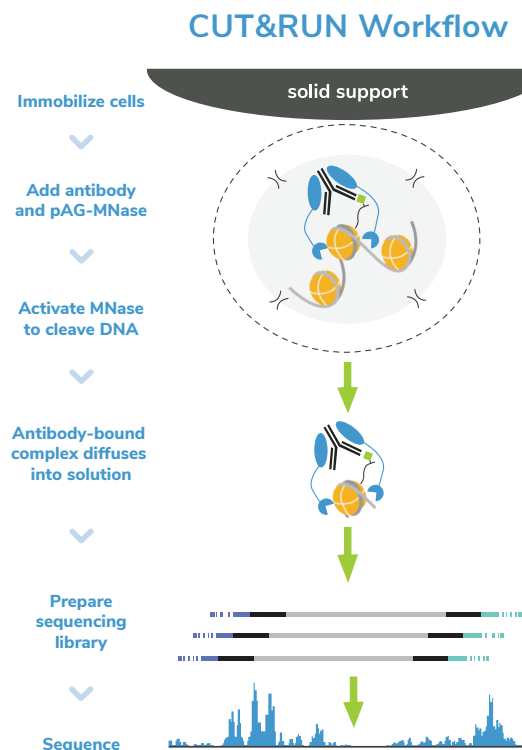


CUTANA™ CUT&RUN Assays for ultra-sensitive genomic mapping

Cleavage Under Targets & Release Using Nuclease (CUT&RUN) is a breakthrough method for genomic mapping of protein-DNA interactions and histone post-translational modifications (PTMs). For new users, our CUTANA™ CUT&RUN platform provides everything you need to get started, including user-friendly kits, protocols, validated antibodies, and more.



Advantages vs. ChIP-seq

- Save 10x in sequencing costs
- Low cell requirements (down to 5k)
- Compatible with diverse targets & sample types
- User-friendly workflow with reproducible results

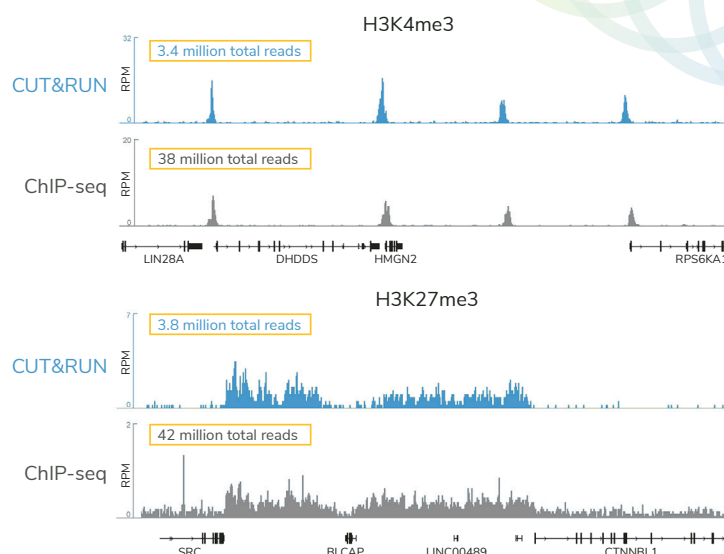


FIGURE 1 CUTANA CUT&RUN outperforms ChIP-seq using fewer sequencing reads.

CUTANA™ CUT&RUN profiles diverse targets

CUT&RUN enables investigation of a wide variety of target classes, including transcription factors, chromatin-interacting proteins, and histone modifications. CUT&RUN also provides access to challenging targets like chromatin remodelers.

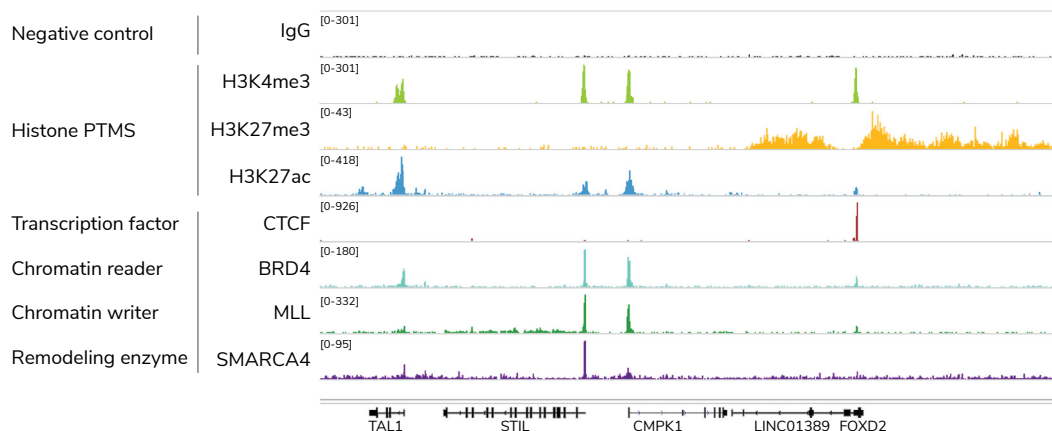
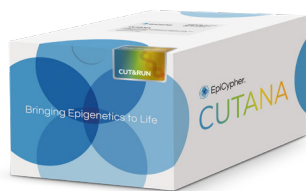


FIGURE 2 Representative genome browser tracks show CUTANA CUT&RUN results using K562 cells. Clear peaks with the expected distribution profile are observed using 3-8 million sequencing reads per sample for a variety of epigenetic targets.

Everything you need to get started!



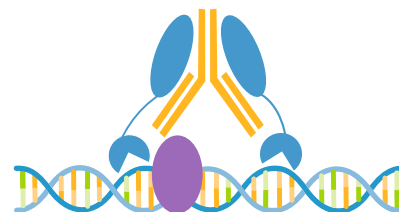
CUTANA CUT&RUN KIT AND LIBRARY PREP KITS

- All-inclusive workflow from cells to sequencing
- Compatible with fresh, frozen, or cross-linked cells or nuclei
- Ideal for setting up CUT&RUN in your lab



CUTANA REAGENTS

- Design and execute custom CUT&RUN experiments
- Include ConA beads, pAG-MNase, DNA purification kit, and spike-in DNA
- Validated using our CUT&RUN protocol

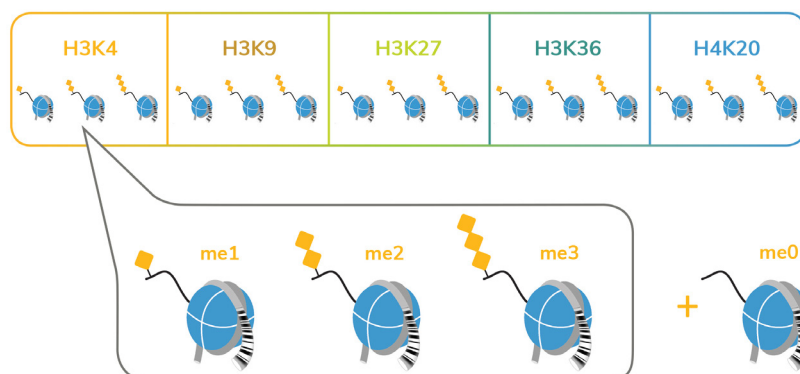


CUTANA CUT&RUN ANTIBODIES

- For histone PTMs, transcription factors, chromatin remodelers, and more
- Lot-validated for robust CUT&RUN mapping assays
- Check our website regularly for new targets

SNAP Spike-ins: Quantitative nucleosome controls for epigenomics

SNAP Spike-ins are the **only** control that provides an accurate readout of CUT&RUN experimental success. Don't question your results - save time and money with SNAP Spike-in Controls.



SNAP Spike-ins are useful for:

- In situ validation of antibody specificity
- Monitoring assay performance
- Quantitative sample comparisons
- Troubleshooting experiments

THE SNAP-CUTANA™ K-METSTAT PANEL comprises 15 DNA-barcoded dNucs carrying disease relevant methyl-lysine PTMs and an unmodified control.

CUTANA™ PRODUCTS AND ORDERING INFO



Let's discuss your project

info@epicypher.com

855.374.2461

epicypher.com

CUT&RUN Kit & Reagents

epicypher.com/cutana-cut-and-run-kit
epicypher.com/cut-and-run-assays

CUTANA CUT&RUN Antibodies

epicypher.com/cut-and-run-antibodies

SNAP-CUTANA Spike-in Controls

epicypher.com/snap-spike-ins