

# PRDM9 Methyltransferase Assay Service

Activity-based PRDM9 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

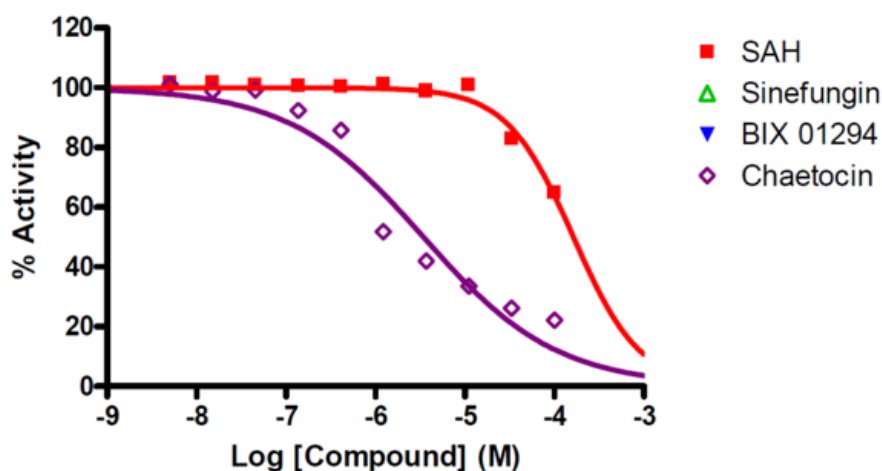
## Target Overview

<b>CAT#</b>	PRDM9
<b>Enzyme</b>	Human PRDM9
<b>Synonym</b>	Histone-lysine N-methyltransferase PRDM9, PR domain zinc finger protein 9, PR domain-containing protein 9, PFM6
<b>Substrates</b>	Histone H3, 5 $\mu$ M SAM, 1 $\mu$ M

## Assay Properties

<b>Reaction</b>	S-adenosyl-L-[methyl- $^3$ H]methionine + histone L-lysine = S-adenosyl-L-homocysteine + histone N-[methyl- $^3$ H]-L-lysine										
<b>Readout</b>	Enzymatic activity										
<b>Reference Compound IC50s</b>	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>SAH</td> <td>160,000</td> </tr> <tr> <td>Sinefungin</td> <td>&gt;200,000</td> </tr> <tr> <td>BIX 01294</td> <td>&gt;200,000</td> </tr> <tr> <td>Chaetocin</td> <td>3,400</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	SAH	160,000	Sinefungin	>200,000	BIX 01294	>200,000	Chaetocin	3,400
Compounds	IC50 (nM)										
SAH	160,000										
Sinefungin	>200,000										
BIX 01294	>200,000										
Chaetocin	3,400										
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/Q9NQV7">Q9NQV7</a> ( <a href="http://www.uniprot.org/uniprot/Q9NQV7">http://www.uniprot.org/uniprot/Q9NQV7</a> )										
<b>Screening Location</b>	Malvern, PA, USA										
<b>Further Information</b>	More information can be found on our website <a href="#">Methyltransferase Assays</a> .										

## Reference compound IC50 for PRDM9



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/prdm9\\_methyl\\_malvern](https://www.reactionbiology.com/datasheet/prdm9_methyl_malvern)

