

DNMT1 Methyltransferase Assay Service

Activity-based DNMT1 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

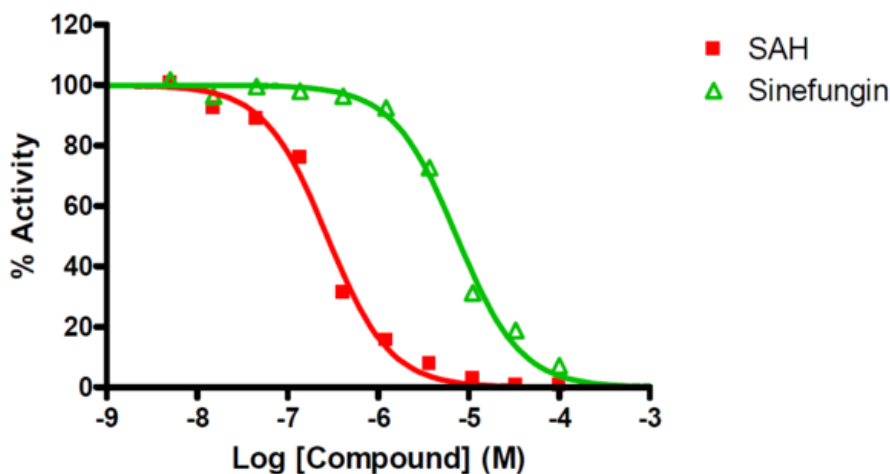
Target Overview

| | |
|-------------------|---|
| CAT# | DNMT1 |
| Enzyme | Human DNMT1 |
| Synonym | DNA (cytosine-5)-methyltransferase 1, CXXC-type zinc finger protein 9, DNA methyltransferase Hsa1, MCMT, AIM, CXXC9, DNMT |
| Substrates | Poly(dI-dC), 0.001 mg/ml SAM, 1 μM |

Assay Properties

| | |
|---------------------------------|--|
| Reaction | S-adenosyl-L-[methyl- ³ H]methionine + DNA = S-adenosyl-L-homocysteine + DNA 5-[methyl- ³ H]-cytosine |
| Readout | Enzymatic activity |
| Reference Compound IC50s | Compounds IC50 (nM) SAH 270 Sinefungin 7,400 |
| Scientific Information | P26358 (http://www.uniprot.org/uniprot/P26358) |
| Screening Location | Malvern, PA, USA |
| Further Information | More information can be found on our website Methyltransferase Assays . |

Reference compound IC50 for DNMT1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/dnmt1_methyl_malvern

