



Acetyltransferase activity kit

Catalog # **ADI-907-026**

Sufficient Reagents for 96 tests
For use with purified in vitro samples

Table of Contents



Reagents require separate storage conditions.



Check our website for additional protocols, technical notes and FAQs.



For proper performance, use the insert provided with each individual kit received.

- 2 Introduction**
- 2 Principle**
- 3 Materials Supplied**
- 4 Storage**
- 4 Materials Needed but Not Supplied**
- 5 Reagent Preparation**
- 7 Sample Handling**
- 7 End point / Kinetic Assay Procedure**
- 8 Example of Endpoint Assay Results**
- 9 Example Kinetic Assay Results**
- 10 Typical Inhibition Assay Procedure**
- 10 Calculation of Results**
- 11 Typical Inhibition Assay Results**
- 12 Interfering Substances**
- 13 References**
- 16 Limited Warranty**

Patent Pending

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

Introduction

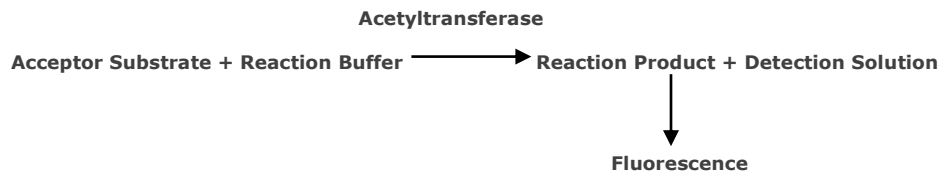
The Acetyltransferase activity kit is a complete kit for the screening of candidate compounds that may alter normal acetyltransferase activity. Please read the complete kit insert before performing this assay.

Acetylation is an important covalent molecular modification. Originally identified as the method by which certain bacteria were able to deactivate anti-microbial compounds, acetylation is now also known as an important partitioning and signaling modification.¹

Acetyltransferases are enzymes that covalently transfer an acetyl group from a donor molecule (Acetyl CoA) to an acceptor. Acetyl CoA serves as a universal donor while the acceptor varies with the acetyltransferase. Acceptors include histones, kinases, transcription factors, receptors, neurotransmitter precursors like choline and serotonin, and anti-microbial agents like chloramphenicol and fluoroquinones.²⁻⁶ Acetylation can signal an increase or decrease in activity based on the context of the message.⁷ Frequently located at critical junctions in metabolic pathways, Acetyltransferases and their regulation have become attractive therapeutic targets to treat everything from insomnia to cancer.⁸⁻¹⁰

Principle

1. Enzyme and reaction mix spiked with acceptor substrate are added to wells of a black 96-well plate. The plate is then incubated.
2. Ice cold isopropyl alcohol is added to stop the reaction. The detection solution is then added and the plate is again incubated.
3. The plate is transferred to a plate reader and fluorescence is measured at 380ex/520em.





Do not mix components from different kit lots or use reagents beyond the kit's expiration date.

Materials Supplied

1. **Black Microtiter Plate**
One plate of 96 wells, Catalog No. 80-1675
The plate is ready to use.
2. **Transferase Assay Buffer Concentrate**
15 mL, Catalog No. 80-1648
A 10X concentrated buffer containing detergent and preservative.
3. **Acetyltransferase Reaction Buffer Concentrate**
0.1 mL, Catalog No. 80-1651
A 50X concentrated buffer containing Acetyl CoA.
4. **Transferase Detection Solution Concentrate.**
0.15 mL, Catalog No. 80-1650
A 100X concentrated solution of fluorescent substrate in DMSO.
5. **Acetyltransferase Positive Control**
0.2 mL, Catalog No. 80-1647
A Stock Solution of buffer and reaction end product.
6. **Foil Plate Sealer**
3 each, Catalog No. 10-3126
7. **Acetyltransferase Assay Layout Sheet**
1 each, Catalog No. 30-0240



Detection Solution contains DMSO. Avoid skin contact. Avoid breathing vapor.



Reagents require separate storage conditions.

Storage

Acetyltransferase Reaction Buffer Concentrate and Transferase Detection Solution concentrate **must** be stored at -20°C , and Acetyltransferase Positive Control **must** be stored at -70°C . Transferase Assay buffer Concentrate can be stored at -20°C or 4°C . All kit components are stable at their recommended storage temperatures until the kit expiration date. Recommended storage temperatures do not necessarily reflect shipping conditions.

Materials Needed but Not Supplied

1. Deionized or distilled water.
2. Acetyl CoA dependent acetyltransferase.
3. Appropriate acceptor substrate.
4. Inhibitor/activator compounds to be screened.
5. Precision pipets for volumes between $5\ \mu\text{L}$ and $1,000\ \mu\text{L}$.
6. Disposable beakers for diluting buffer concentrates.
7. Graduated cylinders.
8. A microplate shaker.
9. Disposable microtubes, 0.5 and $1.5\ \text{mL}$.
10. Microplate reader capable of measuring fluorescence at $380\text{ex}/520\text{em}$.
11. Crushed ice and container.
12. Isopropyl alcohol (ice cold).

Reagent Preparation



Thaw Detection Solution Concentrate at room temperature. Do not place on ice. Thaw all other reagents on ice.



Pre-rinse each pipet tip with reagent.



Store 1X Detection Solution on ice and use within 4 hours of preparation. Discard unused 1X Solution.



A minimum of 2 wells of the positive control are recommended. Adjust the calculations accordingly if more than 2 wells are used.



Detection Solution contains DMSO. Avoid skin contact. Avoid breathing vapor.

1 Transferase Assay Buffer

Prepare the assay buffer by diluting 15mL of the supplied Transferase Assay Buffer Concentrate with 135mL of deionized water. This solution can be stored at 4°C for 3 months, or the kit's expiration, whichever is earlier. The 1X assay buffer is used to prepare dilutions of Acetyltransferase Reaction Buffer, Transferase Detection Solution, enzymes, substrates, and compounds to be screened.

2 Transferase Detection Solution

Count the total number of wells needed for compound screening and add 6 (for the zero, positive control, and blank wells in duplicate). Use the following formula to calculate the volume of 1X Detection Solution required.

A. Total volume required

$$[\text{Total number of wells needed} + 6] \times 100 \mu\text{L} = \text{_____} \mu\text{L}$$

B. Volume of Transferase Detection Solution Concentrate required

$$[\text{Total volume required (from A. above)}] \times 0.01 = \text{_____} \mu\text{L}$$

C. Volume of 1X Transferase Assay Buffer required

$$[\text{Total volume required (from A. above)}] \times 0.99 = \text{_____} \mu\text{L}$$

Prepare 1X Detection Solution by combining the appropriate reagent volumes calculated in B and C above. For example, to prepare 2 mL of 1X Detection Solution, combine the following volumes: 20 μL of the supplied Transferase Detection Solution Concentrate and 1980 μL 1X Transferase Assay Buffer. Diluted Detection Solution should be kept on ice and used within 4 hours of preparation. Any unused 1X Detection Solution should be discarded.

3 Acetyltransferase Reaction Mix

Count the total number of wells needed for compound screening and add 4 (for the zero and blank wells, in duplicate). Use the following formula to calculate the volume of 1X Acetyltransferase Reaction Mix required.

A. Total volume required

$$[\text{Total number of wells needed} + 4] \times 25 \mu\text{L} = \text{_____} \mu\text{L}$$

B. Volume of Acetyltransferase Reaction Buffer Concentrate required

$$[\text{Total volume required (from A. above)}] \times 0.02 = \text{_____} \mu\text{L}$$

C. Volume of acceptor substrate required

$$[\text{Determined empirically based on enzyme used}] = \text{_____} \mu\text{L}$$

D. Volume of 1X Transferase Assay Buffer required

$$[\text{Total volume required (from A. above)}] \times 0.98 - [\text{volume of substrate required}] = \text{_____ } \mu\text{L}$$

Prepare 1X Reaction Mix **on ice** by combining the appropriate reagent volumes calculated in B, C and D above. For example, to prepare 2 mL of Reaction Mix spiked with 22 μL of acceptor substrate, combine the following volumes: 40 μL Acetyltransferase Reaction Buffer Concentrate, 22 μL substrate and 1938 μL 1X Transferase Assay Buffer. Diluted Reaction Mix should be kept on ice and used within 8 hours of preparation. Any unused 1X Reaction Mix should be discarded.



Thaw Acetyltransferase Positive Control on ice.



Store Diluted Positive Control on ice and use within 2 hours of preparation. Discard unused 1X solution.

4 Positive Control

A positive control is included in the kit to verify the activity of the kit components. It should not be used to calculate the concentration of acetyltransferase activity in samples. Store the Positive Control at -70°C and avoid freeze-thaws.

Prepare enough Positive Control to test in duplicate. Use the following formula to calculate the volume of Positive Control required. A minimum of 2 wells of the positive control are recommended per assay.

A. Total volume required

$$[\text{Total number of wells needed}] \times 50\mu\text{L} = \text{_____ } \mu\text{L}$$

B. Dilute Acetyltransferase Positive Control Stock 1:100 by adding 5 μL of Positive Control to 495 μL of 1X Transferase Assay Buffer.

C. Volume of 1:100 Positive Control required

$$[\text{Total volume required (from A. above)}] \times 0.025 = \text{_____ } \mu\text{L}$$

D. Volume of 1X Transferase Assay Buffer required

$$[\text{Total volume required (from A. above)}] \times 0.975 = \text{_____ } \mu\text{L}$$

Prepare Positive Control **on ice** by combining the appropriate reagent volumes calculated in C and D above. For example, to prepare 200 μL of Positive Control, combine the following volumes: 5 μL of (1:100) Positive Control and 195 μL of 1X Transferase Assay Buffer. Diluted Positive Control should be kept on ice and used within 2 hours of preparation. Any unused 1X Positive Control should be discarded.

Sample Handling



If buffers other than those provided are used in the assay, the end-user must determine the appropriate dilution and assay validation.



Before screening colored compounds, the compound should be titrated in the assay to ensure that there is no interference with signal collection.



All samples should be run in duplicate.



Pre-rinse each pipet tip with reagent. Use fresh pipet tips for each sample and reagent.



Pipet the reagents to the sides of the wells to avoid possible contamination.

This assay is suitable for use with all Acetyl CoA-dependent acetyltransferases. It is necessary to titrate each enzyme / substrate system in the assay to determine optimal conditions.

This assay should only be used to screen purified *in vitro* samples in buffer systems without reductants.

It is recommended that an end-point assay be performed to determine the optimal concentration of enzyme/substrate to use prior to screening candidate compounds. Make serial dilutions of the acetyltransferase of interest in the assay buffer. Initial concentrations of 100 nM are recommended. A kinetic assay format is also an available option.

The positive control provided may also be used to test colored compounds for interference in the assay. A suitable protocol follows.

End point / Kinetic Assay Procedure

Refer to the Assay Layout Sheet to determine the number of wells to be used. Cover unused wells tightly with a plate sealer. **DO NOT REUSE WELLS!**

1. Pipet 25 μ L of 1X Transferase Assay Buffer into the blank wells.
2. Pipet 25 μ L of acetyltransferase dilutions into the bottom of the appropriate wells.
3. Pipet 25 μ L of 1X Reaction Mix into each well.
4. Cover plate with foil plate sealer. Incubate for 30 min (for end point format) shaking* at room temperature.

For **Kinetic** measurements, incubate identical reaction wells for the desired periods of time. The duplicate wells for each time point can be stopped as in step #6 below.

5. Pipet 50 μ L of 1X Positive Control into the bottom of the appropriate wells.
6. Pipet 50 μ L of ice cold isopropyl alcohol into each well.
7. Pipet 100 μ L 1X Detection Solution into each well.
8. Cover plate with foil plate sealer. Incubate for 10 min at room temperature, without shaking.
9. Read fluorescence at 380ex/520em.

* Shaking is preferably carried out on a suitable plate or orbital shaker set at a speed to ensure adequate mixing of the contents of the wells. The optimal speed for each shaker will vary and may range from 120-700 rpm.



Make sure to multiply sample concentrations by the dilution factor used during sample preparation.

Example of Endpoint Assay Results

1. Plot the mean of the duplicate relative fluorescence units (RFU) at 380ex/520em versus Enzyme concentration. (Figure 1)
2. Calculate the signal-to-noise ratio : $\frac{\text{Mean RFU for enzyme dilution}}{\text{Mean RFU blank}}$.

Chloramphenicol Acetyltransferase Titration in Acetyltransferase Assay

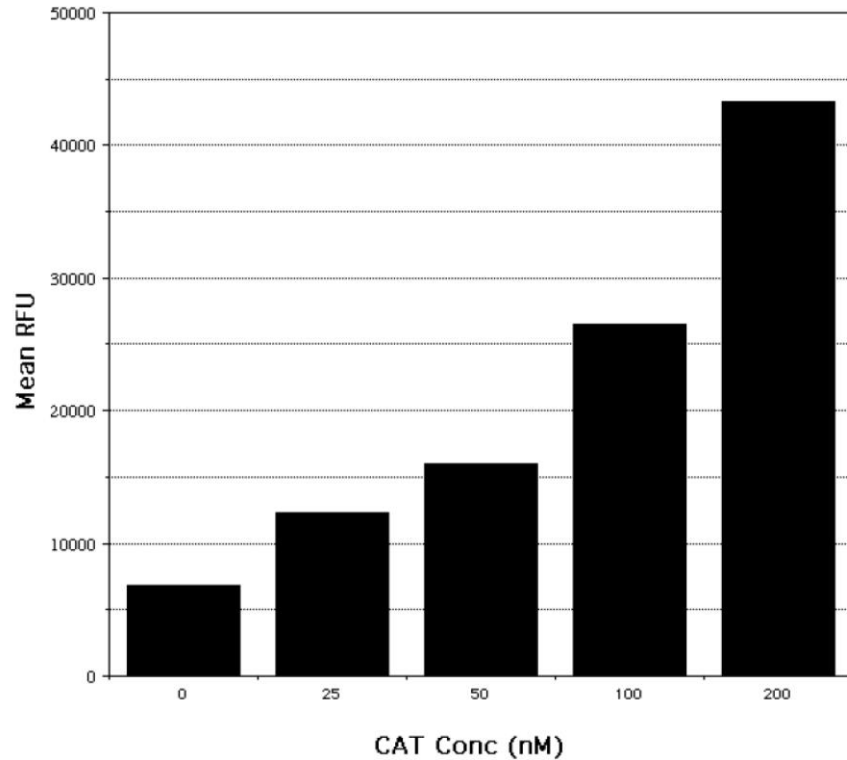


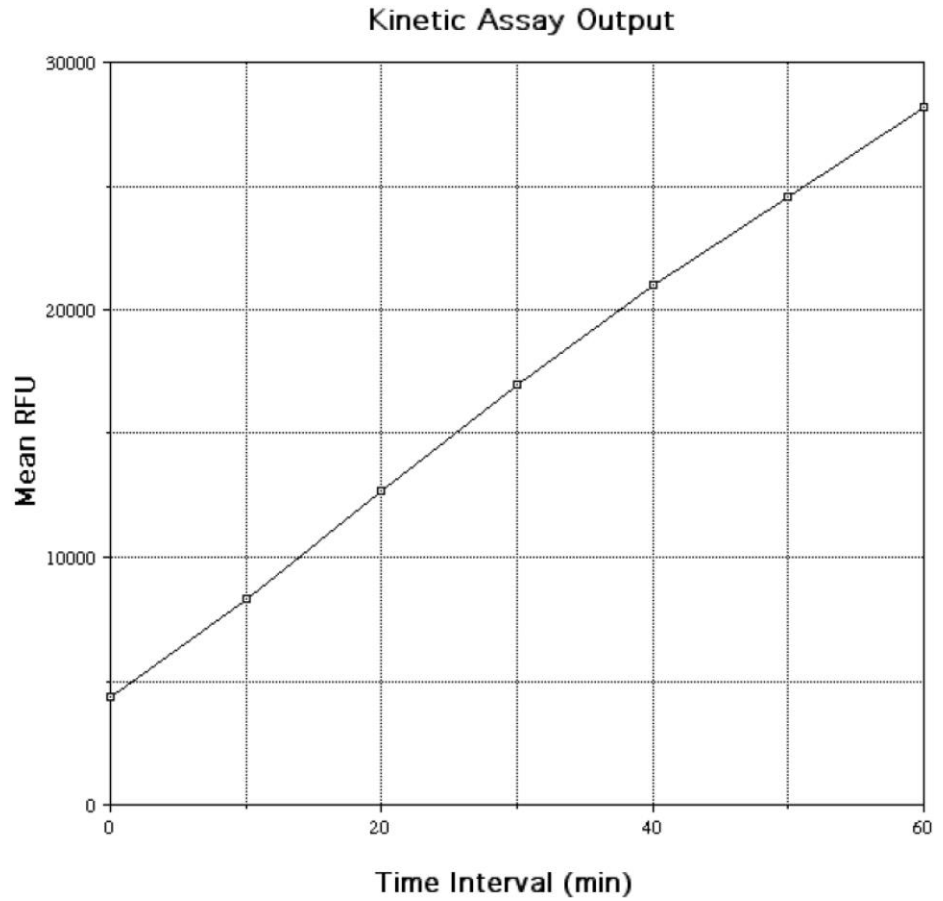
Figure 1:

Chloramphenicol acetyltransferase (CAT) was titrated in the assay using 100 μ M of the Substrate chloramphenicol. Serial dilutions of the enzyme were prepared in Transferase Assay Buffer. Mean relative fluorescence was plotted against CAT concentration to generate the following graph. This is for illustration purposes only. The investigator must titrate their enzyme / substrate system in the assay.

Based on this titration data, the acetyltransferase concentration of 100nM produces a maximum signal within the detection range of the plate reader, with a signal to noise ratio sufficient for easy detection of altered enzyme activity.

Example Kinetic Assay Results

100nM CAT was tested with 100 μ M Chloramphenicol substrate in the kinetic assay format. Mean relative fluorescence was plotted against the stop time interval to generate this graph. This graph is for illustration purposes only. The kinetic assay must be optimized by the investigator, with their enzyme / substrate system.



Typical Inhibition Assay Procedure

Refer to the Assay Layout Sheet to determine the number of wells to be used. Cover unused wells tightly with a plate sealer. **DO NOT REUSE WELLS!**

1. Pipet 25µL of 1X Transferase Assay Buffer into the blank wells.
2. Pipet 10µL of 1X Transferase Assay Buffer into the zero wells.
3. Pipet 10µL of inhibitor dilution into the bottom of the appropriate wells.
4. Pipet 15µL of acetyltransferase at chosen working concentration into appropriate wells.
5. Cover plate with foil plate sealer. Incubate for 10 min at room temperature without shaking.
6. Pipet 25µL of 1X Reaction Mix into each well.
7. Cover plate with foil plate sealer. Incubate for 30 min shaking* at room temperature.
8. Pipet 50µL of ice cold isopropyl alcohol into each well.
9. Pipet 100µL of 1X Detection Solution into each well.
10. Cover plate with foil plate sealer. Incubate for 10 min at room temperature without shaking.
11. Read fluorescence at 380ex/520em.

* Shaking is preferably carried out on a suitable plate or orbital shaker set at a speed to ensure adequate mixing of the contents of the wells. The optimal speed for each shaker will vary and may range from 120-700 rpm.

Calculation of Results

Several options are available for the calculation of the inhibition of acetyltransferase. We recommend that the data be handled by a software package utilizing a suitable curve fitting program to determine the percent inhibition. If data reduction software is not readily available, the data can be calculated as follows:

1. Calculate the mean net RFU for each sample by subtracting the mean blank RFU from the mean RFU for the samples:

$$\text{Mean Net RFU} = \text{Mean Sample RFU} - \text{Mean Blank RFU}$$

2. Percent inhibition should be calculated using the following formula for each inhibitor dilution:

Percent Inhibition =

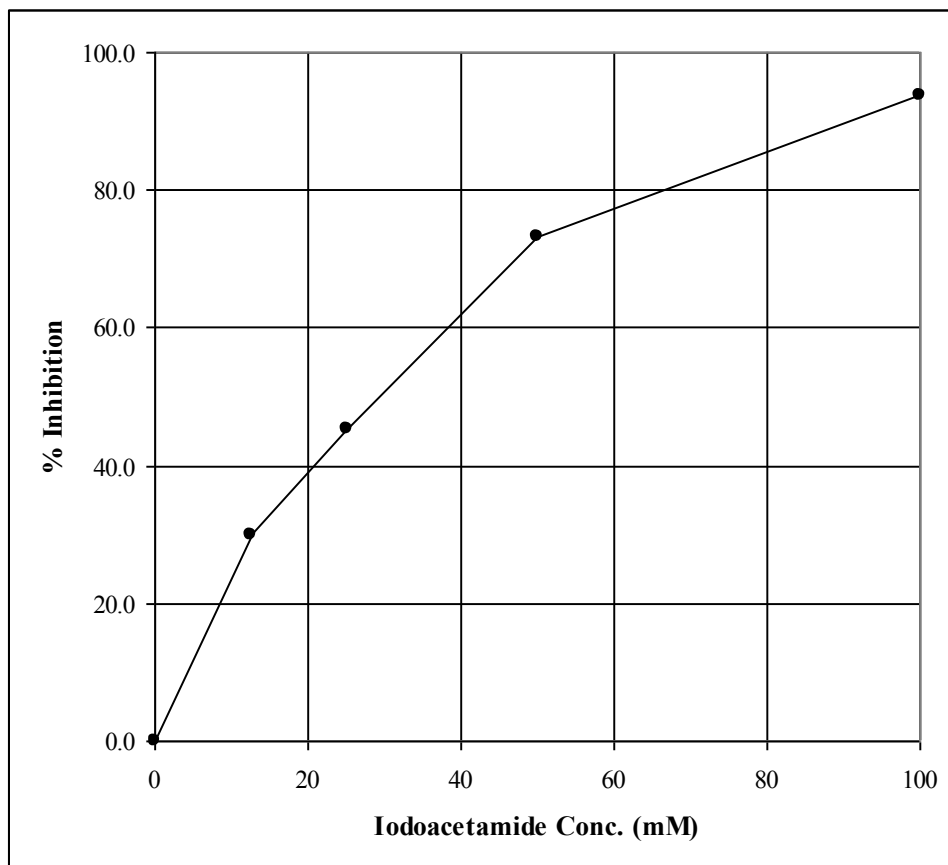
$$\frac{(\text{Mean Net Zero (non-inhibited enzyme) RFU} - \text{Mean Net Inhibited enzyme RFU}) \times 100}{\text{Mean Net Zero (non-inhibited enzyme) RFU}}$$

Typical Inhibition Assay Results

Using Chloramphenicol Acetyltransferase (CAT) and Chloramphenicol as an enzyme substrate system, percent inhibition for dilutions of Iodoacetamide was tested. The results shown below are for illustration only and **should not** be used to calculate results from another assay.

Dilution	Inhib. Conc. (mM)	Mean RFU	Mean Net RFU	% Inhibition
1	100mM	8,707	2,161	93.6%
2	50mM	15,575	9,029	73.1%
3	25mM	24,951	18,405	45.2%
4	12.5mM	30,110	23,564	29.8%
0	0mM	40,127	33,581	0%
blank	-----	6,889	0	-----

Typical Inhibition Curve



Typical Enzyme Interference Assay Procedure

*Prepare dilutions of the colored compound to be screened in 1X Transferase Assay Buffer.

1. Pipet 50 μ L of 1X Transferase Assay Buffer into the blank wells.
2. Pipet 25 μ L of 1X Transferase Assay Buffer into the zero wells.
3. Pipet 25 μ L of colored compound dilution into appropriate wells
4. Pipet 25 μ L of 1X Positive Control into the zero wells, and wells containing colored compound dilutions.
5. Pipet 50 μ L of ice cold isopropyl alcohol into each well.
6. Pipet 100 μ L 1X Detection Solution into each well.
7. Cover plate with foil plate sealer, incubate for 10 min at room temperature without shaking.
8. Read fluorescence at 380ex/520em.

To determine whether or not colored compounds will interfere with the assay, calculate and compare the signal to noise ratio of the colored compound dilutions to the signal to noise ratio of the zero wells.

Interfering Substances

The following solvents were tested for interference with the fluorescent signal generated in the assay. The table lists the percentage of signal in the presence of interferant relative to the zero for each solvent.

% Interferant	DMSO	DMF	Acetonitrile
12.5	108	68	49
6.25	107	82	71
3.12	107	88	86
1.56	112	98	96
0.78	114	104	95
0.39	115	109	100
0	100	100	100

* Percent interferant is relative to a 50 μ L total reaction volume.

Diluents containing bovine serum albumin (BSA), or other Thiol containing reagents, should be treated with N-Ethylmaleimide (NEM) prior to use in the assay. We recommend reacting 10% BSA with 1 mM NEM for 1 hour at room temperature, however this procedure should be optimized for each reagent.

References

1. Walsh, C.T.. Posttranslational Modification of Proteins, Expanding Nature's Inventory. Roberts and Company; Englewood.
2. Schilz, R.L., et. al. Overlapping but Distinct Patterns of Histone Acetylation by the Human Coactivators p300 and PCAF within Nucleosomal Substrates. *J. Biol. Chem.* (1999); 274(3): 1189-1999.
3. Worbly, C.A. & D.E. Dixon. Bacteria Seize Control by Acetylating Host Proteins. *Science* (2006) 312:1150-1151.
4. Kim, M.Y., et. al. A Role for Coactivators and Histone Acetylation in Estrogen Receptor α -mediated Transcription Initiation. *EMBO J* (2001) 20(21):6084-6094.
5. DeAngelis, J., et. al. Kinetic Analysis of the Catalytic Mechanism of Serotonin N-Acetyltransferase (EC 2.3.1.87). *J. Biol. Chem.* (1998) 273(5):3045-3050.
6. Fitton, J.E. & W.V. Shaw. Comparison of Chloramphenicol Acetyltransferase Variants in staphylococci; Purification, Inhibitor Studies and N-terminal Sequences. *Biochem. J.* (1997) 177:575-582.
7. Tai, H.H. et. al. Dynamic Histone Acetylation of Late Embryonic Genes During Seed Germination. *Plant Mol. Biol.* (2005) 59(6):909-925.
8. Khalil, E.M., et. al. Mechanism-based Inhibition of the Melatonin Rhythm Enzyme: Pharmacologic Exploitation of Active Site Functional Plasticity. *PNAS* (1999); (22):12418-12423.
9. Brockton, N. et. al. N-Acetyltransferase Polymorphisms and Colorectal Cancer: A HuGe Review. *J. Epidem.* (2000) 151(9):846-861.
10. Shi, H., et. al. Tripple Analysis of the Cancer Epigenome: An Integrated Microarray System for Assessing Gene Expression, DNA Methylation, and Histone Acetylation. *Cancer Res.* (2003) 63:2164-2171.

Notes

Notes



MSDS (Material Safety Data Sheet) available online



USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

www.enzolifesciences.com
Enabling Discovery in Life Science[®]

Global Headquarters

Enzo Life Sciences Inc.
10 Executive Blvd
Farmingdale, NY 11735
(p) 1-800-942-0430
(f) 1-631-694-7501
(e) info-usa@enzolifesciences.com

Enzo Life Sciences (ELS) AG

Industriestrasse 17, Postfach
CH-4415 Lause / Switzerland
(p) +41/0 61 926 89 89
(f) +41/0 61 926 89 79
(e) info-ch@enzolifesciences.com

Please visit our website at www.enzolifesciences.com for additional contact information.