


Histone H4 (pan-methyl K20) Quantification Kit (Fluorometric)

ab115100

製品の概要

製品名	Histone H4 (pan-methyl K20) Quantification Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Tissue, Adherent cells, Suspension cells
検出感度	1 ng/well
検出範囲	10 ng/well - 2000 ng/well
全工程の試験時間	2h 30m
種交差性	交差種: Mouse, Human 交差が予測される動物種: Mammals 

製品の概要	<p>Methylation of histone H4 at lysine 20 is catalyzed by SET9 and SUV4.20h among other methyl transferases in mammalian cells. H4 (mono-methyl K20) is involved in the maintenance of proper higher order structure of DNA and is consequently essential for chromosome condensation, as well as functioning in gene silencing. H4 (di-methyl K20) has been described as another repressive chromatin domain and is involved in the DNA damage response. H4 (tri-methyl K20) acts as a passive feature or structure determinant for chromatin degradation and release, as well as being an epigenetic marker of early apoptosis. H4 (tri-methyl K20) is also considered as a common hallmark of human cancer.</p> <p>Abcam's Histone H4 (pan-methyl K20) Quantification Kit (Fluorometric) (ab115100) enables the user to measure global mono-, di- and tri-methylation of histone H4K20. The global H4K20 pan-methylation can also be changed by inhibition or activation of HMTs, making quantitative detection of global pan-methyl histone H4K20 a useful tool for better understanding epigenetic regulation of gene activation/repression. ab115100 is suitable for use with a variety of mammalian cells including fresh and frozen tissues, and cultured adherent and suspension cells.</p>
-------	--

試験プラットフォーム	Microplate reader
------------	-------------------

製品の特性

保存方法	Please refer to protocols.
------	----------------------------

内容	ラベル	96 tests
10X Wash Buffer		1 x 20ml

内容	ラベル	96 tests
8 Well Sample Strip (with Frame)		1 x 9 units
8 Well Standard Control Strips	Green Ringed	1 x 3 units
Antibody Buffer		1 x 12ml
Detecting Antibody, 1 mg/mL		1 x 10 μ l
Fluoro Developer		1 x 24 μ l
Fluoro Dilution		1 x 8ml
Fluoro Enhancer		1 x 24 μ l
Signal Enhancer		1 x 240 μ l
Signal Report Solution		1 x 10 μ l
Standard Control, 100 μ g/mL		1 x 20 μ l

機能	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.
配列類似性	Belongs to the histone H4 family.
翻訳後修飾	<p>Acetylation at Lys-6 (H4K5ac), Lys-9 (H4K8ac), Lys-13 (H4K12ac) and Lys-17 (H4K16ac) occurs in coding regions of the genome but not in heterochromatin.</p> <p>Citrullination at Arg-4 (H4R3ci) by PAD4 impairs methylation.</p> <p>Monomethylation and asymmetric dimethylation at Arg-4 (H4R3me1 and H4R3me2a, respectively) by PRMT1 favors acetylation at Lys-9 (H4K8ac) and Lys-13 (H4K12ac).</p> <p>Demethylation is performed by JMJD6. Symmetric dimethylation on Arg-4 (H4R3me2s) by the PRDM1/PRMT5 complex may play a crucial role in the germ-cell lineage.</p> <p>Monomethylated, dimethylated or trimethylated at Lys-21 (H4K20me1, H4K20me2, H4K20me3).</p> <p>Monomethylation is performed by SET8. Trimethylation is performed by SUV420H1 and SUV420H2 and induces gene silencing.</p> <p>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. Monoubiquitinated at Lys-92 of histone H4 (H4K91ub1) in response to DNA damage. The exact role of H4K91ub1 in DNA damage response is still unclear but it may function as a licensing signal for additional histone H4 post-translational modifications such as H4 Lys-21 methylation (H4K20me).</p> <p>Sumoylated, which is associated with transcriptional repression.</p>
細胞内局在	Nucleus. Chromosome.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors