


# Histone H3 Total Quantification Kit (Colorimetric) ab115091

1 References 1 图像

概述

产品名称	Histone H3 Total Quantification试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Tissue, Adherent cells, Suspension cells
灵敏度	10 ng/well
范围	20 ng/well - 1000 ng/well
检测时间	2h 30m
种属反应性	与反应: Mouse, Rat, Human 预测可用于: Mammals 

**产品概述**

Histone H3 Total Quantification Kit (Colorimetric) (ab115091) enables the user to quantify levels of histone H3 proteins independent of their modified states. It can also be used for normalizing the modified histone H3 content of samples when run in parallel with Abcam's histone modification quantification kit series. ab115091 is suitable for specifically measuring total histone H3 on mammalian samples such as human, mouse, and rat, including fresh and frozen tissues and cultured adherent and suspension cells.

Histone H3 can undergo several epigenetic modifications that influence cellular processes such as transcription activation/inactivation, chromosome packaging, and DNA damage/repair. These modifications including acetylation, phosphorylation or methylation occur on the N-terminal tail domains of histone H3 through a variety of histone modifying enzymes. In most species, histone H3 is primarily acetylated at K9, K14, K18, K23 and K56, methylated at K4, K9, K27, K36 and K79, and phosphorylated at S10, S28, T3 and T11, respectively.

平台	Microplate reader
----	-------------------

性能

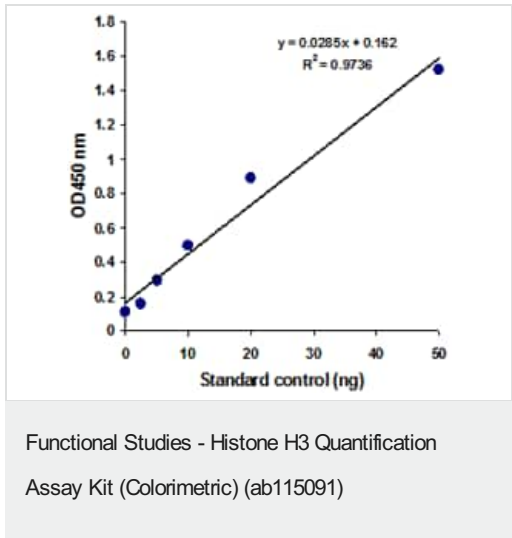
**存放说明** Please refer to protocols.

组件	标识符	48 tests	96 tests
10X Wash Buffer		1 x 10ml	1 x 20ml
8-Well Sample Strip (with Frame)		4 units	9 units

组件	标识符	48 tests	96 tests
8-Well Standard Control Strips	Green Ringed	2 units	3 units
Antibody Buffer		1 x 6ml	1 x 12ml
Color Developer		1 x 5ml	1 x 10ml
Detection Antibody, 1 mg/mL		1 x 5µl	1 x 10µl
Standard Control, 100 µg/mL		1 x 10µl	1 x 20µl
Stop Solution		1 x 3ml	1 x 6ml

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	Nucleus. Chromosome.

图片



Measurement of histone H3 with different concentrations of the included H3 standard in theHistone H3 Total Quantification Kit (Colorimetric) (ab115091).

**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