

PRMT4 Methyltransferase Assay Service

Activity-based PRMT4 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

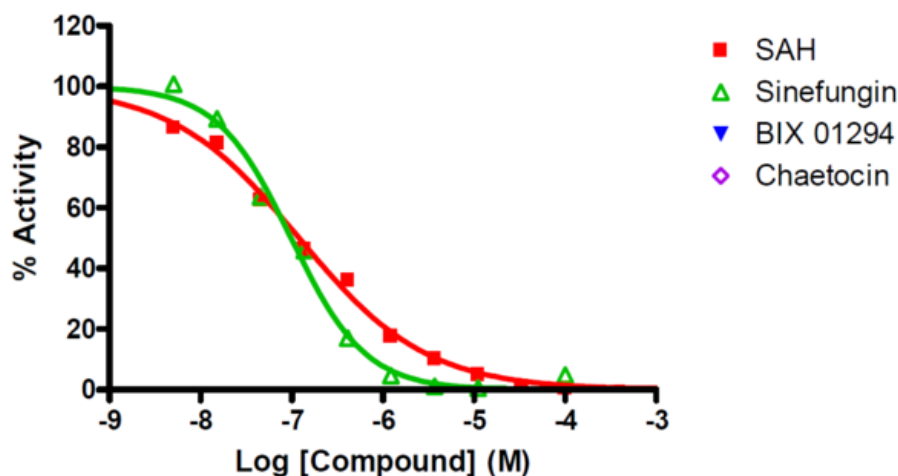
Target Overview

CAT#	PRMT4
Enzyme	Human PRMT4
Synonym	Protein arginine N-methyltransferase 4, Histone-arginine methyltransferase CARM1, Coactivator-associated arginine methyltransferase 1, CARM1
Substrates	Histone H3, 5 μ M SAM, 1 μ M

Assay Properties

Reaction	S-adenosyl-L-[methyl- 3 H]methionine + histone L-arginine = S-adenosyl-L-homocysteine + histone [methyl- 3 H]-L-arginine						
Readout	Enzymatic activity						
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>SAH</td> <td>120</td> </tr> <tr> <td>Sinefungin</td> <td>100</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	SAH	120	Sinefungin	100
Compounds	IC50 (nM)						
SAH	120						
Sinefungin	100						
Scientific Information	Q86X55 (http://www.uniprot.org/uniprot/Q86X55)						
Screening Location	Malvern, PA, USA						
Further Information	More information can be found on our website Methyltransferase Assays .						

Reference compound IC50 for PRMT4



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/prmt4_methyl_malvern

