

Product datasheet

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

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.

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

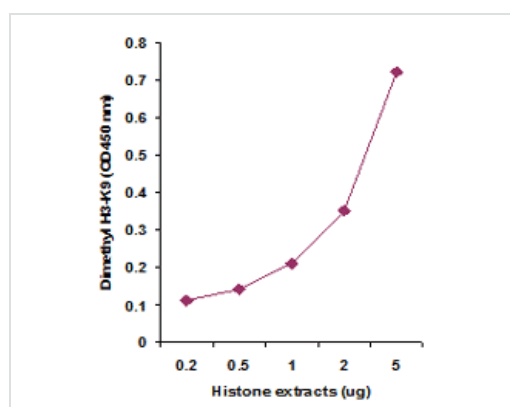
Relevance

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](#).

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.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