


# Histone H3 (K9) Methyltransferase Activity Quantification Assay Kit ab113453

★★★★★ [1 Abreviews](#) [2 References](#) [1 图像](#)

### 概述

产品名称	Histone H3 (K9) Methyltransferase Activity Quantification Assay试剂盒
样品类型	Tissue Extracts, Cell Lysate
检测类型	Quantitative
检测时间	3h 00m
种属反应性	与反应: Mouse, Human 预测可用于: Mammals 

### 产品概述

Methylation of histone H3 at lysine 9 (K9) mediates heterochromatin formation by forming a binding site for HP1 and also participates in silencing expression at euchromatic sites. In mammalian cells, histone lysine methyltransferases such as SuVar39h and ESET among others are responsible for methylation of lysine 9 at histone H3.

Histone H3 (K9) Methyltransferase Activity Quantification Assay Kit (ab113453) allows the user to measure the activity or inhibition of individual histone methyltransferase (HMT) that specifically target histone H3 at lysine 9 (*H3K9*). The kit is ready-to-use and provides all the essential components needed to carry out a successful HMT activity/inhibition quantification experiment without the need for radioactivity or any special equipment.

ab113453 recognizes methylated H3 at lysine site 9.

平台	Microplate reader
----	-------------------

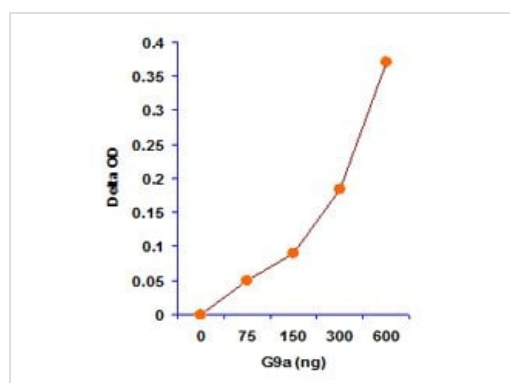
### 性能

存放说明 Please refer to protocols.

组件	48 tests	96 tests
10X Wash Buffer	1 x 11ml	1 x 22ml
8-Well Assay Strips (with Frame)	6 units	12 units
Adomet	1 x 25µl	1 x 50µl
Biotinylated Substrate, 25 µg/mL	1 x 100µl	1 x 200µl

组件	48 tests	96 tests
Capture Antibody, 100 µg/mL	1 x 25µl	1 x 50µl
Control Enzyme (300 µg/mL)	1 x 5µl	1 x 10µl
Detection Antibody, 200 µg/mL	1 x 10µl	1 x 20µl
Developing Solution	1 x 6ml	1 x 12ml
Histone Assay Buffer	1 x 1.5ml	1 x 3ml
HMT Standard, 10 µg/mL	1 x 10µl	1 x 20µl
Stop Solution	1 x 3ml	1 x 6ml

## 图片



Functional Studies - Histone H3 (K9)  
Methyltransferase Activity Quantification Assay Kit  
(ab113453)

Nuclear extracts were prepared from MCF-7 cells using the [ab113474](#) and H3-K9 specific histone methyltransferase activity (G9a) was measured. Purified G9a was spiked into the nuclear extract that has low concentration of nuclear proteins by dilution.

**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.cn/abpromise> or contact our technical team.

## Terms and conditions

---

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