

Anti-Acetyl Lysine Mouse MAb

Cat. # AAC01

Lot: 031

Upon arrival, store at 4°C (desiccated)

See datasheet for storage after reconstitution

Form:	Lyophilized powder
Amount of material:	2 x 100 µl when reconstituted
Validated applications:	WB, IF, IP, ChIP
Species reactivity:	All
Host/Isotype:	Mouse/IgG2b
Clone:	3C6.08.20

Background Information

Acetylation of proteins can occur as a co-translational or post-translational modification (PTM) (1). Co-translational acetylation occurs at the N-terminus of approximately 85% of mammalian proteins, is irreversible, and is thought to be important in protein stability, localization, and synthesis (1). Post-translational acetylation occurs on the epsilon amino group of lysine residues as a reversible and highly dynamic PTM that is known to be a key regulator in multiple cellular events, including chromatin structure, transcription, metabolism, signal transduction, and cytoskeletal regulation (2-3). To date, over 4,000 proteins have been identified as targets for PTM acetylation which is comparable to phosphorylation in cellular prevalence (3). Antibody AAC01 detects acetyl lysine PTMs.

Material

Anti-acetyl lysine antibody is a mouse monoclonal antibody. The antibody was raised against a proprietary mixture of acetylated proteins designed to optimize acetyl lysine recognition in a wide range of sequence contexts. The antibody has been shown to recognize a wide range of acetylated proteins (Fig. 1A), including acetylated tubulin, histones, glutamic dehydrogenase, and chemically acetylated bovine serum albumin (BSA) (Fig. 1B-C). The antibody has also been shown to recognize acetylated actins and Arp 2/3 complex proteins (data not shown). AAC01 is purified by Protein G affinity chromatography and is supplied as a lyophilized white powder.

Storage and Reconstitution

Shipped at ambient temperature. The lyophilized protein can be stored desiccated at 4°C for 6 months. For reconstitution, the product tube should be briefly centrifuged to collect the powder to the bottom of the tube.

Reconstitute each tube in 100 µl of 50% glycerol in water and store at -20°C. Final buffer composition is 250 mM PIPES pH 7.0, 1% sucrose, and 0.5% dextran.

When stored and reconstituted as described, the product is stable for 6 months at -20°C. NOTE: We recommend adding an antibacterial such as sodium azide (0.02% final concentration) to prevent bacterial contamination of the antibody stock.

Control Proteins/Lysates

The following controls are recommended for this antibody.

*For further details and recommendations for control protein/lysates and associated products, visit www.cytoskeleton.com.

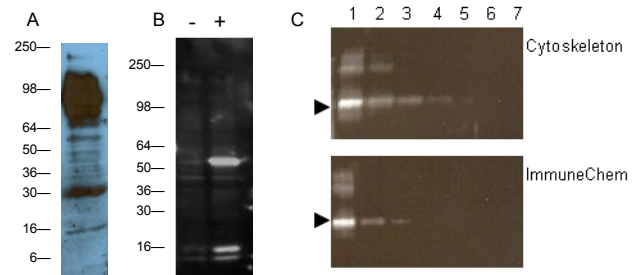
Control Protein/Lysate*	Description	Use	Cat# or Part#
Acetylated BSA	Chemically acetylated BSA	Use at 10-15 µg/ml of Ab final concentration prior to Western blot incubation to inhibit specific membrane bound antigen recognition. Can also be run on a Western blot as a positive control sample.	Cat # AACX1
A431 cell lysate: acetylation control	Trichostatin A (TSA) treated lysate	The HDAC inhibitor TSA stops protein deacetylation and allows accumulation of acetylated proteins. Most notable in A431 cells is the accumulation of tubulin at 55 kDa and histones at 14-16 kDa.	Cat # AACX2

Applications

Western Blot (WB) Applications

Use as indicated below at 1:500-1:1000 dilution, sufficient for 100-200 ml of working strength Ab.

Figure 1 : Western blot applications



Legend: AAC01 was used at a 1:500 dilution following the recommended Western blot protocol (see below). Figure 1A: murine tissue extract, 30 µg liver extract. Figure 1B: 20 µg of A431 cell lysate treated with TSA (+) or untreated (-). Strongly enhanced bands at 55 and 14-16 kDa in TSA-treated lysate correspond to acetylated tubulin and histone proteins, respectively. Figure 1C: Titration of acetylated BSA lanes 1-5 contain 0.5, 0.1, 0.05, 0.01, and 0.005 ng Ac-BSA, lanes 6-7 contain 500 and 1000 ng non-acetylated BSA, respectively. AAC01 recognizes 0.005 ng of chemically acetylated BSA. For comparison ImmuneChem anti-acetyl lysine antibody Cat# ICP0380 was able to detect 0.05 ng of acetylated BSA when used as instructed by manufacturer. Arrowhead indicates acetylated BSA, higher molecular weight bands are acetylated BSA oligomers.

Western Blot Method:

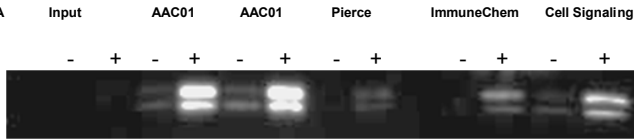
- Run protein samples and control samples in SDS-PAGE.
- We recommend running 20 µg of TSA-treated A431 cell lysate as a control.
- Equilibrate the gel in Western blot buffer (25 mM Tris pH 8.3, 192 mM glycine, and 15% methanol) for 15 min at room temperature prior to electro-blotting.
- Transfer the protein to a PVDF membrane for 60 min at 75 V.
- Wash the membrane with TBST for 10 min. with constant agitation (10 mM Tris-HCl pH 8.0, 150 mM NaCl, 0.05% Tween 20).
- Air dry the membrane at room temperature for 30 min.
- Rehydrate the membrane in TBST for 30 min. at room temperature.
- The membrane may be left in TBST overnight at 4°C if convenient.
- Block the membrane surface with 5% nonfat-dry milk in TBST for 30 min at room temperature with constant agitation.
- Incubate the membrane with a 1:500-1:1000 dilution of anti-acetyl lysine antibody, diluted in TBST/3% non-fat milk, for 1-2 h at room temperature or overnight at 4°C with constant agitation.
- Rinse the membrane three times in 50 ml TBST for 10 min. each at room temperature with constant agitation.
- Incubate the membrane with an appropriate dilution (e.g., 1:20,000) of anti-mouse secondary antibody (e.g., goat anti-mouse HRP conjugated IgG from Jackson Labs., Cat. # 115-035-068) in TBST/3% non-fat milk for 60 min at room temperature with constant agitation.
- Wash the membrane 6 times in TBST for 10 min each with constant agitation.
- Use an enhanced chemiluminescence detection method to detect the signal (e.g., SuperSignal West Dura Extended Duration Substrate; ThermoFisher).

NOTE: This Ab has some cross reactivity with the dyes used for pre-stained molecular weight markers. This does not affect protein specificity for acetyl groups (see Fig. 1B-1C). Also, some molecular weight marker proteins are acetylated, particularly commonly used metabolic enzymes such as glutamic dehydrogenase (m.wt. approx. 55kDa).

Immunoprecipitation (IP) Applications

Use as indicated at 20 μ l per IP reaction (1.5 mg total lysate per IP), sufficient for approximately 10 IP assays.

Figure 2: IP of histone proteins from TSA-treated A431 cells



Legend: A431 cells were either treated (+) or untreated (-) with TSA (0.6 μ M for 6 h). Cell lysates were prepared in a modified RIPA buffer (50 mM Tris pH 7.5, 150 mM NaCl, 1% IGEPAL, 0.1% SDS, 0.5% Na deoxycholate) and 1.5 mg of lysate per reaction was used for IP of acetylated histones. Each IP was performed according to the manufacturer's instructions. IP reactions for AAC01 are shown in duplicate. Western blots of immunoprecipitated proteins were developed using AAC01 at 1:500 dilution. Input signal is from 15 μ g of TSA treated (+) or untreated (-) A431 lysates. Ability of AAC01 to IP histones was compared to other commercially available antibodies, ThermoFisher (Pierce) Cat. # MA1-2021, ImmuneChem Cat. # ICP0388 and Cell Signaling Cat. # 9441. IPs were carried out according to the manufacturer's instructions.

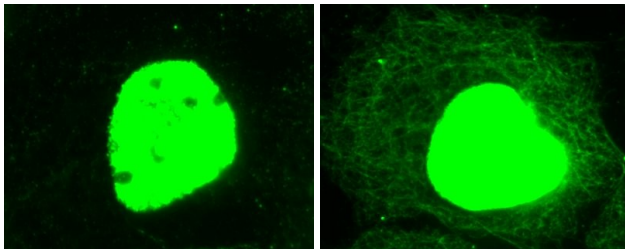
IP Method

1. Add 20 μ l of antibody to 500 μ l of PBS pH 7.4 in a microfuge tube containing 30 μ l of packed Protein G agarose pre-equilibrated in PBS.
2. Gently rotate the reaction for 1 h at 4°C.
3. Add 500 μ l of PBST to the mixture and centrifuge for 1 min at 4°C and 3000 rpm (approx. 960 x g). Addition of the PBST will prevent agarose from sticking to the microfuge tube walls.
4. Discard supernatant and wash beads 3X in PBST.
5. Add 1-1.5 mg of cell lysate (1-1.5 mg/ml protein concentration) to the beads. The lysate must be prepared in an IP compatible buffer. In this assay, the IP protein is in the following buffer: 50 mM Tris pH 7.5, 150 mM NaCl, 1% IGEPAL, 0.05% SDS, 0.25% Na deoxycholate).
6. Gently rotate the reaction at 4°C for 2 h or overnight if convenient.
7. Spin down agarose for 1 min at 4°C at 3000 rpm (approx. 960 x g).
8. Discard supernatant and wash beads with 700 μ l of PBST at 4°C.
9. Repeat wash two more times.
10. Resuspend beads in 30 μ l of 2X Laemmli buffer (125 mM Tris pH 6.8, 20% glycerol, 4% SDS, 0.005% Bromophenol blue, 5% beta-mercaptoethanol) and boil for 5 min prior to loading on SDS-PAGE for subsequent Western blot analysis.

Immunofluorescence (IF) Applications

Use as indicated below at 1:200 dilution, sufficient for 40 ml of working strength Ab, approx. 400 IF slides.

Figure 3: IF of A431 cells untreated and treated with TSA



Legend: Human epidermoid carcinoma A431 cells, untreated (left) or treated (right) with TSA (5 μ M for 16 h), were stained as described in the method. Acetylated cytoplasmic and nuclear proteins were visualized in green fluorescence. Note that in contrast with the untreated control, acetylated microtubule network is clearly visible in TSA-treated sample. The fluorescent nuclear intensities indicate the high abundance of acetylated proteins in the nucleus.

IF Method (TSA treatment of A431 cells)

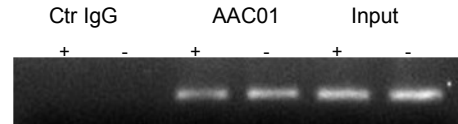
1. Plate A431 cells at 1×10^5 /ml on acid washed coverslips in tissue culture dish with DMEM media containing 10% FBS.
2. Allow cells to grow for 24-48 h, then treat one set of coverslips with TSA (5 μ M for 16 h).
3. Fix cells by dipping coverslips in -20°C methanol for 5 min (it is convenient to use a porcelain coverslip rack for fixing and rinsing samples).
4. Remove coverslips and dip into -20°C acetone for 2 min.

5. Air dry the coverslips. Dip the coverslips into PBS for 5 min. at room temperature.
6. Place coverslips cell side up on parafilm and apply 100 μ l-200 μ l of AAC01 solution to each coverslip (1:200 Ab dilution in PBS/3% BSA).
7. Incubate at room temperature for 30 min.
8. Dip the coverslips into PBS for 5 min. at room temperature.
9. Apply 100 μ l-200 μ l of fluorescently-labeled anti-mouse secondary antibody at manufacturer's recommended dilution. For example, we use fluorescently-labeled goat anti-mouse at 1:200 dilution in PBS.
10. Incubate at room temperature for 30 min.
11. Dip the coverslips into PBS for 5 min. at room temperature.
12. Rinse coverslips for 30 seconds in sterile water.
13. Dip coverslips in absolute ethanol at room temperature for 2 min and transfer to blotting paper, cell side up, allow to air dry.
14. Place coverslips, cell side down, on glass slide with mounting media (e.g., EMS, Cat# 17987-10) and observe cells under fluorescence microscope.

Chromatin Immunoprecipitation (ChIP) Application

Use as indicated below at 1:100 dilution, sufficient for 200 ChIP assays of 100 μ l volume.

Figure 4: ChIP analysis with AAC01 antibody



Legend: Chromatin was prepared from A431 cells, either untreated or TSA-treated (5 μ M, 4 hrs). ChIP was performed according to method see below. Ctr IgG: mouse IgG used for ChIP, AAC01: anti-acetyl antibody used for ChIP, Input: cell lysate prior to ChIP. PCR analysis was carried out using primers for the promoter region of the housekeeping gene GAPDH.

ChIP Method

1. Sheared and cross-linked chromatin from A431 cells was prepared according to published protocols (4).
2. Approximately 25 μ g of chromatin is used per ChIP assay and the reactions are carried out in the following ChIP buffer: 10 mM Tris pH 8.0, 0.5 mM EGTA, 1 mM EDTA, 140 mM NaCl, 1% Triton, 0.15% SDS, 0.1% Na deoxycholate, 1 mM PMSF, 2 μ g/ml pepstatin/leupeptin, and aprotinin.
3. Make the final volume to 100 μ l in ChIP reaction buffer and add 1 μ l of AAC01 per ChIP. NOTE: it is also recommended to run mouse IgG (5 μ g) as a non-specific control reaction (see Ctr IgG lanes above) and a no-antibody as input control (see Input lanes above).
4. Incubate on a rotator at 4°C for 1-2 h or overnight.
5. Add 20 μ l of a 50% slurry of Protein G beads to all reactions and rotate at 4°C for 2-4 h.
6. Pellet beads by centrifugation at 3000 rpm (approx. 960 x g) for 1 min.
7. Transfer the no-antibody control supernatant to a fresh tube on ice, this will be processed later and serve as the total input control.
8. Discard all other supernatants.
9. Wash beads five times with 500 μ l of RIPA buffer (10 mM Tris pH 8.0, 0.5 mM EGTA, 1 mM EDTA, 140 mM NaCl, 1% Triton, 0.1% SDS, 0.1% Na deoxycholate, and 1 mM PMSF).
10. Wash beads once with 500 μ l of LiCl buffer (0.25 M LiCl, 0.5% NP-40, 0.5% Na deoxycholate, 1 mM EDTA, 10 mM Tris pH 8.0).
11. Wash beads once with 500 μ l of TE buffer (10 mM Tris pH 8.0, 1 mM EDTA) and resuspend beads in 100 μ l of TE buffer.
12. The immunoprecipitated chromatins bound to the bead samples (and chromatins in the no-antibody control supernatant) were processed for DNA isolation and PCR analysis according to published protocols (4).

References

1. Bogdan P. and Sherman F. 2002. The diversity of acetylated proteins. *Genome Biol.* 3 (5): reviews 0006.
2. Lundby A. et al. 2012. Proteomic analysis of lysine acetylation sites in rat tissues reveals organ specificity and cellular patterns. *Cell Reports* 2:419-431.
3. Sadouli K. et al. 2010. The tale of protein lysine acetylation in the cytoplasm. *J. Biomed. Biotech.* 2011:1-15.
4. Golemis EA et. Al, Protein-Protein Interactions, *CSHLP*, 2005, p67.

Product Citations/Related Products

For the latest citations and related products please visit www.cytoskeleton.com