

DNMT3B/3L Methyltransferase Assay Service

Activity-based DNMT3B/3L Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

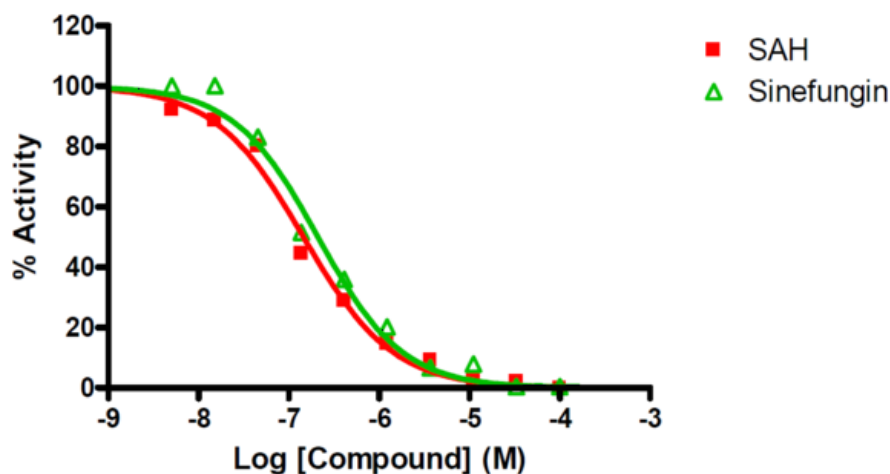
Target Overview

CAT#	DNMT3B/3L
Enzyme	Co-expressed Human DNMT3B/3L catalytic domain
Synonym	DNMT3B: DNA (cytosine-5)-methyltransferase 3B, DNA methyltransferase HsaIIIB, DNMT3B; DNMT3L: DNA (cytosine-5)-methyltransferase 3-like, DNMT3L
Substrates	Lambda DNA, 0.0075 mg/mlSAM, 1 μM

Assay Properties

Reaction	S-adenosyl-L-[methyl- ³ H]methionine + DNA = S-adenosyl-L-homocysteine + DNA 5-[methyl- ³ H]-cytosine	
Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (nM)
	SAH	146
	Sinefungin	210
Scientific Information	DNMT3B: Q9UBC3 (http://www.uniprot.org/uniprot/Q9UBC3)DNMT3L: Q9UJW3 (http://www.uniprot.org/uniprot/Q9UJW3)	
Screening Location	Malvern, USA	

Reference compound IC50 for DNMT3B/3L



Let's Discover Together.

