

Histone H4 Total Quantification Kit (Colorimetric) ab156909

1 References 5 图像

概述	
产品名称	Histone H4 Total Quantification试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Cell culture extracts, Tissue Extracts, Nuclear Extracts
检测类型	Quantitative
灵敏度	>= 0.5 ng/well
范围	50 ng - 1000 ng
检测时间	3h 30m
种属反应性	与反应: Mouse, Rat, Human
产品概述	<p>Abcam's Histone H4 Total Quantification Kit (Colorimetric) is a complete set of essential components which enables an experimenter to colorimetrically quantify levels of histone H4 proteins independent of their modified states. It can also be used for normalizing the modified histone H4 content of samples when run in parallel with Abcam's histone modification quantification kit series.</p> <p>The kit is suitable for specifically measuring total histone H4 using human, mouse, and rat samples, including fresh and frozen tissues and cultured adherent and suspension cells.</p>
平台	Microplate reader

性能

存放说明 Please refer to protocols.

组件	48 tests	96 tests
1000X Capture Antibody	1 x 5µl	1 x 10µl
10X Wash Buffer	1 x 14ml	1 x 28ml
2000X Signal Reporter	1 x 6µl	1 x 12µl
8-Well Assay Strips (with Frame)	1 x 6 units	1 x 12 units
Color Developer	1 x 5ml	1 x 10ml

组件	48 tests	96 tests
Enhancer Solution	1 x 120µl	1 x 240µl
Histone Assay Buffer	1 x 4ml	1 x 8ml
Standard Control, 100 µg/mL	1 x 10µl	1 x 20µl
Stop Solution	1 x 5ml	1 x 10ml

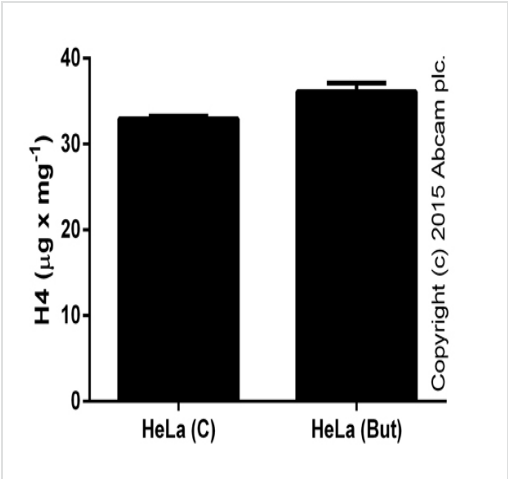
相关性

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.

细胞定位

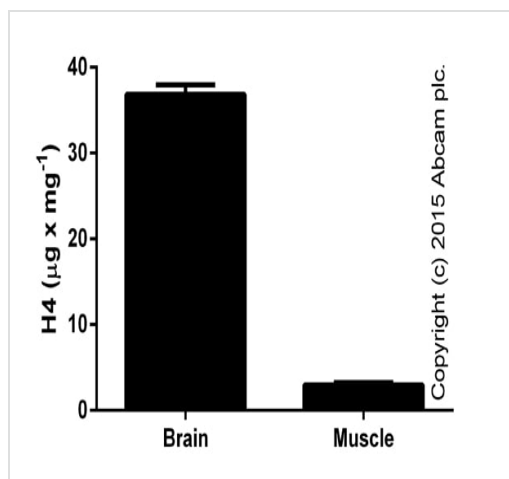
Nuclear

图片



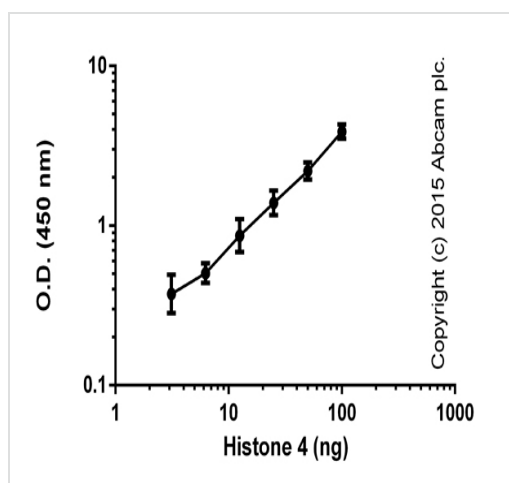
Total H4 per mg of extracted histones from HeLa control cells (C) or cells treated with 10 micromolar sodium butyrate (But) (extracted using [ab113476](#); 50-500 ng per well) (duplicates +/- SD).

Functional Assay: ab156909 Histone H4 Total Quantification Kit (Colorimetric)



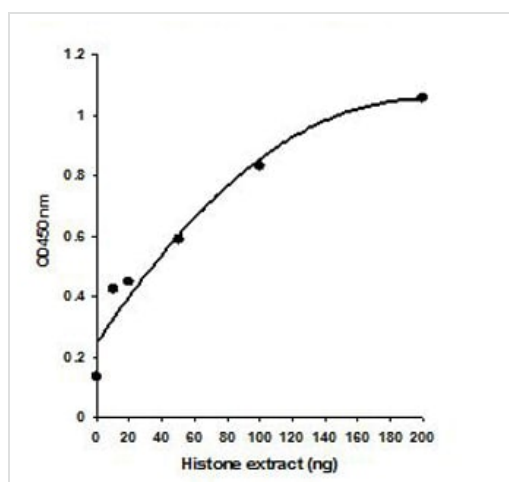
Functional Assay: ab156909 Histone H4 Total  
Quantification Kit (Colorimetric)

Total H4 per mg of extracted histones from murine tissue preparations (extracted using [ab113476](#); 50-500 ng per well) (duplicates +/- SD).



Functional Assay: ab156909 Histone H4 Total  
Quantification Kit (Colorimetric)

Standard curve with background signal subtracted (duplicates; +/- SD).



Histone H4 Quantification

Histone nuclear extracts were prepared from MDA-231 cells using the [Histone Extraction Kit \(ab113476\)](#) and the ODs generated from Histone H4 are measured.

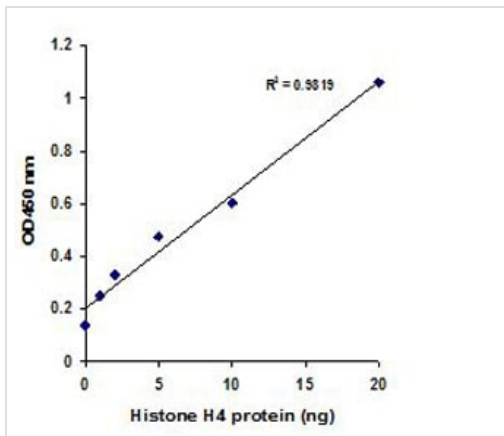


Figure 1. Illustrated standard curve generated with H4 standard.

Illustrated Standard Curve

**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