

Rabbit IgG Antibody, CUTANA™ CUT&RUN Negative Control

Catalog No. 13-0042
Lot No. 20036001-52
Pack Size 100 µg

Type Isotype control

Target Size N/A

Reactivity Negative Control

Host Rabbit

Format Aff. Pur. IgG

Appl. CUT&RUN

Product Description:

This Rabbit IgG Antibody can be used to control for non-specific background signal in CUT&RUN and other applications. A positive control antibody is also available (H3K4me3, EpiCypher Catalog No. 13-0041).

Immunogen:

None

Formulation:

Affinity-purified antibody (11.5 mg/mL) in PBS, pH 7.6

Storage and Stability:

Stable for 1 year at 4°C from date of receipt

Application Notes:

Prior to use, prepare a 1:10 working stock dilution in CUT&RUN Antibody Buffer (20 mM HEPES pH 7.5, 150 mM NaCl, 0.5 mM Spermidine, 0.01% Digitonin, 2 mM EDTA). Store at 4°C for 1 month. For each CUT&RUN assay, dilute 1:100 from the working stock. A typical yield for a CUT&RUN assay using 500,000 K562 cells is 2-5 ng DNA.

References:

Applications Key: ChIP: Chromatin immunoprecipitation; ChIP-seq: ChIP-sequencing; E: ELISA; FACS: Flow cytometry; IF: Immunofluorescence; IHC: Immunohistochemistry; ICC: Immunocytochemistry; IP: Immunoprecipitation; WB: Western Blotting; L: Luminex
Reactivity Key: B: Bovine; Ce: *C. elegans*; Ch: Chicken; Dm: *Drosophila*; Eu: Eukaryote; H: Human; M: Mouse; Ma: Mammal; R: Rat; Sc: *S. cerevisiae*; Sp: *S. pombe*; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish



EpiCypher®

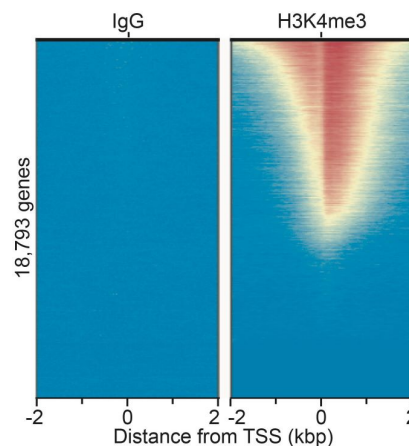


Figure 1. CUT&RUN data was generated using the EpiCypher CUTANA™ CUT&RUN Protocol (EpiCypher.com/cutana-protocol). Shown is a heatmap for Rabbit IgG CUT&RUN negative control antibody (left) compared to H3K4me3 CUT&RUN positive control antibody (right, EpiCypher Catalog No. 13-0041). Signal from 2.2 and 3.9 million paired-end reads, respectively, are aligned to the transcription start site (TSS, +/- 2kb) of 18,793 genes. High and low signal are ranked by intensity (top to bottom) and reflected by red and blue colors, respectively.

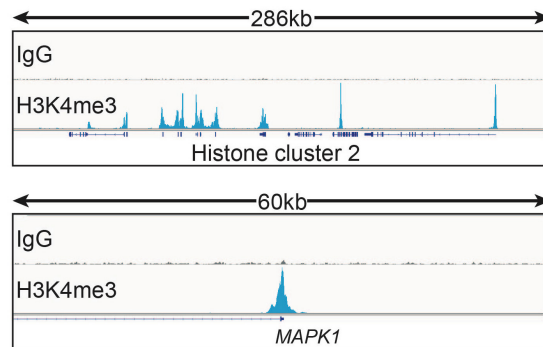


Figure 2. Rabbit IgG (top track) and H3K4me3 (bottom track) CUT&RUN data was generated as described above. Two representative loci are shown in the Integrative Genomics Viewer (IGV, Broad Institute). Window size is denoted at the top of each panel.