

# Histone H3 Modification Multiplex Assay Kit (Colorimetric) ab185910

★★★★★ [1 Abreviews](#) [3 References](#) [1 图像](#)

### 概述

产品名称	Histone H3 Modification Multiplex Assay试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Tissue, Adherent cells, Suspension cells
检测类型	Quantitative
灵敏度	0.5 ng/well
范围	20 ng/well - 500 ng/well
检测时间	2h 30m
种属反应性	与反应: Mouse, Rat, Human
产品概述	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.
平台	Microplate reader

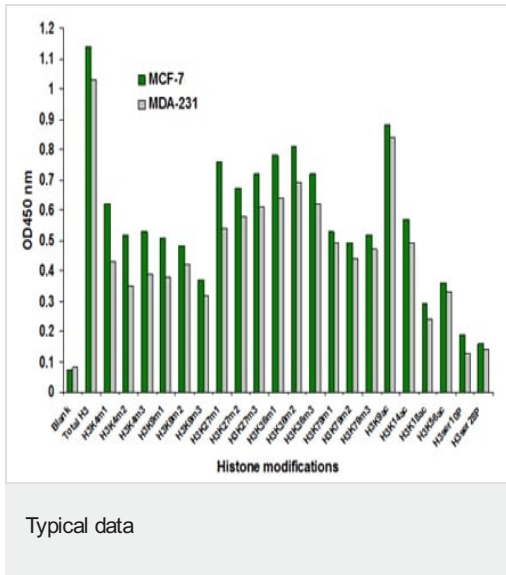
### 性能

**存放说明** Please refer to protocols.

组件	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

组件	96 tests
Detection Antibody	1 x 12 $\mu$ l
Developer Solution	1 x 12ml
Stop Solution	1 x 12ml

图片



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.cn/abpromise> or contact our technical team.

**Terms and conditions**

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