

Anti-Mouse Secondary Antibody for CUTANA™ CUT&Tag



EpiCypher®

Catalog No. 13-0048

Lot No. 20302002-10

Pack Size 50 reactions

Type Recombinant Polyclonal

Host Goat

Reactivity Anti-Mouse

Mol. Wgt. N/A

Format Aff. Pur. IgG

Applications CUT&Tag

Product Description:

Anti-Mouse Secondary Antibody for CUTANA™ CUT&Tag is affinity-purified and specific for mouse immunoglobulins. Minimal cross reactivity with rabbit, 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. For best results, use with CUTANA™ pAG-Tn5 (EpiCypher Catalog No. 15-1017) and CUTANA Compatible Antibodies.

Immunogen:

Recombinant full-length mouse immunoglobulin protein.

Formulation:

Protein G affinity-purified antibody (1 mg/mL) in PBS, pH 7.2, with 0.09% sodium azide.

Storage and Stability:

Store at 4°C short term. For long term storage, store at -20°C. Avoid freeze/thaw cycles.

Application Notes:

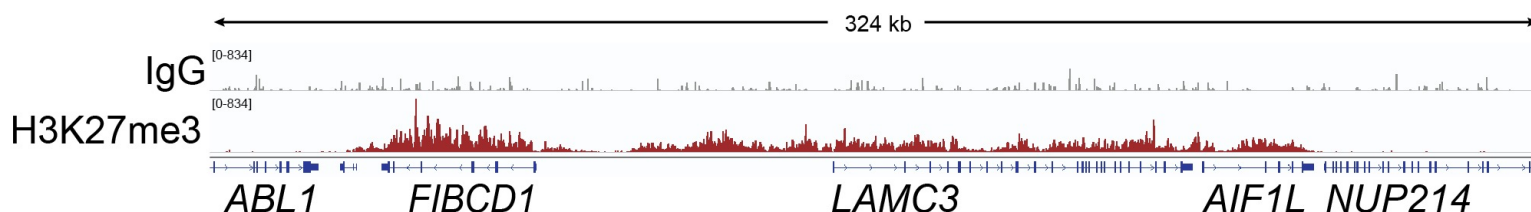
Recommended usage amounts:

CUT&Tag: 0.5 µg per sample

For detailed instructions for use, see EpiCypher CUTANA Direct-to-PCR CUT&Tag protocol:

EpiCypher.com/cutana-protocol

References Using this Product:



CUT&Tag Data: Representative sequencing track obtained using Anti-Mouse Secondary Antibody for CUTANA™ CUT&Tag (0.5 µg) is shown at a 324 kb window centered at the LAMC3 gene. CUT&Tag was performed using the EpiCypher CUTANA Direct-to-PCR CUT&Tag protocol (EpiCypher.com/cutana-protocol) using 100,000 K-562 nuclei with mouse anti-H3K27me3 antibody. Image was generated using the Integrative Genomics Viewer (IGV, Broad Institute). Anti-Mouse Secondary Antibody for CUTANA™ CUT&Tag produced clear peaks with a genomic distribution pattern consistent with H3K27me3.

Applications Key: ChIP: Chromatin Immunoprecipitation; ChIC: Chromatin Immunocleavage; CUT&RUN: Cleavage Under Targets and Release Using Nuclease; CUT&Tag: Cleavage Under Targets and Tagmentation; 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.cerevesiae; Sp: S. pombe; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish

This product is for *in vitro* research use only and is not intended for use in humans or animals.

EpiCypher, Inc • Phone: 855-374-2461 • Fax: 855-420-6111 • www.epicypher.com