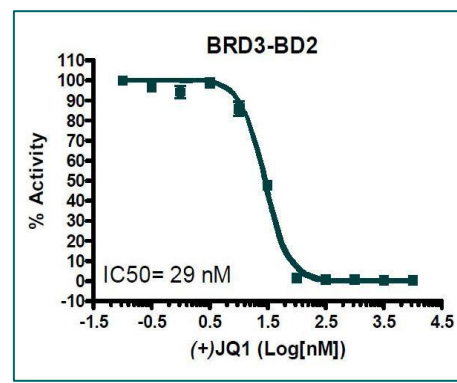
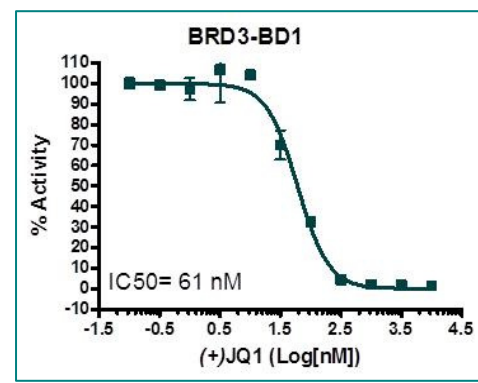
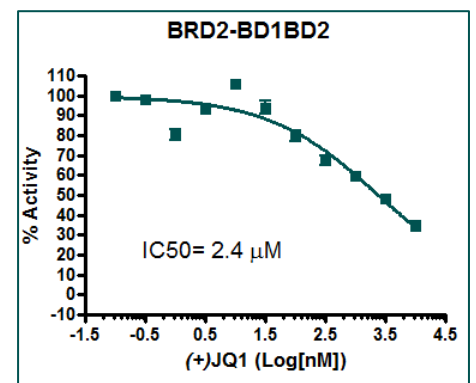
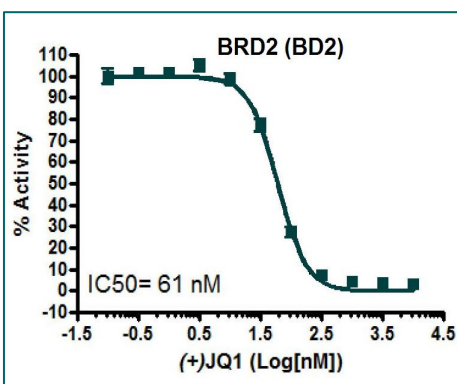
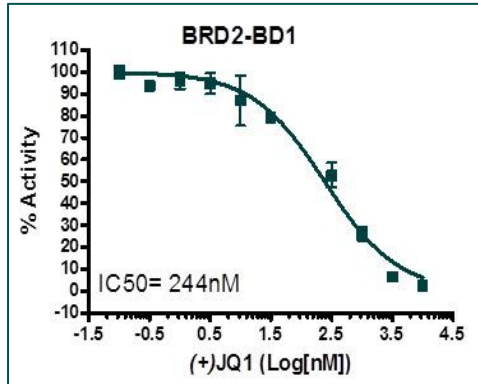
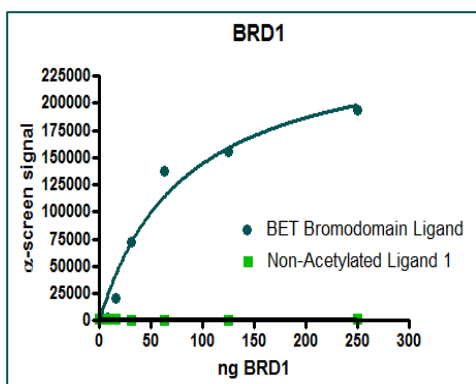
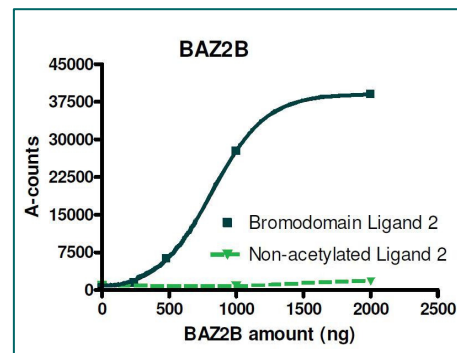
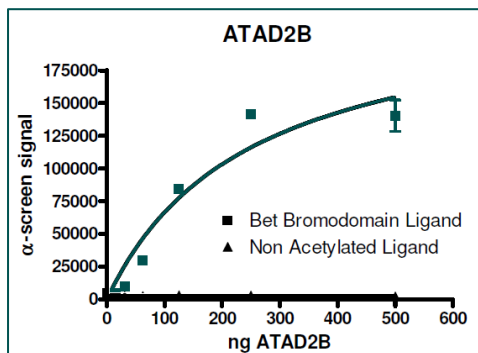
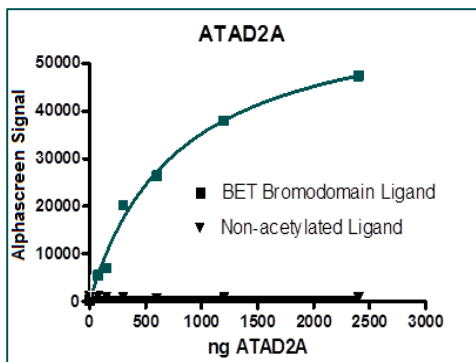


Bromodomain Profiling Services

BPS can provide full IC50 assays, or screen your compounds against a panel of bromodomains.



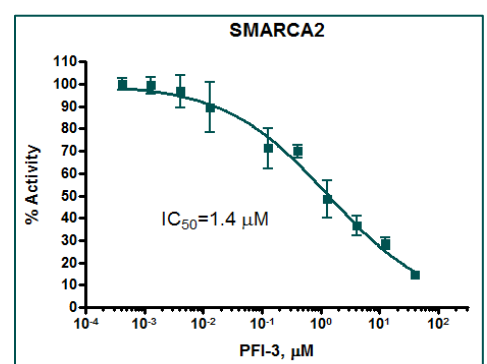
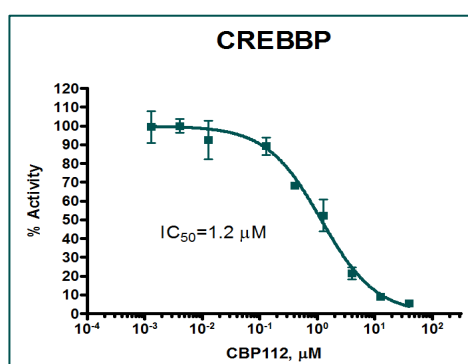
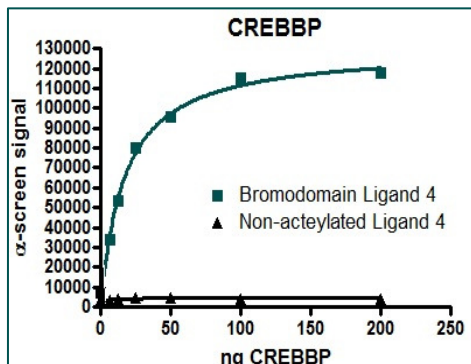
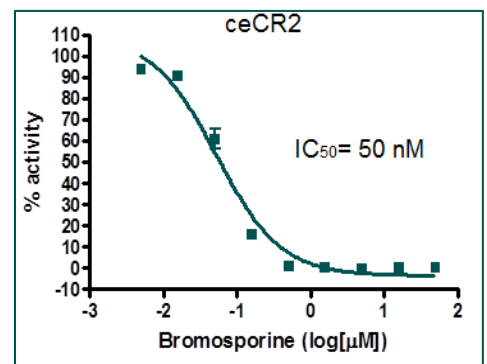
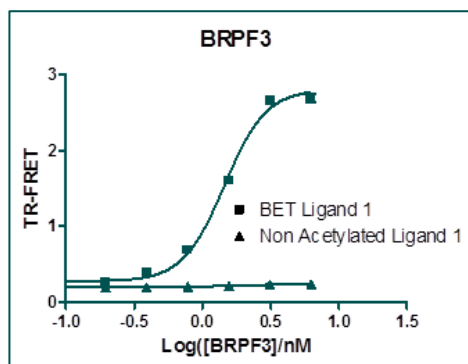
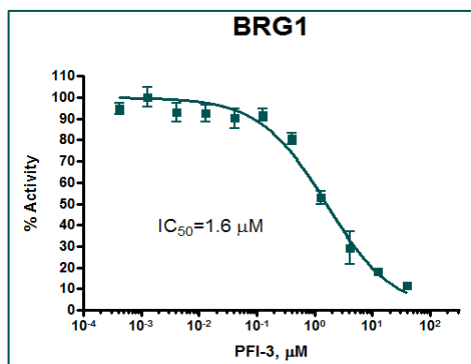
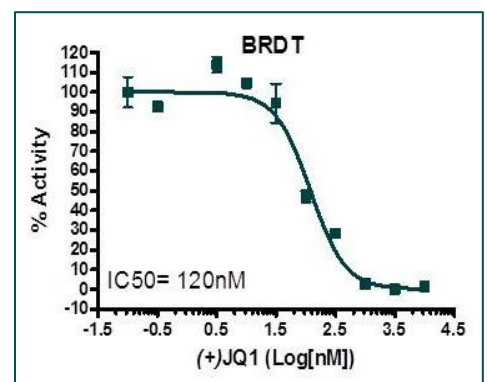
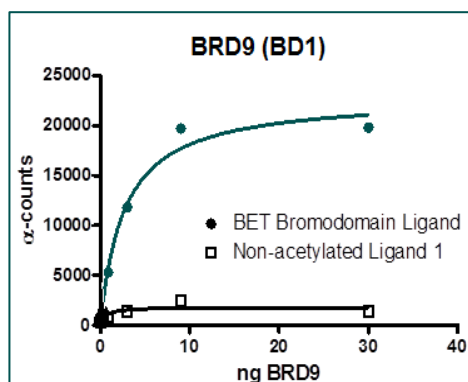
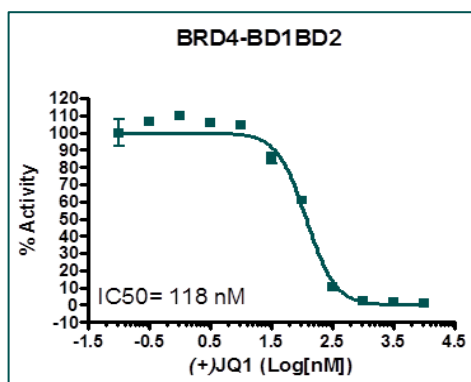
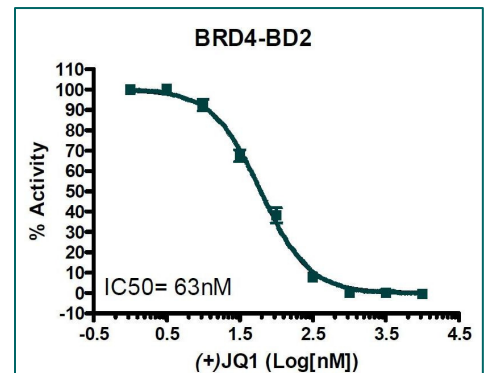
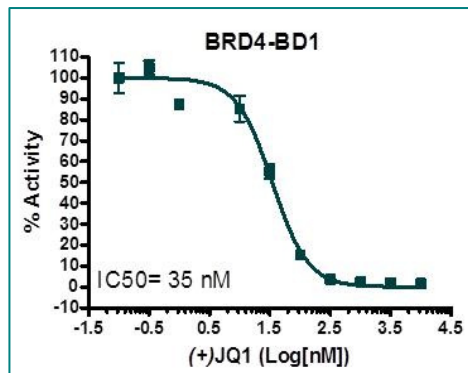
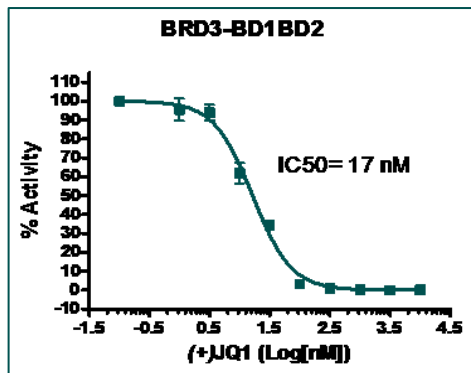
Other Screening/Profiling Services: HDACs, Histone Demethylases, Histone Methyltransferases, DNMTs, SIRT, and PARPs

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

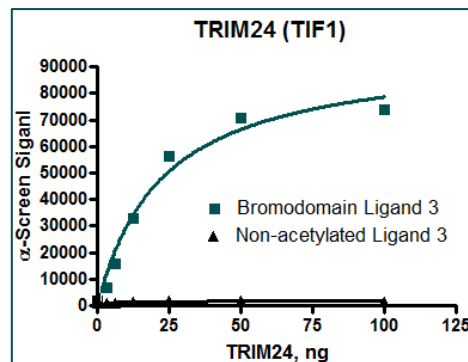
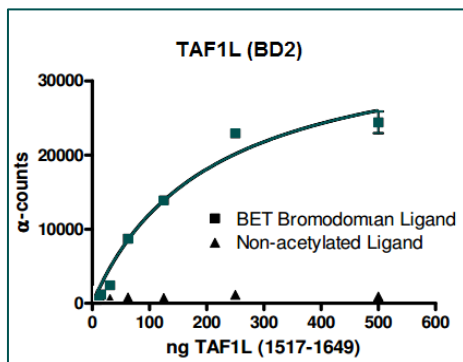
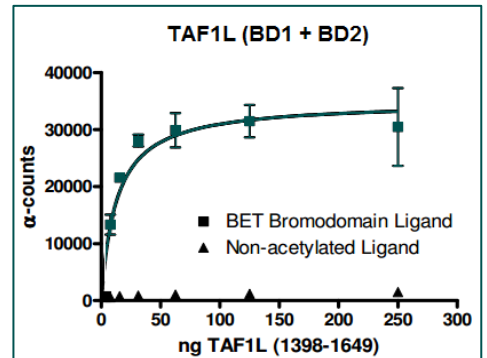
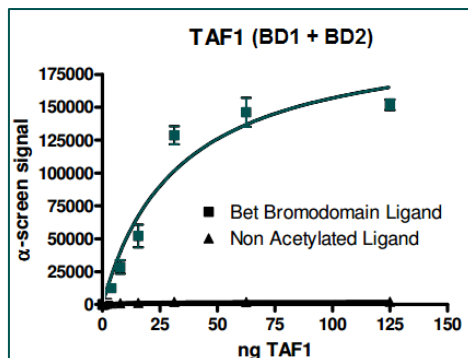
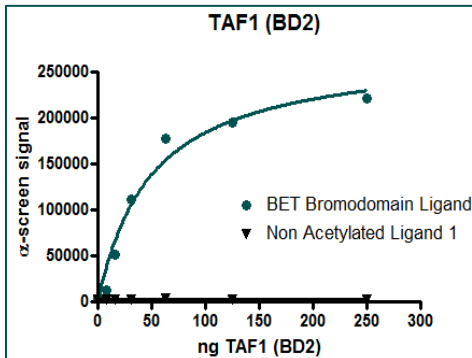
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



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



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