

SMYD2 Methyltransferase Assay Service

Activity-based SMYD2 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

Target Overview

CAT#	SMYD2
Enzyme	Human SMYD2
Synonym	SET and MYND domain-containing protein 2, HSKM-B Lysine N-methyltransferase 3C, KMT3C
Substrates	Histone H3, 5 μ M SAM, 1 μ M

Assay Properties

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	Compounds IC50 (nM) SAH 526 Sinefungin 15,000
Scientific Information	Q9NRG4 (http://www.uniprot.org/uniprot/Q9NRG4)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website Methyltransferase Assays .

Reference compound IC50 for SMYD2

