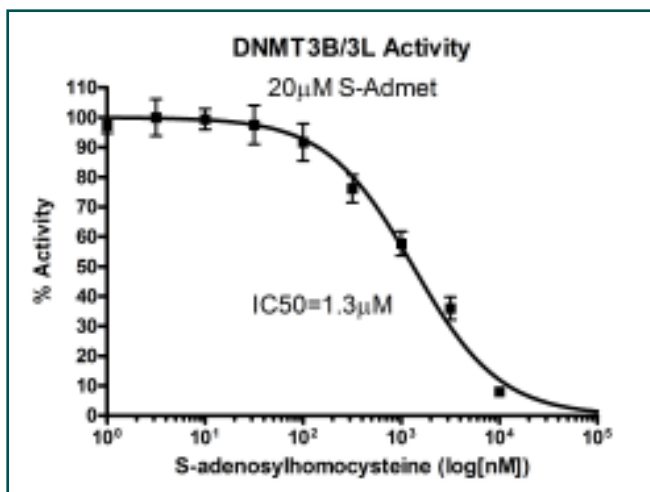
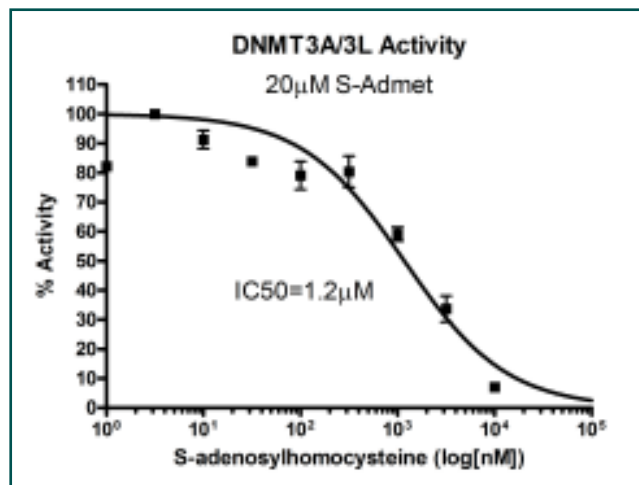
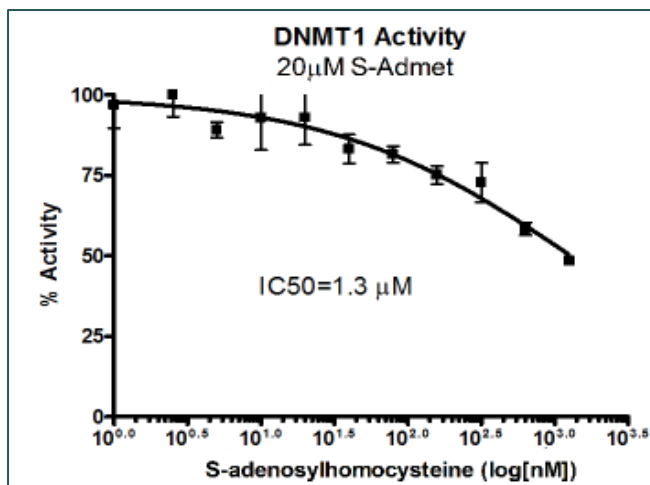


DNA Methyltransferase Profiling



Other Screening/Profiling Services:

Histone Methyltransferases (13 available)

EZH1, EZH2, G9a, PRMT1, PRMT3, PRMT4 (CARM1), PRMT5, PRMT6, PRMT8, SET7/9, SET8, SUV39H1, and SUV39H2

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