

Anti-Rabbit Secondary Antibody for CUTANA™ CUT&Tag



EpiCypher®

Catalog No 13-0047

Lot No 21175003-61

Pack Size 50 Reactions

Type Mixed Monoclonal*

Host Goat

Reactivity Anti-Rabbit

Mol. Wgt. N/A

Format Aff. Pur. IgG

Applications CUT&Tag

Product 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. For best results, use with CUTANA pAG-Tn5 (EpiCypher 15-1017 & 15-1117) and CUTANA™ validated antibodies.

*Mixed Monoclonal: a pool of multiple monoclonal antibodies.

Immunogen:

Recombinant full-length rabbit 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 dilution:

CUT&Tag: 0.5 µg per reaction

For detailed instructions for use, see EpiCypher's CUTANA™ CUT&Tag protocol: epicypher.com/protocols

References:

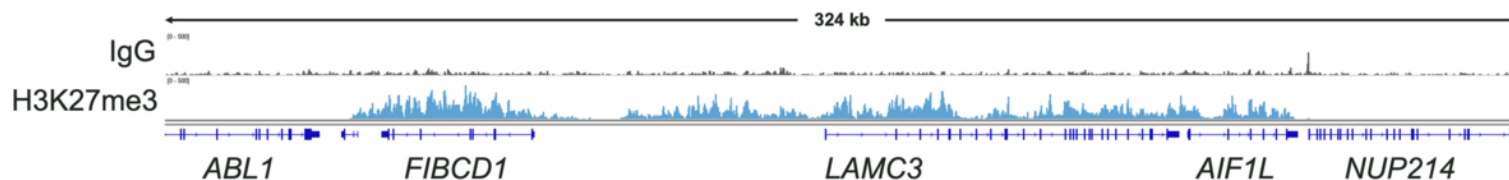


Figure 1: CUT&Tag data. CUT&Tag was performed on 100k K562 nuclei with rabbit anti-H3K27me3 antibody (EpiCypher 13-0030) and 0.5 µg of anti-rabbit secondary antibody for CUTANA™ CUT&Tag using the EpiCypher CUTANA™ Direct-to-PCR CUT&Tag protocol (epicypher.com/protocols). Gene browser shots generated using the Integrative Genomics Viewer (IGV, Broad Institute) show a representative 324 kb window centered at the LAMC3 gene. The genomic distribution pattern was consistent with that of H3K27me3.

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