


Histone H3 (pan-methyl K9) Quantification Kit (Fluorometric) ab115067

1 Image

Overview

|                    |   |
|--------------------|---|
| Product name       | Histone H3 (pan-methyl K9) Quantification Kit (Fluorometric)  |
| Detection method   | Fluorescent   |
| Sample type        | Tissue, Adherent cells, Suspension cells  |
| Assay type         | Quantitative  |
| Sensitivity        | 0.4 ng/well   |
| Range              | 5 ng/well - 2000 ng/well  |
| Assay time         | 3h 00m  |
| Species reactivity | <b>Reacts with:</b> Mouse, Human<br><b>Predicted to work with:</b> Mammals   |
| Product overview   | <p>Methylation of histone H3 at lysine 9 by ESET, G9a among other methyl transferases is involved in euchromatin gene silencing as well as heterochromatin formation. H3 (mono-methyl K9) is enriched in certain euchromatic domains, which have been postulated to be transcriptionally silent. H3 (di-methyl K9) and H3 (tri-methyl K9) mediate heterochromatin formation by forming a binding site for HP1 and also participate in silencing gene expression at euchromatic sites.</p> <p>Histone H3 (pan-methyl K9) Quantification Kit (Fluorometric) (ab115067) allows the user to specifically measure global mono-, di-, and tri-methylation of histone H3K9 fluorometrically using a variety of mammalian cells including fresh and frozen tissues, cultured adherent and suspension cells.</p> |
| Platform           | Microplate reader   |

Properties

Storage instructions Please refer to protocols.

| Components                        | Identifier | 96 tests    |
|-----------------------------------|------------|-------------|
| 10X Wash Buffer                   |            | 1 x 20ml    |
| 8-Well Sample Strips (with Frame) |            | 1 x 9 units |
|                                   |            |             |

| Components                     | Identifier   | 96 tests    |
|--------------------------------|--------------|-------------|
| 8-Well Standard Control Strips | Green Ringed | 1 x 3 units |
| Antibody Buffer                |              | 1 x 12ml    |
| Detection Antibody, 1 mg/mL    |              | 1 x 10µl    |
| Fluoro Developer               |              | 1 x 24µl    |
| Fluoro Dilution                |              | 1 x 8ml     |
| Fluoro Enhancer                |              | 1 x 24µl    |
| Standard Control, 100 µg/mL    |              | 1 x 20µl    |

### Function

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.

### Sequence similarities

Belongs to the histone H3 family.

### Developmental stage

Expressed during S phase, then expression strongly decreases as cell division slows down during the process of differentiation.

### Post-translational modifications

Acetylation is generally linked to gene activation. Acetylation on Lys-10 (H3K9ac) impairs methylation at Arg-9 (H3R8me2s). Acetylation on Lys-19 (H3K18ac) and Lys-24 (H3K24ac) favors methylation at Arg-18 (H3R17me).

Citrullination at Arg-9 (H3R8ci) and/or Arg-18 (H3R17ci) by PAD4 impairs methylation and represses transcription.

Asymmetric dimethylation at Arg-18 (H3R17me2a) by CARM1 is linked to gene activation. Symmetric dimethylation at Arg-9 (H3R8me2s) by PRMT5 is linked to gene repression. Asymmetric dimethylation at Arg-3 (H3R2me2a) by PRMT6 is linked to gene repression and is mutually exclusive with H3 Lys-5 methylation (H3K4me2 and H3K4me3). H3R2me2a is present at the 3' of genes regardless of their transcription state and is enriched on inactive promoters, while it is absent on active promoters.

Methylation at Lys-5 (H3K4me), Lys-37 (H3K36me) and Lys-80 (H3K79me) are linked to gene activation. Methylation at Lys-5 (H3K4me) facilitates subsequent acetylation of H3 and H4. Methylation at Lys-80 (H3K79me) is associated with DNA double-strand break (DSB) responses and is a specific target for TP53BP1. Methylation at Lys-10 (H3K9me) and Lys-28 (H3K27me) are linked to gene repression. Methylation at Lys-10 (H3K9me) is a specific target for HP1 proteins (CBX1, CBX3 and CBX5) and prevents subsequent phosphorylation at Ser-11 (H3S10ph) and acetylation of H3 and H4. Methylation at Lys-5 (H3K4me) and Lys-80 (H3K79me) require preliminary monoubiquitination of H2B at 'Lys-120'. Methylation at Lys-10 (H3K9me) and Lys-28 (H3K27me) are enriched in inactive X chromosome chromatin.

Phosphorylated at Thr-4 (H3T3ph) by GSG2/haspin during prophase and dephosphorylated during anaphase. Phosphorylation at Ser-11 (H3S10ph) by AURKB is crucial for chromosome condensation and cell-cycle progression during mitosis and meiosis. In addition phosphorylation at Ser-11 (H3S10ph) by RPS6KA4 and RPS6KA5 is important during interphase because it enables the transcription of genes following external stimulation, like mitogens, stress, growth factors or UV irradiation and result in the activation of genes, such as c-fos and c-jun.

Phosphorylation at Ser-11 (H3S10ph), which is linked to gene activation, prevents methylation at

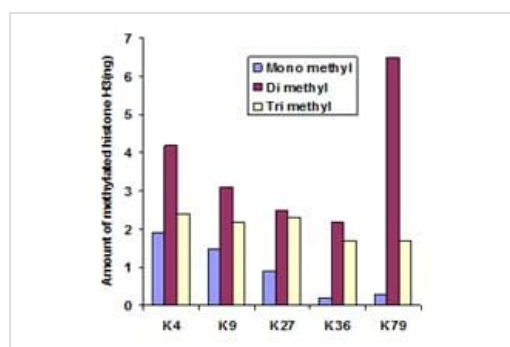
Lys-10 (H3K9me) but facilitates acetylation of H3 and H4. Phosphorylation at Ser-11 (H3S10ph) by AURKB mediates the dissociation of HP1 proteins (CBX1, CBX3 and CBX5) from heterochromatin. Phosphorylation at Ser-11 (H3S10ph) is also an essential regulatory mechanism for neoplastic cell transformation. Phosphorylated at Ser-29 (H3S28ph) by MLTK isoform 1, RPS6KA5 or AURKB during mitosis or upon ultraviolet B irradiation. Phosphorylation at Thr-7 (H3T6ph) by PRKCBB is a specific tag for epigenetic transcriptional activation that prevents demethylation of Lys-5 (H3K4me) by LSD1/KDM1A. At centromeres, specifically phosphorylated at Thr-12 (H3T11ph) from prophase to early anaphase, by DAPK3 and PKN1. Phosphorylation at Thr-12 (H3T11ph) by PKN1 is a specific tag for epigenetic transcriptional activation that promotes demethylation of Lys-10 (H3K9me) by KDM4C/JMJD2C. Phosphorylation at Tyr-42 (H3Y41ph) by JAK2 promotes exclusion of CBX5 (HP1 alpha) from chromatin.

Monoubiquitinated by RAG1 in lymphoid cells, monoubiquitination is required for V(D)J recombination (By similarity). Ubiquitinated by the CUL4-DDB-RBX1 complex in response to ultraviolet irradiation. This may weaken the interaction between histones and DNA and facilitate DNA accessibility to repair proteins.

## Cellular localization

Nucleus. Chromosome.

## Images



Functional Studies - Histone H3 (pan-methyl K9)  
Quantification Kit (Fluorometric) (ab115067)

ab115067 used to specifically measure global histone H3K9 mono-, di-, and tri-methylation. The ratio of mono-, di-, and tri-methylated H3K9 is proportional to the intensity of fluorescence. The absolute amount of the methylated H3K9 can be quantitated by comparing to the standard control.

**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