

# MLL2 Methyltransferase Assay Service

Activity-based MLL2 Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

## Target Overview

<b>CAT#</b>	MLL2
<b>Enzyme</b>	Human MLL2; 5 component complex
<b>Synonym</b>	Complex of MLL2: Histone-lysine N-methyltransferase MLL2, ALL1-related protein, ALR, KMT2B 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
<b>Substrates</b>	Core Histone, 0.05 mg/ml SAM, 1 $\mu$ M

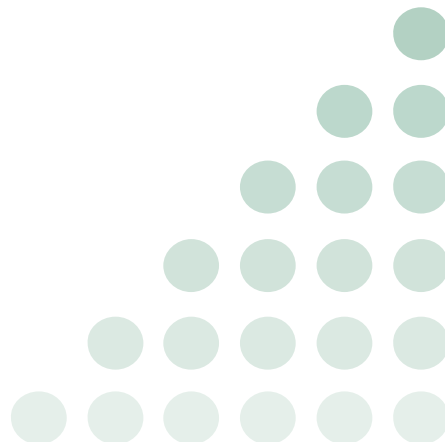
## Assay Properties

<b>Reaction</b>	S-adenosyl-L-[methyl- <sup>3</sup> H]methionine + histone L-lysine = S-adenosyl-L-homocysteine + histone [methyl- <sup>3</sup> H]-L-lysine										
<b>Readout</b>	Enzymatic activity										
<b>Reference Compound IC50s</b>	<table border="1"> <thead> <tr> <th colspan="2">Compounds IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>SAH</td> <td>22,500</td> </tr> <tr> <td>Sinefugin</td> <td>38,000</td> </tr> <tr> <td>BIX01294</td> <td>&gt;500,000</td> </tr> <tr> <td>Chaetocin</td> <td>400</td> </tr> </tbody> </table>	Compounds IC50 (nM)		SAH	22,500	Sinefugin	38,000	BIX01294	>500,000	Chaetocin	400
Compounds IC50 (nM)											
SAH	22,500										
Sinefugin	38,000										
BIX01294	>500,000										
Chaetocin	400										
<b>Scientific Information</b>	MLL2 <a href="http://www.uniprot.org/uniprot/O14686">O14686</a> ( <a href="http://www.uniprot.org/uniprot/O14686">http://www.uniprot.org/uniprot/O14686</a> ) WDR5 <a href="http://www.uniprot.org/uniprot/P61964">P61964</a> ( <a href="http://www.uniprot.org/uniprot/P61964">http://www.uniprot.org/uniprot/P61964</a> ) RBBP5 <a href="http://www.uniprot.org/uniprot/Q15291">Q15291</a> ( <a href="http://www.uniprot.org/uniprot/Q15291">http://www.uniprot.org/uniprot/Q15291</a> ) ASH2L <a href="http://www.uniprot.org/uniprot/Q9UBL3">Q9UBL3</a> ( <a href="http://www.uniprot.org/uniprot/Q9UBL3">http://www.uniprot.org/uniprot/Q9UBL3</a> ) DPY30 <a href="http://www.uniprot.org/uniprot/Q9C005">Q9C005</a> ( <a href="http://www.uniprot.org/uniprot/Q9C005">http://www.uniprot.org/uniprot/Q9C005</a> )										
<b>Screening Location</b>	Malvern, PA, USA										
<b>Further Information</b>	More information can be found on our website <a href="#">Methyltransferase Assays</a> .										

Let's Discover Together.

© Oct 2022 Reaction Biology. All rights reserved.

[https://www.reactionbiology.com/datasheet/ml2\\_complex\\_methyl\\_malvern](https://www.reactionbiology.com/datasheet/ml2_complex_methyl_malvern)



Reference compound IC50 for MLL2

