

Histone H4 Total Acetylation Assay Kit (Colorimetric)

ab115125

4 References

Overview

| | |
|---------------------------|---|
| Product name | Histone H4 Total Acetylation Assay Kit (Colorimetric) |
| Detection method | Colorimetric |
| Sample type | Tissue, Adherent cells, Suspension cells |
| Sensitivity | 2 ng/well |
| Range | 5 ng/well - 2000 ng/well |
| Assay time | 2h 30m |
| Species reactivity | Reacts with: Mouse, Human |

Predicted to work with: Mammals 

Product overview

Acetylation of histones such histone H4 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 H4 acetylation. Reversible acetylation of nucleosomal histone H4 is believed to correlate with potential transcriptional activity of eukaryotic chromatin domains. The reversible lysine acetylation of histone H4 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 H4 Total Acetylation Detection Fast Kit (Colorimetric) enables the user to colorimetrically detect and quantify if histone H4 is acetylated. The kit is ready-to-use and provides all the essential components needed to carry out a successful assay experiment. Histone H4 Total Acetylation Detection Fast Kit (Colorimetric) is suitable for specifically measuring total histone H4 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 | 48 tests | 96 tests |
|-----------------------------------|-----------|-----------|
| 10X Wash Buffer | 1 x 10ml | 1 x 20ml |
| 8-Well Sample Strips (with Frame) | 4 units | 9 units |
| 8-Well Standard Control Strips | 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 |
| Signal Enhancer | 1 x 120µl | 1 x 240µl |
| Signal Report Solution | 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 |

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