

Histone H3K36ac Antibody, SNAP-ChIP[®] Certified



EpiCypher[®]

Catalog No 13-0035
Lot No 20336002-40
Pack Size 100 µg

Type Monoclonal **Host** Rabbit
Reactivity Human, Mouse, Wide Range

Target Size 15 kDa **Format** Aff. Pur. IgG
Applications ChIP, WB, ICC, Luminex

Product Description:

This antibody meets EpiCypher's "SNAP-ChIP[®] Certified" criteria for specificity and efficient target enrichment in a ChIP experiment (<20% cross-reactivity across the panel, >5% recovery of target input) based on technology originating from Grzybowski et al. [1] and profiling standards from Shah et al. [2]. This antibody reacts to H3K36ac and exhibits no cross reactivity to other lysine acylations in the EpiCypher SNAP-ChIP K-AcylStat panel (EpiCypher 19-3001) is detected.

Immunogen:

A synthetic peptide corresponding to histone H3 acetylated at lysine 36.

Formulation:

Protein A affinity-purified recombinant monoclonal antibody (1 mg/mL) in PBS, with 0.09% sodium azide, 1% BSA, and 50% glycerol.

Storage and Stability:

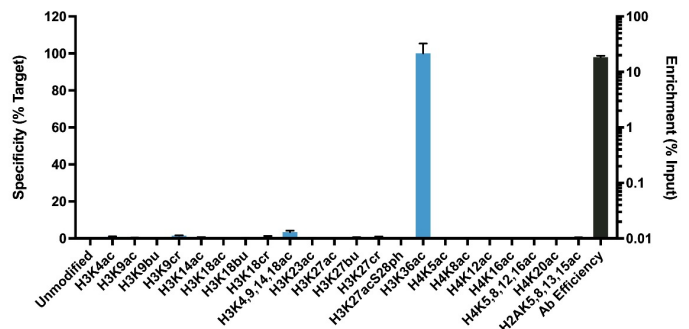
Stable for 1 year at -20°C from date of receipt.

Recommended Dilution:

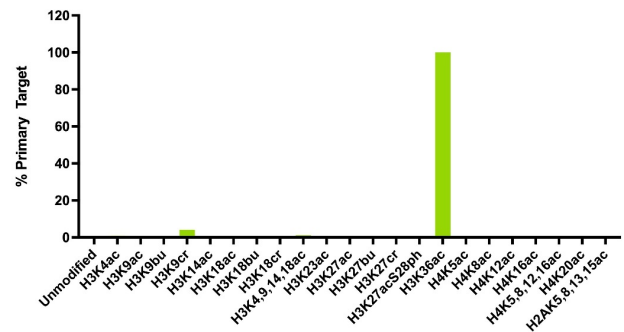
ChIP: 2 - 5 µg per 10⁶ cells **WB:** 0.5 - 2 µg/mL
ICC: 0.5 - 2 µg/mL **Luminex:** 0.25 - 4 µg/mL

References:

- [1] Grzybowski et al (2015) *Mol Cell* 58:886
 [2] Shah et al (2018) *Mol Cell* 72:162

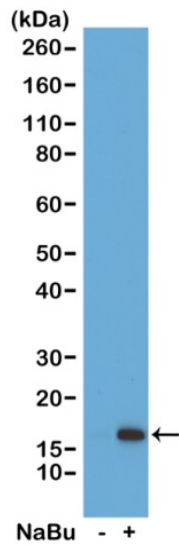


SNAP-ChIP-qPCR: Histone H3K36ac antibody (3 µg) was tested in a native ChIP experiment using chromatin from K-562 cells (3 µg) with the SNAP-ChIP K-AcylStat Panel (EpiCypher 19-3001) spiked-in prior to micrococcal nuclease digestion. Specificity (left y-axis) was determined by qPCR for the DNA barcodes corresponding to modified nucleosomes in the SNAP-ChIP panel (x-axis). Black bar represents antibody efficiency (right y-axis; log scale) and indicates percentage of the target immunoprecipitated relative to input.

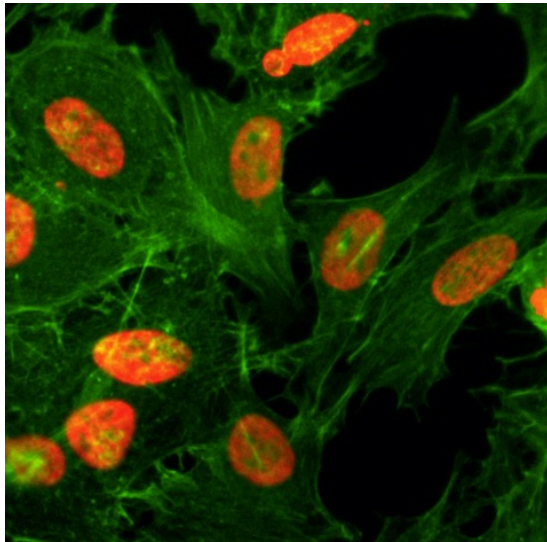


Luminex Data: Histone H3K36ac antibody was assessed using a Luminex[®] based approach employing dCypher[™] Nucleosome K-AcylStat Panel (EpiCypher 16-9003). The panel comprises biotinylated designer nucleosomes (x-axis) individually coupled to uniquely identifiable Luminex MagPlex[®] beads. Antibody binding to nucleosomes was tested in multiplex (24-plex) at a 1:1,000 dilution, and detected with an anti-IgG*PE secondary. Data was generated using a Luminex FlexMAP3D[®] and is shown normalized to on-target signal (H3K36ac; set to 100).

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



Western Blot Data: Western blot of acid extracts from HeLa cells untreated (-) or treated with sodium butyrate (+), using H3K36ac antibody at 1 μ g/mL.



Immunocytochemistry Data: Immunocytochemistry of HeLa cells treated with sodium butyrate, using H3K36ac antibody (red). Actin filaments have been labeled with fluorescein phalloidin (green).

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