

Product datasheet

Histone Extraction Kit ab113476

★★★★☆ 4 Abreviews 16 References 3 Images

Overview

Product name	Histone Extraction Kit
Sample type	Tissue, Adherent cells, Suspension cells
Assay time	1h 00m
Species reactivity	Reacts with: Mouse, Human, Drosophila melanogaster Predicted to work with: Mammal

Product overview

Histone Extraction Kit (ab113476) is suitable for a quick preparation of total histone extracts from mammalian cells and tissue samples. The minimal amount of starting materials can be as low as 10^5 cells or 1 mg of tissue. For the best results, the cell number should be greater than 10^6 cells or the tissue amount should be greater than 10 mg. A total of 100 standard extractions (use 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.

Notes

Shipped at +4°C. Store refrigerated at +4°C before preparation.

Tested applications

Suitable for: Functional Studies

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

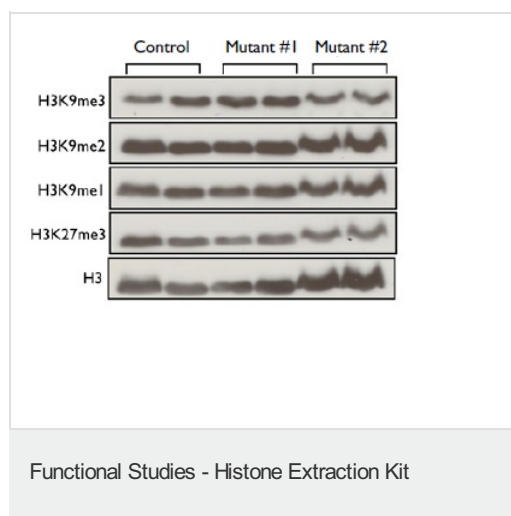
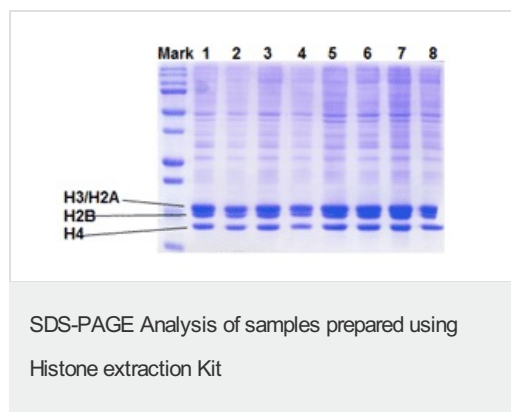
Applications

Our [Abpromise guarantee](#) covers the use of **ab113476** in the following tested applications.

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

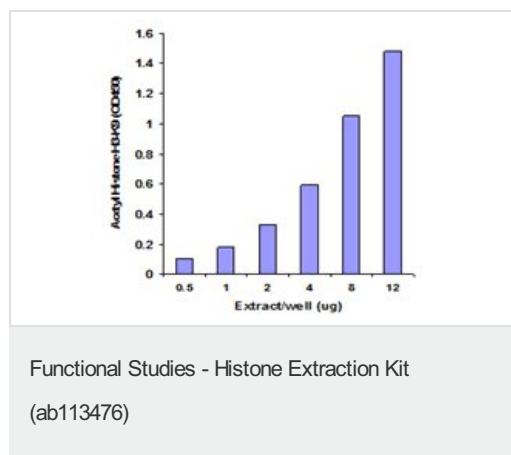
Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration. Total histone protein yield = $0.4\text{mg}/10^7$ or 100mg tissue (yield may vary depending on cell or tissue type).

Images



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.

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.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