





The Protein  
Experts

Cytoskeleton, Inc.

# Datasheet

## Western Blot Method:

1. Run protein samples and control samples in SDS-PAGE.
2. 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 electroblotting.
3. Transfer the protein to a PVDF membrane for 60 min at 75 V.
4. 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).
5. Air dry the membrane at room temperature for 30 min.
6. Rehydrate the membrane in TBST for 30 min. at room temperature.
7. The membrane may be left in TBST overnight at 4° C if convenient.
8. Block the membrane surface with 5% nonfat-dry milk in TBST for 30 min at room temperature with constant agitation.
9. Incubate the membrane with a 1:500-1:1000 dilution of anti-acetyl lysine antibody, diluted in TBST, for 1-2 h at room temperature or overnight at 4°C with constant agitation. If Ac-BSA is to be used as a competitor, then include this at 10 µg/ml.
10. Rinse the membrane three times in 50 ml TBST for 10 min. each at room temperature with constant agitation.
11. 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/0.5% non-fat milk for 60 min at room temperature with constant agitation.
12. Wash the membrane 6 times in TBST for 10 min each with constant agitation.
13. 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. 1 & 2). Also, some molecular weight marker proteins are acetylated, particularly commonly used metabolic enzymes such as glutamic dehydrogenase (m.wt. approx. 55kDa).

## References

1. Yang X-J. 2004. The diverse superfamily of lysine acetyltransferases and their roles in leukemia and other diseases. *Nucleic Acids Res.* **32**, 959-976.
2. Komatsu Y. et al. 2003. Four different clones of mouse anti-acetyllysine monoclonal antibodies having different recognition properties share a common immunoglobulin framework structure. *J. Immunol. Methods* **272**, 161-175.

## Product Citations / Related Products

For the latest citations and related products please visit [www.cytoskeleton.com](http://www.cytoskeleton.com).