

## H3K36me3 Antibody, SNAP-Certified™ for CUT&RUN

<b>Catalog No</b>	13-0058	<b>Type</b>	Monoclonal [2091-1E2]
<b>Lot No</b>	23072002-89	<b>Host</b>	Rabbit
<b>Pack Size</b>	100 µg	<b>Concentration</b>	0.5 mg/mL
<b>Applications</b>	CUT&RUN	<b>Reactivity</b>	Human, Wide Range (Predicted)

### DESCRIPTION

This H3K36me3 (histone H3 lysine 36 trimethyl) antibody meets EpiCypher's lot-specific SNAP-Certified™ criteria for specificity and efficient target enrichment in CUT&RUN. This requires <20% cross-reactivity to related histone PTMs determined using the SNAP-CUTANA™ K-MetStat Panel of spike-in controls (EpiCypher 19-1002, **Figure 1**). High target efficiency is confirmed by consistent genomic enrichment at 500k and 50k starting cells (**Figures 2-3**). This antibody targets histone H3 trimethylated at lysine 36, which is enriched in promoters and gene bodies of active genes [1].

### TECHNICAL INFORMATION

<b>Immunogen</b>	A synthetic peptide corresponding to histone H3 trimethylated at lysine 36
<b>Storage</b>	Stable for 1 year at 4°C from date of receipt
<b>Formulation</b>	Protein A affinity-purified recombinant monoclonal antibody in borate buffered saline pH 8.0, 0.09% sodium azide
<b>Target Size</b>	15 kDa

### RECOMMENDED DILUTION

**CUT&RUN:** 0.5 µg per reaction

### GENE & PROTEIN INFORMATION

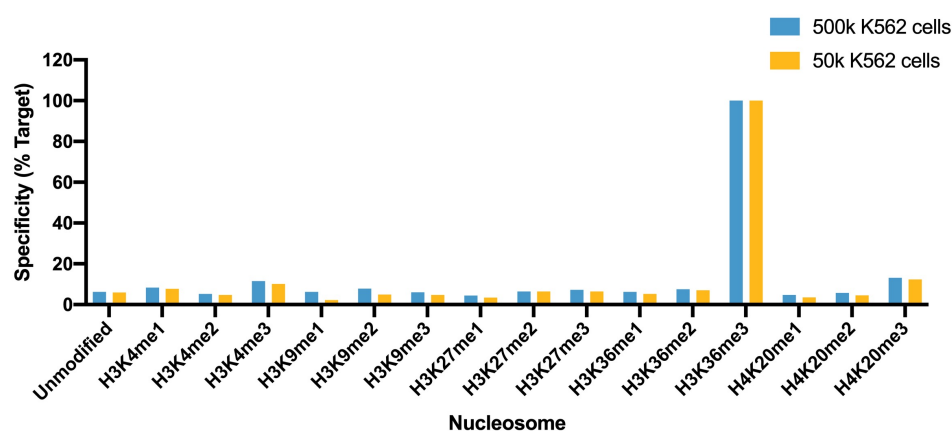
**UniProt ID** H3.1 - P68431

### REFERENCES

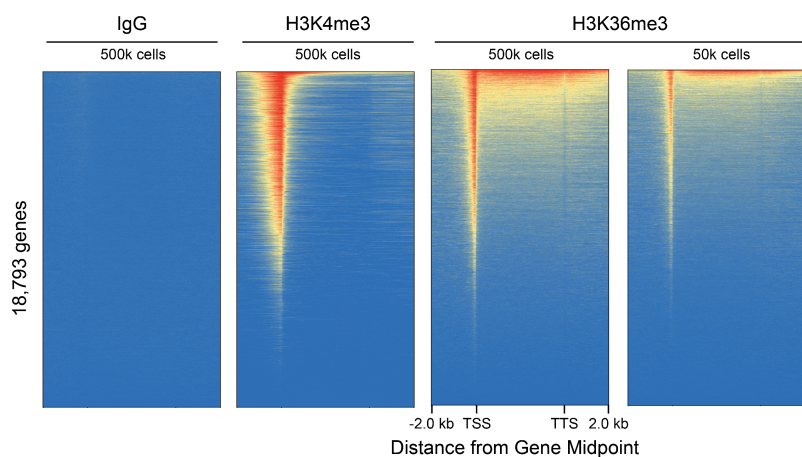
[1] Zhang et al. Nature Communications (2022). PMID: 35680905

**CUT&RUN Methods**

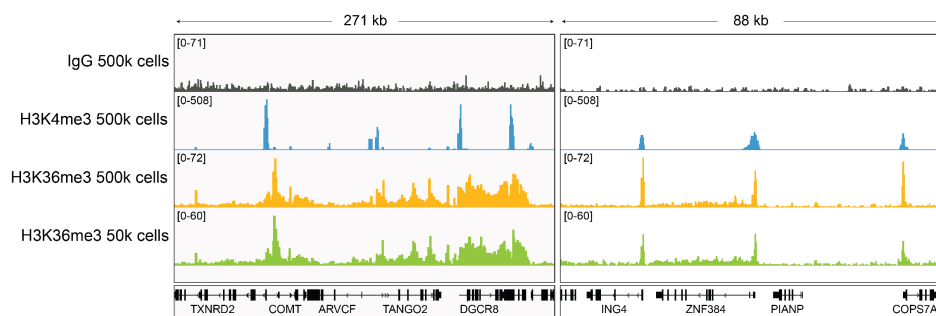
CUT&RUN was performed on 500k and 50k K562 cells with the SNAP-CUTANA™ K-MetStat Panel (EpiCypher 19-1002) spiked-in prior to the addition of 0.5 µg of either IgG negative control (EpiCypher 13-0042), H3K4me3 positive control (EpiCypher 13-0041), or H3K36me3 antibodies. The experiment was performed using the CUTANA™ ChIC/CUT&RUN Kit v3.0 (EpiCypher 14-1048). Library preparation was performed with 5 ng of CUT&RUN enriched DNA (or the total amount recovered if less than 5 ng) using the CUTANA™ CUT&RUN Library Prep Kit (EpiCypher 14-1001/14-1002). Both kit protocols were adapted for high throughput Tecan liquid handling. Libraries were run on an Illumina NextSeq2000 with paired-end sequencing (2x50 bp). Sample sequencing depth was 16.8 million reads (IgG 500k cell input), 14.4 million reads (H3K4me3 500k cell input), 24.4 million reads (H3K36me3 500k cell input) and 16.4 million reads (H3K36me3 50k cell input). Data were aligned to the hg19 genome using Bowtie2. Data were filtered to remove duplicates, multi-aligned reads, and ENCODE DAC Exclusion List regions.



**FIGURE 1 SNAP specificity analysis in CUT&RUN.** CUT&RUN was performed as described above. CUT&RUN sequencing reads were aligned to the unique DNA barcodes corresponding to each nucleosome in the K-MetStat panel (x-axis). Data are expressed as a percent relative to on-target recovery (H3K36me3 set to 100%).



**FIGURE 2 CUT&RUN genome-wide enrichment.** CUT&RUN was performed as described above. Sequence reads were aligned to 18,793 annotated transcription start sites (TSSs ± 2 kbp). Signal enrichment was sorted from highest to lowest (top to bottom) relative to the H3K36me3 - 500k cells sample (all gene rows aligned). High, medium, and low intensity are shown in red, yellow, and blue, respectively. H3K4me3 positive control and H3K36me3 antibodies produced the expected enrichment pattern, which was consistent between 500k and 50k cells and greater than the IgG negative control.



**FIGURE 3 H3K36me3 CUT&RUN representative browser tracks.** CUT&RUN was performed as described above. Gene browser shots were generated using the Integrative Genomics Viewer (IGV, Broad Institute). H3K36me3 antibody tracks display the characteristic enrichment known to be consistent with the function of this PTM [1]. Similar results in peak structure and location were observed for both 500k and 50k cell inputs.