

PRMT5 (C449S)/MEP50 Methyltransferase Assay

Service

Activity-based PRMT5 (C449S)/MEP50 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™](#) and [FlashPlate methyltransferase assay](#).

Target Overview

CAT#	PRMT5 (C449S)/MEP50
Enzyme	Human PRMT5 (C449S)/MEP50
Synonym	PRMT5: Protein arginine N-methyltransferase 5, Shk1 kinase-binding protein 1 homolog, Jak-binding protein 1, 72 kDa IκBn-binding protein, HRMT1L5, IBP72, JBP1, SKB1, ANM5 MEP50: Methylosome protein 50, Androgen receptor cofactor p44 WD repeat-containing protein 77, p44/Mep50, WDR77, HKMT1069
Substrates HotSpot™ assay:	Histone H2A, 5 μM SAM, 1 μM
Substrates FlashPlate assay:	H4 (1-15)-Biotin, 40 nM SAM, 1 μM

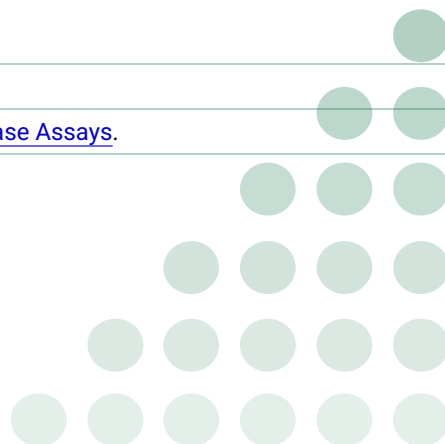
Assay Properties

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 HotSpot™ assay	Compounds IC50 (nM) SAH 1,000 Sinefugin 640 MTA 87 EPZ015666 2,600
Reference Compound IC50s FlashPlate Assay	Compounds IC50 (nM) SAH 1,100 Sinefugin 440 MTA 1,400 EPZ015666 14
Scientific Information	PRMT5; O14744 (http://www.uniprot.org/uniprot/O14744) MEP50; Q9BQA1 (http://www.uniprot.org/uniprot/Q9BQA1)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website Methyltransferase Assays .

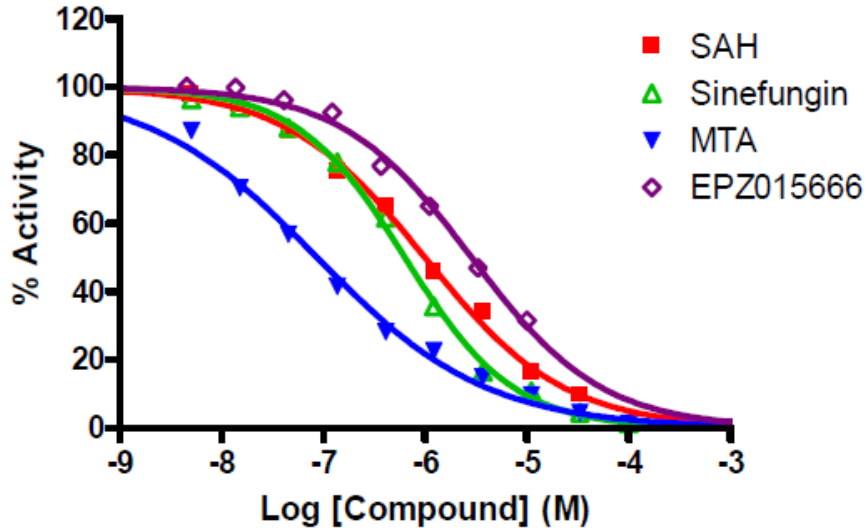
Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/prmt5_c449s_mep50_methyl_malvern



Reference compound IC50 for PRMT5(C449S)/MEP50 - HotSpot



Reference compound IC50 for PRMT5(C449S)/MEP50 - FlashPlate

