


Product datasheet

Histone H3 Total Acetylation Detection Fast Kit (Colorimetric) ab115124

★★★★★ 1 Abreviews 7 References 3 Images

Overview

| | |
|--------------------|--|
| Product name | Histone H3 Total Acetylation Detection Fast Kit (Colorimetric) |
| Detection method | Colorimetric |
| Sample type | Tissue, Adherent cells, Suspension cells |
| Sensitivity | 2 ng |
| Range | 2 ng - 100 ng |
| Assay time | 2h 30m |
| Species reactivity | Reacts with: Mouse, Human Predicted to work with: Mammals  |

Product overview

Acetylation of histones such histone H3 has been involved in the regulation of chromatin structure and the recruitment of transcription factors to gene promoters. HATs (histone acetyltransferases) and HDACs (histone deacetylases) play a critical role in controlling histone H3 acetylation. Histone acetylation is tightly involved in cell cycle regulation, cell proliferation and apoptosis. The reversible lysine acetylation of histone H3 may play a vital role in the regulation of many cellular processes including chromatin dynamics and transcription, gene silencing, cell cycle progression, apoptosis, differentiation, DNA replication and repair, nuclear import and neuronal repression. Histone H3 Total Acetylation Detection Fast Kit (Colorimetric) allows the user to colorimetrically detect and quantify if histone H3 is acetylated. The kit is ready-to-use and provides all the essential components needed to carry out a successful assay experiment. ab115124 is suitable for specifically measuring total histone H3 acetylation using a variety of mammalian cells including fresh and frozen tissues, and cultured adherent and suspension cells.

| | |
|----------|-------------------|
| Platform | Microplate reader |
|----------|-------------------|

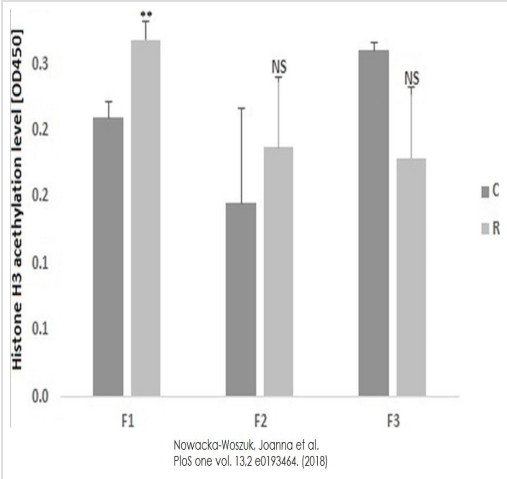
Properties

| | |
|----------------------|----------------------------|
| Storage instructions | Please refer to protocols. |
|----------------------|----------------------------|

| Components | Identifier | 48 tests | 96 tests |
|-----------------|------------|----------|----------|
| 10X Wash Buffer | | 1 x 10ml | 1 x 20ml |
| | | | |

| Components | Identifier | 48 tests | 96 tests |
|-----------------------------------|--------------|----------|----------|
| 8-Well Sample Strips (with Frame) | | 4 units | 9 units |
| 8-Well Standard Control Strips | Green Ringed | 2 units | 3 units |
| Antibody Buffer | | 1 x 6ml | 1 x 12ml |
| Color Developer | | 1 x 5ml | 1 x 10ml |
| Detection Antibody, 1 mg/ml | | 1 x 5µl | 1 x 10µl |
| Standard Control (100 µg/ml) | | 1 x 10µl | 1 x 20µl |
| Stop Solution | | 1 x 3ml | 1 x 6ml |

Images



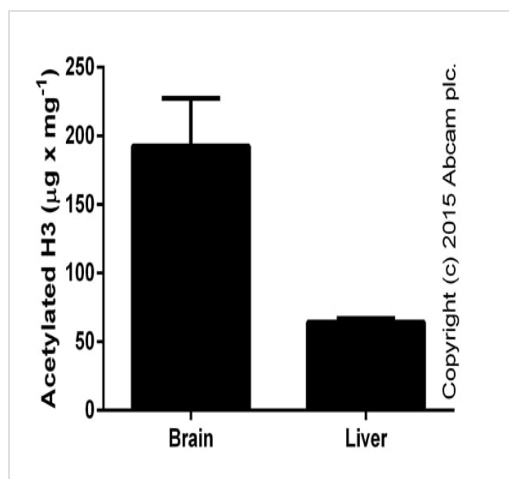
The differences in H3 acetylation level in fetal livers (female fetuses aged 19 dpc) from F1, F2, and F3 generations. measured with **ab113476**. Histones were prepared using ab115124

Asterisk indicate a significant difference between the respective C and R groups: □ P < 0.01; NS not significant. Data presented as the means ± SE (n = 6 per group).

Functional Studies - Histone H3 Total Acetylation

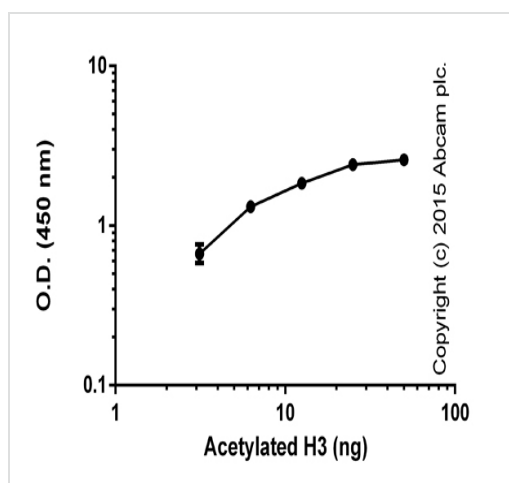
Detection Fast Kit (Colorimetric) (ab115124)

Nowacka-Wozuk, Joanna et al., PloS one?vol. 13,2 e0193464., Fig 4, doi:10.1371/journal.pone.0193464



Acetylation Histone 3 per mg of extracted histones from murine tissue preparations (extracted using **ab113476**; 20-500 ng per well) (duplicates +/- SD).

Functional Assay: ab115124 Histone H3 Total
Acetylation Detection Fast Kit (Colorimetric)



Standard curve with background signal subtracted (duplicates; +/- SD).

Functional Assay: ab115124 Histone H3 Total
Acetylation Detection Fast Kit (Colorimetric)

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