

Product datasheet

Histone H3 (acetyl K18) Quantification Kit (Fluorometric) ab115109

概述

| | |
|-------|---|
| 产品名称 | Histone H3 (acetyl K18) Quantification试剂盒(Fluorometric) |
| 样品类型 | Tissue, Adherent cells, Suspension cells |
| 灵敏度 | 0.4 ng/well |
| 范围 | 5 ng/well - 2000 ng/well |
| 检测时间 | 2h 30m |
| 种属反应性 | 与反应: Mouse, Human 预测可用于: Mammal |

产品概述

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. Lysine 18 at histone H3, along with K9 and K14, are primary acetylated sites of histone H3. Histone H3 (acetyl K18) is tightly involved in cell cycle regulation, cell proliferation and apoptosis. It also correlates with transcription activation.

Abcam's Histone H3 (acetyl K18) Quantification Kit (Fluorometric) (ab115109) is a convenient package of tools that allows the experimenter to measure global acetylation of histone H3K18 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 H3K18 acetylation using a variety of mammalian cells including fresh and frozen tissues, and cultured adherent and suspension cells.

说明

经测试应用 **适用于: Functional Studies**

性能

存放说明 Please refer to protocols.

| 组件 | 标识符 | 48 tests | 96 tests |
|-----------------|-----|----------|----------|
| 10X Wash Buffer | | 1 x 10ml | 1 x 20ml |

| 组件 | 标识符 | 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 |
| 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 |

应用

Our [Abpromise guarantee](#) covers the use of **ab115109** in the following tested applications.

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

| 应用 | Ab评论 | 说明 |
|--------------------|------|--|
| 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.cn/abpromise> or contact our technical team.

Terms and conditions

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