

GCN5 Histone Acetyltransferase Assay Service

Activity-based GCN5 Histone Acetyltransferase Assay for compound screening and profiling via [radiometric HAT assay](#).

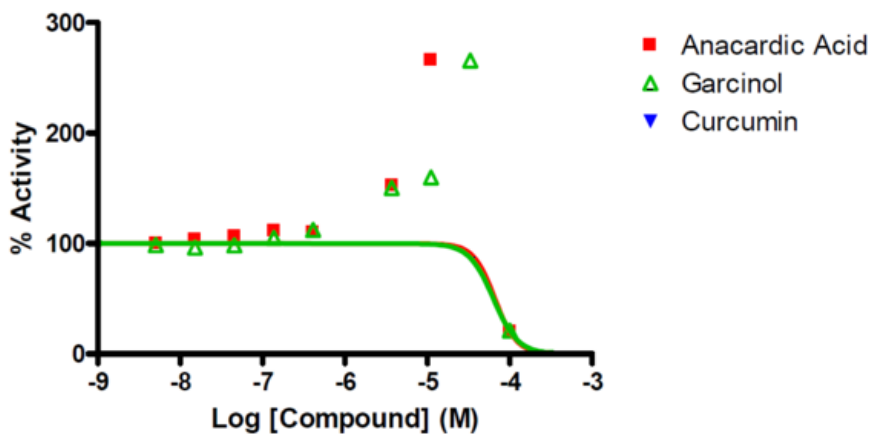
Target Overview

CAT#	GCN5
Enzyme	Human GCN5
Synonyms	Histone acetyltransferase KAT2A, Lysine acetyltransferase 2A Histone acetyltransferase GCN5, General control of amino acid synthesis protein 5-like 2, STAF97, GCN5L2, HGCN5, KAT2A
Substrates	Histone H3, 5 μ M Acetyl CoA, 3 μ M

Assay Properties

Reaction	[Acetyl- ³ H]-Acetyl Coenzyme A + histone L-lysine = Coenzyme A + [Acetyl- ³ H]-L-lysine						
Readout	Enzymatic activity						
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Anacardic Acid</td> <td>66,000</td> </tr> <tr> <td>Garcinol</td> <td>64,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Anacardic Acid	66,000	Garcinol	64,000
Compounds	IC50 (nM)						
Anacardic Acid	66,000						
Garcinol	64,000						
Scientific Information	Q92830 (http://www.uniprot.org/uniprot/Q92830)						
Screening Location	Malvern, USA						

Reference Compound IC50 for GCN5



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/hgcn5_hat_malvern

