

NRMT1 Methyltransferase Assay Service

Activity-based NRMT1 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

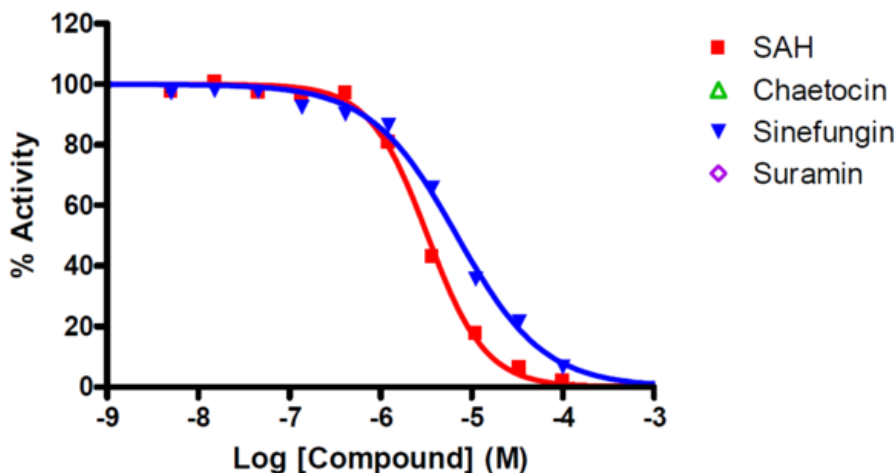
Target Overview

CAT#	NRMT1
Enzyme	Human NRMT1
Synonym	N-terminal Xaa-Pro-Lys N-methyltransferase 1, NTMT1, N-terminal RCC1 methyltransferase, Methyltransferase-like protein 11A, METTL11A
Substrates	RCC1, 5 μ M SAM, 1 μ M

Assay Properties

Reaction	S-adenosyl-L-[methyl- ³ H]methionine + N-terminal-A/S/PPK-protein = S-adenosyl-L-homocysteine + protein [methyl- ³ H]-L- lysine										
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>3,300</td> </tr> <tr> <td>Chaetocin</td> <td>>100,000</td> </tr> <tr> <td>Sinefungin</td> <td>6,900</td> </tr> <tr> <td>Suramin</td> <td>>100,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	SAH	3,300	Chaetocin	>100,000	Sinefungin	6,900	Suramin	>100,000
Compounds	IC50 (nM)										
SAH	3,300										
Chaetocin	>100,000										
Sinefungin	6,900										
Suramin	>100,000										
Scientific Information	Q9BV86 (http://www.uniprot.org/uniprot/Q9BV86)										
Screening Location	Malvern, USA										

Reference compound IC50 for NRMT1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/nrmt1_methyl_malvern

