

Histone H3 (di-methyl K9) Quantification Kit (Colorimetric) ab115062

画像数 1

製品の概要

製品名	Histone H3 (di-methyl K9) Quantification Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Tissue, Adherent cells, Suspension cells
アッセイタイプ	Quantitative
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

Methylation of histone H3 at lysine 9 by ESET, G9a among other methyl transferases is involved in euchromatin gene silencing as well as in heterochromatin formation. Specifically, *H3 (di-methyl K9)* mediates heterochromatin formation by forming a binding site for HP1.

Histone H3 (di-methyl K9) Quantification Kit (Colorimetric) (ab115062) allows the user to specifically measure global di-methylation of histone H3K9 colorimetrically using a variety of mammalian cells (human, mouse, etc.) including fresh and frozen tissues, cultured adherent and suspension cells.

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.

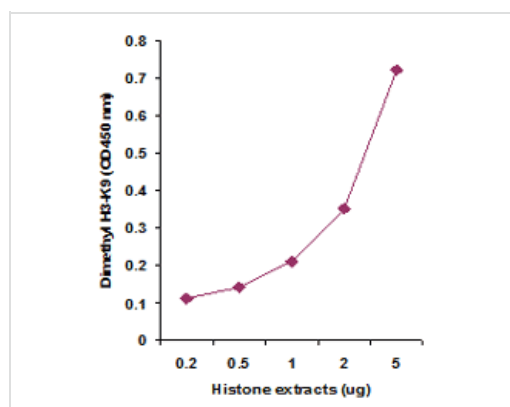
内容	ラベル	48 tests	96 tests
10X Wash Buffer		1 x 10ml	1 x 20ml
8-Well Sample Strips (with frame)		4 units	9 units
8-Well Standard Control Strips	Green Ringed	2 units	3 units
Antibody Buffer		1 x 6ml	1 x 2ml
Color Developer		1 x 5ml	1 x 10ml

内容	ラベル	48 tests	96 tests
Detection Antibody, 1 mg/mL		1 x 5 μ l	1 x 10 μ l
Standard Control (100 μ g/mL)		1 x 10 μ l	1 x 20 μ l
Stop Solution		1 x 3ml	1 x 6ml

関連性 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.

細胞内局在 Nucleus. Chromosome.

画像



Histone extracts were prepared from MDA-MB-231 cells using the Histone Extraction Kit ([ab113476](#)) and the amount of dimethyl-H3-K9 was measured using ab115062.

Functional Studies - Histone H3 (di-methyl K9)
Quantification Kit (Colorimetric) (ab115062)

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