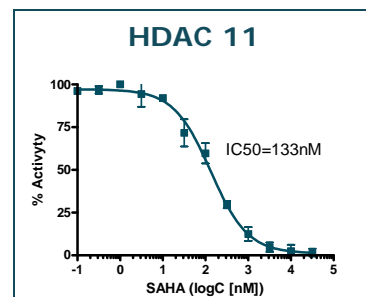
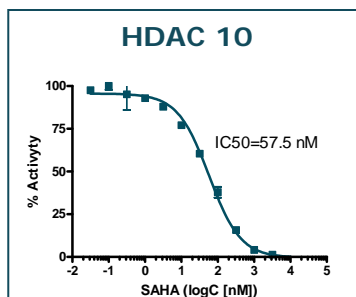
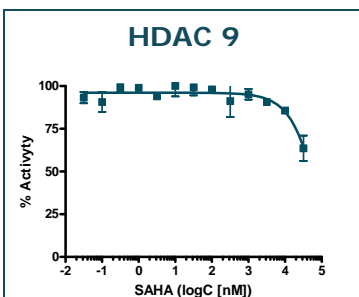
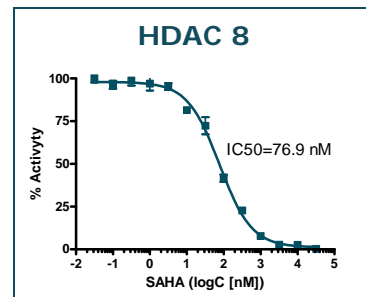
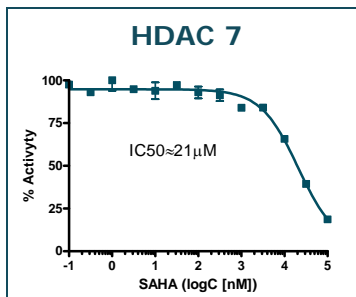
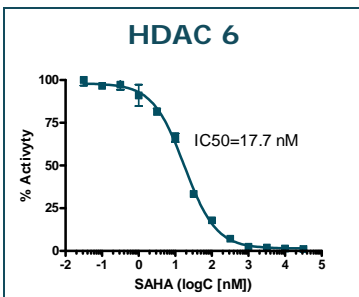
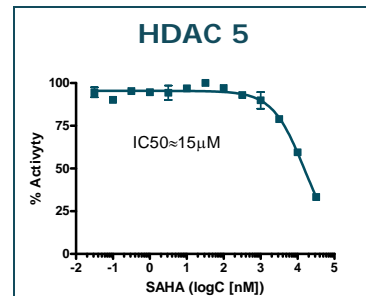
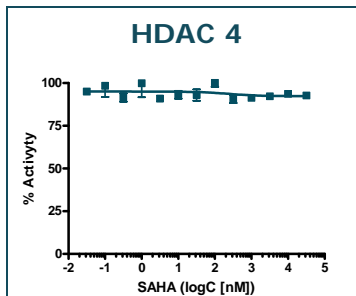
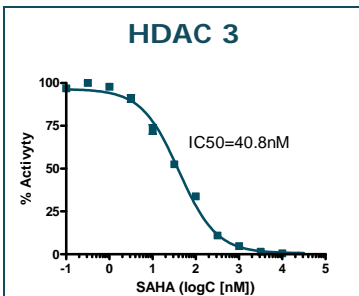
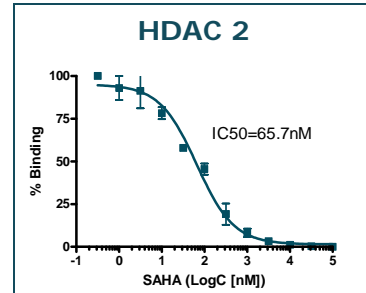
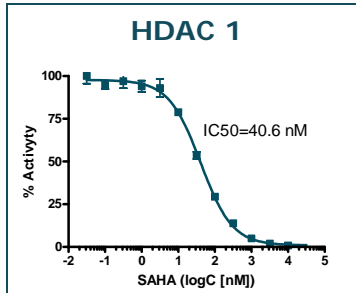




**SAHA Screening**  
 Assay Buffer: BPS#50031.  
 Assay Developer: BPS#50030.  
 Substrate BPS#50037 for HDAC 1, 2, 3, 6, 8, 10 and 11.  
 Substrate BPS#50040 for HDAC 4, 5, 7 and 9.



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

BPS Bioscience Inc. 6044 Cornerstone Court West, Suite E, San Diego, CA 92121.

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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)