

Product datasheet

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

[1 Image](#)

Overview

Product name	Histone H3 (pan-methyl K9) Quantification Kit (Fluorometric)
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	Reacts with: Mouse, Human Predicted to work with: Mammal

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 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.

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.

Tested applications

Suitable for: Functional Studies

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
8-Well Standard Control Strips	Green Ringed	1 x 3 units

Components	Identifier	96 tests
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

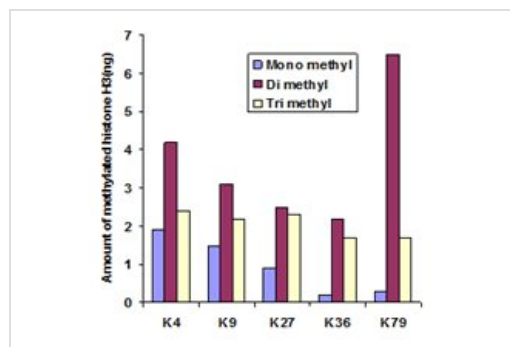
Applications

Our [Abpromise guarantee](#) covers the use of **ab115067** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration.

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 AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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