

DOT1L Methyltransferase Assay Service

Activity-based DOT1L Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

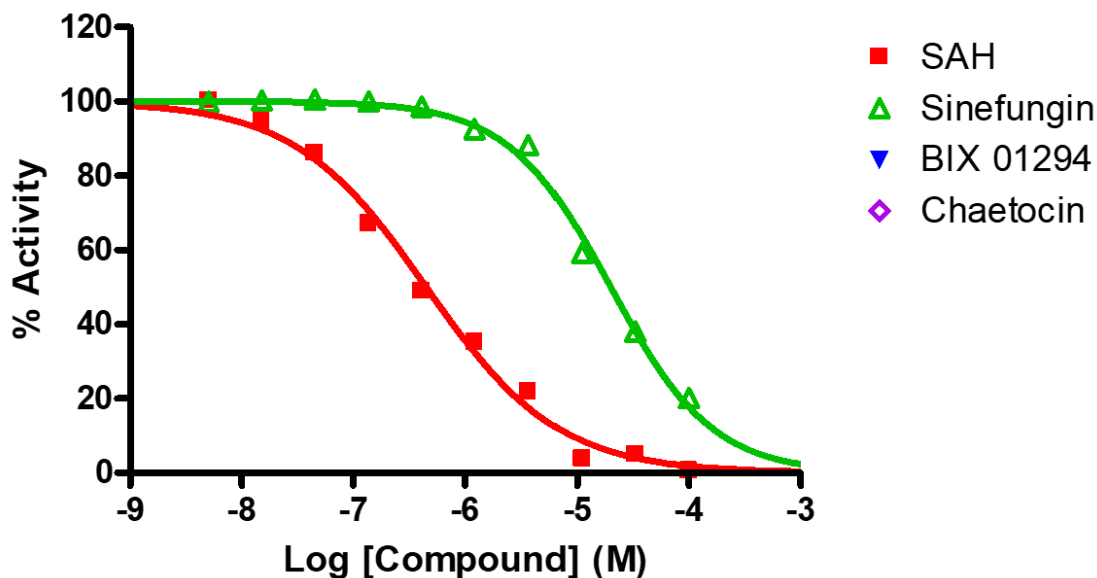
Target Overview

CAT#	DOT1L
Enzyme	Human DOT1L
Synonym	Histone-lysine N-methyltransferase, H3 lysine-79 specific, Histone H3-K79 methyltransferase, H3-K79-HMTase, DOT1-like protein, Lysine N-methyltransferase 4, KIAA1814, KMT4
Scientific Information	Q8TEK3 (http://www.uniprot.org/uniprot/Q8TEK3)

Assay Properties

Substrates	Nucleosomes, 0.05 mg/mL SAM, 1 μM						
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	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>SAH</td> <td>450</td> </tr> <tr> <td>Sinefungin</td> <td>20,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	SAH	450	Sinefungin	20,000
Compounds	IC50 (nM)						
SAH	450						
Sinefungin	20,000						
Screening Location	Malvern, PA, USA						
Further Information	More information can be found on our website Methyltransferase Assays .						

Reference compound IC50 for DOT1



Let's Discover Together.

© Sep 2023 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/dot1L_methyl_malvern

