

MLL4 Methyltransferase Assay Service

Activity-based MLL4 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

Target Overview

CAT#	MLL4
Enzyme	Human MLL4; 5 component complex
Synonym	Complex of MLL4: Histone-lysine N-methyltransferase MLL4, WW domain-binding protein 7, HRX2, KIAA0304, KMT2D, TRX2, WBP7 WDR5: WD repeat-containing protein 5, BMP2-induced 3-kb gene protein RBBP5: Retinoblastoma-binding protein 5, Retinoblastoma-binding protein RBQ-3 ASH2L: Set1/Ash2 histone methyltransferase complex subunit ASH2, ASH2-like protein DPY30: Protein dpy-30 homolog, Dpy-30-like protein
Substrates	Core Histone, 0.05 mg/ml SAM, 1 μM

Assay Properties

Reaction	S-adenosyl-L-[methyl- ³ H]methionine + histone L-lysine = S-adenosyl-L-homocysteine + histone [methyl- ³ H]-L-lysine										
Readout	Enzymatic activity										
Reference Compound IC50s	<table border="1"> <thead> <tr> <th colspan="2">Compounds IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>SAH</td> <td>8,300</td> </tr> <tr> <td>Sinefugin</td> <td>42,000</td> </tr> <tr> <td>BIX01294</td> <td>>500,000</td> </tr> <tr> <td>Chaetocin</td> <td>310</td> </tr> </tbody> </table>	Compounds IC50 (nM)		SAH	8,300	Sinefugin	42,000	BIX01294	>500,000	Chaetocin	310
Compounds IC50 (nM)											
SAH	8,300										
Sinefugin	42,000										
BIX01294	>500,000										
Chaetocin	310										
Scientific Information	MLL4 Q9UMN6 (http://www.uniprot.org/uniprot/Q9UMN6) WDR5 P61964 (http://www.uniprot.org/uniprot/P61964) RBBP5 Q15291 (http://www.uniprot.org/uniprot/Q15291) ASH2L Q9UBL3 (http://www.uniprot.org/uniprot/Q9UBL3) DPY30 Q9C005 (http://www.uniprot.org/uniprot/Q9C005)										
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/ml4_complex_methyl_malvern



Reference compound IC50 for MLL4

