

## Product datasheet

# Histone H3 Total Acetylation Assay Kit (Fluorometric) ab131561

★★★★☆ [3 Abreviews](#) [6 References](#)

### Overview

---

<b>Product name</b>	Histone H3 Total Acetylation Assay Kit (Fluorometric)
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Tissue, Adherent cells, Suspension cells
<b>Assay type</b>	Quantitative
<b>Sensitivity</b>	> 0.4 ng
<b>Range</b>	0.4 ng - 50 ng
<b>Assay time</b>	2h 30m
<b>Species reactivity</b>	<b>Reacts with:</b> 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. 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.

ab131561 is designed for measuring total histone H3 acetylation in a fast format. In an assay with this kit, the acetyl histone H3 is captured to the strip wells coated with an anti-acetyl histone H3 antibody. The captured acetyl histone H3 can then be detected with a labeled detection antibody followed by a fluorescent development reagent. The ratio of acetyl histone H3 is proportional to the intensity of fluorescence. The absolute amount of acetyl histone H3 can be quantified by comparing to the standard control.

**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
8-Well Assay 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
Detection Antibody, 1 mg/mL		1 x 5µl	1 x 10µl
Fluoro Developer		1 x 12µl	1 x 24µl
Fluoro Dilution		1 x 4ml	1 x 8ml
Fluoro Enhancer		1 x 12µl	1 x 24µl
Standard Control (100 µg/mL)		1 x 10µl	1 x 20µl

**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