

CUTANA™ ChIC/CUT&RUN Kit

Catalog No	14-1048-24	Pack Size	24 Reactions
Lot No	24326001-81	Kit Version	v5

DESCRIPTION

The CUTANA™ ChIC/CUT&RUN Kit enables streamlined chromatin profiling of histone post-translational modifications (PTMs) and chromatin associated proteins. The CUT&RUN Kit Version 5 (v5) now includes an additional control antibody (H3K27me3). Positive (H3K4me3 and H3K27me3) and negative (IgG) control antibodies pair with SNAP-CUTANA™ spike-in controls for assay optimization and continuous assay monitoring (**Figure 2**). *E. coli* DNA is included for data normalization. SPRI magnetic beads are used for DNA purification, enabling seamless multi-channel pipetting throughout the workflow to maximize throughput and reproducibility. The kit is compatible with a variety of inputs including cells or nuclei derived from native, cryopreserved, or cross-linked samples. While it is recommended to start with 500,000 cells, comparable data can be generated using as few as 5,000 cells. The inclusion of controls, as well as compatibility with diverse target types, sample inputs, and low cell numbers, make this kit the go-to solution for chromatin mapping experiments.

KIT CONTENTS

<u>Item</u>	<u>CAT</u>	<u>Item</u>	<u>CAT</u>
8-strip Tubes	10-0009-01	Stop Buffer	21-1003-01
0.5 M EDTA	21-1006-01	5% Digitonin	21-1004-01
100 mM Calcium Chloride	21-1007-01	1 M Spermidine	21-1005-01
SPRIselect Reagent from Beckman Coulter, Inc. ¹	21-1405-01	SNAP-CUTANA K-MetStat Panel	19-1002-02
0.1X TE Buffer	21-1025-01	H3K4me3 Positive Control Antibody	13-0060-02
ConA Beads	21-1401-01	H3K27me3 Positive Control Antibody	13-0055-02
<i>E. coli</i> Spike-in DNA	18-1401-01	Rabbit IgG Negative Control Antibody	13-0042-02
Bead Activation Buffer	21-1001-01	pAG-MNase	15-1016-01
Pre-Wash Buffer	21-1002-01		

TECHNICAL INFORMATION

Storage	OPEN KIT IMMEDIATELY and store components at room temperature, 4°C, and -20°C as indicated (see User Manual corresponding to Kit Version 5). Stable for 12 months upon date of receipt.
Instructions for Use	See User Manual corresponding to Kit Version 5. This kit is not compatible with previous user manuals.

VALIDATION DATA

CUT&RUN Methods	CUT&RUN was performed using the CUTANA™ ChIC/CUT&RUN Kit starting with 500k K562 cells with 0.5 µg of IgG (EpiCypher 13-0042), H3K4me3 (EpiCypher 13-0060), H3K27me3 (EpiCypher 13-0055), or 0.125 µg of CTCF (EpiCypher 13-2014) antibodies in duplicate. Library preparation was performed with 5 ng of DNA (or the total amount recovered if less than 5 ng) using the CUTANA™ CUT&RUN Library Prep Kit (EpiCypher 14-1001/14-1002). Libraries were run on an Illumina NextSeq2000 with paired-end sequencing (2x50 bp). Sample sequencing depth was 5.5/18.8 million reads (IgG Rep 1/Rep 2), 14.2/17.0 million reads (H3K4me3 Rep 1/Rep 2), 24.7/18.1 million reads (H3K27me3 Rep 1/Rep 2), and 8.6/5.5 million reads (CTCF Rep 1/Rep 2). Data were aligned to the T2T-CHM13v2.0 genome using Bowtie2. Data were filtered to remove duplicates, multi-aligned reads, and ENCODE DAC Exclusion List regions.
----------------------------	--

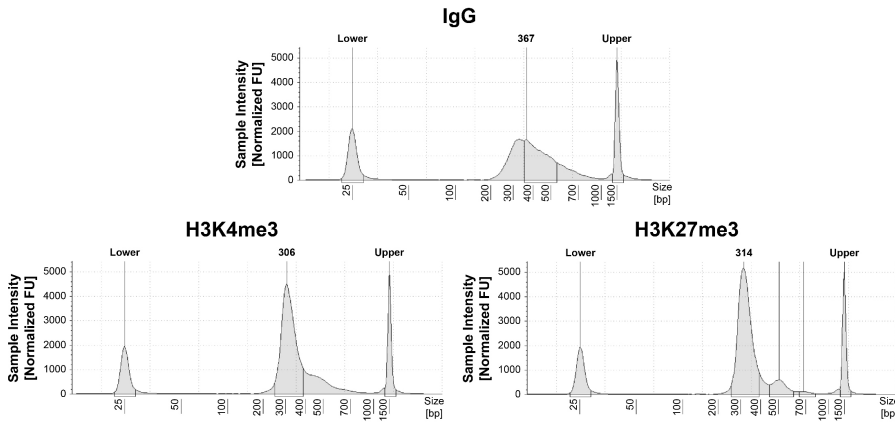


FIGURE 1 CUT&RUN DNA fragment size distribution analysis. CUT&RUN was performed as described above. Library DNA was analyzed by Agilent TapeStation®. This analysis confirmed that mononucleosomes were predominantly enriched in CUT&RUN (~300 bp peaks represent 150 bp nucleosomes + sequencing adapters).

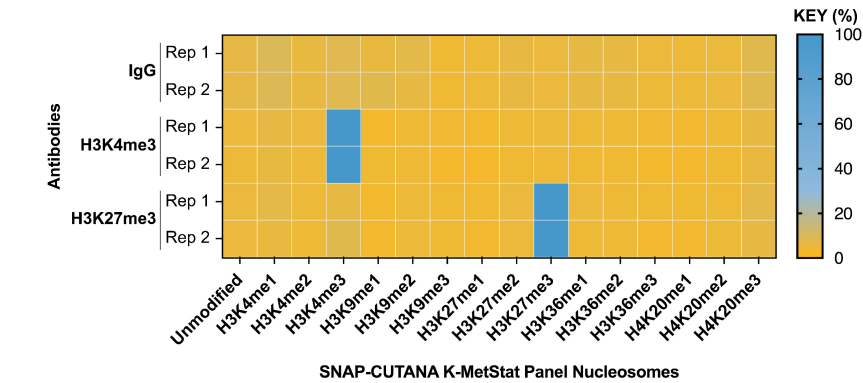


FIGURE 2 SNAP-CUTANA™ Spike-in controls. DNA-barcoded designer nucleosomes (dNucs) harboring distinct K-methyl PTMs were spiked into CUT&RUN reactions prior to antibody addition. Spike-in barcodes were analyzed using the shell script at epicypher.com/19-1002. Barcodes for IgG (normalized to total reads), H3K4me3 and H3K27me3 (normalized to on-target) antibodies are shown. The spike-ins confirmed H3K4me3 and H3K27me3 antibodies specifically recovered the target dNucs, while IgG showed no preferential enrichment.

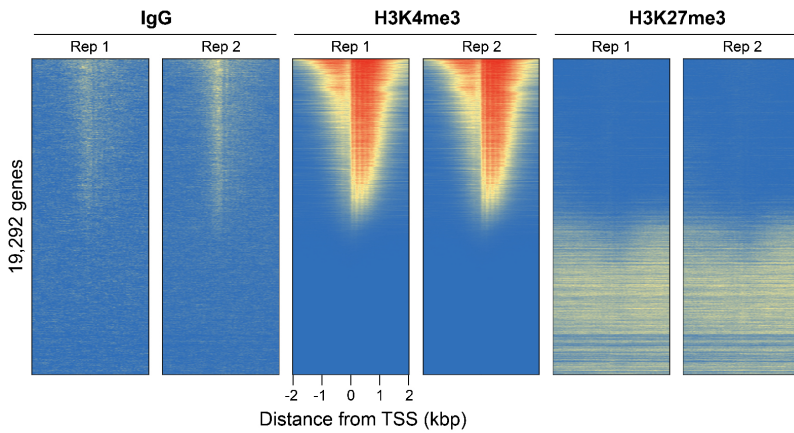


FIGURE 3 CUT&RUN genome-wide heatmaps. CUT&RUN was performed as described above. Heatmaps show two replicates (“Rep”) of IgG, H3K4me3, and H3K27me3 kit control antibodies in aligned rows ranked by intensity (top to bottom) relative to the H3K4me3 Rep 1 reaction and colored such that red indicates high localized enrichment and blue denotes background signal.

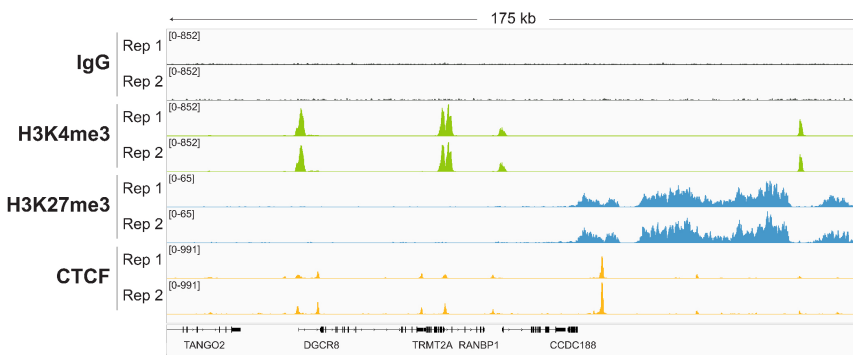


FIGURE 4 Representative gene browser tracks. CUT&RUN was performed as described above. A representative 175 kb window at the TRMT2A gene is shown for two replicates (“Rep”) of IgG, H3K4me3, and H3K27me3 kit control antibodies. Representative tracks are also shown for the transcription factor CTCF. The CUT&RUN kit produced the expected genomic distribution for each target. Images were generated using the Integrative Genomics Viewer (IGV, Broad Institute).

¹ Beckman Coulter, the stylized logo, and SPRIselect are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

US Pat. No. 7790379, 11885814 and related patents and pending applications