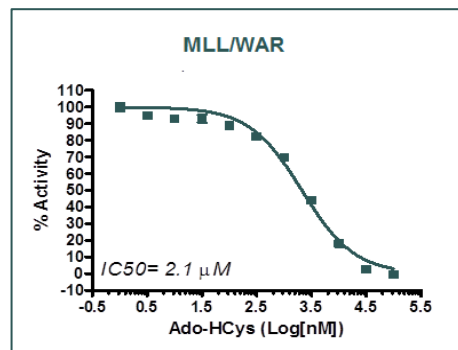
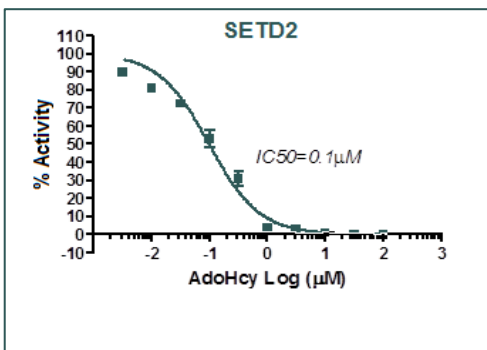
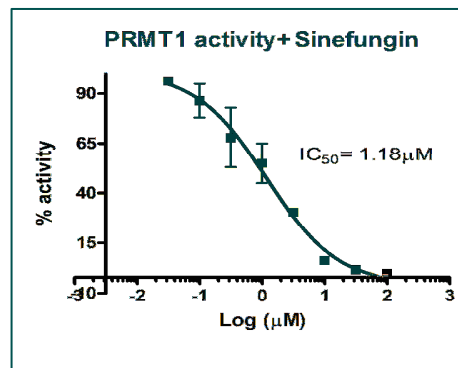
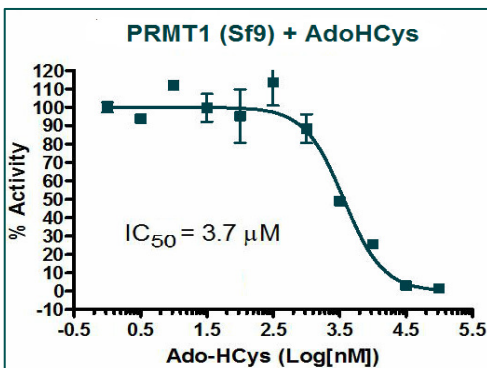
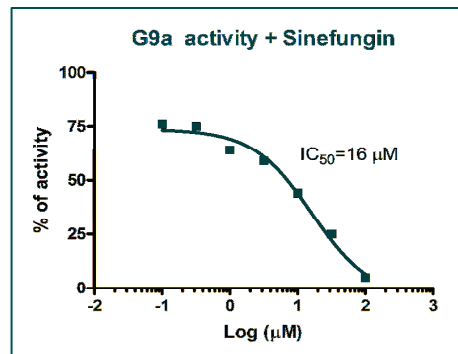
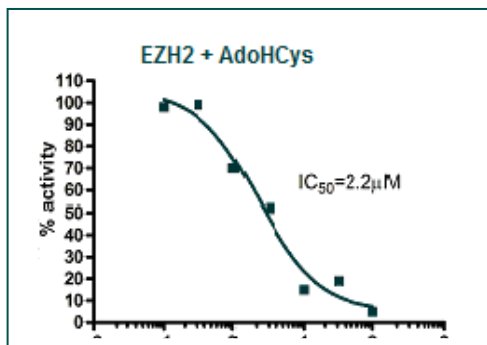
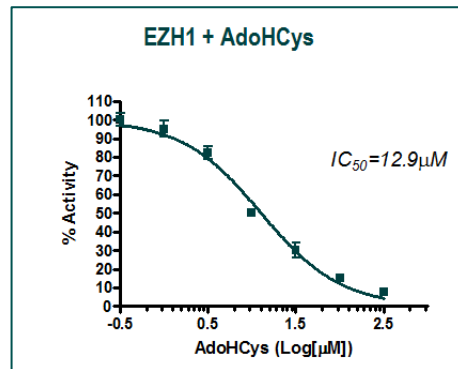


Methyltransferase Profiling

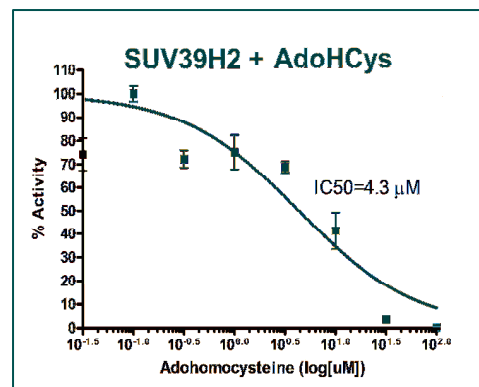
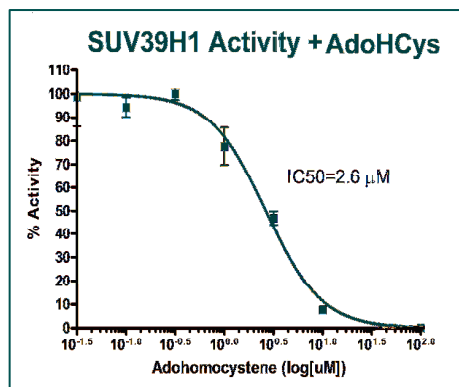
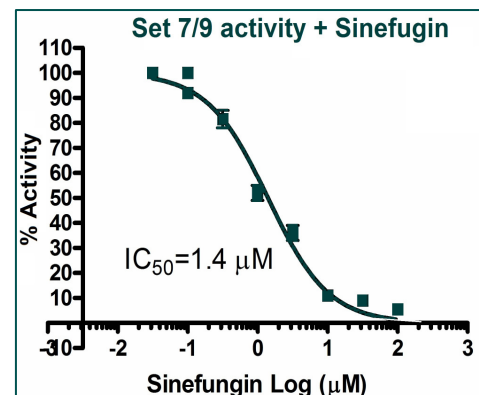
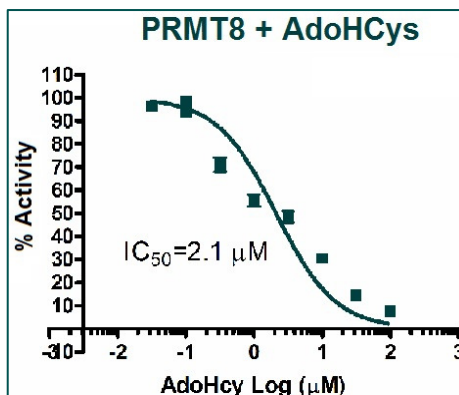
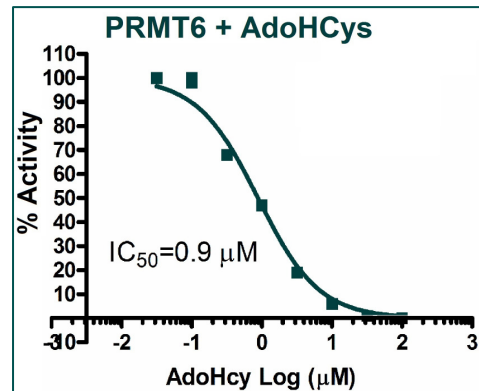
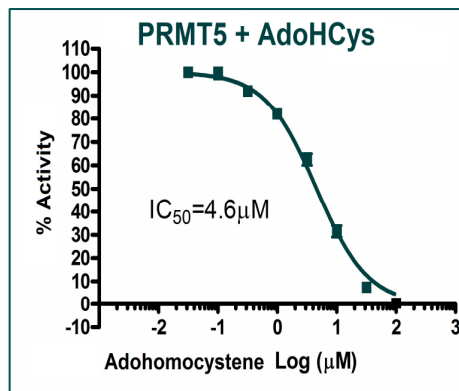
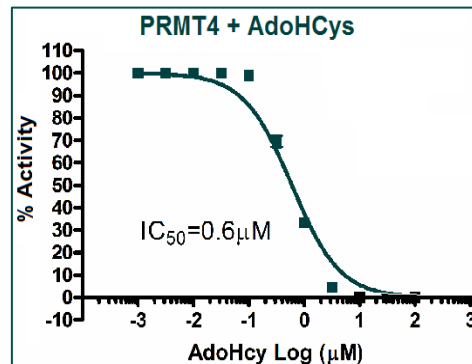
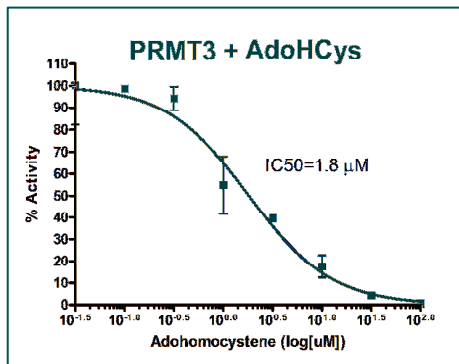
15 Different HMTs

EZH1, EZH2, G9a, PRMT1,
 PRMT3, PRMT4, PRMT5,
 PRMT6, PRMT8,
 SETD2, SET7/9, SET8,
 SUV39H1, SUV39H2, and
 MLL Complex



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: HDAC, 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