

Histone H3 Modification Multiplex Assay Kit (Colorimetric)
ab185910

★★★★★ [1 Abreviews](#) [3 References](#) [画像数 1](#)

製品の概要

製品名	Histone H3 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
製品の概要	The Histone H3 Modification Multiplex Assay Kit (Colorimetric) (ab185910) is a complete set of optimized reagents to detect and quantify up to 21 modified Histone H3 patterns simultaneously in a simple, ELISA-like format with use of a standard microplate reader.
試験プラットフォーム	Microplate reader

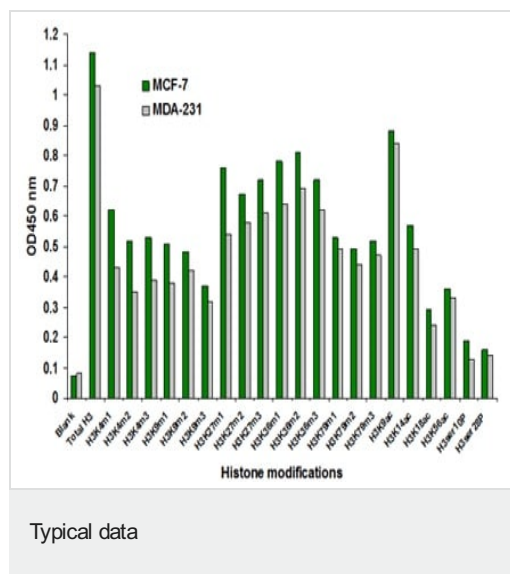
製品の特性

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 20µl

内容	96 tests
Detection Antibody	1 x 12 μ l
Developer Solution	1 x 12ml
Stop Solution	1 x 12ml

画像



Histone extracts were prepared from MCF-7 and MDA-231 cells and multiple Histone H3 modifications were screened and measured using the Histone H3 Modification Multiplex Assay Kit (Colorimetric). 100 ng of total histone proteins 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