

Histone Extraction Kit ab113476

★★★★★ 6 Abreviews 86 References 4 图像

概述

产品名称	Histone Extraction试剂盒
样品类型	Tissue, Adherent cells, Suspension cells
检测时间	1h 00m
产品概述	<p>Histone Extraction Kit (ab113476) is suitable for a quick isolation of total histone extracts from cells and tissue samples from mammals and other species.</p> <p>The amount of starting material can be as low as 10^5 cells or 1 mg of tissue. For best results, the cell number should be $>10^6$ cells or the tissue amount should be >10 mg.</p> <p>A total of 100 standard extractions (using 10^7 cells or 100 mg of tissue per extraction) can be performed with this kit. Yield of the total histone proteins can be up to 0.4 mg per 10^7 cells or per 100 mg of tissue. The yield may vary depending on the cell or tissue type.</p>

说明

Histone purification is most often performed before the analysis of the global levels within a sample of post-translational histone modifications in western blot.

Chromatin immunoprecipitation (ChIP) is the most common technique used for the analysis of histone modifications at specific sites within the genome. Use a chromatin extraction kit instead of a histone purification kit for the preparation of samples for ChIP.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

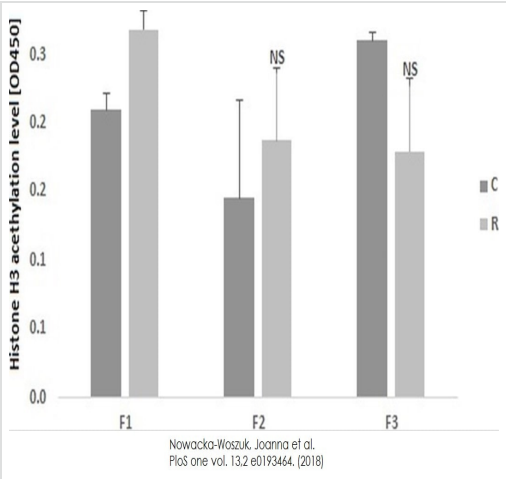
性能

存放说明 Store at +4°C. Please refer to protocols.

组件	100 tests
10X Pre-Lysis Buffer	1 x 10ml

组件	100 tests
Balance Buffer	1 x 8ml
DTT Solution	1 x 20μl
Lysis Buffer	1 x 20ml

图片



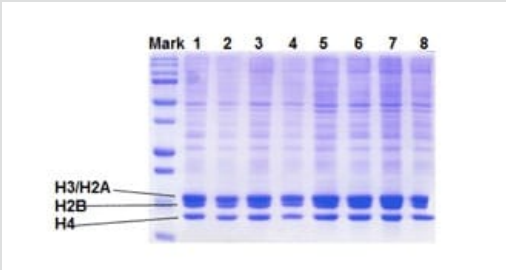
The differences in H3 acetylation level in fetal livers (female fetuses aged 19 dpc) from F1, F2, and F3 generations. measured with ab113476. Histones were prepared using **ab115124**

Asterisk indicate a significant difference between the respective C and R groups: ** P < 0.01; NS not significant. Data presented as the means ± SE (n = 6 per group).

Functional Studies - Histone Extraction Kit

(ab113476)

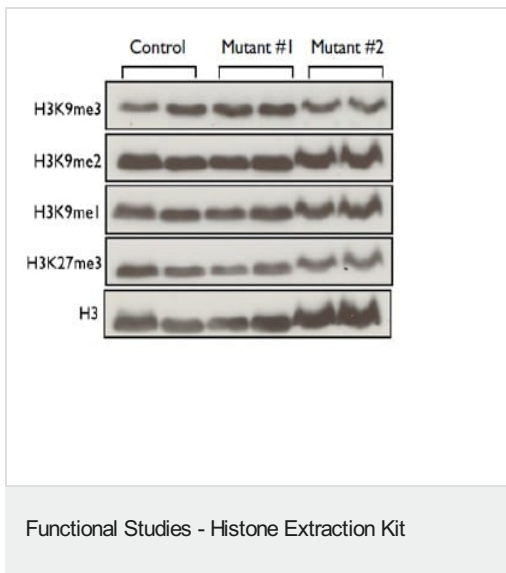
Nowacka-Woszuk, Joanna et al., PloS one?vol. 13,2 e0193464., Fig 4, doi:10.1371/journal.pone.0193464



SDS-PAGE Analysis of samples prepared using

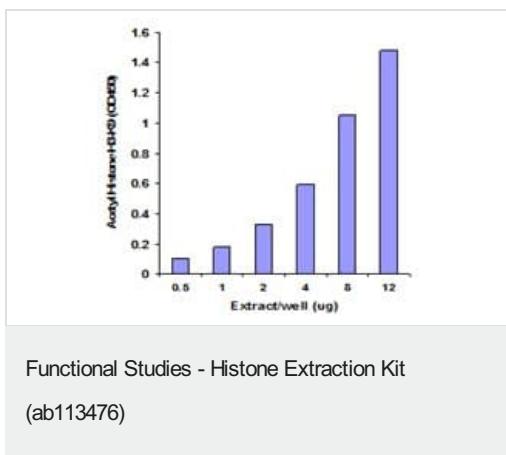
Histone extraction Kit

SDS-PAGE Analysis of samples prepared using Histone extraction Kit.



Histones were purified from mouse liver cells using ab113476. Each lane was loaded with 15µg of purified histone.

Image courtesy of Kevin Dalgaard, MPI for Immunobiology & Epigenetics, Freiburg, DE



Histone extracts were prepared from MCF-7 cells using the ab113476 and acetyl histone H3-K9 was quantified.

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.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