

Anti-Rabbit Secondary Antibody for CUTANA™ CUT&Tag

Catalog No	13-1047	Type	Mixed Monoclonal*
Lot No	22250005-81	Host	Goat
Pack Size	250 Reactions	Concentration	1 mg/mL
Applications	CUT&Tag	Reactivity	Anti-Rabbit

DESCRIPTION

Anti-Rabbit Secondary Antibody for CUTANA™ CUT&Tag is affinity purified and specific for rabbit immunoglobulins. Minimal cross reactivity with mouse, rat, human, bovine, guinea pig, and donkey IgG is observed. Use of a secondary antibody in CUT&Tag aids in signal amplification by increasing bound IgG per target epitope [1]. For best results, use with CUTANA™ pAG-Tn5 (EpiCypher 15-1017 & 15-1117). For detailed instructions for use, see EpiCypher's CUTANA™ CUT&Tag protocol: epicypher.com/protocols.

*Mixed Monoclonal: a pool of multiple monoclonal antibodies.

TECHNICAL INFORMATION

Immunogen	Recombinant full-length rabbit immunoglobulin protein
Storage	Store at 4°C short term. For long term storage, store at -20°C. Avoid freeze/thaw cycles.
Formulation	Protein G affinity-purified antibody in PBS pH 7.2, 0.09% sodium azide

RECOMMENDED DILUTION

CUT&Tag	0.5 µg per reaction
--------------------	---------------------

REFERENCES

[1] Kaya-Okur et al. Nat. Commun. (2019). PMID: 31036827

VALIDATION DATA

CUT&Tag Methods

CUT&Tag was performed on 100k K562 nuclei with 0.5 µg of either IgG (EpiCypher 13-0042), H3K4me3 (EpiCypher 13-0041), or H3K27me3 (ThermoFisher MA5-11198) antibodies, followed by 0.5 µg of anti-rabbit secondary antibody for CUTANA™ CUT&Tag. The EpiCypher Direct-to-PCR CUT&Tag protocol (epicypher.com/protocols) was used. Libraries were run on an Illumina NextSeq2000 with paired-end sequencing (2x50 bp). Sample sequencing depth was 8.5 million reads (IgG), 6.0 million reads (H3K4me3), and 12.5 million reads (H3K27me3). Data were aligned to the hg19 genome using Bowtie2. Data were filtered to remove duplicates, multi-aligned reads, and ENCODE DAC Exclusion List regions.

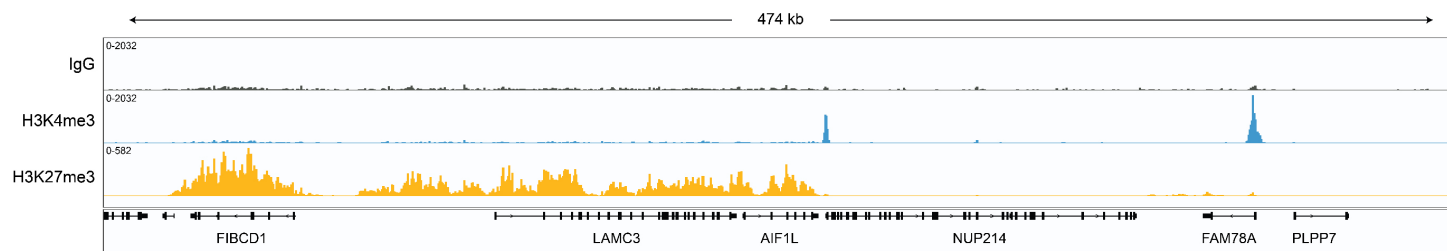


FIGURE 1 CUT&Tag data. CUT&Tag was performed as described above. Gene browser shots generated using the Integrative Genomics Viewer (IGV, Broad Institute) show a representative 474 kb window centered at the *LAMC3* gene. The genomic distribution pattern was consistent with that expected for H3K4me3 and H3K27me3.