


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

1 Image

Overview

Product name	Histone H3 (di-methyl K9) Quantification Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Tissue, Adherent cells, Suspension cells
Assay type	Quantitative
Assay time	2h 30m
Species reactivity	Reacts with: Mouse, Human Predicted to work with: 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 in heterochromatin formation. Specifically, <i>H3 (di-methyl K9)</i> mediates heterochromatin formation by forming a binding site for HP1.</p> <p>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.</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
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

Components	Identifier	48 tests	96 tests
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

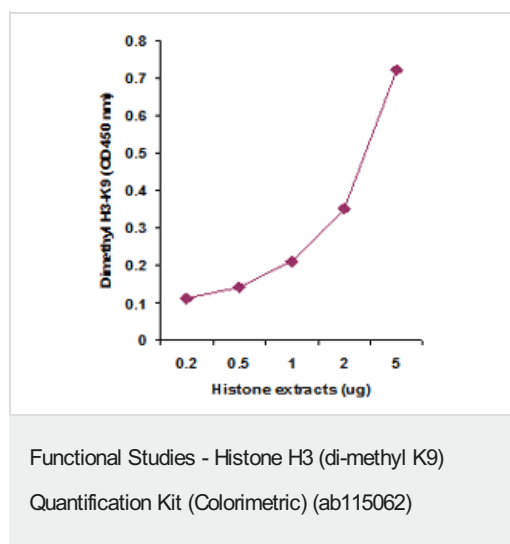
Relevance

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



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.

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