

# MYST4 Histone Acetyltransferase Assay Service

Activity-based MYST4 Histone Acetyltransferase Assay for compound screening and profiling via [radiometric HAT assay](#).

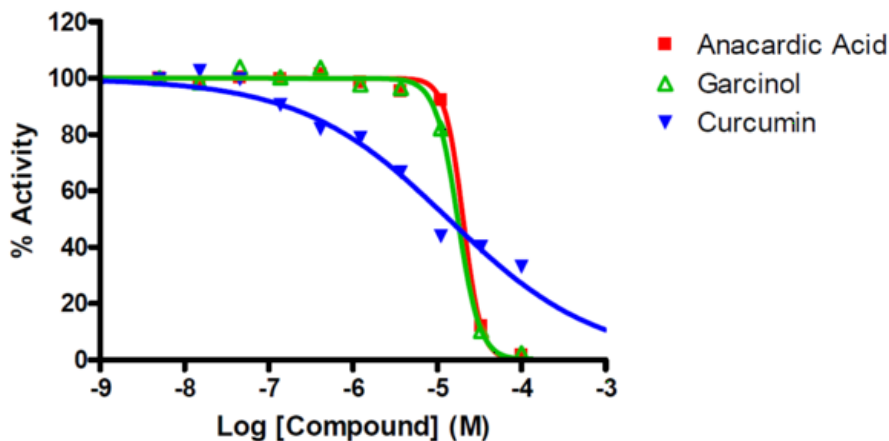
## Target Overview

<b>CAT#</b>	MYST4
<b>Enzyme</b>	Human MYST4
<b>Synonyms</b>	Histone acetyltransferase KAT6B, MOZ, YBF2/SAS3, SAS2 and TIP60 protein 4, Monocytic leukemia zinc finger protein-related factor, KIAA0383, MORF, MOZ2
<b>Substrates</b>	Histone H4, 5 $\mu$ M Acetyl CoA, 3 $\mu$ M

## Assay Properties

<b>Reaction</b>	[Acetyl- $^3$ H]-Acetyl Coenzyme A + histone L-lysine = Coenzyme A + [Acetyl- $^3$ H]-L-lysine								
<b>Readout</b>	Enzymatic activity								
<b>Reference Compound IC50s</b>	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Anacardic Acid</td> <td>20,000</td> </tr> <tr> <td>Garcinol</td> <td>17,500</td> </tr> <tr> <td>Curcumin</td> <td>14,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Anacardic Acid	20,000	Garcinol	17,500	Curcumin	14,000
Compounds	IC50 (nM)								
Anacardic Acid	20,000								
Garcinol	17,500								
Curcumin	14,000								
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/Q8WYB5">Q8WYB5</a> ( <a href="http://www.uniprot.org/uniprot/Q8WYB5">http://www.uniprot.org/uniprot/Q8WYB5</a> )								
<b>Screening Location</b>	Malvern, USA								

## Reference Compound IC50 for MYST4



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/myst4\\_hat\\_malvern](https://www.reactionbiology.com/datasheet/myst4_hat_malvern)

