



SensoLyte[®] HAT (p300) Assay Kit **Fluorimetric**

Catalog #	72172
Kit Size	100 Assays (96-well plate)

- **Optimized Performance:** This kit is optimized to detect p300 activity
- **Enhanced Value:** It provides ample reagents to perform 100 assays in a 96-well plate format
- **High Speed:** The entire process can be completed in one hour
- **Assured Reliability:** Detailed protocol and references are provided

Kit Components, Storage and Handling

Component	Description	Quantity
Component A	Acetyl CoA	200 µL
Component B	Histone H3 peptide (1-21)	1 mM, 200 µL
Component C	p53 peptide (368-386)	4 mM, 200 µL
Component D	HAT, Recombinant p300 enzyme	100 µL, 0.1 mg/mL
Component E	Assay Buffer	40 mL
Component F	p300 Inhibitor	10 mM, 10 µL
Component G	p300 Developer	400 µL
Component H	Stop Solution	10 mL
Component I	CoASH standard	5 mM, 10 µL

Other Materials Required (but not provided)

- Microplate: Black, flat-bottom 96-well plate with non-binding surface.
- Fluorescence microplate reader: Capable of detecting emission at 513 nm with excitation at 389 nm.

Storage and Handling

- Store all kit components, except Components A and D, at -20°C
- Store Components A and D at -80°C. Aliquot as needed to avoid multiple freeze-thaw cycles.
- Protect Component G from light and moisture.
- Components E and H can be stored at room temperature for convenience.

Introduction

Histone acetyltransferases (HATs) enzymes regulate the acetylation of histones and non-histone proteins.^{1, 2} The acetylation of the ϵ -amino groups of lysine residues present at histone tails correlates largely with transcriptional activation, but it is also involved in DNA replication, DNA repair and protein–protein interactions.³ HATs play major roles in the control of cell fate and misregulation is implicated in the development of some human tumors.^{4,5} HAT p300 is a transcriptional coactivator that acetylates core histones facilitating chromatin decondensation and recruiting basic RNA polymerase machinery.⁶ Many non-histone proteins, such as p53, STATs, and alpha interferon receptor, serve as substrates for p300.⁷

The SensoLyte[®] HAT (p300) Assay Kit provides a convenient assay for screening of enzyme inhibitors and for continuous measurement of p300 activity. For added convenience, this kit includes two peptide substrates: histone H3 (1-21) peptide and non-histone p53 peptide (368-386). After incubation with acetyl CoA and the substrate, the p300 enzyme generates acetylated H3 or p53 peptide and CoASH. The thiol groups of CoASH can be detected with fluorogenic reagent (ABD-F) at excitation/emission=389nm/513nm.

Protocol

Note: For standard curve, please refer to [Appendix I](#) (optional).

1. Prepare working solutions.

Note: Thaw all kit components to room temperature before starting the experiments.

1.1 Acetyl CoA solution: Dilute Acetyl CoA (Component A) 10-fold in assay buffer (Component D) for one experiment. Refer to Table 1. For each experiment, prepare fresh solution of Acetyl CoA.

Table 1. Acetyl CoA solution for one 96-well plate (100 assays)

Components	Volume
Acetyl CoA (Component A)	100 μ L
Assay Buffer (Component E)	900 μ L
Total volume	1 mL

1.2 H3 peptide or p53 peptide solution: Dilute peptide (Component B or C) 10-fold in assay buffer (Component E). Refer to Table 2.

Table 2. Substrate peptide solutions for one 96-well plate (100 assays)

Components	Volume
H3 or p53 peptide (Component B or C)	200 μ L
Assay Buffer (Component E)	1800 μ L
Total volume	2 mL

1.3 p300 diluent: Dilute the p300 enzyme (Component D) 10-fold in assay buffer (Component E). Refer to Table 3. This amount of enzyme is enough for a full 96-well plate. If not using the entire plate, adjust the amount of enzyme to be diluted accordingly.

Table 3. p300 enzyme solution for one 96-well plate (100 assays)

Components	Volume
p300 enzyme (Component D)	100 μ L
Assay buffer (Component E)	900 μ L
Total volume	1 mL

1.4 p300 inhibitor (Anacardic acid): Dilute the 10 mM inhibitor solution (Component F) 100-fold in assay buffer (Component E) to 100 μ M.

1.5 p300 Developer solution: Dilute the p300 Developer (Component G) 50-fold in assay buffer (Component E). Refer to Table 4.

Table 4. Developer solution for one 96-well plate (100 assays)

Components	Volume
p300 Developer (Component G)	200 μ L
Assay buffer (Component E)	9800 μ L
Total volume	10 mL

2. Set up the enzymatic reaction.

2.1 Add test compounds and diluted p300 enzyme solution to the microplate wells. For one well of a 96-well plate, the suggested volume of p300 enzyme solution is 10 μ L and 10 μ L for test compound.

2.2 Simultaneously establish the following control wells, as deemed necessary:

- Positive control contains the p300 without test compound.
- Inhibitor control contains p300 and inhibitor Anacardic acid from Step 1.4.
- Vehicle control contains p300 and vehicle used in delivering test compound (e.g. DMSO, concentration not to exceed 1%).
- Test compound control contains assay buffer (Component E) and test compound. Some test compounds have strong autofluorescence and may give false results.
- Background control contains the p300 without test compound.

2.3 Using the assay buffer (Component E), bring the total volume of all controls to 20 μ L.

2.4 Pre-incubate the plate for 10 min at room temperature.

3. Run the enzymatic reaction.

3.1 Add 10 μ L of Acetyl CoA solution from Step 1.1 into each well.

3.2 Add 20 μ L of substrate peptide solution from Step 1.2 into each well, except background control wells. Mix the reagents completely by shaking the plate gently for 30 sec.

3.3 Incubate the mixture for 15 min at 37 °C and add 50 μ L of Stop Solution (Component H).

Note: For best accuracy, it is advisable to have the solutions equilibrated to the assay temperature.

3.4 Add 20 μ L of substrate peptide solution from Step 1.2 into background control wells.

3.5 Add 100 μ L of Developer solution from Step 1.5 into each well. Incubate the reaction for 30 min at room temperature.

3.6 Measure fluorescence signal: Keep plate from direct light. Measure fluorescence intensity at Ex/Em=389/513nm. Fluorescence signal is stable at room temperature for at least 2 hours.

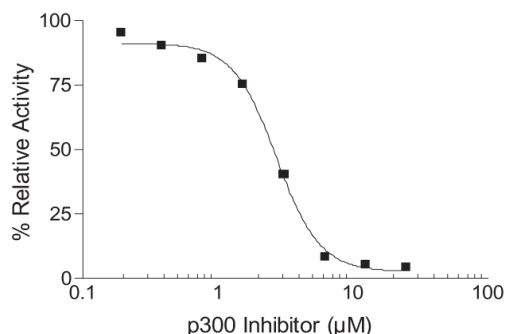


Figure 1. Inhibition of p300 activity (H3 peptide) was measured using Sensolyte[®] HAT(p300) Assay Kit.

Appendix I. Standard Curve

- **CoASH standard curve:** Dilute 5 mM CoASH (Component I) 100-fold to 50 μM with assay buffer (Component E). Perform 2-fold serial dilutions to get concentrations of 25, 12.5, 6.25, 3.13, 1.56 and 0.78 μM, and include an assay buffer blank. Add 50 μL/well of these serially diluted CoASH standard solutions.
- Add 50 μL/well of the Stop Solution (Component H).
- Add 100 μL/well of the Developer solution from Step 1.5. Incubate the reaction for 30 min at room temperature.
- Measure the fluorescence of the CoASH standard at Ex/Em=389/513 nm. Use the same setting of sensitivity as used in the enzyme reaction.
- Plot the CoASH standard curve as RFU (relative fluorescent units) versus concentration as shown in Figure 2.
- The final concentrations of CoASH standard are 12.5, 6.25, 3.13, 1.56, 0.78, 0.39, 0.20, and 0 μM. This reference standard is used to calibrate the variation of different instruments and different experiments. It is also an indicator of the amount of final product of the enzymatic reaction.

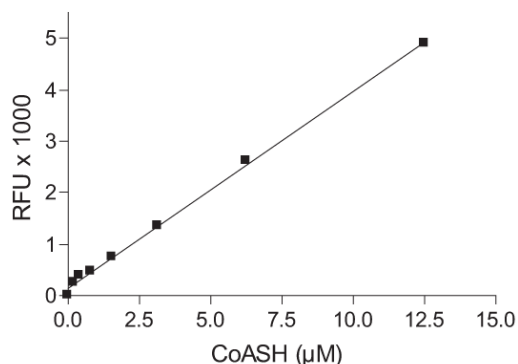


Figure 2. CoASH standard. CoASH was serially diluted in assay buffer and after 30 min incubation with Developer solution, fluorescence was recorded at Ex/Em=389/513 nm (Flexstation 384II, Molecular Devices).

References :

1. Roth, SY. et al. *Annu Rev Biochem.* **70**, 81 (2001).
2. Glozak, MA. et al. *Gene.* **363**, 15 (2005).
3. Turner, BM. *Bioessays* **22**, 836 (2000).
4. Timmermann, S. et al. *Cell Mol Life Sci.* **58**, 728 (2001).
5. Iyer, NG. et al. *Oncogene.* **23**, 4225 (2004).
6. Chan, HM. et al. *J Cell Sci.* **114**, 2363 (2001).
7. Grossman, SR. *Eur J Biochem.* **268**, 2773 (2001).

Revised: May 24, 2011