

Product datasheet

Histone H3 (acetyl K14) Quantification Kit (Colorimetric) ab115106

Overview

Product name	Histone H3 (acetyl K14) Quantification Kit (Colorimetric)
Sample type	Tissue, Adherent cells, Suspension cells
Sensitivity	> 2 ng/well
Range	20 ng/well - 5000 ng/well
Assay time	2h 30m
Species reactivity	Reacts with: Mouse, Human Predicted to work with: Mammal

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. Lysine 14 at histone H3, along with K9 and K18, are primary acetylated sites of histone H3. Histone H3 (acetyl K14) is tightly involved in cell cycle regulation, cell proliferation and apoptosis. It also correlates with transcription activation.

Abcam's Histone H3 (acetyl K14) Quantification Kit (Colorimetric) (ab115106) allows the user to measure global histone H3K14 acetylation. 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 H3K14 acetylation using a variety of mammalian cells including fresh and frozen tissues, and cultured adherent and suspension cells.

Notes

Tested applications **Suitable for:** Functional Studies

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

Applications

Our [Abpromise guarantee](#) covers the use of **ab115106** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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