

## Product datasheet

# Histone H4 (acetyl K5) Quantification Kit (Colorimetric) ab115116

### 1 References

#### 製品の概要

製品名	Histone H4 (acetyl K5) Quantification Kit (Colorimetric)
サンプルの種類	Tissue, Adherent cells, Suspension cells
検出感度	2 ng/well
検出範囲	20 ng/well - 5000 ng/well
全工程の試験時間	2h 30m
種交差性	交差種: Mouse, Human 交差が予測される動物種: Mammal

#### 製品の概要

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. Histone H4 (acetyl K5) reflects the hyperacetylated state in histone H4 and strongly correlates with active state of genes. H4 (acetyl K5) is also considered to be a marker of dexamethasone transactivation and a change in H4 (acetyl K5) is observed in cancer and inflammatory diseases. Abcam's Histone H4 (acetyl K5) Quantification Kit (Colorimetric) (ab115116) allows the experimenter to measure global acetylation of histone H4K5 quickly and consistently. The kit is ready-to-use and provides all the essential components needed to carry out a successful assay experiment and it is suitable for specifically measuring global histone H4K5 acetylation using a variety of mammalian cells including fresh and frozen tissues, and cultured adherent and suspension cells.

#### 特記事項

アプリケーション **適用あり:** Functional Studies

#### 製品の特性

保存方法 Please refer to protocols.

内容	ラベル	48 tests	96 tests
10X Wash Buffer		1 x 10ml	1 x 20ml
8-Well Sample Strips (with Frame)		4 units	9 units

内容	ラベル	48 tests	96 tests
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 5 $\mu$ l	1 x 10 $\mu$ l
Signal Enhancer		1 x 120 $\mu$ l	1 x 240 $\mu$ l
Signal Report Solution		1 x 5 $\mu$ l	1 x 10 $\mu$ l
Standard Control, 100 $\mu$ g/mL		1 x 10 $\mu$ l	1 x 20 $\mu$ l
Stop Solution		1 x 3ml	1 x 6ml

## アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab115116** in the following tested applications.

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

アプリケーション	Abreviews	特記事項
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.co.jp/abpromise> or contact our technical team.

## Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors