

Product datasheet

Histone Extraction Kit ab113476

★★★★☆ 4 Abreviews 26 References 3 Images

Overview

Product name	Histone Extraction Kit
Sample type	Tissue, Adherent cells, Suspension cells
Assay time	1h 00m
Product overview	<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>

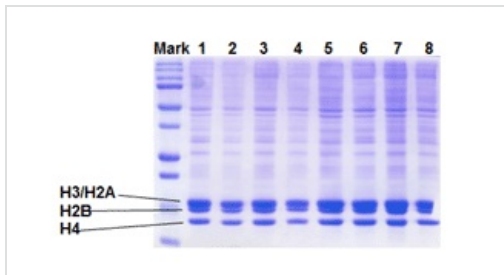
Notes	<p>Histone purification is most often performed before the analysis of the global levels within a sample of post-translational histone modifications in western blot.</p> <p>Chromatin immunoprecipitation (ChIP) is the most common technique used for the analysis of histone modifications at specific sites within the genome. Use a chromatin extraction kit instead of a histone purification kit for the preparation of samples for ChIP.</p>
--------------	------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------

Properties

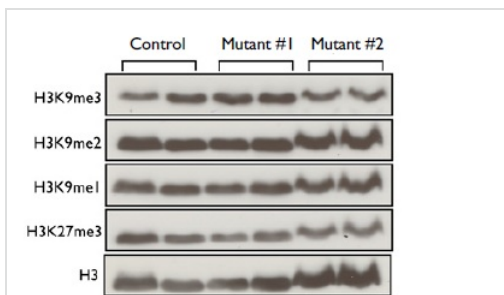
Storage instructions	Store at +4°C. Please refer to protocols.
-----------------------------	-------------------------------------------

Components	100 tests
10X Pre-Lysis Buffer	1 x 10ml
Balance Buffer	1 x 8ml
DTT Solution	1 x 20µl
Lysis Buffer	1 x 20ml

Images



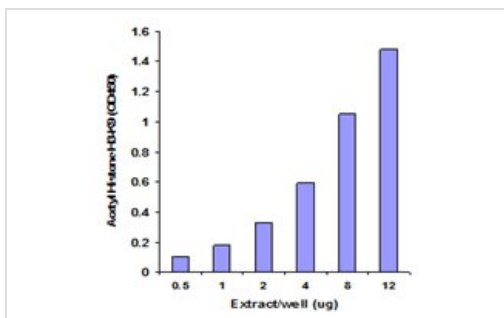
SDS-PAGE Analysis of samples prepared using Histone extraction Kit



Functional Studies - 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



Functional Studies - Histone Extraction Kit (ab113476)

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 AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

Our Promise 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