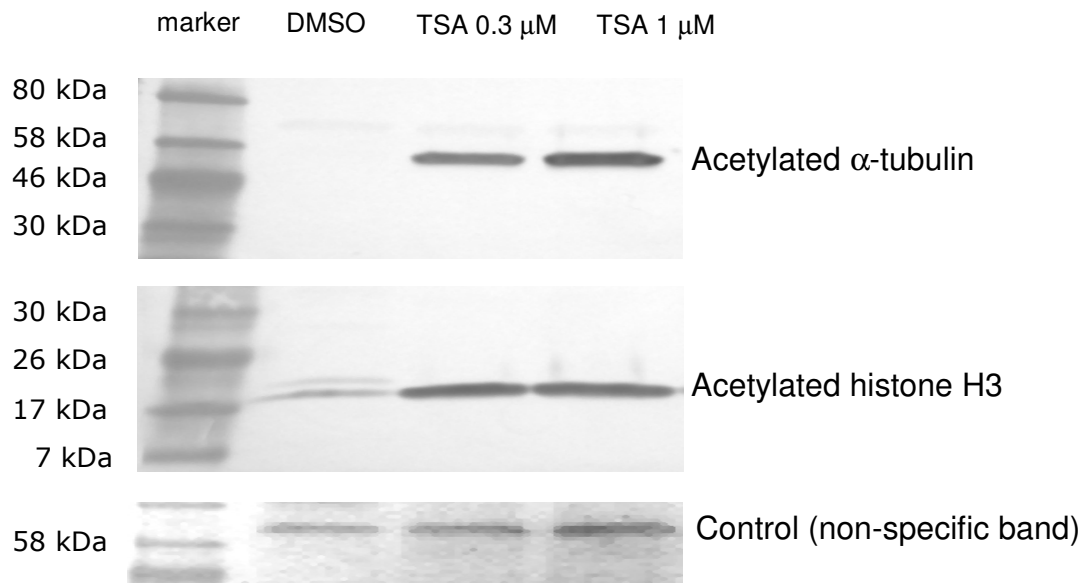


Histone Deacetylase (HDAC) Cell-Based Screening Services

BPS Bioscience now provides services to determine IC₅₀'s or screen inhibitors of HDACs using our **cell-based assays!** We can perform assays in tumor cell lines such as A549, PC-12, or other cell lines of your choice.

Contact us today so we can help you accelerate your drug discovery research!

Western blot-based assay to determine inhibition of HDAC activity in whole cells



PC-12 cells were treated with TSA (in DMSO stock solution) or DMSO alone. Following treatment, cells were harvested and cell lysates were prepared. Cell lysates were run on SDS-PAGE gel and transferred to nitrocellulose membrane. Western blotting was performed with anti-acetylated tubulin or anti-acetylated histone H3 antibodies. A non-specific band reactive with acetylated histone H3 antibody was used as a control to confirm protein normalization.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com