

DNMT3A Methyltransferase Assay Service

Activity-based DNMT3A Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

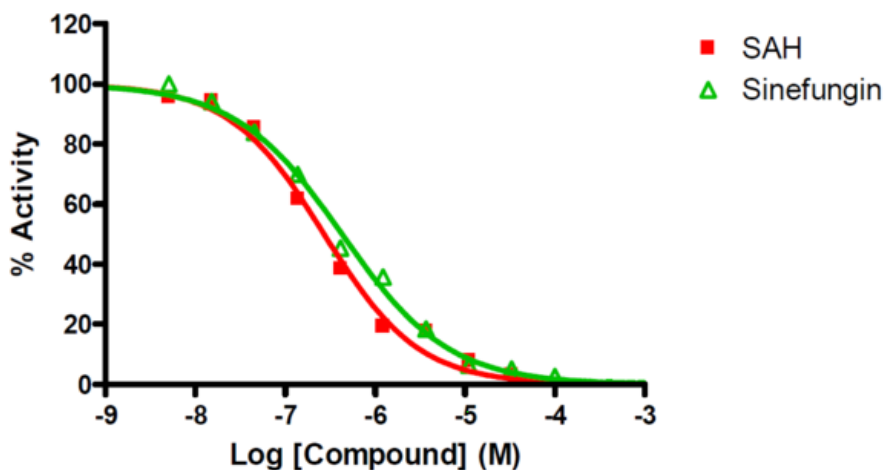
Target Overview

CAT#	DNMT3A
Enzyme	Human DNMT3A
Synonym	DNA (cytosine-5)-methyltransferase 3A, DNA methyltransferase HsaIIIa, DNMT3A
Substrates	Lambda DNA, 0.0075 mg/ml SAM, 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 271 Sinefungin 430
Scientific Information	Q9Y6K1 (http://www.uniprot.org/uniprot/Q9Y6K1)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website Methyltransferase Assays .

Reference compound IC50 for DNMT3A



Let's Discover Together.

© Oct 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/dnmt3a_methyl_malvern

