


Histone H4 Modification Multiplex Assay Kit (Colorimetric) ab185914

1 References [画像数 1](#)

製品の概要

製品名	Histone H4 Modification Multiplex Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Tissue, Adherent cells, Suspension cells
アッセイタイプ	Quantitative
検出感度	0.5 ng/well
検出範囲	20 ng/well - 500 ng/well
全工程の試験時間	2h 30m
種交差性	交差種: Mouse, Rat, Human 交差が予測される動物種: Plants, Fungi 

製品の概要

Abcam's Histone H4 Modification Multiplex Assay Kit (Colorimetric) (ab185914) is a complete set of optimized reagents to detect and quantify nearly all histone H4 modifications (10 different types) simultaneously in a simple, ELISA-like format with use of a standard microplate reader.

ab185914 is designed for measuring multiple histone H4 modifications simultaneously. In an assay with this kit, each histone H4 modified at specific sites will be captured by an antibody that is coated on the strip wells and specifically targets the appropriate histone H4 modification pattern. The captured histone modified at specific sites will be detected with a detection antibody, followed by a color development reagent. The ratio of modified histone is proportional to the intensity of absorbance measured by a microplate reader at a wavelength of 450 nm.

Starting Materials

Input materials can be histone extracts or purified histone H4 proteins obtained from human, mouse, rat, as well as a broad range of species including most plants, fungi, and bacteria, based on high sequence homology of histone H4. The amount of histone extracts for each assay can be 20 ng to 500 ng with an optimal range of 50 to 100 ng depending on the purity of histone extracts. The amount of purified histone H4 proteins for each assay can be 1 ng to 25 ng with an optimal range of 4 to 5 ng.

試験プラットフォーム	Microplate reader
------------	-------------------

製品の特性

保存方法

Please refer to protocols.

内容	96 tests
10X Wash Buffer	1 x 28ml
8-Well Strips	2 units
96-Well Strip Plate (with Frame)	1 unit
Adhesive Covering Film	1 unit
Antibody Buffer	1 x 8ml
Assay Control Protein	1 x 20 μ l
Detection Antibody	1 x 12 μ l
Developer Solution	1 x 12ml
Stop Solution	1 x 12ml

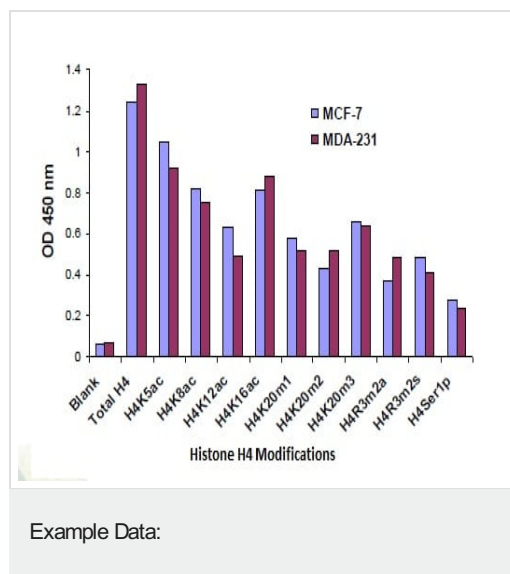
関連性

Core component of nucleosome. Nucleosomes wrap and compact DNA into chromatin, limiting DNA accessibility to the cellular machineries which require DNA as a template. Histones thereby play a central role in transcription regulation, DNA repair, DNA replication and chromosomal stability. DNA accessibility is regulated via a complex set of post-translational modifications of histones, also called histone code, and nucleosome remodeling.

細胞内局在

Nuclear

画像



Histone extracts were prepared from MCF-7 and MDA-231 cells using the Histone Extraction Kit ([ab113476](#)). 10 histone H4 modifications were measured using the Histone H4 Modification Multiplex Assay Kit (Colorimetric) ([ab185914](#)). 100 ng of total histone proteins per well were used.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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