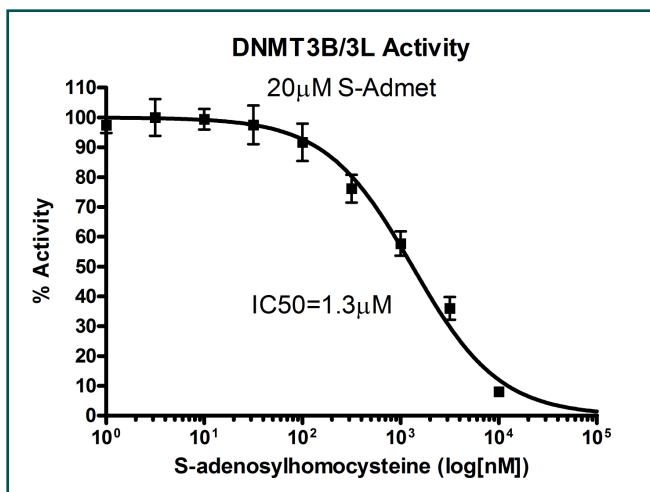
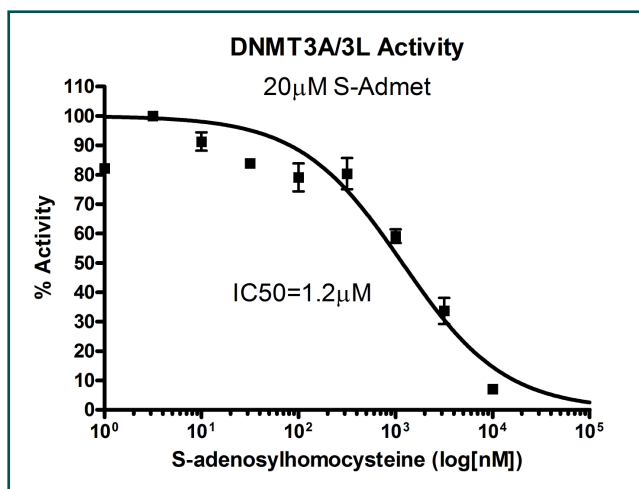
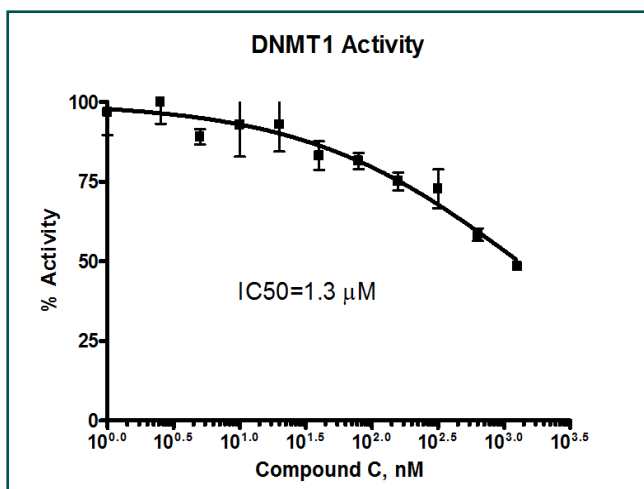


DNA Methyltransferase Profiling



Other Screening/Profiling Services:

Histone Methyltransferases

(14 available)

EZH1, EZH2, G9a, MLL, PRDM2, PRMT1, PRMT3, PRMT5, SET7/9, SET8, SUV39H1, SUV39H2, SUV4-20H1, and SUV4-20H2

Histone Demethylases

LSD1, JMJD2A, JMJD2B, JMJD2C

HDACs (all 11 HDACs)

HDACs 1-11, SIRT1-7

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