

CBP Histone Acetyltransferase Assay Service

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

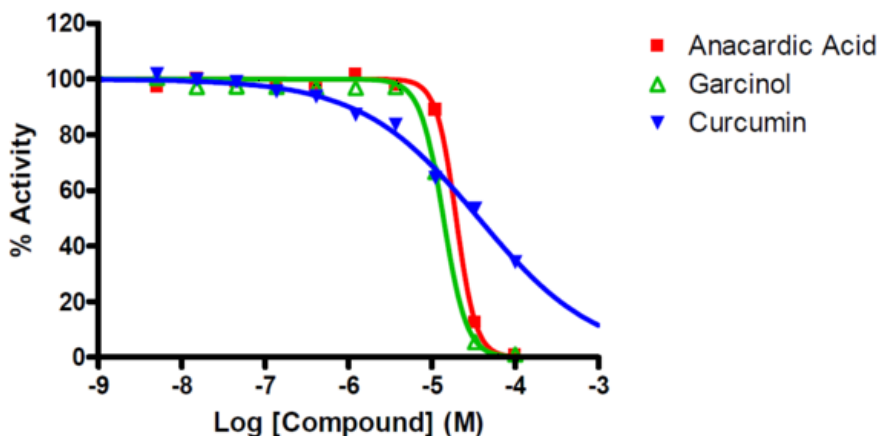
Target Overview

CAT#	CBP
Enzyme	Human CBP
Synonyms	CREB-binding protein, CREBBP
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>20,000</td> </tr> <tr> <td>Garcinol</td> <td>14,000</td> </tr> <tr> <td>Curcumin</td> <td>37,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Anacardic Acid	20,000	Garcinol	14,000	Curcumin	37,000
Compounds	IC50 (nM)								
Anacardic Acid	20,000								
Garcinol	14,000								
Curcumin	37,000								
Scientific Information	Q92793 (http://www.uniprot.org/uniprot/Q92793)								
Screening Location	Malvern, USA								

Reference Compound IC50 for CBP



Let's Discover Together.

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

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

