# Datasheet

Anti-Acetyl Lysine Mouse Monoclonal Antibody Horseraddish peroxidase conjugates

Cat. # AAC03-HRP

Upon arrival, store at 4°C (desiccated) See datasheet for storage after reconstitution

Online Datasheet Contains AAC03-HRP (V1.0) and AAC03-HRP-S (V1.0)

# **Background Information**

Acetylation of proteins can occur as a co-translational or post-translational modification (PTM) (1). Co-translational acetylation occurs at the N-terminal of approximately 85% of mammalian proteins, it 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 (3).

# Material

AAC03-HRP 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 broad range of acetylated proteins, including acetylated tubulin, histones, and chemically acetylated bovine serum albumin (Fig. 1). AAC03-HRP is supplied as a lyophilized white powder. Each Lot of antibody is quality controlled to provide high batch-to-batch consistency. The Lot specific µg per tube can be found in the Lot specific COA documents at www.cytoskeleton.com.

# Storage and Reconstitution

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

Reconstitute each tube with 100  $\mu$ l of 50% glycerol (room temperature). We do not recommend using 50% glycerol at 4°C as this can cause the lyophilized antibody to stick to the pipet tip during resuspension. Store reconstituted antibody at –20°C. Final buffer composition is 200 mM PIPES, 50% glycerol, 5% sucrose,1% dextran and 10mg/ml BSA.

When stored and reconstituted as described, the product is stable for 12 months at -20°C. NOTE: Sodium azide is an irreversible inhibitor of HRP. Do not add sodium azide to APY03-HRP antibody.

## Applications

### Western Blot (WB) Applications

Use as indicated below at 1:3000-1:6000 dilution, sufficient for 300-600 ml of working strength Ab.

### Western Blot Method:

- 1. Run protein samples and control samples in SDS-PAGE.
- We recommend running 30 µg of TSA/nicotinamide-treated Cos-7 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.
- 4. Transfer the protein to a PVDF membrane for 60 min at 70 V.
- Wash the membrane once with TBST (10 mM Tris-HCl pH 8.0, 150 mM NaCl, 0.05% Tween 20).
- 6. The membrane may be left in TBST overnight at 4°C if convenient.
- 7. Block the membrane surface with 3% nonfat-dry milk in TBST for 60 min at room temperature with constant agitation.
- Incubate the membrane with a 1:3000-1:6000 dilution of anti-acetyl lysine antibody, diluted in 3% nonfat-dry milk in TBST, for 1-2 h at room temperature or overnight at 4°C with constant agitation.
- 9. Rinse the membrane five times in 50 ml TBST for 10 min. each at room temperature with constant agitation.
- 10. Use an enhanced chemiluminescence detection method to detect the signal (e.g., SuperSignal West Dura Extended Duration Substrate; ThermoFisher).

Form:	Lyophilized powder
Amount of material:	1 x 100 $\mu$ l when reconstituted
Validated applications:	WB
Species reactivity:	All
Host/Isotype:	Mouse/IgG1
Clone:	19C4B2.1

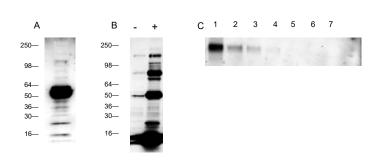


Fig 1: Utilization of AAC03-HRP for western blotting. A: Murine tissue extract, 30 μg brain extract. B: 30 μg of Cos-7 cell lysate treated with TSA and nicotinamide (+) or untreated (-). Strongly enhanced bands at 55 and 14-16 kDa in TSA-treated lysate correspond to acetylated tubulin and histone proteins, respectively. C: 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. AAC03-HRP was used at a 1:3000 dilution following the recommended Western blot protocol.

### 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 Sadoul 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

cytoskeleton.com

Phone: (303) 322.2254 Fax: (303) 322.2257 Customer Service: cserve@cytoskeleton.com Technical Support: tservice@cytoskeleton.com

V.1.0

# Datasheet

# Anti-Acetyl Lysine Mouse Monoclonal Antibody Horseraddish peroxidase conjugates

Cat. # AAC03-HRP-S

Upon arrival, store at 4°C (desiccated) See datasheet for storage after reconstitution

## **Background Information**

Acetylation of proteins can occur as a co-translational or post-translational modification (PTM) (1). Co-translational acetylation occurs at the N-terminal of approximately 85% of mammalian proteins, it 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 (3).

### Material

AAC03-HRP 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 broad range of acetylated proteins, including acetylated tubulin, histones, and chemically acetylated bovine serum albumin (Fig. 1). AAC03-HRP is supplied as a lyophilized white powder. Each Lot of antibody is quality controlled to provide high batch-to-batch consistency. The Lot specific µg per tube can be found in the Lot specific COA documents at www.cytoskeleton.com.

### Storage and Reconstitution

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

Reconstitute each tube with 25  $\mu$ I of 50% glycerol (room temperature). We do not recommend using 50% glycerol at 4°C as this can cause the lyophilized antibody to stick to the pipet tip during resuspension. Store reconstituted antibody at –20°C. Final buffer composition is 200 mM PIPES, 50% glycerol, 5% sucrose,1% dextran and 10mg/mI BSA.

When stored and reconstituted as described, the product is stable for 12 months at -20°C. NOTE: Sodium azide is an irreversible inhibitor of HRP. Do not add sodium azide to APY03-HRP antibody.

## Applications

### Western Blot (WB) Applications

Use as indicated below at 1:3000-1:6000 dilution, sufficient for 75-150 ml of working strength Ab.

### Western Blot Method:

- 1. Run protein samples and control samples in SDS-PAGE.
- We recommend running 30 µg of TSA/nicotinamide-treated Cos-7 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.
- 4. Transfer the protein to a PVDF membrane for 60 min at 70 V.
- Wash the membrane once with TBST (10 mM Tris-HCl pH 8.0, 150 mM NaCl, 0.05% Tween 20).
- 6. The membrane may be left in TBST overnight at 4°C if convenient.
- 7. Block the membrane surface with 3% nonfat-dry milk in TBST for 60 min at room temperature with constant agitation.
- Incubate the membrane with a 1:3000-1:6000 dilution of anti-acetyl lysine antibody, diluted in 3% nonfat-dry milk in TBST, for 1-2 h at room temperature or overnight at 4°C with constant agitation.
- 9. Rinse the membrane five times in 50 ml TBST for 10 min. each at room temperature with constant agitation.
- 10. Use an enhanced chemiluminescence detection method to detect the signal (e.g., SuperSignal West Dura Extended Duration Substrate; ThermoFisher).

Lyophilized powder
1 x 25 µl when reconstituted
WB
All
Mouse/IgG1
19C4B2.1

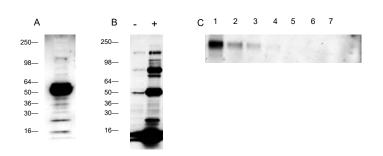


Fig 1: Utilization of AAC03-HRP for western blotting. A: Murine tissue extract, 30 µg brain extract. B: 30 µg of Cos-7 cell lysate treated with TSA and nicotinamide (+) or untreated (-). Strongly enhanced bands at 55 and 14-16 kDa in TSA-treated lysate correspond to acetylated tubulin and histone proteins, respectively. C: 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. AAC03-HRP was used at a 1:3000 dilution following the recommended Western blot protocol.

#### 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 Sadoul 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

cytoskeleton.com

# Phone: (303) 322.2254 Fax: (303) 322.2257 Customer Service: cserve@cytoskeleton.com Technical Support: tservice@cytoskeleton.com

V.1.0