

## Product datasheet

# Histone Extraction Kit ab113476

★★★★☆ 4 Abreviews 16 References 3 Images

### Overview

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

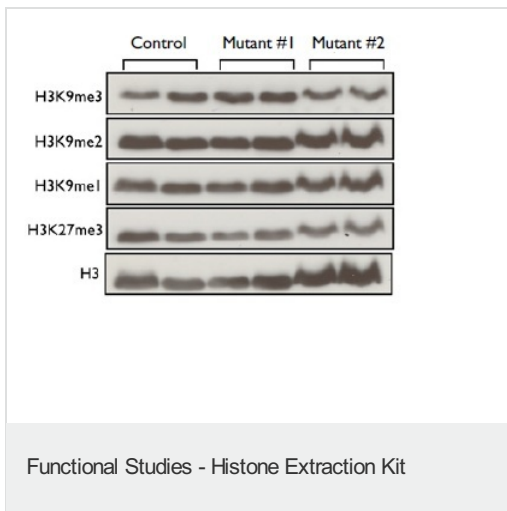
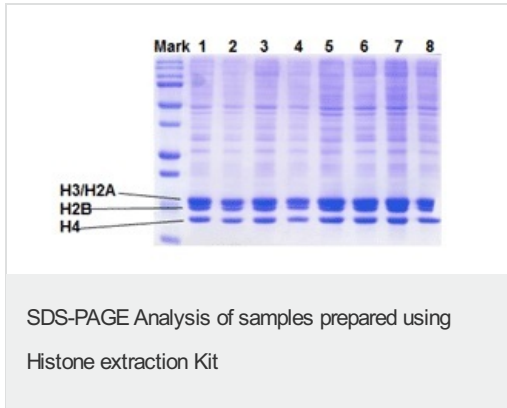
### Applications

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.

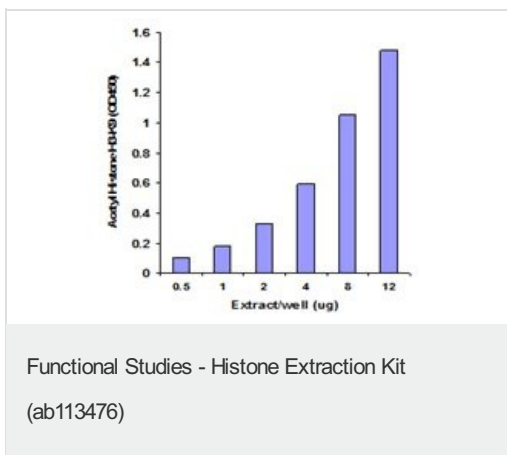
Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration. Total histone protein yield = 0.4mg/10 <sup>7</sup> or 100mg tissue (yield may vary depending on cell or tissue type).

## Histone Extraction Kit 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.

**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.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