

Histone H4 Total Quantification Kit (Colorimetric)

ab156909

[1 References](#) [5 Images](#)

Overview

Product name	Histone H4 Total Quantification Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Cell culture extracts, Tissue Extracts, Nuclear Extracts
Assay type	Quantitative
Sensitivity	>= 0.5 ng/well
Range	50 ng - 1000 ng
Assay time	3h 30m
Species reactivity	Reacts with: Mouse, Rat, Human
Product overview	<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>
Platform	Microplate reader

Properties

Storage instructions Please refer to protocols.

Components	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

Components	48 tests	96 tests
Color Developer	1 x 5ml	1 x 10ml
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

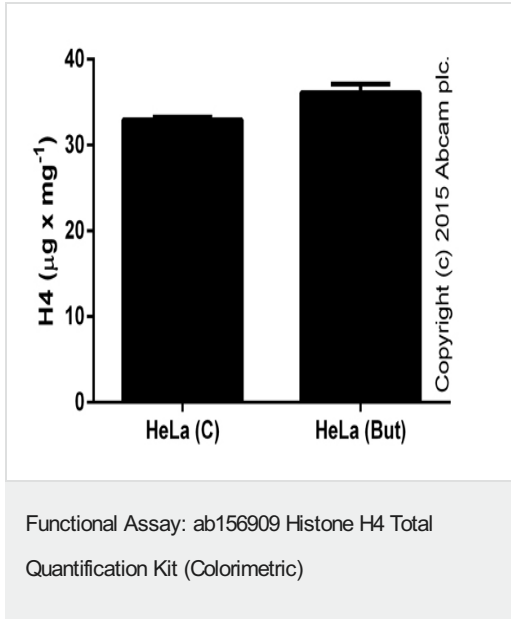
Relevance

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.

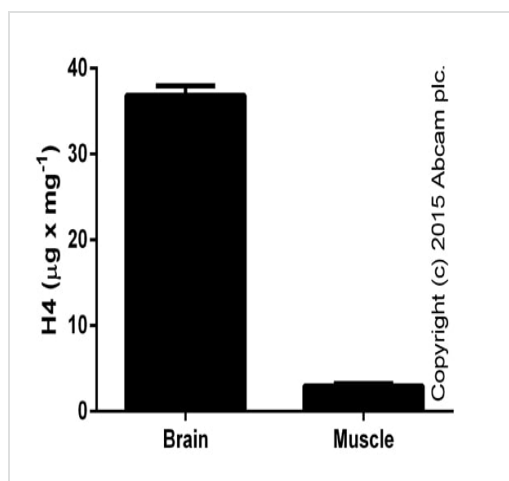
Cellular localization

Nuclear

Images

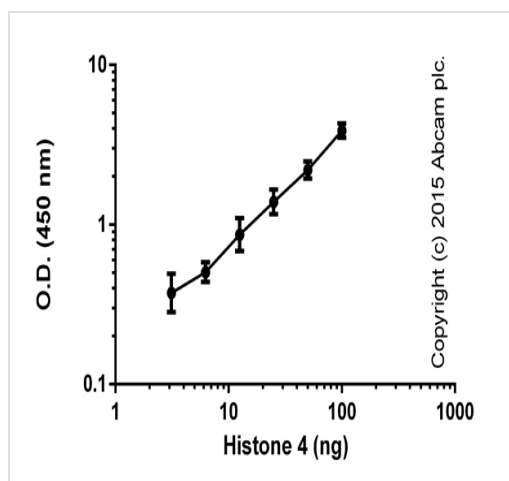


Total H4 per mg of extracted histones from HeLa control cells (C) or cells treated with 10 micromolar sodium butyrate (But) (extracted using [ab113476](#); 50-500 ng per well) (duplicates +/- SD).



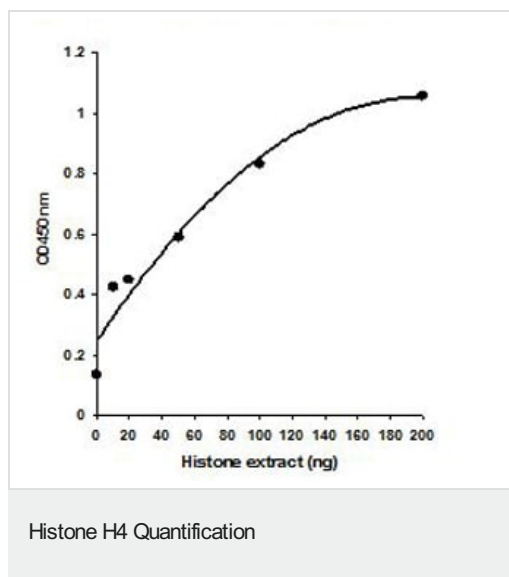
Total H4 per mg of extracted histones from murine tissue preparations (extracted using [ab113476](#); 50-500 ng per well) (duplicates \pm SD).

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



Standard curve with background signal subtracted (duplicates; \pm SD).

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



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

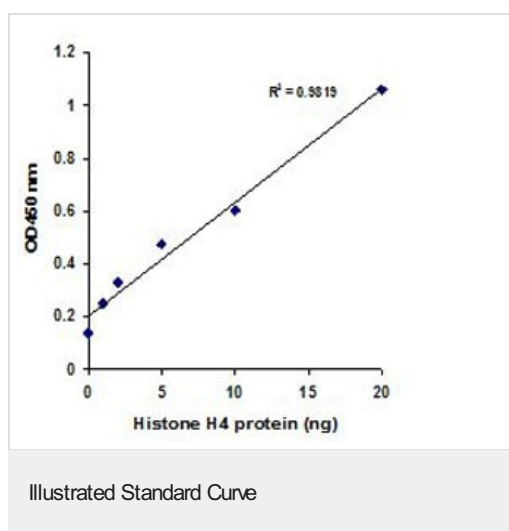


Figure 1. Illustrated standard curve generated with H4 standard.

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.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