

HDAC Screening and Profiling

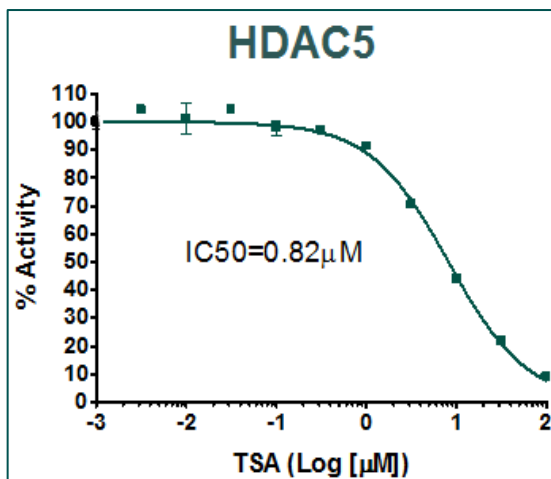
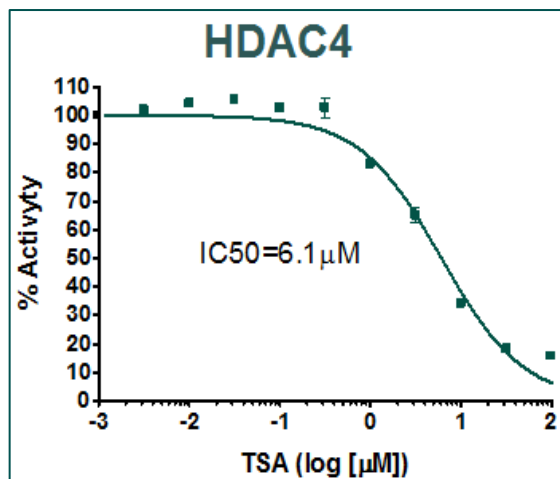
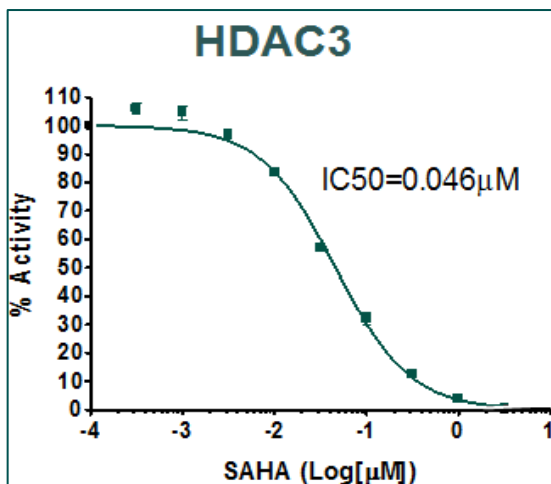
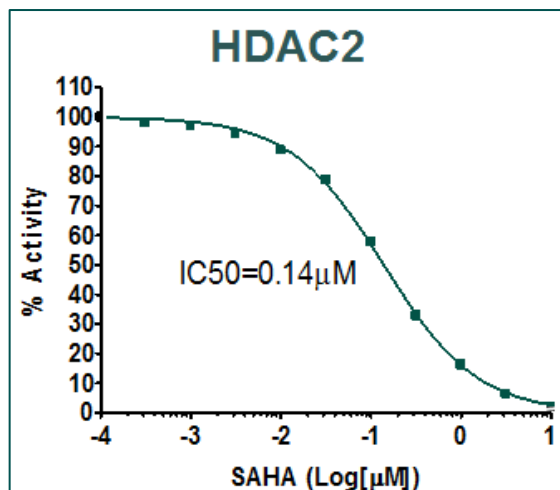
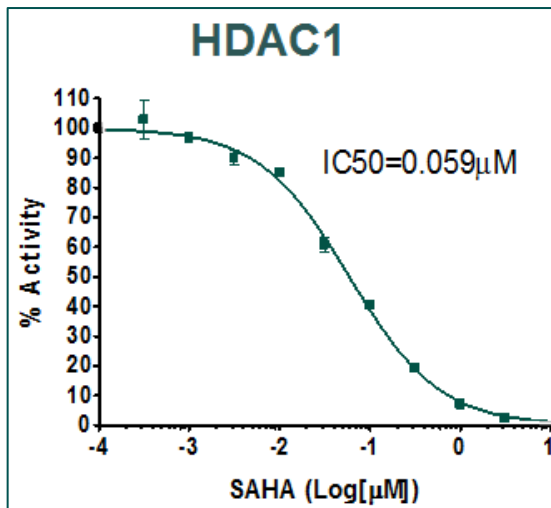
SAHA & TSA Screening

Assay Buffer: BPS #50031

Assay Developer: BPS #50030

Substrate: BPS #50037 for
 HDACs 1, 2, 3, 6, 10

Substrate: BPS #50040 for
 HDACs 4, 5, 7, 8, 9, 11

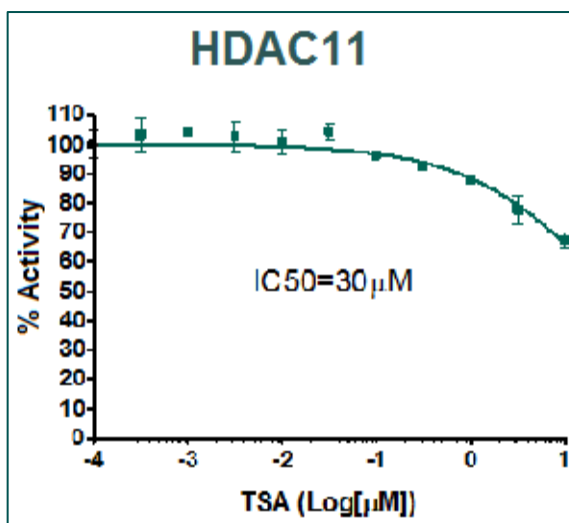
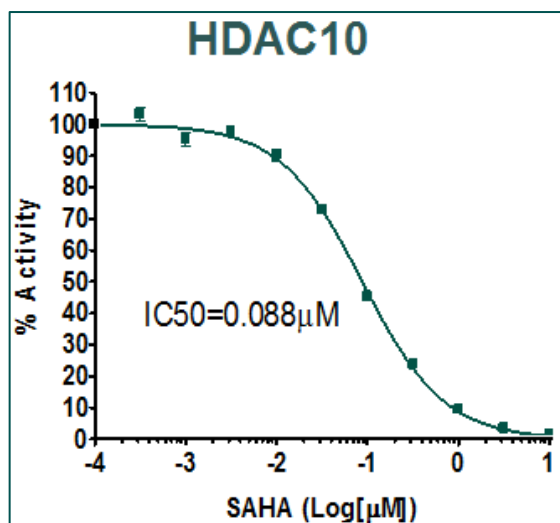
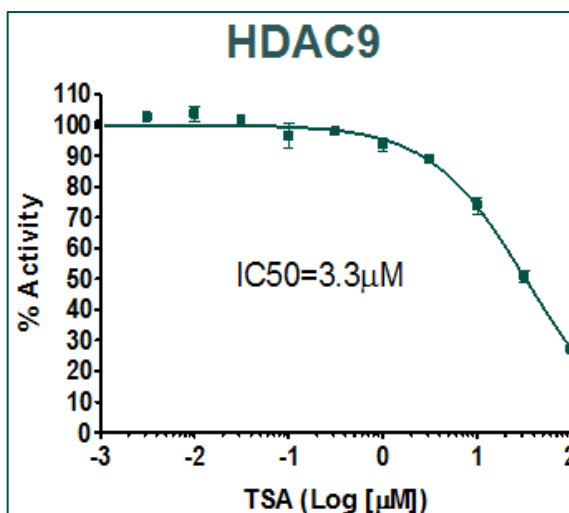
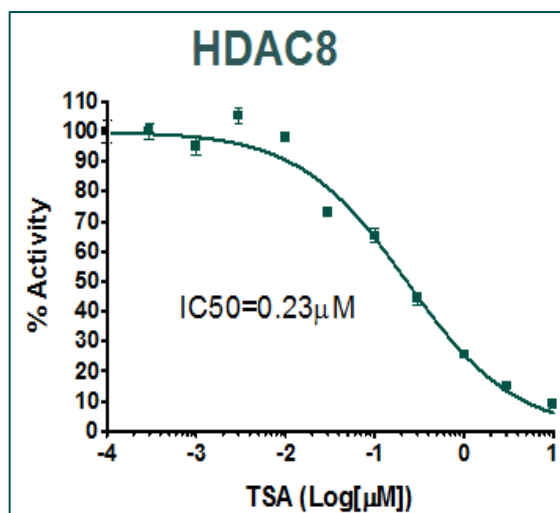
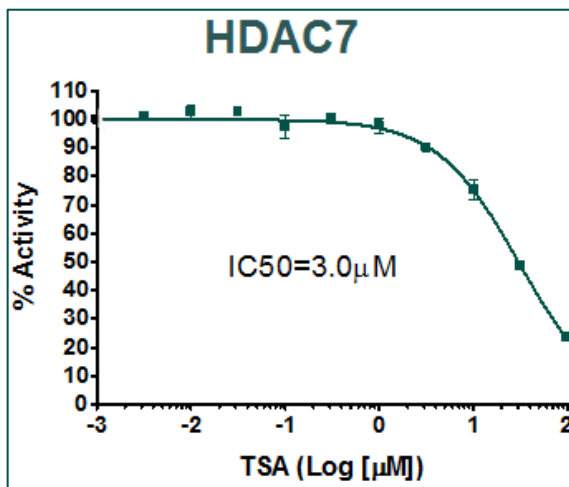
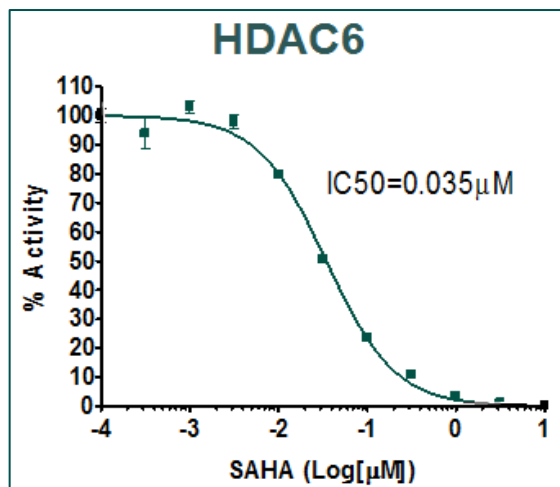


OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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



Also Available: HMT, DNMT, PARP, and Demethylase Profiling

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