

SET8 Methyltransferase Assay Service

Activity-based SET8 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

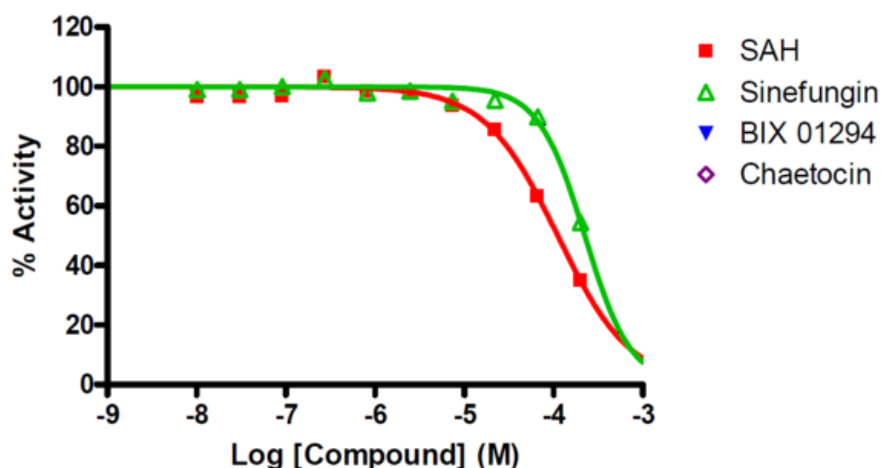
Target Overview

CAT#	SET8
Enzyme	Human SET8
Synonym	Histone-lysine N-methyltransferase SETD8, H4-K20-HMTase SETD8, SET domain-containing protein 8, PR/SET domain-containing protein 07, Lysine N-methyltransferase 5A, KMT5A, PRSET7, SET07
Substrates	Nucleosomes, 0.05 mg/ml SAM, 1 μM

Assay Properties

Reaction	S-adenosyl-L-[methyl- ³ H]methionine + histone L-lysine = S-adenosyl-L-homocysteine + histone N - [methyl- ³ H]-L-lysine
Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (nM) SAH 110,000 Sinefungin 220,000
Scientific Information	Q9NQR1 (http://www.uniprot.org/uniprot/Q9NQR1)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website Methyltransferase Assays .

Reference compound IC50 for SET8



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/set8_methyl_malvern

