

# PRODUCT INFORMATION



## Albumin (bovine) Acetylated

Item No. 10010566

### Overview and Properties

<b>Contents:</b>	This vial contains 1 mg of acetylated bovine serum albumin (BSA), lyophilized.
<b>Synonym:</b>	(Ac)BSA
<b>Form:</b>	Solid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Applications:</b>	ELISA or Western blot (WB); reconstitute the lyophilized protein with double distilled water. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.
<b>Special Conditions:</b>	Use within one month after reconstitution or store reconstituted aliquots at -20°C to avoid freeze-thaw cycles.

### Description

Acetylated BSA may be used as a positive control for ELISA or WB when employing an acetyl-lysine monoclonal or polyclonal antibody. Cayman's Acetyl Lysine Monoclonal Antibody (Clone 7F8) (Item No. 10010567) can detect less than 1 ng of this positive control by immunoblotting. Histone subunit modifications such as lysine acetylation are regulated by the activity of histone acetyltransferases (HATs) and histone deacetylases (HDACs).<sup>1,2</sup> Epigenetic modifications such as protein acetylation directly influence cellular genetic programs including those contributing to cancer cell viability.<sup>1,3,4</sup>

### References

1. Loidl, P. Histone acetylation: Facts and questions. *Chromosoma* **103(7)**, 441-449 (1994).
2. Cheung, W.L., Briggs, D.B., and Allis, C.D. Acetylation and chromosomal functions. *Curr. Opin. Cell Biol.* **12(3)**, 326-333 (2000).
3. Marks, P.A. and Breslow, R. Dimethyl sulfoxide to vorinostat: Development of this histone deacetylase inhibitor as an anticancer drug. *Nat. Biotechnol.* **25(1)**, 84-90 (2007).
4. Bolden, J.E., Peart, M.J., and Johnstone, R.W. Anticancer activities of histone deacetylase inhibitors. *Nat. Rev. Drug Discov.* **5(9)**, 769-784 (2006).

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/14/2023

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM