

Product datasheet

Histone Acetyltransferase Activity Assay Kit (Colorimetric) ab65352

★★★★★ [1 Abreviews](#) [44 References](#) [3 Images](#)

Overview

Product name	Histone Acetyltransferase Activity Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Nuclear Extracts, Purified protein, Tissue Lysate
Assay type	Enzyme activity
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Abcam's Histone Acetyltransferase Activity Assay Kit offers a convenient, nonradioactive system for a rapid and sensitive detection of HAT activity in mammalian samples. The kit utilizes active Nuclear Extract (NE) as a positive control and acetyl-CoA as a cofactor. Acetylation of peptide substrate by active HAT releases the free form of CoA which then serves as an essential coenzyme for producing NADH. NADH can easily be detected spectrophotometrically upon reacting with a soluble tetrazolium dye. The detection can be continuous and suitable for kinetic studies. The kit provides a simple, straightforward protocol for a complete assay. Visit our FAQs page for tips and troubleshooting.</p> <p>HAT assay protocol summary:</p> <ul style="list-style-type: none">- add samples to wells- add HAT assay mix- incubate for 1-4 hrs- analyze with microplate reader
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K332 HAT Activity Colorimetric Assay Kit. K332-100 is the same size as the 100 test size of ab65352.</p> <p>Histone acetyltransferases (HATs) have been implicated to play a crucial role in various cellular functions, such as gene transcription, differentiation, and proliferation.</p> <p>Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.</p>
Platform	Microplate reader

Properties

Storage instructions

Store at -80°C. Please refer to protocols.

Components	Identifier	100 tests
2X HAT Assay Buffer	Amber	1 x 7.5ml
HAT Reconstitution Buffer	Clear	1 x 1.8ml
HAT Substrate I	Blue	1 vial
HAT Substrate II	Red	1 vial
NADH Generating Enzyme	Green	1 x 800µl
Nuclear Extract (4 mg/ml)	Violet	1 x 50µl

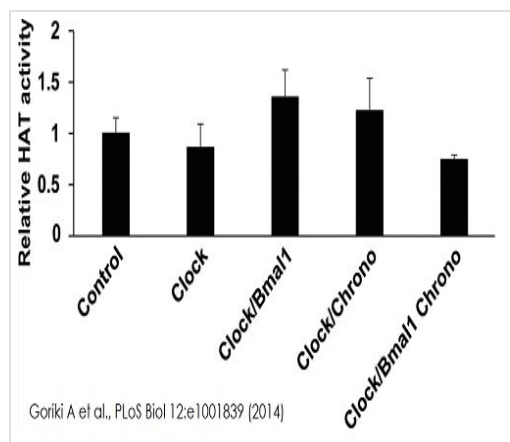
Relevance

Histone acetyltransferases (HAT) are enzymes that acetylate conserved lysine amino acids on histone proteins. They have been implicated to play a crucial role in various cellular functions, such as gene transcription, differentiation, and proliferation.

Cellular localization

Cytoplasmic and Nuclear. Note=Nuclear in S-phase cells and cytoplasmic.

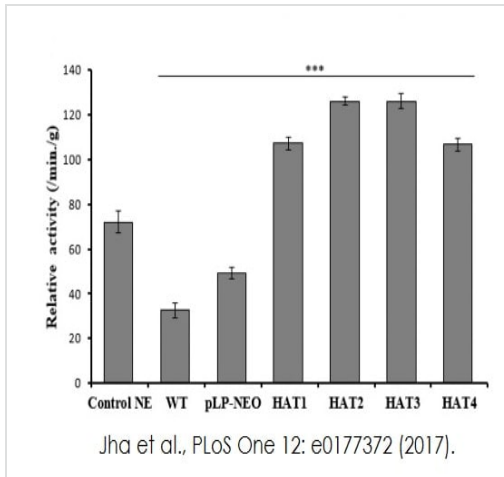
Images



HAT activity in NIH3T3 cells, which were transfected with the desired plasmids by using Lipofectamine 2000. After 24 hours from transfection, the nuclear protein was extracted and the Histone acetyltransferase activity was measured by using ab65352.

Functional studies - ab65352

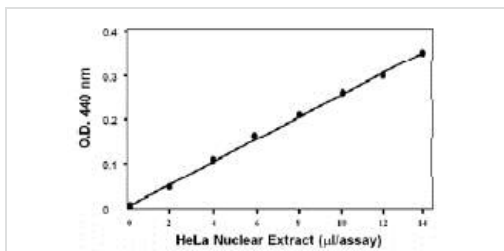
Image from Goriki A et al., PLoS Biol 12(4), Figure S11. doi:10.1371/journal.pbio.1001839. Reproduced under the Creative Commons license <http://creativecommons.org/licenses/by/4.0/>



Histone acetylation activity assay for *L. donovani* cells over-expressing HAT1, HAT2, HAT3 and HAT4. The activities were compared to wild-type (WT) cells and pLPneo2 (vector) only transfected cells. The relative activity was determined using Histone Acetyltransferase activity assay kit (ab65352).

Functional studies - ab65352

Image from Jha et al., PLoS One 12(5), fig 2C. doi: 10.1371/journal.pone.0177372. Reproduced under the Creative Commons license <http://creativecommons.org/licenses/by/4.0/>



Analysis of Histone Acetyltransferase Activity in HeLa Nuclear Extract using ab65352. HeLa nuclear extract in various amounts was incubated with HAT substrate. Activity was analyzed in a micro plate reader at 440 nm according to the kit instructions.

Functional Studies - Histone Acetyltransferase Activity Assay Kit (Colorimetric) (ab65352)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you

- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors