

CUTANA<sup>™</sup> CUT&RUN Services For scaled epigenomic analyses

#### Chromatin mapping made simple

## The challenge

Biomedical and pharmaceutical researchers are increasingly turning to epigenomic analyses for identification of new cell types, biomarkers, and drug targets. Mapping the genomic enrichment of transcription factors, chromatin modifying enzymes, and histone post-translational modifications (PTMs) is essential to drive these efforts.

To fully achieve these goals, scientists need access to genomics expertise and supporting infrastructure, including customized high-throughput workflows and bioinformatics pipelines.

### Our solution

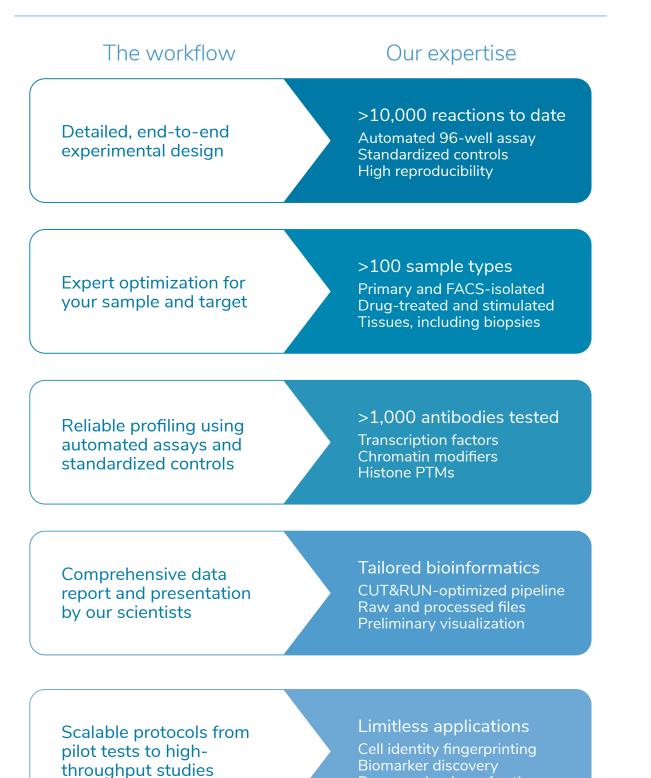
EpiCypher is pioneering the development of CUT&RUN chromatin mapping assays, which provide massive improvements in throughput, data quality, and sensitivity compared to ChIP-based methods.

To meet the needs of our research partners, we developed automated CUTANA CUT&RUN Services. Our services offer exclusive access to EpiCypher's genomics experts, guaranteeing strong support from experimental design to data analysis. This turnkey solution delivers rapid, high-resolution chromatin mapping at scale – unlocking the epigenome to advance biological discovery.



The perfect partnership: your project & our experts

At a glance

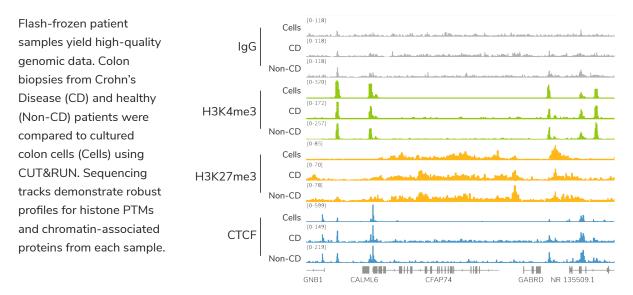


Drug mechanism of action

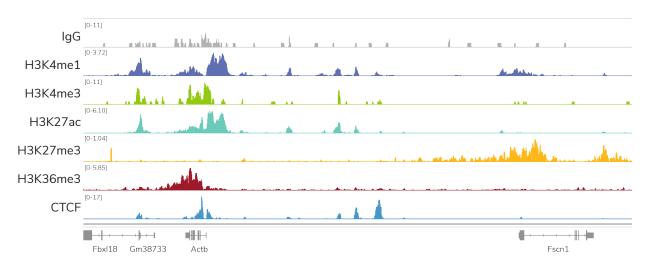
#### Validated applications

Our team provides detailed, end-to-end experimental design and expert optimization for targets and cell types, including high-value clinical samples.

## High-resolution chromatin profiling from biopsies



### Map diverse targets to illuminate the chromatin landscape

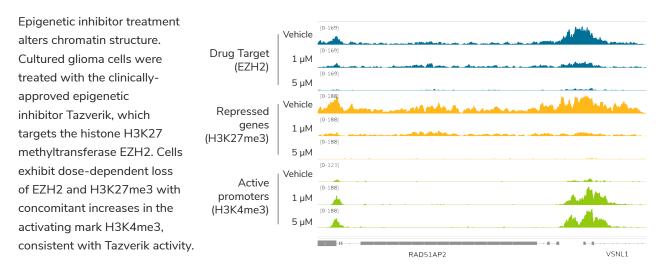


High-resolution chromatin profiling using 10,000 FACS-isolated innate lymphoid cells identifies unique genomic compartments, including active regulatory elements (H3K4me1, H3K27ac), promoters (H3K4me3), gene bodies (H3K36me3), repressed genes (H3K27me3), and transcription factor binding sites (CTCF).

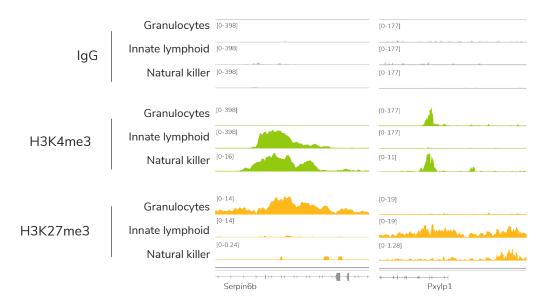
#### Validated applications

Our scaled workflows have limitless applications, ranging from pilot studies aimed at defining gene regulatory mechanisms to large-scale drug screening.

## Uncover drug mechanism of action



## Reveal cell-specific epigenomic profiles



Investigate unique cell types and uncover novel regulatory pathways using CUTANA CUT&RUN Services. Here, FACS-isolated immune cells from a multi-site consortium reveal distinct H3K4me3 (promoters) and H3K27me3 (repressed genes) profiles across granulocytes, innate lymphoid cells, and natural killer cells.

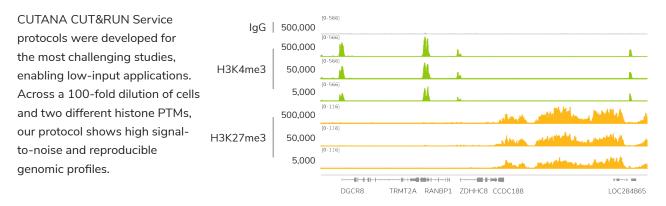
#### Our approach

Our automated workflows include multiple quality control steps. Together we review tailored bioinformatic analyses and quality control metrics to ensure you receive data you can trust.

## Standardized controls and stringent quality metrics

Designated controls create checkpoints for ongoing project risk evaluation. Sample preparation: Viability staining to assess cell and nuclei integrity Spike-in controls: Proprietary approach to confirm experimental success Pre-sequencing: DNA yield, sequencing library size and concentration Post-sequencing: Sequencing and control metrics, genomic enrichment

### Robust, proven protocols



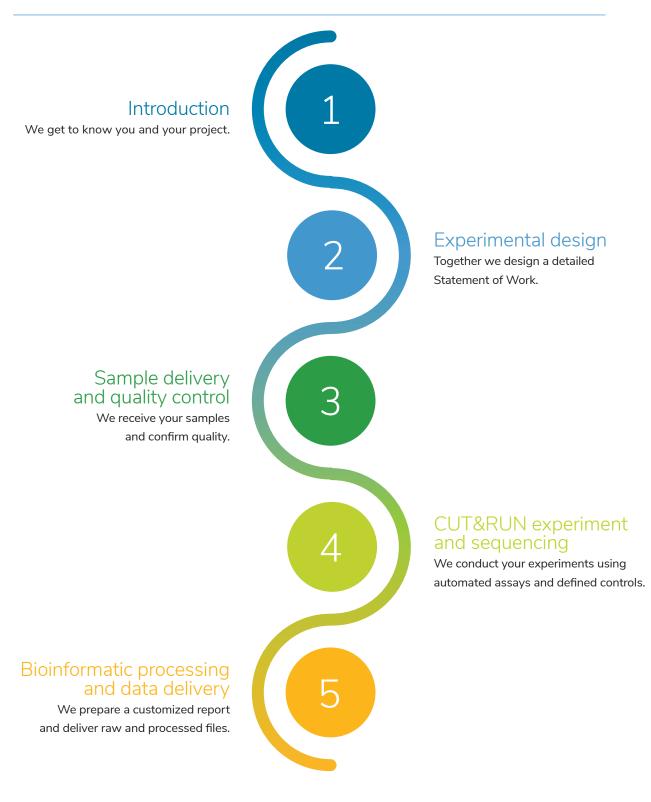
## Expert bioinformatic processing

Our bioinformaticians developed a CUT&RUN-optimized pipeline that we use to analyze sequencing data. You receive raw and processed files that can immediately be used to perform downstream analyses and derive biological insights in your lab.

To see a sample data report, email services@epicypher.com



#### Project roadmap



Ready to get started? Email services@epicypher.com