

# EZH2 (Y641F) Methyltransferase Assay Service

Activity-based EZH2 (Y641F) Methyltransferase Assay for compound screening and profiling via [radiometric HotSpot™ methyltransferase assay](#).

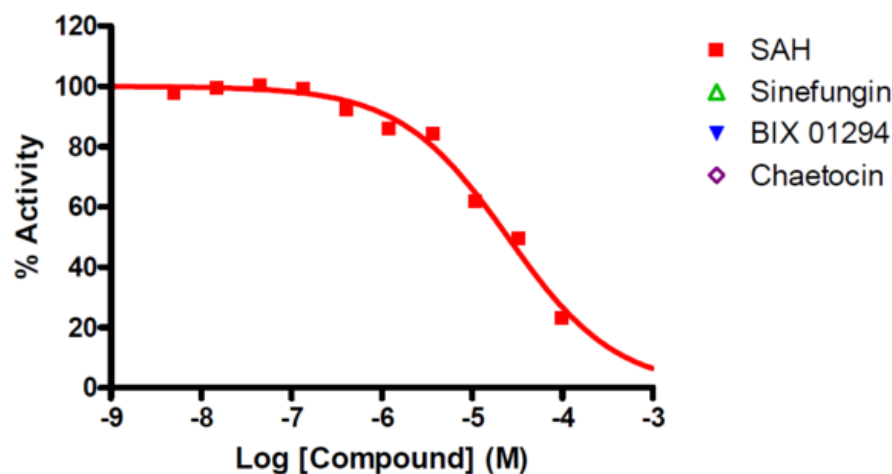
## Target Overview

<b>CAT#</b>	EZH2 (Y641F)
<b>Enzyme</b>	Human EZH2 (Y641F); 5 component complex
<b>Synonym</b>	EZH2: Polycomb protein EZH2, Enhancer of zeste homolog 2, ENX-1, KMT6, EED: WD protein associating with integrin cytoplasmic tails 1, SUZ12: Suppressor of zeste 12 protein homolog, RBAP48: Histone-binding protein RBBP4, AEBP2: Zinc finger protein AEBP2
<b>Substrates</b>	Core Histone, 0.05 mg/mlSAM, 1 μM

## Assay Properties

<b>Reaction</b>	S-adenosyl-L-[methyl- <sup>3</sup> H]methionine + histone L-lysine = S-adenosyl-L-homocysteine + histone N -[methyl- <sup>3</sup> H]-L-lysine	
<b>Readout</b>	Enzymatic activity	
<b>Reference Compound IC50s</b>	<b>Compounds</b>	<b>IC50 (nM)</b>
	SAH	26,000
	Sinefungin	>200,000
	BIX01294	>200,000
	Chaetocin	>200,000
<b>Scientific Information</b>	EZH2 <a href="http://www.uniprot.org/uniprot/Q15910">Q15910</a> ( <a href="http://www.uniprot.org/uniprot/Q15910">http://www.uniprot.org/uniprot/Q15910</a> )EED <a href="http://www.uniprot.org/uniprot/O75530">O75530</a> ( <a href="http://www.uniprot.org/uniprot/O75530">http://www.uniprot.org/uniprot/O75530</a> )SUZ12 <a href="http://www.uniprot.org/uniprot/Q15022">Q15022</a> ( <a href="http://www.uniprot.org/uniprot/Q15022">http://www.uniprot.org/uniprot/Q15022</a> )RBBP4 <a href="http://www.uniprot.org/uniprot/Q09028">Q09028</a> ( <a href="http://www.uniprot.org/uniprot/Q09028">http://www.uniprot.org/uniprot/Q09028</a> )AEBP2 <a href="http://www.uniprot.org/uniprot/Q6ZN18">Q6ZN18</a> ( <a href="http://www.uniprot.org/uniprot/Q6ZN18">http://www.uniprot.org/uniprot/Q6ZN18</a> )	
<b>Screening Location</b>	Malvern, USA	

## Reference compound IC50 for EZH2 (Y641F)



Let's Discover Together.

