

PRMT1 Methyltransferase Assay Service

Activity-based PRMT1 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

Target Overview

| | |
|-------------------------------------|---|
| CAT# | PRMT1 |
| Enzyme | Human PRMT1 |
| Synonym | Protein arginine N-methyltransferase 1, Interferon receptor 1-bound protein 4, HMT2, HRMT1L2, IR1B4, ANM1 |
| Substrates HotSpot assay: | Histone H4, 5 μM; SAM, 1 μM |
| Substrates FlashPlate assay: | H4 (1-15)-Biotin, 40 nM; SAM, 1 μM |

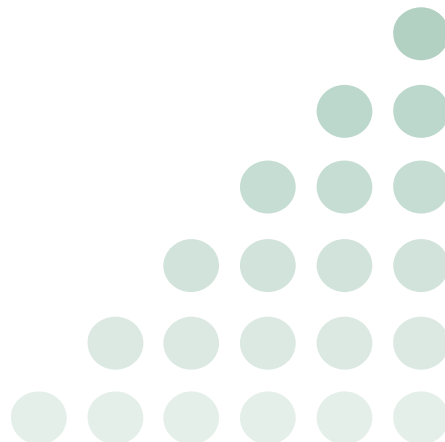
Assay Properties

| | | |
|--|--|------------------|
| 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 | 97 |
| | Sinefugin | 110 |
| | GSK 3368715 | 65 |
| Reference Compound IC50s FlashPlate Assay | Compounds | IC50 (nM) |
| | SAH | 5,500 |
| | Sinefugin | 720 |
| | GSK 3368715 | 2.3 |
| Scientific Information | Q99873 (http://www.uniprot.org/uniprot/Q99873) | |
| Screening Location | Malvern, USA | |

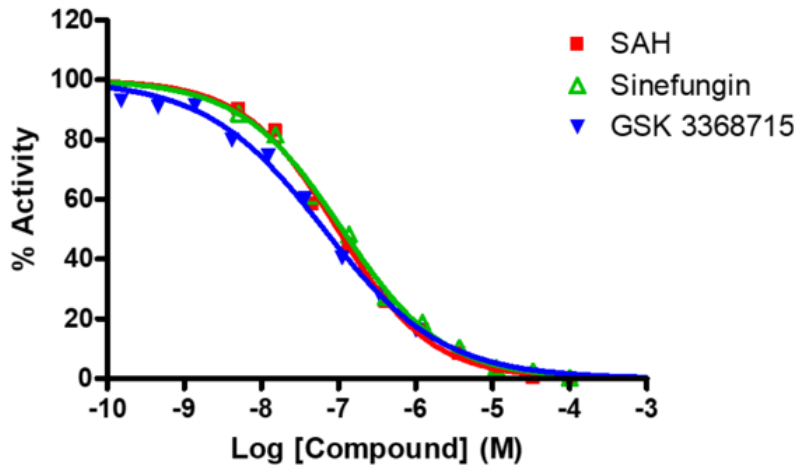
Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/prmt1_methyl_malvern



Reference Compound IC50 for PRMT1 - HotSpot -



Reference Compound IC50 for PRMT1 - FlashPlate -

