

Product datasheet

Histone H3 Acetylation Assay Kit ab115102

★★★★☆ 2 Abreviews 3 References

Overview

Product name	Histone H3 Acetylation Assay Kit
Sample type	Tissue, Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Sensitivity	1 ng
Range	1 ng - 100 ng
Assay time	5h 00m
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 acylation. Histone acetylation is tightly involved in cell cycle regulation, cell proliferation and apoptosis. An imbalance in the equilibrium of histone acetylation has been associated with tumorigenesis and cancer progression.

Histone H3 Acetylation Assay Kit (ab115102) allows the user to measure global acetylation of histone H3 at tremendously fast speeds and consistency, superior and safer than all other current methods. The kit is ready-to-use and provides all the essential components needed to carry out a successful assay experiment. The kit is suitable for specifically measuring global 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	48 tests	96 tests
10X Lysis Buffer	1 x 5ml	1 x 10ml
10X Wash Buffer	1 x 14ml	1 x 28ml

Components	48 tests	96 tests
8-Well Assay Strips (with Frame)	6 units	12 units
Acetylated Histone H3 Control (60 µg/mL)	1 x 10µl	1 x 20µl
Antibody Buffer	1 x 6ml	1 x 12ml
Blocking Buffer	1 x 10ml	1 x 20ml
Capture Antibody, 100 µg/mL	1 x 25µl	1 x 50µl
Detection Antibody, 400 µg/mL	1 x 10µl	1 x 20µl
Developing Solution	1 x 5ml	1 x 10ml
Extraction Buffer	1 x 8ml	1 x 16ml
Histone Buffer	1 x 0.5ml	1 x 1ml
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