

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

Cleavage Under Targets and Release Using Nuclease (CUT&RUN) is an advanced method for chromatin profiling with improved sensitivity and reduced background compared to ChIP-seq. CUTANA™ CUT&RUN offers optimized kits, expert validated antibodies, and complete workflows for reliable results across target classes.

What can CUT&RUN do for you?

- Reveal chromatin changes driving disease and therapeutic response
- Detect transient and low abundance DNA-binding
- Generate high-quality data from limited primary cells and tissues

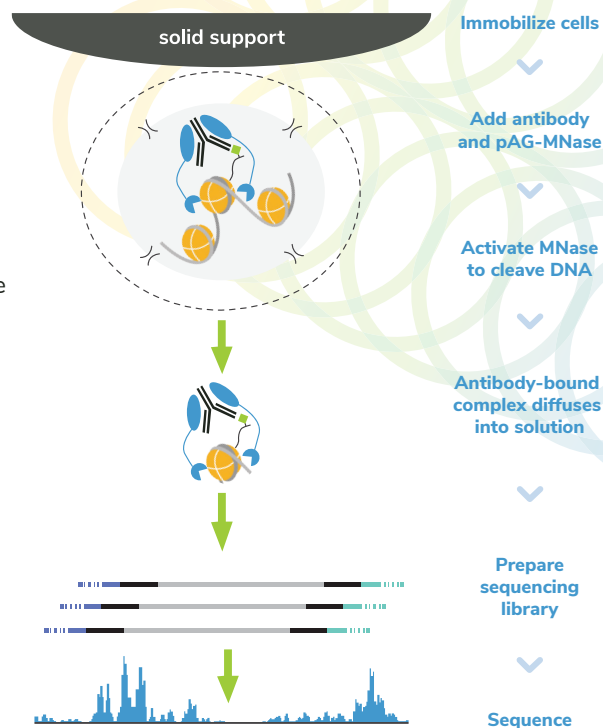
Advantages vs. ChIP-seq

- High signal-to-noise reduces sequencing requirements
- Effective with inputs as low as 5,000 cells
- Native profiling without crosslinking or sonication
- Streamlined workflow with improved reproducibility

For high-quality profiling, choose CUTANA™ assays

FEATURES	ChIP-seq	CUTANA™ CUT&RUN	CUTANA™ CUT&Tag
Cells required	>1 Million	5,000 - 500,000	1* - 100,000
Compatible targets	Histone PTMs, TFs	Histone PTMs, TFs, and chromatin remodelers	Histone PTMs
Sequencing depth (reads)	>30 Million	3-8 Million	5-8 Million
Signal-to-noise	Low	High	High
Assay automation	Difficult	Yes	Yes

*Single-cell CUT&Tag has been described in the literature; see Kaya-Okur et al. Nat Commun. 10 (2019) and others.



CUTANA™ CUT&RUN profiles diverse targets

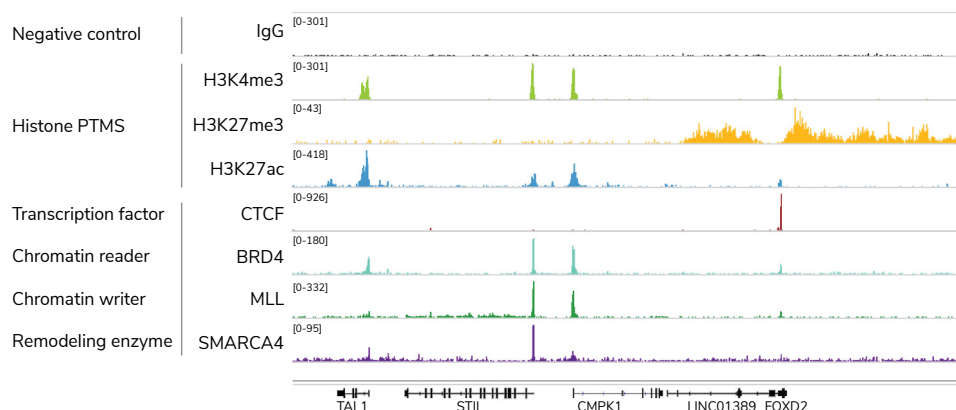


FIGURE 1 CUT&RUN supports native chromatin profiling across transcription factors, chromatin remodelers, and histone modifications, enabling detection of low abundance and transient targets without crosslinking.

Everything you need to get started



CUT&RUN KITS & LIBRARY PREP KITS

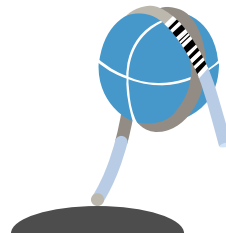
- All-inclusive kits with built in controls and validated protocols
- Compatible with nuclei, cells, and tissue samples
- Pair both kits for a seamless end-to-end workflow

CUTANA CLOUD DATA ANALYSIS

- Biologist-friendly no-code platform with rigorously optimized pipelines
- Transparent pay-as-you-go pricing, with no platform license required
- Secure cloud infrastructure powered by DNAnexus

SERVICES

- Scalable support for biomarker discovery and translational studies
- Fully managed start-to-finish studies led by CUT&RUN experts
- Standardized automated assays
- Custom data reporting



CUT&RUN ANTIBODIES

- Over 1,000 antibodies screened to identify top performers
- Lot-validated directly in CUT&RUN for reliable enrichment
- Available for histone PTMs, transcription factors, remodelers, and more

SPIKE-IN CONTROLS

- SNAP-CUTANA controls confirm antibody specificity & performance
- Epitope-tagged standards troubleshoot antibody and expression failures
- *E. coli* spike-in supports cross-sample normalization

REAGENTS & ACCESSORIES

- Build custom workflows with pAG-MNase and core reagents
- Validated accessories including ConA beads & SPRI reagent
- Precision-engineered magnetic racks for consistent sample recovery

CUTANA™ PRODUCTS AND ORDERING INFO



Let's discuss your project!

info@epicypher.com
855.374.2461
epicypher.com

CUT&RUN Kits, Reagents & Services

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

CUTANA CUT&RUN Antibodies

epicypher.com/cut-and-run-antibodies

SNAP-CUTANA Spike-in Controls

epicypher.com/snap-spike-ins