


Histone H3 Total Quantification Kit (Fluorometric)

ab115092

1 Image

Overview

| | |
|--------------------|----------------------------------------------------------------------------------------------------------------------------|
| Product name | Histone H3 Total Quantification Kit (Fluorometric) |
| Detection method | Fluorescent |
| Sample type | Tissue, Adherent cells, Suspension cells |
| Sensitivity | 2 ng/well |
| Range | 10 ng/well - 1000 ng/well |
| Assay time | 2h 30m |
| Species reactivity | Reacts with: Mouse, Rat, Human |
| | Predicted to work with: Mammals  |

| | |
|------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Product overview | <p>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.</p> <p>Histone H3 Total Quantification Kit (Fluorometric) (ab115092) 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.</p> |
| Platform | Microplate reader |

Properties

| | |
|----------------------|----------------------------|
| Storage instructions | Please refer to protocols. |
|----------------------|----------------------------|

| Components | Identifier | 48 tests | 96 tests |
|-----------------|------------|----------|----------|
| 10X Wash Buffer | | 1 x 10ml | 1 x 20ml |
| | | | |

| Components | Identifier | 48 tests | 96 tests |
|-----------------------------------|--------------|----------|----------|
| 8 well Sample Strips (with frame) | | 4 units | 9 units |
| 8 well Standard Control Strips | Green Ringed | 2 units | 3 units |
| Antibody Buffer | | 1 x 6ml | 1 x 12ml |
| Detection Antibody, 1 mg/mL | | 1 x 5µl | 1 x 10µl |
| Fluoro Developer | | 1 x 12µl | 1 x 24µl |
| Fluoro Dilution | | 1 x 4ml | 1 x 8ml |
| Fluoro Enhancer | | 1 x 12µl | 1 x 24µl |
| Standard Control (100 µg/mL) | | 1 x 10µl | 1 x 20µl |

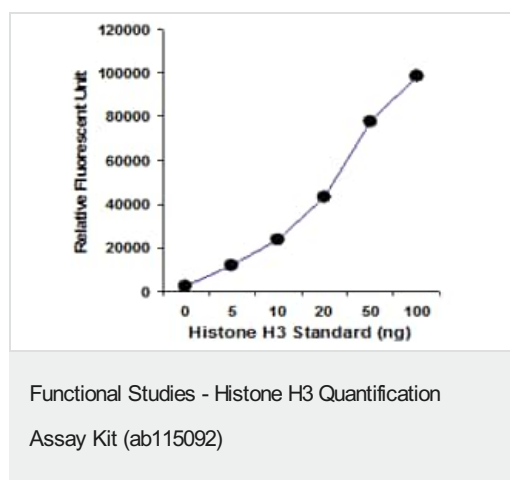
Relevance

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

Nucleus. Chromosome.

Images



A histone H3 standard curve was generated as described in the protocol of the Histone H3 Total Quantification Kit (Fluorometric) (ab115092).

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