

HAT Assay Kit

Fluorescent

(version B1)

Catalog No. 56100

Active Motif North America

1914 Palomar Oaks Way, Suite 150

Carlsbad, California 92008, USA

Toll free: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

Active Motif Europe

104 Avenue Franklin Roosevelt

B-1330 Rixensart, Belgium

UK Free Phone: 0800 169 31 47

France Free Phone: 0800 90 99 79

Germany Free Phone: 0800 181 99 10

Telephone: +32 (0)2 653 0001

Fax: +32 (0)2 653 0050

Active Motif Japan

Azuma Bldg, 7th Floor

2-21 Ageba-Cho, Shinjuku-Ku

Tokyo, 162-0824, Japan

Telephone: +81 3 5225 3638

Fax: +81 3 5261 8733

Copyright 2009 Active Motif, Inc.

Information in this manual is subject to change without notice and does not constitute a commitment on the part of Active Motif, Inc. It is supplied on an “as is” basis without any warranty of any kind, either explicit or implied. Information may be changed or updated in this manual at any time.

This documentation may not be copied, transferred, reproduced, disclosed, or duplicated, in whole or in part, without the prior written consent of Active Motif, Inc. This documentation is proprietary information and protected by the copyright laws of the United States and international treaties.

The manufacturer of this documentation is Active Motif, Inc.

© 2009 Active Motif, Inc., 1914 Palomar Oaks Way, Suite 150; Carlsbad, CA 92008. All rights reserved.

All trademarks, trade names, service marks or logos referenced herein belong to their respective companies.

TABLE OF CONTENTS	Page
Overview	1
Introduction	2
Kit Performance and Benefits	3
Kit Components and Storage	
Additional Materials Required	4
Reagent Preparation	
Assay Buffer AM1	5
Acetyl-CoA	5
Recombinant p300, catalytic domain	5
Histone H3 Peptide	5
Histone H4 Peptide	5
Anacardic Acid	5
DMSO	5
Stop Solution (Fluor)	5
Developer (Fluor)	5
Protocols	
Preparation of Standard Curve	6
Assay Protocol	7
Data Analysis	8
References	8
Appendix	
Section A. Troubleshooting Guide	9
Section B. Related Products	10
Technical Services	12

Overview

Active Motif's fluorescent HAT Assay Kit is an easy and sensitive assay to either determine the activity of your own source of purified histone acetyltransferases (HATs) or to screen potential inhibitors of HAT activity.

The HAT Assay Kit contains a purified recombinant active p300 catalytic domain protein, histone H3 and H4 N-terminal substrate peptides, anacardic acid as a control inhibitor, and all the essential buffers for a full 96-well plate assay.

This fluorescent HAT Assay works in three easy steps: first, the HAT protein is incubated with acetyl-CoA and either a histone H3 or histone H4 substrate peptide. During this reaction, the HAT enzyme catalyzes the enzymatic transfer of acetyl groups from the acetyl-CoA to the histone peptide thereby generating the acetylated peptide and CoA-SH. The reaction is stopped by adding stop solution, and after adding the complete developing solution, the developer reacts with the free sulfhydryl groups on the CoA-SH to give a fluorescent reading of acetyltransferase activity.

product	format	catalog no.
HAT Assay Kit (Fluorescent)	1 x 96 rxns	56100

Introduction

Histone Acetylation

The basic structural unit of chromatin is the nucleosome, which consists of 146 base pairs (bp) of DNA wrapped around a histone octamer. The histone octamer consists of two copies each of the core histone H2A-H2B dimers and H3-H4 dimers, and the N-terminal tails of the histones are free from the bound core of the nucleosome structure where they can be modified by protein interactions in the nucleus. Histones, especially the N-terminal tails, are highly conserved from yeast to humans, and the modifications are commonly acetylation, phosphorylation and methylation. The modifications that occur at specific amino acid residues on the histone tails are very important for downstream signaling events and are of much interest in recent studies¹.

Histone acetylation is a modification that effects the nucleosome structure and therefore the ability of transcription factors to access the DNA and regulate gene expression. Recent studies have found that hyperacetylation of histones can be associated with relaxed chromatin structure and active gene transcription, while hypoacetylated histones can lead to transcriptional repression by condensing the structure of chromatin and restricting the access of transcription factors to the DNA².

Histone acetylation is a reversible process. Two superfamilies of enzymes are involved in histone acetylation: histone acetyltransferases (HAT) and histone deacetylases (HDAC), which catalyze the addition and removal of acetyl groups on histones. Histone acetyltransferases (HAT) are enzymes that catalyze the acetylation of lysine residues in the N-terminal tails of histone proteins.

The first nuclear HATs discovered were the general control of nuclear-5 (GCN5), and its paralog p300/CBP accessory factor (PCAF). Both are conserved in organisms from yeast to humans, and GCN5/PCAF HATs have served as paradigms for the biochemical analysis of HAT structure, mechanism and function³. In addition to histones, GCN5/PCAF HATs target non-histone protein substrates, such as transcription factors (*i.e.*, p53) and structural proteins (*i.e.* α -tubulin). More than 20 HATs have been identified and these have been classified into one of five families (GNAT1, MYST (MOZ, YBF2/Sas3, Sas2, Tip60), TAFII250, nuclear receptor co-activators, and p300/cAMP response element-binding protein (CBP)). To date, a small number of HAT inhibitors have been reported. A few natural small molecules such as anacardic acid, garcinol and curcumin have been described as potent p300 and PCAF inhibitors⁴.

Kit Performance and Benefits

The HAT Assay Kit (Fluorescent) is for research use only. Not for use in diagnostic procedures.

Recombinant p300 catalytic domain protein, active

p300 catalytic domain is a purified recombinant human protein expressed in insect cells. p300 catalytic domain is supplied at 0.1 mg/ml. This recombinant protein corresponds to amino acids 965-1810 from the GenBank sequence accession number U01877.

Anacardic Acid

Anacardic acid is supplied at 28.7 mM in DMSO. Anacardic acid is a small molecule compound from cashew nut shell liquid. 15 μM anacardic acid has been shown to quench HAT activity more than 90%. The IC_{50} value of anacardic acid for p300 using histones as substrates was found to be 8.5 μM and the IC_{50} value of anacardic acid for p300 using H3 peptide as the substrate was found to be 500 nM⁴.

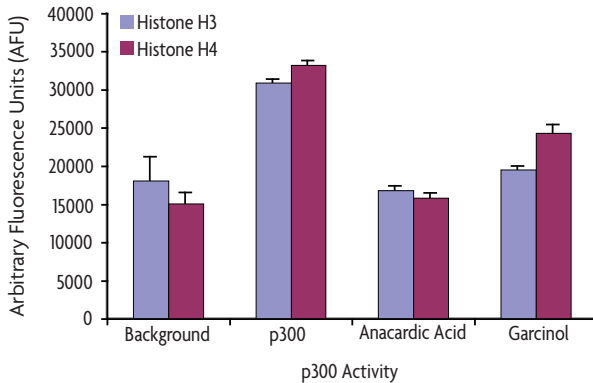


Figure 1: Effect of inhibitors on p300 HAT activity.

Fluorescent HAT Assay Kit results using 50 ng p300 catalytic domain per well with 50 μM Histone H3 peptide, 50 μM Acetyl-CoA, 15 μM Anacardic Acid, 25 μM Garcinol and 30 minute incubation at room temperature. The background fluorescence was subtracted from the data.

Kit Components and Storage

The HAT Assay Kit (Fluorescent) is for research use only. Not for use in diagnostic procedures. Store components at the temperatures indicated below. All components are guaranteed stable for six months when stored at the indicated temperature.

Reagent	Quantity	Storage / Stability
5X Assay Buffer AMI	1.8 ml	-20°C to room temp.
Acetyl-CoA	105 µl	-80°C
Recombinant p300 catalytic domain	50 µl (0.1 mg/ml)	-80°C
Histone H3 Peptide (MW=2068 Da)	0.275 µmoles (568 µg)	-20°C
Histone H4 Peptide (MW=1905 Da)	0.275 µmoles (524 µg)	-20°C
Stop Solution (Fluor)	5 ml	room temp.
Developer (Fluor), 120X	84 µl	-20°C
Developing Solution (Fluor)	10 ml	-20°C to room temp.
96-well plate	1 plate	room temp.
Anacardic Acid	4 µl (10 mg/ml)	-80°C to -20°C
Plate sealer	1	room temp.

Additional materials required

- β-mercaptoethanol or Coenzyme A
- DMSO
- multi-channel pipette
- multi-channel pipette reservoir
- Fluorescent plate reader capable of exciting at 360-390 nm and reading emission at 460-490 nm

Reagent Preparation

5X Assay Buffer AM1

Prior to the assay, dilute to 1X with sterile water. Allow solution to warm to room temperature before use. This assay buffer is used in the assay and for diluting the Acetyl-CoA, Histone peptide substrates, and inhibitors. Prepare the quantity required for your assay including what you need to dilute the components. The diluted 1X Assay Buffer AM1 is stable for 3 months.

Acetyl-CoA

Acetyl-CoA is provided as a 6 mM stock solution. Prior to the assay, dilute 1:12 with the 1X Assay Buffer to a final concentration of 0.5 mM. 5 μ l are used per assay well. Use freshly diluted Acetyl-CoA for each assay.

Recombinant p300 catalytic domain

Avoid multiple freeze-thaws of the protein. Please aliquot and store frozen at -80°C if only a portion of the enzyme is needed for your assays.

Histone H3 Peptide and Histone H4 Peptide

Peptides are supplied lyophilized. Rehydrate each peptide with 2.2 ml of the 1X Assay Buffer to obtain a 125 μ M stock solution (125 pmol/ μ l). The peptide is supplied in a tube with a 1.5 ml volume capacity. Start by adding a portion of the 2.2 ml 1X Assay Buffer to transfer the peptide to a new tube to resuspend in a 2.2 ml final volume. After the peptide is in solution, store at -20°C.

Use 20 μ l per assay well for a 50 μ M final concentration in the reaction. This concentration may be reduced or increased with the 1X Assay Buffer if another final concentration is preferable.

Anacardic Acid

Anacardic Acid is supplied at 28.7 mM in DMSO. Dilute in 1X Assay Buffer AM1 to make a 150 μ M solution. For example, dilute 1 μ l of Anacardic Acid stock solution in 190 μ l 1X Assay Buffer. For a 15 μ M Anacardic Acid treatment, add 5 μ l of the 150 μ M solution per assay well.

DMSO

Prepare a DMSO dilution equivalent to the Anacardic Acid dilution to serve as a background control within the assay. Add 1 μ l of DMSO to 190 μ l of 1X Assay Buffer. Use 5 μ l of diluted DMSO per assay well.

Stop Solution (Fluor) Ready to use as supplied.

Developer (Fluor)

Developer stock is 3 mM in DMSO. Prior to use, dilute the developer 1:120 with Developing Solution to obtain a final developing solution concentration of 25 μ M. Protect the Developer and final Developing Solution from light. The final diluted developing solution is stable for 6 hours at room temperature.

Note: The assay will not work if the developer is diluted in assay buffer instead of developing solution.

Protocol

PLEASE READ THE ENTIRE PROTOCOL BEFORE STARTING!

Minimize the use of sample buffers containing DTT or β -mercaptoethanol for your HAT source. Buffers containing these reducing agents or other sources of reactive thiol groups will result in high background.

This protocol has been optimized for use with recombinant proteins. If you wish to calculate enzymatic activity, a standard curve must be prepared.

It is recommended to run test samples, controls, standards and background wells at least in duplicate.

Preparation of Standard Curve (Optional)

If you want to calculate enzymatic activity, you will need to run a standard curve alongside your samples in the assay. A standard curve may be prepared with Coenzyme A or β -mercaptoethanol.

For example, to prepare a standard curve with β -mercaptoethanol, make dilutions in duplicate ranging from 0 to 14 μ M in 2 μ M increments in a 50 μ l volume to result in a 0 to 500 pmol range for the X-axis (Figure 1). The volumes below are enough for duplicate wells.

	0 μ M	2 μ M	4 μ M	6 μ M	8 μ M	10 μ M	12 μ M	14 μ M
1X Assay Buffer	120 μ l	117.6 μ l	115.2 μ l	112.8 μ l	110.4 μ l	108 μ l	105.6 μ l	103.2 μ l
β -mercaptoethanol	0 μ l	2.4 μ l	4.8 μ l	7.2 μ l	9.6 μ l	12 μ l	14.4 μ l	16.8 μ l

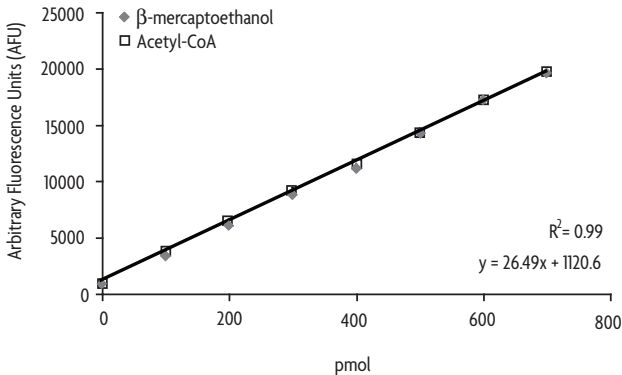


Figure 1: Standard Curve Data.

Example standard curve data of both acetyl-CoA and β -mercaptoethanol diluted from 0 to 14 μ M. pmol of β -mercaptoethanol = μ M of β -mercaptoethanol X 50 μ l per well

Run standard curve wells by adding 50 μ l of each dilute per well in duplicate and continue with step 4 in the following protocol.

Assay Protocol

1. Design your assay to test all samples in duplicate. Prepare the quantity required of 1X Assay Buffer, diluted Acetyl-CoA and desired inhibitors. If using the Anacardic acid inhibitor provided in the kit, prepare a DMSO dilution for background comparison. Prepare the amount of final developing solution required and keep protected from light. The HAT enzyme and Acetyl-CoA should be kept on ice.
2. Using any desired microcentrifuge tubes for samples including background, positive control and inhibition wells, gently mix the following components together in the quantities indicated. 30 μ l of each mix is needed per assay well. Keep the reactions at room temperature. The volumes listed in the following table are enough for duplicate wells.

Reagent	Background Wells	Positive Control Wells	Inhibitor Wells	Standard Curve Wells
1X Assay Buffer	48.75 μ l	48.75 μ l	48.75 μ l	
0.5 mM Acetyl-CoA	12.5 μ l	12.5 μ l	12.5 μ l	
150 μ M Anacardic Acid			12.5 μ l	
Diluted DMSO	12.5 μ l	12.5 μ l		
p300 catalytic domain (0.1 mg/ml)	1.25 μ l	1.25 μ l	1.25 μ l	
β -mercaptoethanol standard				120 μ l
Total volume per well	30 μl	30 μl	30 μl	50 μl

3. Add 20 μ l of Histone H3 or Histone H4 peptide to each assay well. Do not add peptide substrate to background wells so that autoacetylation activity of the acetyltransferase is included in the background measurement. Mix solution by pipetting contents up-and-down with multichannel pipettor. Cover plate with plate sealer and incubate for 10 to 30 minutes at room temperature.

Note: If you are determining the activity of a new HAT protein, we recommend running a time course to establish the correct incubation time for your protein.

4. Add 50 μ l of Stop Solution per well. To background wells, add 20 μ l peptide substrate after adding the stop solution.
5. Add 100 μ l final Developing Solution to each well and incubate 15 minutes in the dark at room temperature.
6. Read fluorescence with excitation at 360-390 nm and emission at 450 to 470 nm.

Data Analysis

Plot the standard curve data with fluorescence units on the Y-axis and pmol of β -mercaptoethanol on the X-axis.

Determine slope of the line through the linear part of the curve as arbitrary fluorescence units per pmol. Subtract the fluorescence of background from that of the assay samples and divide the difference by the incubation time (AFU/minute). Divide the AFU/minute by the slope of the standard curve (AFU/pmol) to obtain the rate in pmol/minute.

References

1. Trievel, R. C., et al., (2000) *Analytical Biochemistry* 287: 319-328.
2. Wang, L., et al., (1997) *Mol. and Cell. Biol.* 17(1): 519-527.
3. Marmorstein, R. (2001) *Cell. Mol. Life Sci.* 58: 693-703.
4. Balasubramanyam K, et al., (2003) *J. Biol. Chem.* 278(21):19134-19140.

Appendix

Section A. Troubleshooting Guide

Problem/question	Recommendation
Must the assay be incubated at room temperature?	The assay may be performed at 30 °C or room temperature without effecting results.
Is it possible to use diluted Acetyl-CoA?	No. Fresh Acetyl-CoA should be prepared for each assay.
Is it necessary to run a standard curve?	No. The standard curve is only necessary if the specific activity needs to be calculated (pmol/min/ μ g).
Why was no HAT activity detected?	The recombinant p300 supplied in this kit may lose activity after 3 freeze-thaws. Please aliquot at first use to store and reduce the number of freeze-thaws.
	Check that the developer has been diluted into developing solution. Do not use Assay Buffer AM1 to dilute!
Why is there high background?	The background is due to the free sulfhydryl groups in the formulation buffer. Some background may be due to the protein's auto-acetylation activity. Run controls with and without acetyl-CoA to determine background and auto-acetylation levels.

Section B. Related Products

Histone ELISAs	Format	Catalog No.
Histone H3 monomethyl Lys4 ELISA	1 x 96 rxns	53101
Histone H3 dimethyl Lys4 ELISA	1 x 96 rxns	53102
Histone H3 trimethyl Lys4 ELISA	1 x 96 rxns	53103

Recombinant Methylated Histones	Format	Catalog No.
Recombinant Histone H3 (C110A)	50 µg	31207
Recombinant Histone H3 monomethyl Lys4	50 µg	31208
Recombinant Histone H3 dimethyl Lys4	50 µg	31209
Recombinant Histone H3 trimethyl Lys4	50 µg	31210
Recombinant Histone H3 monomethyl Lys9	50 µg	31211
Recombinant Histone H3 dimethyl Lys9	50 µg	31212
Recombinant Histone H3 trimethyl Lys9	50 µg	31213
Recombinant Histone H3 monomethyl Lys27	50 µg	31214
Recombinant Histone H3 dimethyl Lys27	50 µg	31215
Recombinant Histone H3 trimethyl Lys27	50 µg	31216
Recombinant Histone H4	50 µg	31223
Recombinant Histone H4 monomethyl Lys20	50 µg	31224
Recombinant Histone H4 dimethyl Lys20	50 µg	31225
Recombinant Histone H4 trimethyl Lys20	50 µg	31226

Control Acid Extracts	Format	Catalog No.
HeLa acid extract	100 µg	36200
HeLa acid extract (Paclitaxel treated)	100 µg	36201
HeLa acid extract (Sodium Butyrate treated)	100 µg	36202
HeLa acid extract (Etoposide treated)	100 µg	36203
HeLa acid extract (Anacardic acid treated)	100 µg	36204

DNA Methylation	Format	Catalog No.
MethylDetector™	50 rxns	55001
MethylCollector™	25 rxns	55002
Fully Methylated Jurkat DNA	10 µg	55003

Histone Purification	Format	Catalog No.
Histone Purification Kit	10 rxns	40025
Histone Purification Mini Kit	20 rxns	40026

Chromatin Assembly	Format	Catalog No.
Chromatin Assembly Kit	10 rxns	53500

Histone Acetyltransferase and Deacetylase Activity	Format	Catalog No.
HAT Assay Kit (Fluorescent)	1 x 96 rxns	56100
Recombinant p300 protein, catalytic domain	5 µg	31205
HDAC Assay Kit (Fluorescent)	1 x 96 rxns	56200
HDAC Assay Kit (Colorimetric)	1 x 96 rxns	56210

Co-Immunoprecipitation	Format	Catalog No.
Nuclear Complex Co-IP Kit	50 rxns	54001
Universal Magnetic Co-IP Kit	25 rxns	54002

SUMOylation	Format	Catalog No.
SUMOLink™ SUMO-1 Kit	20 rxns	40120
SUMOLink™ SUMO-2/3 Kit	20 rxns	40220

ChIP-IT™ Kits	Format	Catalog No.
ChIP-IT™ Express	25 rxns	53008
ChIP-IT™ Express Enzymatic	25 rxns	53009
ChIP-IT™ Express HT	96 rxns	53018
ChIP-IT™ Protein G Magnetic Beads	25 rxns	53014
Re-ChIP-IT™	25 rxns	53016
ChIP-IT™	25 rxns	53001
ChIP-IT™ w/o controls	25 rxns	53004
ChIP-IT™ Shearing Kit	10 rxns	53002
ChIP-IT™ Enzymatic	25 rxns	53006
ChIP-IT™ Enzymatic w/o controls	25 rxns	53007
Enzymatic Shearing Kit	10 rxns	53005
Salmon Sperm DNA/Protein G agarose	25 rxns	53003
ChIP-IT™ Control Kit – Human	5 rxns	53010
ChIP-IT™ Control Kit – Mouse	5 rxns	53011
ChIP-IT™ Control Kit – Rat	5 rxns	53012
Ready-to-ChIP HeLa Chromatin	10 rxns	53015
Ready-to-ChIP Hep G2 Chromatin	10 rxns	53019
Ready-to-ChIP K-562 Chromatin	10 rxns	53020
Ready-to-ChIP NIH/3T3 Chromatin	10 rxns	53021

Transcription Factor ELISAs	Format	Catalog No.
TransAM™ AP-1 Family	2 x 96-well plates	44296
TransAM™ AP-1 c-Jun	1 x 96-well plate	46096
TransAM™ GR	1 x 96-well plate	45496
TransAM™ HIF-1	1 x 96-well plate	47096
TransAM™ IRF-3 (Human)	1 x 96-well plate	48396
TransAM™ IRF-7	1 x 96-well plate	50196
TransAM™ NFATc1	1 x 96-well plate	40296
TransAM™ NFκB Family	2 x 96-well plates	43296
TransAM™ NFκB p50	1 x 96-well plate	41096
TransAM™ NFκB p52	1 x 96-well plate	48196
TransAM™ NFκB p65	1 x 96-well plate	40096
TransAM™ p53	1 x 96-well plate	41196
TransAM™ STAT Family	2 x 96-well plates	42296

For a complete list of the over 40 TransAM Kits available, please visit www.activemotif.com/transam.

Technical Services

If you need assistance at any time, please call Active Motif Technical Service at one of the numbers listed below.

Active Motif North America

1914 Palomar Oaks Way, Suite 150

Carlsbad, CA 92008

USA

Toll Free: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

E-mail: tech_service@activemotif.com

Active Motif Europe

104 Avenue Franklin Roosevelt

B-1330 Rixensart, Belgium

UK Free Phone: 0800 169 31 47

France Free Phone: 0800 90 99 79

Germany Free Phone: 0800 181 99 10

Telephone: +32 (0)2 653 0001

Fax: +32 (0)2 653 0050

E-mail: eurotech@activemotif.com

Active Motif Japan

Azuma Bldg, 7th Floor

2-21 Ageba-Cho, Shinjuku-Ku

Tokyo, 162-0824, Japan

Telephone: +81 3 5225 3638

Fax: +81 3 5261 8733

E-mail: japantech@activemotif.com

Visit Active Motif on the worldwide web at <http://www.activemotif.com>

At this site:

- Read about who we are, where we are, and what we do
- Review data supporting our products and the latest updates
- Enter your name into our mailing list to receive our catalog, MotifVations newsletter and notification of our upcoming products
- Share your ideas and results with us
- View our job opportunities

Don't forget to bookmark our site for easy reference!