

# NSD2 (T1150A) Methyltransferase Assay Service

Activity-based NSD2 (T1150A) Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

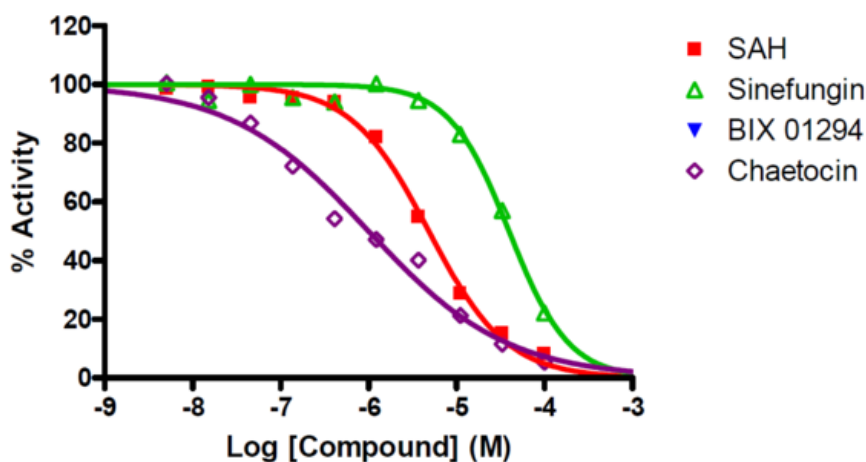
## Target Overview

<b>CAT#</b>	NSD2 (T1150A)
<b>Enzyme</b>	Human NSD2 (T1150A)
<b>Synonym</b>	Probable histone-lysine N-methyltransferase NSD2, Multiple myeloma SET domain-containing protein Nuclear SET domain-containing protein 2, Protein trithorax-5, Wolf-Hirschhorn syndrome candidate 1 protein, WHSC1, KIAA1090, MMSET, TRX5
<b>Substrates</b>	Nucleosomes, 0.05 mg/ml SAM, 1 μM

## Assay Properties

<b>Reaction</b>	S-adenosyl-L-[methyl- <sup>3</sup> H]methionine + histone L-lysine = S-adenosyl-L-homocysteine + histone [methyl- <sup>3</sup> 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>4,900</td> </tr> <tr> <td>Sinefungin</td> <td>40,000</td> </tr> <tr> <td>BIX01294</td> <td>&gt;100,000</td> </tr> <tr> <td>Chaetocin</td> <td>1,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	SAH	4,900	Sinefungin	40,000	BIX01294	>100,000	Chaetocin	1,000
Compounds	IC50 (nM)										
SAH	4,900										
Sinefungin	40,000										
BIX01294	>100,000										
Chaetocin	1,000										
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/O96028">O96028</a> ( <a href="http://www.uniprot.org/uniprot/O96028">http://www.uniprot.org/uniprot/O96028</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 NSD2 (T1150A)



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/nsd2\\_t1150a\\_methyl\\_malvern](https://www.reactionbiology.com/datasheet/nsd2_t1150a_methyl_malvern)

